

Active Motif Epigenetic Services Publications

The papers below contain data generated by Active Motif's Epigenetic Services team. To learn more about our services, please give us a call or visit us at [activemotif.com/services](https://www.activemotif.com/services).

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K27me3	Nat Commun.	2024	Zoe Veneti <i>et al.</i> Polycomb-mediated silencing of miR-8 is required for maintenance of intestinal stemness in <i>Drosophila melanogaster</i> . DOI: 10.1038/s41467-024-46119-9
ModSpec		Sci Adv.	2024	Anthony KN Chan <i>et al.</i> Therapeutic targeting Tudor domains in leukemia via CRISPR-Scan Assisted Drug Discovery. DOI: 10.1126/sciadv.adk3127
ATAC-Seq		Nature	2024	Violetta Karwacki-Neisius <i>et al.</i> WNT signalling control by KDM5C during development affects cognition. DOI: 10.1038/s41586-024-07067-y
ChIP-Seq	FOXO1	Mol Cell Endocrinol.	2024	Caitlin E Stallings <i>et al.</i> FOXO1 regulates expression of Neurod4 in the pituitary gland. DOI: 10.1016/j.mce.2023.112128
ATAC-Seq		Life Sci Alliance	2024	Shivakshi Sulekh <i>et al.</i> A feedback loop that drives cell death and proliferation and its defect in intestinal stem cells. DOI: 10.26508/lsa.202302238
ChIP-Seq	H3K27me3, EZH2	Genome Biol.	2024	Georgette Tanner <i>et al.</i> IDHwt glioblastomas can be stratified by their transcriptional response to standard treatment, with implications for targeted therapy. DOI: 10.1186/s13059-024-03172-3
ChIP-Seq	Hoxa5, Pbx1	bioRxiv	2024	KC Ritesh <i>et al.</i> Multimodal Hox5 activity generates motor neuron diversity. DOI: 10.1101/2024.02.08.579338
ATAC-Seq, RNA-Seq		EMBO Rep.	2024	Panagiotis Papoutsoglou <i>et al.</i> TGF β -induced long non-coding RNA LINC00313 activates Wnt signaling and promotes cholangiocarcinoma. DOI: 10.1038/s44319-024-00075-z
ChIP-Seq	H3K27ac	Cancer Res.	2024	Tathiane M Malta <i>et al.</i> The Epigenetic Evolution of Glioma Is Determined by the IDH1 Mutation Status and Treatment Regimen. DOI: 10.1158/0008-5472.CAN-23-2093
snRNA-Seq		Aging Cell	2024	Spiros Palikyras <i>et al.</i> Rapid and synchronous chemical induction of replicative-like senescence via a small molecule inhibitor. doi: 10.1111/ace1.14083
RNA-Seq		Cell Rep.	2024	Clement Maghe <i>et al.</i> The paracaspase MALT1 controls cholesterol homeostasis in glioblastoma stem-like cells through lysosome proteome shaping. DOI: 10.1016/j.celrep.2023.113631
RNA-Seq		iScience	2023	Yufei Wei <i>et al.</i> Sirt6 regulates the proliferation of neural precursor cells and cortical neurogenesis in mice. DOI: 10.1016/j.isci.2023.108706
ChIP-Seq	Vitamin D Receptor	JCI Insight	2023	Yuan Wen <i>et al.</i> Vitamin D status associates with skeletal muscle loss after anterior cruciate ligament reconstruction. DOI: 10.1172/jci.insight.170518
CUT&Tag	H3K9me3	Genes (Basel)	2023	Takahiro Sano <i>et al.</i> Impaired Repopulating Ability of Uhrf2-/- Hematopoietic Progenitor Cells in Mice. DOI: 10.3390/genes14081531

Technique	Target	Journal	Year	Reference
ChIP	H3K4me1	Front Endocrinol.	2023	Yeongseik Oh <i>et al.</i> The NR2F2-HAND2 signaling axis regulates progesterone actions in the uterus at early pregnancy. Doi: 10.3389/fendo.2023.1229033
RRBS		Cancers	2023	Yomoki Yamamoto <i>et al.</i> DNA Methylation Aberrations in Dimethylarsinic Acid-Induced Bladder Carcinogenesis. Doi: 10.3390/cancers15215274
ChIP-Seq	H3K79me2	Nat Immunol.	2023	Liangyu Lin <i>et al.</i> Oleic acid availability impacts thymocyte preprogramming and subsequent peripheral Treg cell differentiation. Doi: 0.1038/s41590-023-01672-1
ChIP-Seq	C/EBPα	Nat Commun.	2023	Dongmei Wang <i>et al.</i> Homodimer-mediated phosphorylation of C/EBPα-p42 S16 modulates acute myeloid leukaemia differentiation through liquid-liquid phase separation. Doi: 10.1038/s41467-023-42650-3
CUT&Tag, RNA-Seq	H3K9Ac	Cells	2023	Shah Alam <i>et al.</i> S1P Lyase Deficiency in the Brain Promotes Astroglia and NLRP3 Inflammasome Activation via Purinergic Signaling. DOI: 10.3390/cells12141844
ChIP-Seq	CTCF	FASEB J.	2023	Sylvia C Hewitt <i>et al.</i> Chromatin architectural factor CTCF is essential for progesterone-dependent uterine maturation. doi: 10.1096/fj.202300862R
ChIP-Seq	AhR (aryl hydrocarbon receptor)	Toxicol Sci.	2023	David Filipovic <i>et al.</i> Interpretable predictive models of genome-wide aryl hydrocarbon receptor-DNA binding reveal tissue-specific binding determinants. DOI: 10.1093/toxsci/kfad094
ChIP-Seq, RIME	H3K27ac	Life Sci Alliance	2023	Victor Bessenneau_Gaborit <i>et al.</i> Exploring the impact of dexamethasone on gene regulation in myeloma cells. doi: 10.26508/lsa.202302195
ChIP-Seq	Klf9	Cell Signal.	2023	Chandni Thakkar <i>et al.</i> Klf9 plays a critical role in GR – dependent metabolic adaptations in cardiomyocytes. DOI: 10.1016/j.cellsig.2023.110886
RIME		J Virol.	2023	Lauren McKenzie Spires <i>et al.</i> KSHV RTA utilizes the host E3 ubiquitin ligase complex RNF20/40 to drive lytic reactivation. DOI: 10.1128/jvi.01389-23
ChIP-qPCR	Myc	Mol Oncol.	2023	Marta Kazimierska <i>et al.</i> CRISPR/Cas9 screen for genome-wide interrogation of essential MYC-bound E-boxes in cancer cells. DOI: 10.1002/1878-0261.13493
ChIP	H3K23Pr	J Clin Invest.	2023	Zhi Yang <i>et al.</i> Reducing branched-chain amino acids improves cardiac stress response in mice by decreasing histone H3K23 propionylation. doi: 10.1172/JCI169399
ChIP-Seq	PHF8	Sci Adv.	2023	Song Peng <i>et al.</i> PHF8-GLUL axis in lipid deposition and tumor growth of clear cell renal cell carcinoma. DOI: 10.1126/sciadv.adf3566
ATAC-Seq		Sci Adv.	2023	Raisa A Reyes-Castro <i>et al.</i> Phosphorylated nuclear DICER1 promotes open chromatin state and lineage plasticity of AT2 tumor cells in lung adenocarcinomas. DOI: 10.1126/sciadv.adf6210
ATAC-Seq, hMeDIP-Seq, RNA-Seq		Int J Dev Biol.	2022	Rosa Guzzo <i>et al.</i> Single-cell transcriptomics defines Dot1L interacting partners and downstream target genes in the mouse molar dental pulp. DOI: 10.1387/ijdb.220141db
ChIP-Seq	H3K27ac, PLAGL1, PLAGL2	Acta Neuropathol.	2023	Michaela-Kristina Keck <i>et al.</i> Amplification of the PLAG-family genes—PLAGL1 and PLAGL2—is a key feature of the novel tumor type CNS embryonal tumor with PLAGL amplification. 10.1007/s00401-022-02516-2

Technique	Target	Journal	Year	Reference
RNA-Seq, hMeDIP-Seq		Cleft Palate Craniofac J.	2023	Badam Enkhmandakh <i>et al.</i> 2023. Single-cell Transcriptome Landscape of DNA Methylome Regulators Associated with Orofacial Clefts in the Mouse Dental Pulp. doi: 10.1177/10556656231172296
RNA-Seq		bioRxiv	2023	Marissa Saenz <i>et al.</i> Longitudinal analysis of a dietary mouse model of non-alcoholic fatty liver disease (NAFLD) and non-alcoholic steatohepatitis (NASH). doi: 10.1101/2023.05.19.540989
Mod Spec		Cells	2023	Monika krosel <i>et al.</i> Bromodomain Protein Inhibitors Reorganize the Chromatin of Synovial Fibroblasts. doi: 10.3390/cells12081149
ChIP	Myc Tag	J Mol Biol.	2023	Christalle Chow <i>et al.</i> ZBTB2 is Recruited to a Specific Subset of HIF-1 Target Loci to Facilitate Full Gene Expression Under Hypoxia. doi: 10.1016/j.jmb.2023.168162
ATAC-Seq		Proc Natl Acad Sci USA	2023	Mateus Mota <i>et al.</i> Targeting SWI/SNF ATPases in H3.3K27M diffuse intrinsic pontine gliomas. 10.1073/pnas.2221175120
ChIP-Seq	Eralpha	Nat Commun.	2023	Jose M Ramon-Pittol <i>et al.</i> Dax1 modulates ER α -dependent hypothalamic estrogen sensing in female mice. doi: 10.1038/s41467-023-38618-y
Mod Spec		Adv Sci.	2023	Mingli Li <i>et al.</i> Epigenetic Control of Translation Checkpoint and Tumor Progression via RUVBL1-EEF1A1 Axis. doi: 10.1002/advs.202206584
RNA-Seq		Int Immunopharmacol.	2023	Arijita Subuddhi <i>et al.</i> Comparative transcriptome profile of mouse macrophages treated with the RhoA/Rock pathway inhibitors Y27632, Fingolimod (Gilenya), and Rezerock (Belumosudil, SLx-2119). doi: 10.1016/j.intimp.2023.110017
RNA-Seq		Cell Rep.	2023	Marion Dejosez <i>et al.</i> Regulatory architecture of housekeeping genes is driven by promoter assemblies. doi: 10.1016/j.celrep.2023.112505
Hi-C			2023	Robbin A Nameki <i>et al.</i> Rewiring of master transcription factor cisomes during high-grade serous ovarian cancer development. doi: 10.1101/2023.04.11.536378
ChIP-Seq	H3K27ac	Epigenetics	2023	Tanner B Jefferson <i>et al.</i> Multiple tissue-specific epigenetic alterations regulate persistent gene expression changes following developmental DES exposure in mouse reproductive tissues.
ATAC-Seq		Nature	2023	Hideko Isozaki <i>et al.</i> Therapy-induced APOBEC3A drives evolution of persistent cancer cells. doi: 10.1038/s41586-023-06303-1
ChIP-Seq	H3K27me3, H3K27ac	Nat Commun.	2023	Thaddeus J Kunkel <i>et al.</i> The cholesterol transporter NPC1 is essential for epigenetic regulation and maturation of oligodendrocyte lineage cells. doi: 10.1038/s41467-023-39733-6
ChIP-Seq		Cancer Res.	2023	Dipika R Mohan <i>et al.</i> β -Catenin-Driven Differentiation Is a Tissue-Specific Epigenetic Vulnerability in Adrenal Cancer. doi: 10.1158/0008-5472
RNA-Seq		Nat Commun.	2023	Daniel Schator <i>et al.</i> Legionella para-effectors target chromatin and promote bacterial replication. doi: 10.1038/s41467-023-37885-z
CUT&Tag	H3K9me3	eLife	2023	Suhas Sureshchandra <i>et al.</i> Maternal obesity blunts antimicrobial responses in fetal monocytes. doi: 10.7554/eLife.81320

Technique	Target	Journal	Year	Reference
ChIP-qPCR	RelA, RUNX2	eLife	2023	Daniel Richard <i>et al.</i> Lineage-specific differences and regulatory networks governing human chondrocyte development. doi: 10.7554/eLife.79925
ChIP-Seq	TEAD1	Cell Rep.	2023	Feng Li <i>et al.</i> VGLL4 and MENIN function as TEAD1 corepressors to block pancreatic β cell proliferation. doi: 10.1016/j.celrep.2022.111904
ChIP-Seq	HDAC1, H3K27ac	Nat Commun.	2023	Dasol Han <i>et al.</i> ZBTB12 is a molecular barrier to dedifferentiation in human pluripotent stem cells. doi: 10.1038/s41467-023-36178-9
ChIP-Seq	EZH2, H3K27me3	Nat Commun.	2023	Fan Chen <i>et al.</i> Polycomb deficiency drives a FOXP2-high aggressive state targetable by epigenetic inhibitors. doi: 10.1038/s41467-023-35784-x
CUT&Tag, RRBS	H3K4me3	Adv Sci.	2023	David A McBride <i>et al.</i> Immunomodulatory Microparticles Epigenetically Modulate T Cells and Systemically Ameliorate Autoimmune Arthritis. doi: 10.1002/adv.202202720
RIME	Basonuclin 1 (BNC1)	bioRxiv	2023	Rafik Boudra <i>et al.</i> PRMT1 Inhibition Selectively Targets BNC1-Dependent Proliferation, but not Migration in Squamous Cell Carcinoma. doi: 10.1101/2023.03.27.533164
ATAC-Seq		Nat Commun.	2023	Kuniyuji Aso <i>et al.</i> Itaconate ameliorates autoimmunity by modulating T cell imbalance via metabolic and epigenetic reprogramming. doi: 10.1038/s41467-023-36594-x
ATAC-Seq		NAR Cancer	2023	Anke Koeniger <i>et al.</i> Tumor-suppressive disruption of cancer subtype-associated super enhancer circuits by small molecule treatment. doi: 10.1093/narcan/zcad007
ChIP-Seq	H3K27me3	bioRxiv	2023	Yingqin Hou <i>et al.</i> Transient EZH2 suppression by Tazemetostat during in vitro expansion maintains T cell stemness and improves adoptive T cell therapy. doi: 10.1101/2023.02.07.527459
ChIP-Seq	DUX4	Sci Adv.	2023	Kaoru Inoue <i>et al.</i> DUX4 double whammy: The transcription factor that causes a rare muscular dystrophy also kills the precursors of the human nose. doi: 10.1126/sciadv.abq7744
ChIP-qPCR	H3K9Ac	Cell Rep.	2023	Alison Iroz <i>et al.</i> A Specific ChREBP and PPAR α Cross-Talk Is Required for the Glucose-Mediated FGF21 Response. doi: 10.1016/j.celrep.2017.09.065
ChIP-Seq	IRF-4	Sci Immunol.	2023	Oriol Fornes <i>et al.</i> A multimorphic mutation in IRF4 causes human autosomal dominant combined immunodeficiency. doi: 10.1126/sciimmunol.ade7953
Mod Spec, RRBS		Cell Rep.	2023	Tetsuya Hirabayashi <i>et al.</i> Hepatic phosphatidylcholine catabolism driven by PNPLA7 and PNPLA8 supplies endogenous choline to replenish the methionine cycle with methyl groups. doi: 10.1016/j.celrep.2022.111940
ATAC-Seq		Nat Commun.	2023	Bogang Wu <i>et al.</i> RNA polymerase II pausing factor NELF in CD8 $^{+}$ T cells promotes antitumor immunity. doi: 10.1038/s41467-022-29869-2
ChIP-Seq	H3K27ac	Int J Mol Sci.	2022	Tomoaki Saito <i>et al.</i> Genome-Wide Super-Enhancer-Based Analysis: Identification of Prognostic Genes in Oral Squamous Cell Carcinoma. Int. J. Mol. Sci. doi: 10.3390/ijms23169154.
ATAC-Seq		Cell Rep.	2022	Yayun Gu <i>et al.</i> Multi-omics profiling visualizes dynamics of cardiac development and functions. doi: 10.1016/j.celrep.2022.111891

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K4me1, H3K-4me3, H3K27ac, RNA pol II	Cell Rep.	2022	Yayun Gu <i>et al.</i> Multi-omics profiling visualizes dynamics of cardiac development and functions. doi: 10.1016/j.celrep.2022.111891
ATAC-Seq		Nucleic Acids Res.	2022	Sebastian Gregoricchio <i>et al.</i> HDAC1 and PRC2 mediate combinatorial control in SPI1/PU.1-dependent gene repression in murine erythroleukaemia. doi: 10.1093/nar/gkac613
ChIP-Seq	H3K4me3, H3K-9me3	Mol. Ther.	2022	Britta Handyside <i>et al.</i> Mol. Ther. Vector genome loss and epigenetic modifications mediate decline in transgene expression of AAV5 vectors produced in mammalian and insect cells. doi: 10.1016/j.jymthe.2022.11.001
Hi-C		Cells	2022	Sylvia C. Hesitt <i>et al.</i> Progesterone Signaling in Endometrial Epithelial Organoids. doi: 10.3390/cells11111760
ChIP-Seq	H3K27ac	NPJ Precis Oncol.	2022	Steven W. Criscione <i>et al.</i> The landscape of therapeutic vulnerabilities in EGFR inhibitor osimertinib drug tolerant persister cells. doi: 10.1038/s41698-022-00337-w
scATAC-Seq		Cell Rep Med.	2022	Yimiing Zhang <i>et al.</i> Pegylated arginine deiminase drives arginine turnover and systemic autophagy to dictate energy metabolism. doi: 10.1016/j.xcrm.2021.100498
RNA-Seq, ATAC-Seq		Inflamm Regen.	2022	Sakurako Kobayashi <i>et al.</i> Collagen type I-mediated mechanotransduction controls epithelial cell fate conversion during intestinal inflammation. doi: 10.1186/s41232-022-00237-3
ATAC-Seq		Cell Death Differ.	2022	Jeehyun Yoon <i>et al.</i> E2F and STAT3 provide transcriptional synergy for histone variant H2AZ activation to sustain glioblastoma chromatin accessibility and tumorigenicity. doi: 10.1038/s41418-021-00926-5
ChIP-Seq	p300	Nucleic Acids Res.	2022	Avital Sarusi Portuguese <i>et al.</i> Ep300 sequestration to functionally distinct glucocorticoid receptor binding loci underlie rapid gene activation and repression. doi: 10.1093/nar/gkac488
ChIP-Seq	KLF4	J. Invest Dermatol.	2022	Yao Yuan <i>et al.</i> Oncogenic Hedgehog-Smoothed Signaling Depends on YAP1-TAZ/TEAD Transcription to Restrain Differentiation in Basal Cell Carcinoma. doi: 10.1016/j.jid.2021.06.020
ChIP-Seq	H3K27ac, H3K-27me3, BRG1	Cancer Discov.	2022	Eshini Padnitharatna <i>et al.</i> BAF Complex Maintains Glioma Stem Cells in Pediatric H3K27M Glioma. doi: 10.1158/2159-8290.CD-21-1491
RNA-Seq		Nat Metab.	2022	Ana P. Gomes <i>et al.</i> Altered propionate metabolism contributes to tumour progression and aggressiveness. doi: 10.1038/s42255-022-00553-5
ChIP-Seq		eLife	2022	Antigoni Gogolou <i>et al.</i> Early anteroposterior regionalisation of human neural crest is shaped by a pro-mesodermal factor. doi: 10.7554/eLife.74263
ChIP-Seq	TEAD1	Nucleic Acids Res.	2022	Feng Li <i>et al.</i> TEAD1 regulates cell proliferation through a pocket-independent transcription repression mechanism. doi: 10.1093/nar/gkac1063
ChIP-qPCR, ATAC-Seq	H3R2me1, H3R2me2s	Nat Commun.	2022	Gizem Gunes Günsel <i>et al.</i> The arginine methyltransferase PRMT7 promotes extravasation of monocytes resulting in tissue injury in COPD. doi: 10.1038/s41467-022-28809-4
ATAC-Seq		Blood Adv.	2022	Genki Yamato <i>et al.</i> Genome-wide DNA methylation analysis in pediatric acute myeloid leukemia. doi: 10.1182/bloodadvances.2021005381.

Technique	Target	Journal	Year	Reference
ATAC-Seq, RNA-Seq		Cell Death Differ.	2022	Kristina Seiler <i>et al.</i> Hexokinase 3 enhances myeloid cell survival via non-glycolytic functions. doi: 10.1038/s41419-022-04891-w.
ATAC-Seq		Mucosal Immunol.	2022	Soh Yamazaki <i>et al.</i> I κ B ζ controls IL-17-triggered gene expression program in intestinal epithelial cells that restricts colonization of SFB and prevents Th17-associated pathologies. doi: 10.1038/s41385-022-00554-3.
ATAC-Seq		Sci Adv.	2022	Jean Personnaz <i>et al.</i> Nuclear HMGB1 protects from nonalcoholic fatty liver disease through negative regulation of liver X receptor. doi: 10.1126/sciadv.abg9055.
ATAC-Seq		Front Med (Lausanne)	2022	Jehan Alam <i>et al.</i> IL-17 Producing Lymphocytes Cause Dry Eye and Corneal Disease With Aging in RXRa Mutant Mouse. doi: 10.3389/fmed.2022.849990.
RNA-Seq		Mol Oncol.	2022	Carlotta Zampieri <i>et al.</i> p53 mutations define the chromatin landscape to confer drug tolerance in pancreatic cancer. doi: 10.1002/1878-0261.13161.
ChIP-Seq	Androgen Receptor	Oncogene	2022	Oh-Joon Kwon <i>et al.</i> Elevated expression of the colony-stimulating factor 1 (CSF1) induces prostatic intraepithelial neoplasia dependent of epithelial-Gp130. doi: 10.1038/s41388-021-02169-7.
ChIP-Seq	Estrogen Receptor alpha	Endocrinology	2022	Sylvia C. Hewitt <i>et al.</i> The Estrogen Receptor α Cistrome in Human Endometrium and Epithelial Organoids. doi: 10.1210/endocr/bqac116.
ATAC-Seq		Mol Cancer Ther.	2022	Aaron Burmeister <i>et al.</i> Establishment and Evaluation of Dual HDAC/BET Inhibitors as Therapeutic Options for Germ Cell Tumors and Other Urological Malignancies. doi: 10.1158/1535-7163.MCT-22-0207.
ChIP-Seq	H3K36me3	Nat Commun.	2022	Zhaoyun Ding <i>et al.</i> Setd2 supports GATA3+ST2+ thymic-derived Treg cells and suppresses intestinal inflammation. doi: 10.1038/s41467-022-35250-0.
ChIP-Seq	p53, H3K4me3, H3K9me3	Cells	2022	Takatsune Shimizu <i>et al.</i> Depletion of R270C Mutant p53 in Osteosarcoma Attenuates Cell Growth but Does Not Prevent Invasion and Metastasis In Vivo. doi: 10.3390/cells11223614.
ChIP-Seq	Progesterone Receptor	Reprod Sci.	2022	Ariel J. Dotts <i>et al.</i> In Vivo Genome-Wide PGR Binding in Pregnant Human Myometrium Identifies Potential Regulators of Labor. doi: 10.1007/s43032-022-01002-0.
ATAC-Seq, RNA-Seq		Acta Biochim Pol.	2022	Pujan Joshi <i>et al.</i> The chromatin accessibility landscape in the dental pulp of mouse molars and incisors. doi: 10.18388/abp.2020_5771.
ChIP-qPCR	FOXM1	Int J Mol Sci.	2022	Erkut Ilaslan <i>et al.</i> Distinct Roles of NANOS1 and NANOS3 in the Cell Cycle and NANOS3-PUM1-FOXM1 Axis to Control G2/M Phase in a Human Primordial Germ Cell Model. doi: 10.3390/ijms23126592.
ChIP-Seq	Estrogen Receptor beta	J Biomed Sci.	2022	Yuri Park <i>et al.</i> Oleuropein suppresses endometriosis progression and improves the fertility of mice with endometriosis. doi: 10.1186/s12929-022-00883-2.
ChIP-Seq	CTCF	Int J Mol Sci.	2022	Clarissa Boschiero <i>et al.</i> Differentially CTCF-Binding Sites in Cattle Rumen Tissue during Weaning. doi: 10.3390/ijms23169070.
ChIP-Seq	CTCF	Biomolecules	2022	Clarissa Boschiero <i>et al.</i> Butyrate Induces Modifications of the CTCF-Binding Landscape in Cattle Cells. doi: 10.3390/biom12091177.

Technique	Target	Journal	Year	Reference
ATAC-Seq, RNA-Seq		Autophagy	2022	Eutteum Jeong <i>et al.</i> The FACT complex facilitates expression of lysosomal and antioxidant genes through binding to TFEB and TFE3. <i>Autophagy</i> . doi: 10.1080/15548627.2022.2029671.
ChIP-Seq	BRD4	Biomed Pharmacother.	2022	Laura M. Tsujikawa <i>et al.</i> Breaking boundaries: Pan BETi disrupt 3D chromatin structure, BD2-selective BETi are strictly epigenetic transcriptional regulators. <i>Biomed. Pharmacother.</i> doi: 10.1016/j.biopha.2022.113230.
ATAC-Seq		Cancer Immunol Res.	2022	Andrew Fedoriw <i>et al.</i> Inhibiting Type I arginine methyltransferase activity promotes the T cell mediated antitumor immune response. <i>Cancer Immunol. Res.</i> doi: 10.1158/2326-6066.CIR-21-0614.
ATAC-Seq		Cell Rep Med.	2022	Sally Mortlock <i>et al.</i> A multi-level investigation of the genetic relationship between endometriosis and ovarian cancer histotypes. <i>Cell Rep Med.</i> doi: 10.1016/j.xcrm.2022.100542.
ChIP-Seq	RNA Pol II	Cell Rep.	2022	Shin-Ichiro Ohno <i>et al.</i> Nuclear microRNAs release paused Pol II via the DDX21-CDK9 complex. <i>Cell Rep.</i> doi: 10.1016/j.celrep.2022.110673.
ChIP-Seq	MAX	Cell Tissue Res.	2022	Weixu Ma <i>et al.</i> MAX deficiency impairs human endometrial decidualization through down-regulating OSR2 in women with recurrent spontaneous abortio. <i>Cell Tissue Res.</i> doi: 10.1007/s00441-022-03579-z.
ChIP-Seq	MAX	Cell Tissue Res.	2022	Weixu Ma <i>et al.</i> MAX deficiency impairs human endometrial decidualization through down-regulating OSR2 in women with recurrent spontaneous abortion. <i>Cell Tissue Res.</i> doi: 10.1007/s00441-022-03579-z.
ChIP-Seq	SMARCA4, H3K- 27me3	Cells	2022	Amelie Alfert <i>et al.</i> Smarcb1 Loss Results in a Deregulation of esBAF Binding and Impacts the Expression of Neurodevelopmental Genes. <i>Cells.</i> doi: 10.3390/cells11081354.
ATAC-Seq		Commun Biol.	2022	Hidenori Machino <i>et al.</i> The metabolic stress-activated checkpoint LKB1-MARK3 axis acts as a tumor suppressor in high-grade serous ovarian carcinoma. <i>Commun. Biol.</i> doi: 10.1038/s42003-021-02992-4.
ATAC-Seq		Commun Biol.	2022	Hidenori Machino <i>et al.</i> The metabolic stress-activated checkpoint LKB1-MARK3 axis acts as a tumor suppressor in high-grade serous ovarian carcinoma. <i>Commun. Biol.</i> doi: 10.1038/s42003-021-02992-4.
ChIP-Seq	H3K27ac	CRISPR J.	2022	Antonia A. Dominguez <i>et al.</i> CRISPR-Mediated Synergistic Epigenetic and Transcriptional Control. <i>CRISPR J.</i> doi: 10.1089/crispr.2021.0099.
ChIP-Seq	H3K27ac	CRISPR J.	2022	Antonia A. Dominuez <i>et al.</i> CRISPR-mediated Synergistic Epigenetic and Transcriptional Control. <i>CRISPR J.</i> doi: 10.1089/crispr.2021.0099.
MeDIP-Seq		EMBO Rep.	2022	Antonia Piazzesi <i>et al.</i> CEST-2.2 overexpression alters lipid metabolism and extends longevity of mitochondrial mutants. <i>EMBO Rep.</i> doi: 10.15252/embr.202152606.
ChIP-Seq	H3K4me1, H3K- 4me3	FASEB J.	2022	Matthew Baxter <i>et al.</i> Circadian clock function does not require the histone methyltransferase MLL3. <i>FASEB J.</i> doi: 10.1096/fj.202200368R.

Technique	Target	Journal	Year	Reference
ChIP-Seq	BRD4, H3K27ac	Front Cell Dev Biol.	2022	Cho-Hao Lin <i>et al.</i> AZD5153, a Bivalent BRD4 Inhibitor, Suppresses Hepatocarcinogenesis by Altering BRD4 Chromosomal Landscape and Modulating the Transcriptome of HCC Cells. <i>Fron Cell Dev Biol.</i> doi: 10.3389/fcell.2022.853652.
ChIP-Seq	DNMT1	Gastroenterology	2022	Shikai Hu <i>et al.</i> NOTCH-YAP1/TEAD-DNMT1 Axis Drives Hepatocyte Reprogramming Into Intrahepatic Cholangiocarcinoma. <i>Gastroenterology.</i> doi: 10.1053/j.gastro.2022.05.007.
ATAC-Seq		Genes (Basel)	2022	Clarissa Boschiero <i>et al.</i> Characterization of Accessible Chromatin Regions in Cattle Rumen Epithelial Tissue during Weaning. <i>Genes (Basel).</i> doi: 10.3390/genes13030535.
ATAC-Seq		Genes (Basel)	2022	Clarissa Boschiero <i>et al.</i> Characterization of Accessible Chromatin Regions in Cattle Rumen Epithelial Tissue during Weaning. <i>Genes (Basel)</i> doi: 10.3390/genes13030535.
ChIP-Seq	CTCF, H3K4me3, H3K27ac	Genome Res.	2022	Blair W. Perry <i>et al.</i> Snake venom gene expression is coordinated by novel regulatory architecture and the integration of multiple co-opted vertebrate pathways. <i>Genome Res.</i> doi: 10.1101/gr.276251.121.
ATAC-Seq, ChIP-Seq	H3K27ac, H3K-27me3, H3K4me1, H3K4me3, CTCF	Genomics	2022	Yahui Gae <i>et al.</i> Functional annotation of regulatory elements in cattle genome reveals the roles of extracellular interaction and dynamic change of chromatin states in rumen development during weaning. <i>Genomics.</i> doi: 10.1016/j.ygeno.2022.110296
ChIP-Seq	H3K4me3	Hepatology Commun.	2022	Michael Schonfeld <i>et al.</i> Alcohol-associated fibrosis in females is mediated by female-specific activation of lysine demethylases KDM5B and KDM5C. <i>Hepatology Commun.</i> doi: 10.1002/hep4.1967.
ChIP-Seq	H3K4me3	Hepatology Commun.	2022	Michael Schonfeld <i>et al.</i> Male-Specific Activation of Lysine Demethylases 5B and 5C Mediates Alcohol-Induced Liver Injury and Hepatocyte Dedifferentiation. <i>Hepatology Commun.</i> doi: 10.1002/hep4.1895.
ChIP-Seq	H3K4me3	Hepatology Commun.	2022	Michael Schonfeld <i>et al.</i> Male-Specific Activation of Lysine Demethylases 5B and 5C Mediates Alcohol-Induced Liver Injury and Hepatocyte Dedifferentiation. <i>Hepatology Commun.</i> doi: 10.1002/hep4.1895.
RIME	EOMES	iScience	2022	James A. Heslop <i>et al.</i> Chromatin remodeling is restricted by transient GATA6 binding during iPSC differentiation to definitive endoderm. <i>iScience.</i> doi: 10.1016/j.isci.2022.104300.
ChIP-Seq	YAP	J Clin Invest.	2022	Toshihide Kashihara <i>et al.</i> YAP mediates compensatory cardiac hypertrophy through aerobic glycolysis in response to pressure overload. <i>J. Clin. Invest.</i> doi: 10.1172/JCI150595.
ChIP-Seq	H3K9me3	J Immunol.	2022	Jun Hyung Sin <i>et al.</i> ATF7ip Targets Transposable Elements for H3K9me3 Deposition to Modify CD8+ T Cell Effector and Memory Responses. <i>J. Immunol.</i> doi: 10.4049/jimmunol.2100996.
ChIP-Seq	H3K9me3	J Immunol.	2022	Jun Hyung Sin <i>et al.</i> ATF7ip Targets Transposable Elements for H3K9me3 Deposition to Modify CD8 + T Cell Effector and Memory Responses. <i>J. Immunol.</i> doi: 10.4049/jimmunol.2100996.

Technique	Target	Journal	Year	Reference
ATAC-Seq		Mol Cancer Res.	2022	Aditya Wirawan <i>et al.</i> A Novel Therapeutic Strategy Targeting the Mesenchymal Phenotype of Malignant Pleural Mesothelioma by Suppressing LSD1. Mol. Cancer Res. doi: 10.1158/1541-7786.MCR-21-0230.
ChIP-qPCR	WT1	Mol Oncol.	2022	Alain Chebly <i>et al.</i> Exploring hTERT promoter methylation in cutaneous T-cell lymphomas. Mol Oncol. doi: 10.1002/1878-0261.12946
ChIP-qPCR	H3K27ac, H3K-4me2	Nat Commun.	2022	Emily V. Dutrow <i>et al.</i> Modeling uniquely human gene regulatory function via targeted humanization of the mouse genome. Nat. Commun. doi: 10.1038/s41467-021-27899-w.
ChIP-Seq	H3K27ac, RNA pol II, ΔNp63	Nat Commun.	2022	Marco Napoli <i>et al.</i> ΔNp63 regulates a common landscape of enhancer associated genes in non-small cell lung cancer. Nat. Commun. doi: 10.1038/s41467-022-28202-1.
ChIP-qPCR, Antibody Validation	H3K27ac, H3K-4me2	Nat Commun.	2022	Emily V. Dutrow <i>et al.</i> Modeling uniquely human gene regulatory function via targeted humanization of the mouse genome. Nat. Commun. doi: 10.1038/s41467-021-27899-w.
ChIP-qPCR, Antibody Validation	KLF5	Sci Adv.	2022	Juliane Mooz <i>et al.</i> ARAF suppresses ERBB3 expression and metastasis in a subset of lung cancers. Sci. Adv. doi: 10.1126/sciadv.abk1538.
ChIP-Seq	H3K9ac	Theranostics	2022	Grace Huang <i>et al.</i> Diabetes impairs cardioprotective function of endothelial progenitor cell-derived extracellular vesicles via H3K9Ac inhibition. Theranostics. doi: 10.7150/thno.70821.
ChIP-Seq	H3K27ac, RelA	Acta Neuropathol.	2021	Daniela Lotsch <i>et al.</i> Targeting fibroblast growth factor receptors to combat aggressive ependymoma. Acta Neuropathol. doi: 10.1007/s00401-021-02327-x.
ChIP	H3K27ac, RelA	Acta Neuropathol.	2021	Daniela Lotsch <i>et al.</i> Targeting fibroblast growth factor receptors to combat aggressive ependymoma. Acta Neuro. doi: 10.1007/s00401-021-02327
RNA-Seq		Am J Physiol Heart Circ Physiol.	2021	Satvik Mareedu <i>et al.</i> Sarcolipin haploinsufficiency prevents dystrophic cardiomyopathy in mdx mice. Am J Physiol Heart Circ Physiol. doi: 10.1152/ajpheart.00601.2020.
ChIP-Seq	Sp6	Biochem Biophys Res Commun.	2021	Craig S. Rhodes <i>et al.</i> Sp6/Epiprofin is a master regulator in the developing tooth. Biochem. Biophys. Res. Commun. doi: 10.1016/j.bbrc.2021.10.017.
Mod Spec		Blood	2021	Zachary C. Murphy <i>et al.</i> Regulation of RNA polymerase II activity is essential for terminal erythroid maturation. Blood. doi: 10.1182/blood.2020009903.
ATAC-Seq		Blood	2021	Julie Agopian <i>et al.</i> GlcNAc is a mast-cell chromatin-remodeling oncometabolite that promotes systemic mastocytosis aggressiveness. Blood. doi: 10.1182/blood.2020008948.
ChIP-qPCR	LXRalpha	Br J Pharmacol.	2021	Louise Menegaut <i>et al.</i> Regulation of glycolytic genes in human macrophages by oxysterols: a potential role for liver X receptors. Br. J. Pharmacol. doi: 10.1111/bph.15358.
ChIP-Seq	Multi-Histones	Cancer Cell	2021	Daniel Haag <i>et al.</i> H3.3-K27M drives neural stem cell-specific gliomagenesis in a human iPSC-derived model. Cancer Cell. doi: 10.1016/j.ccell.2021.01.005.

Technique	Target	Journal	Year	Reference
ATAC-Seq		Cancer Discov.	2021	Sonia Iyer <i>et al.</i> Genetically Defined Syngeneic Mouse Models of Ovarian Cancer as Tools for the Discovery of Combination Immunotherapy. <i>Cancer Discov.</i> doi: 10.1158/2159-8290
ATAC-Seq, RNA-Seq		Cancer Immunol Res.	2021	Justin C. Boucher <i>et al.</i> CD28 Costimulatory Domain-Targeted Mutations Enhance Chimeric Antigen Receptor T-cell Function. <i>Cancer Immunol. Res.</i> doi: 10.1158/2326-6066.CIR-20-0253.
ATAC-Seq		Cancers (Basel)	2021	Dongho Kim <i>et al.</i> Metformin Reduces Histone H3K4me3 at the Promoter Regions of Positive Cell Cycle Regulatory Genes in Lung Cancer Cells. <i>Cancers (Basel)</i> . doi: 10.3390/cancers13040739.
Mod Spec		Cancers (Basel)	2021	Anke Koeniger <i>et al.</i> Activation of Cilia-Independent Hedgehog/GLI1 Signaling as a Novel Concept for Neuroblastoma Therapy. <i>Cancers (Basel)</i> . doi: 10.3390/cancers13081908
ATAC-Seq		Cell	2021	Florian Wimmers <i>et al.</i> The single-cell epigenomic and transcriptional landscape of immunity to influenza vaccination. <i>Cell</i> . doi: 10.1016/j.cell.2021.05.039.
ChIP-Seq	TDF4	Cell Chem. Biol.	2021	Jiongjia Cheng <i>et al.</i> Small-molecule probe reveals a kinase cascade that links stress signaling to TCF/LEF and Wnt responsiveness. <i>Cell Chem. Biol.</i> doi: 10.1016/j.chembiol.2021.01.001
ATAC-Seq		Cell Rep.	2021	Ho Sun Jung <i>et al.</i> SOX17 integrates HOXA and arterial programs in hemogenic endothelium to drive definitive lympho-myeloid hematopoiesis. <i>Cell Rep.</i> doi: 10.1016/j.celrep.2021.108758.
RIME	RelA	Cell Rep.	2021	James A. Heslop <i>et al.</i> GATA6 defines endoderm fate by controlling chromatin accessibility during differentiation of human-induced pluripotent stem cells. <i>Cell Rep.</i> doi: 10.1016/j.celrep.2021.109145
RNA-Seq		Cells	2021	Melanie Lavaud <i>et al.</i> Overexpression of the Ubiquitin Specific Proteases USP43, USP41, USP27x and USP6 in Osteosarcoma Cell Lines: Inhibition of Osteosarcoma Tumor Growth and Lung Metastasis Development by the USP Antagonist PR619. <i>Cells</i> . doi: 10.33990/cells10092268
ChIP-Seq	p300	Clin Transl Med.	2021	Chuanjie Zhang <i>et al.</i> Epigenome screening highlights that JMJD6 confers an epigenetic vulnerability and mediates sunitinib sensitivity in renal cell carcinoma. <i>Clin. Transl. Med.</i> doi: 10.1002/ctm2.328.
ChIP-Seq	FANCD2	Commun Biol.	2021	Philippe Fernandes <i>et al.</i> FANCD2 modulates the mitochondrial stress response to prevent common fragile site instability. <i>Commun Biol.</i> doi: 10.1038/s42003-021-01647-8
Histone PTM Analysis		Dev Cell	2021	Igor L. Bado <i>et al.</i> The bone microenvironment increases phenotypic plasticity of ER + breast cancer cells. <i>Dev. Cell.</i> doi: 10.1016/j.devcel.2021.03.008
RNA-Seq		Development	2021	Matthias Godart <i>et al.</i> The murine intestinal stem cells are highly sensitive to the modulation of the T3/TRa1-dependent pathway. <i>Development</i> . doi: 10.1242/dev.194357
ChIP-Seq	H3K79me2	Diabetes	2021	Lin Shuai <i>et al.</i> DOT1L Regulates Thermogenic Adipocyte Differentiation and Function via Modulating H3K79 Methylation. <i>Diabetes</i> . doi: 10.2337/db20-1110

Technique	Target	Journal	Year	Reference
PTM Quantitation		eLife	2021	Helen M. Tauc <i>et al.</i> Age-related changes in polycomb gene regulation disrupt lineage fidelity in intestinal stem cells. eLife. doi: 10.7554/eLife.62250
ATAC-Seq		EMBO Rep.	2021	Raquel Sales-Gil <i>et al.</i> Non-redundant functions of H2A.Z.1 and H2A.Z.2 in chromosome segregation and cell cycle progression. EMBO Rep. doi: 10.15252/embr.202052061.
ATAC-Seq		EMBO Rep.	2021	Raquel Sales-Gil <i>et al.</i> Non-redundant functions of H2A.Z.1 and H2A.Z.2 in chromosome segregation and cell cycle progression. EMBO Reports. doi: 10.15252/embr.202052061
ChIP-Seq	ZNF263	ERJ Open Res.	2021	Sai Sneha Priya Nemani <i>et al.</i> COL4A3 expression in asthmatic epithelium depends on intronic methylation and ZNF263 binding. ERJ Open Res. doi: 10.1183/23120541.00802-2020
ATAC-Seq		Int J Mol Sci.	2021	Lilas Courtot <i>et al.</i> Low Replicative Stress Triggers Cell-Type Specific Inheritable Advanced Replication Timing. Int. J. Mol. Sci. doi.org/10.3390/ijms22094959
ChIP-qPCR	FOXO3	J Cell Physiol.	2021	Siyun Wang <i>et al.</i> cMET promotes metastasis and epithelial-mesenchymal transition in colorectal carcinoma by repressing RKIP. J. Cell Physiol. doi: 10.1002/jcp.30142.
ChIP-Seq	BRD2, BRD4, RNA Pol II	J Immunol.	2021	Vinod Krishna <i>et al.</i> Integration of the Transcriptome and Genome-Wide Landscape of BRD2 and BRD4 Binding Motifs Identifies Key Superenhancer Genes and Reveals the Mechanism of Bet Inhibitor Action in Rheumatoid Arthritis Synovial Fibroblasts. J. Immunol. doi: doi: 10.4049/jimmunol.2000286.
ATAC-Seq		J Mol Cell Cardiol.	2021	Douglas J. Chapski <i>et al.</i> Early adaptive chromatin remodeling events precede pathologic phenotypes and are reinforced in the failing heart. J. Mol. Cell Cardiol. doi: 10.1016/j.jmcc.2021.07.002.
RIME	Estrogen Receptor- α	Mol Cancer Res.	2021	Joseph L. Sottnik <i>et al.</i> Mediator of DNA Damage Checkpoint 1 (MDC1) Is a Novel Estrogen Receptor Coregulator in Invasive Lobular Carcinoma of the Breast. Mol. Cancer Res. doi: 10.1158/1541-7786.MCR-21-0025.
ChIP-qPCR	H3K27ac	Mol Cell	2021	Liang Ma <i>et al.</i> Co-condensation between transcription factor and coactivator p300 modulates transcriptional bursting kinetics. Mol Cell. doi: 10.1016/j.molcel.2021.01.031
ChIP-qPCR	WT1	Mol Oncol.	2021	Alain Chebly <i>et al.</i> Exploring hTERT promoter methylation in cutaneous T-cell lymphomas. Mol. Oncol. doi: 10.1002/1878-0261.12946.
ChIP-Seq	BRD4, H3K27ac, H3K4me1	Mol Ther Methods Clin Dev.	2021	Qingwei Wang <i>et al.</i> A hierarchical and collaborative BRD4/CEBPD partnership governs vascular smooth muscle cell inflammation. Mol. Ther. Methods Clin. Dev. doi: 10.1016/j.omtm.2021.02.021
ATAC-Seq		Nat Cancer	2021	Katherine R. Morel <i>et al.</i> EZH2 inhibition activates a dsRNA–STING–interferon stress axis that potentiates response to PD-1 checkpoint blockade in prostate cancer. Nat Cancer. doi: 0.1038/s43018-021-00185-w
ChIP-Seq	BRD4	Nat Chem Biol.	2021	Gabi Schutzius <i>et al.</i> BET bromodomain inhibitors regulate keratinocyte plasticity. Nat. Chem. Biol. doi: 10.1038/s41589-020-00716-z.
ChIP-Seq	H2AK119Ub, Ring1b	Nat Chem Biol.	2021	Shirish Shukla <i>et al.</i> Small-molecule inhibitors targeting Polycomb repressive complex 1 RING domain. Nature Chemical Biology. doi: 10.1038/s41589-021-00815-5

Technique	Target	Journal	Year	Reference
ATAC-Seq		Nat Commun.	2021	Joana Esteves de Lima <i>et al.</i> HIRA stabilizes skeletal muscle lineage identity. Nat. Commun. doi: 10.1038/s41467-021-23775-9.
ChIP-Seq	BMAL1	Nat Commun.	2021	Eleonore Maury <i>et al.</i> Circadian clock dysfunction in human omental fat links obesity to metabolic inflammation. Nat. Commun. doi: 10.1038/s41467-021-22571-9.
ATAC-Seq		Nat Commun.	2021	Joana Esteves de Lima <i>et al.</i> HIRA stabilizes skeletal muscle lineage identity. Nature. doi: 10.1038/s41467-021-23775-9
RRBS		Nucleic Acids Res.	2021	Miriam Recalde <i>et al.</i> The splicing regulator SLU7 is required to preserve DNMT1 protein stability and DNA methylation. Nuc. Acids Res. doi: 10.1093/narlgkab649
ATAC-Seq		PLoS Pathog.	2021	Naveen Parmar <i>et al.</i> Intestinal-epithelial LSD1 controls goblet cell maturation and effector responses required for gut immunity to bacterial and helminth infection. PLoS Pathog. doi: 10.1371/journal.ppat.1009476.
ChIP-qPCR	H3K27ac	RNA	2021	Ryan L. Setten <i>et al.</i> CRED9: A differentially expressed elncRNA regulates expression of transcription factor CEBPA. RNA. doi: 10.1261/rna.078752.121.
RRBS		Sci Adv.	2021	Rina Baba <i>et al.</i> LSD1 enzyme inhibitor TAK-418 unlocks aberrant epigenetic machinery and improves autism symptoms in neurodevelopmental disorder models. Sci Adv. doi: 10.1126/sciadv.aba1187
ChIP-Seq	NR2F2	Sex Dev.	2021	Fei Zhao <i>et al.</i> Molecular Actions Underlying Wolffian Duct Regression in Sexual Differentiation of Murine Reproductive Tracts. Sex Dev. doi: 10.1159/000513878
ChIP-Seq	Histone PTMs	Toxicol Sci.	2021	Jo Jongpyo Lim <i>et al.</i> Neonatal Exposure to BPA, BDE-99, and PCB Produces Persistent Changes in Hepatic Transcriptome Associated With Gut Dysbiosis in Adult Mouse Livers. Toxicol. Sci. doi: 10.1093/toxsci/kfab104.
ChIP-Seq	H3K9me3, H3K-27me3	Arch Toxicol.	2020	Masaki Fujioka <i>et al.</i> Dimethylarsinic acid (DMA) enhanced lung carcinogenesis via histone H3K9 modification in a transplacental mouse model. Arch. Toxicol. doi: 10.1007/s00204-020-02665-x.
RNA-Seq		Blood Adv.	2020	Iosifina P. Foskolou <i>et al.</i> The S enantiomer of 2-hydroxyglutarate increases central memory CD8 populations and improves CAR-T therapy outcome. Blood Adv. doi: 10.1182/bloodadvances.2020002309.
Mod Spec		Blood Adv.	2020	Tiziana Bruno <i>et al.</i> Che-1/AATF-induced transcriptionally active chromatin promotes cell proliferation in multiple myeloma. Blood Adv. doi: 10.1182/bloodadvances.2020002566.
ChIP-Seq	H3K27me3	Cancer Cell	2020	Huairui Yuan <i>et al.</i> SETD2 Restricts Prostate Cancer Metastasis by Integrating EZH2 and AMPK Signaling Pathways. Cancer Cell. doi: 10.1016/j.ccell.2020.05.022
ATAC-Seq		Cancer Discov.	2020	Di Zhao <i>et al.</i> Chromatin Regulator, CHD1, Remodels the Immunosuppressive Tumor Microenvironment in PTEN-Deficient Prostate Cancer. Cancer Discov. 2020 doi: 10.1158/2159-8290.CD-19-1352
RNA-Seq		Cancers (Basel)	2020	Mathilde Mullard <i>et al.</i> Sonic Hedgehog Signature in Pediatric Primary Bone Tumors: Effects of the GLI Antagonist GANT61 on Ewing's Sarcoma Tumor Growth. Cancers (Basel). doi: 10.3390/cancers12113438.

Technique	Target	Journal	Year	Reference
RNA-Seq		Cancers (Basel)	2020	Sarah Morice <i>et al.</i> The YAP/TEAD Axis as a New Therapeutic Target in Osteosarcoma: Effect of Verteporfin and CA3 on Primary Tumor Growth. <i>Cancers (Basel)</i> . doi: 10.3390/cancers12123847.
ChIP-qPCR	BRD2, BRD4	Cardiovasc Ther.	2020	Sylvia Wasiak <i>et al.</i> Epigenetic Modulation by Apabetalone Counters Cytokine-Driven Acute Phase Response In Vitro, in Mice and in Patients with Cardiovascular Disease. <i>Cardiovasc. Ther.</i> doi: 10.1155/2020/9397109.
RIME	Bcl-6	Cell	2020	Leandro Venturutti <i>et al.</i> TBL1XR1 Mutations Drive Extranodal Lymphoma by Inducing a Pro-tumorigenic Memory Fate. <i>Cell</i> . doi: 10.1016/j.cell.2020.05.049.
ChIP-Seq	CEBP, H3K4me3, H3K27ac	Cell	2020	Ada Ndoja <i>et al.</i> Ubiquitin Ligase COP1 Suppresses Neuroinflammation by Degrading c/EBP β in Microglia. <i>Cell</i> . doi: 10.1016/j.cell.2020.07.011.
ChIP-qPCR	HIF1alpha	Cell Rep.	2020	Louise Ménégaud <i>et al.</i> Interplay between Liver X Receptor and Hypoxia Inducible Factor 1a Potentiates Interleukin-1b Production in Human Macrophages. <i>Cell Reports</i> . doi: 10.1016/j.celrep.2020.107665
ChIP-Seq	H3K27ac, AFF4	Commun Biol.	2020	Yujing Gao <i>et al.</i> Acetylation of histone H3K27 signals the transcriptional elongation for estrogen receptor alpha. <i>Commun Biol</i> . 2020 doi: 10.1038/s42003-020-0898-0
ATAC-Seq		Diabetes	2020	Marion de Toledo <i>et al.</i> Lamin C counteracts glucose intolerance in aging, obesity and diabetes. <i>Diabetes</i> . 2020 doi: 10.2337/db19-0377.
ChIP-Seq	HOXB13, H3K27ac, H3K27me3, H3K-4me3, RNA Pol II	Exp Cell Res.	2020	Paul-Joseph Aspuria <i>et al.</i> HOXB13 controls cell state through super-enhancers. <i>Exp Cell Res</i> . 2020 doi: 10.1016/j.yexcr.2020.112039
ChIP-Seq	b-catenin	Exp Mol Med.	2020	Jung-Yoon Yoo <i>et al.</i> β -catenin activates TGF- β -induced epithelial-mesenchymal transition in adenomyosis. <i>Exp. Mol. Med.</i> doi: 10.1038/s12276-020-00514-6
ChIP-Seq	Estrogen Receptor	FASEB J.	2020	Yukitomo Arao <i>et al.</i> The genomic regulatory elements for estrogen receptor alpha transactivation-function-1 regulated genes. <i>FASEB J.</i> doi: 10.1096/fj.202001435R.
ChIP-Seq	N-MYC, C-MYC	Front Oncol.	2020	Robyn T. Sussman <i>et al.</i> CAMKV Is a Candidate Immunotherapeutic Target in MYCN Amplified Neuroblastoma. <i>Front. Oncol.</i> 2020 doi: 10.3389/fonc.2020.00302
ChIP-Seq	H3K4me1, H3K-4me3, H3K9me3, H3K27me3, H3K27ac, H3K-36me3, BRD4, MED1, MYC	Haematologica	2020	"Aneta Mikulasova <i>et al.</i> Microhomology-mediated end joining drives complex rearrangements and over expression of MYC and PVT1 in multiple myeloma. <i>Haematologica</i> .doi:10.3324/haematol.2019.217927"
RNA-Seq		iScience	2020	Aarti Mishra <i>et al.</i> Dynamic Neuroimmune Profile during Mid-life Aging in the Female Brain and Implications for Alzheimer Risk. <i>iScience</i> . doi: 10.1016/j.isci.2020.101829.
Hi-C		J Biol Chem.	2020	Sylvia Hewitt <i>et al.</i> Estrogen receptor α (ER α)-binding super enhancers drive key mediators that control uterine estrogen responses in mice. <i>JBC</i> . 2020 doi: 10.1074/jbc.RA120.013666
ChIP-Seq	HA-tagged AHR1	mBio.	2020	Sofia Ruben <i>et al.</i> Ahr1 and Tup1 Contribute to the Transcriptional Control of Virulence-Associated Genes in <i>Candida albicans</i> . <i>mBio</i> . 2020 doi: 10.1128/mBio.00206-20

Technique	Target	Journal	Year	Reference
ChIP-qPCR	BRG1, MITF	Mol Cancer Ther.	2020	Florencia Rago <i>et al.</i> The Discovery of SWI/SNF Chromatin Remodeling Activity as a Novel and Targetable Dependency in Uveal Melanoma. Mol. Cancer Ther. doi: 10.1158/1535-7163
ChIP-Seq	H3K27ac	Mol Neurobiol.	2020	Sunil Bhattarai <i>et al.</i> Modulation of Brain Pathology by Enhancer RNAs in Cerebral Ischemia. Mol. Neurobiol. Mol. Neurobiol. doi: 10.1007/s12035-020-02194-9
ChIP-Seq	IRF-2	Nat Cell Biol.	2020	Taku Sato <i>et al.</i> Regulated IFN signalling preserves the stemness of intestinal stem cells by restricting differentiation into secretory-cell lineages. Nat. Cell Biol. doi: 10.1038/s41556-020-0545-5.
ChIP-Seq		Nat Commun.	2020	Gulfem D. Guler <i>et al.</i> Detection of early stage pancreatic cancer using 5-hydroxymethylcytosine signatures in circulating cell free DNA. Nat. Commun. doi: 10.1038/s41467-020-18965-w.
ChIP-Seq	SMARCA2, SMARCA4	Nat Commun.	2020	Tharu M. Fernando <i>et al.</i> Functional characterization of SMARCA4 variants identified by targeted exome-sequencing of 131,668 cancer patients. Nat. Commun. doi: 10.1038/s41467-020-19402-8.
ChIP-Seq	H3K27ac, H3K-27me3, H3K3me1, H3K9me3, H3K-36me3	Nat Commun.	2020	Yonghe Wu <i>et al.</i> Glioblastoma epigenome profiling identifies SOX10 as a master regulator of molecular tumour subtype. Nat. Commun. doi: 10.1038/s41467-020-20225-w.
ChIP-Seq	N-MYC, BRD4, H3K9me3, RNA Pol II	Nat Commun.	2020	"Maged Zeineldin <i>et al.</i> MYCN amplification and ATRX mutations are incompatible in neuroblastoma. Nat. Commun. doi: 10.1038/s41467-020-14682-6"
ATAC-Seq		Nat Genet.	2020	Mark M. Pomerantz <i>et al.</i> Prostate cancer reactivates developmental epigenomic programs during metastatic progression. Nat. Genet. doi: 10.1038/s41588-020-0664-8.
ChIP-Seq	FLAG-tagged MRG15	Nat Metab.	2020	Yuda Wei <i>et al.</i> MRG15 orchestrates rhythmic epigenomic remodelling and controls hepatic lipid metabolism. Nat Metab. 2020 doi: 10.1038/s42255-020-0203-z
RNA-Seq		Nature	2020	Ana P. Gomes <i>et al.</i> Age-induced accumulation of methylmalonic acid promotes tumour progression. Nature. doi: 10.1038/s41586-020-2630-0.
ATAC-Seq		Nature	2020	"Biraj Mahato <i>et al.</i> Pharmacologic fibroblast reprogramming into photoreceptors restores vision. Nature. 2020 doi: 10.1038/s41586-020-2201-4"
ChIP-Seq	N-MYC, H3K27ac	Neoplasia	2020	Dongdong Chen <i>et al.</i> LIN28B promotes neuroblastoma metastasis and regulates PDZ binding kinase. Neoplasia. 2020 doi: 10.1016/j.neo.2020.04.001
ChIP-Seq	RNA Pol II	PLoS Genet.	2020	Katherine I. Farley-Barnes <i>et al.</i> Paired Box 9 (PAX9), the RNA polymerase II transcription factor, regulates human ribosome biogenesis and craniofacial development. PLoS Genet. doi: 10.1371/journal.pgen.1008967.
ATAC-Seq		Proc Natl Acad Sci USA	2020	Krutika Bhat <i>et al.</i> The dopamine receptor antagonist trifluoperazine prevents phenotype conversion and improves survival in mouse models of glioblastoma. PNAS. 2020 doi: 10.1073/pnas.1920154117

Technique	Target	Journal	Year	Reference
ATAC-Seq, ChIP-Seq	N-MYC, C-MYC, H3K4me1, H3K-4me3, H3K27ac, H3K27me3	Sci Data	2020	Kristen Upton <i>et al.</i> Epigenomic profiling of neuroblastoma cell lines. Sci Data. 2020 doi: 10.6084/m9.figshare.11993883
RIME	Estrogen Receptor- α	Sci Signal.	2020	Yin Li <i>et al.</i> A mutant form of ER α associated with estrogen insensitivity affects the coupling between ligand binding and coactivator recruitment. Sci. Signal. doi: 10.1126/scisignal.aaw4653.
ChIP-Seq	Progesterone Receptor	Sci Signal.	2020	Margeaux Wetendorf <i>et al.</i> Constitutive expression of progesterone receptor isoforms promotes the development of hormone-dependent ovarian neoplasms. Sci. Signal. doi: 10.1126/scisignal.aaz9646.
ChIP-Seq	H3K4me3, H3K-9me3, H3K27me3, H3R8me2a, H4K20me3	Theranostics	2020	Qian Zhang <i>et al.</i> Mdig promotes oncogenic gene expression through antagonizing repressive histone methylation markers. Theranostics 2020; 10(2):602-614.
ChIP-Seq	H3K27ac	Transl Oncol.	2020	Jie Cui <i>et al.</i> A zinc finger family protein, ZNF263, promotes hepatocellular carcinoma resistance to apoptosis via activation of ER stress-dependent autophagy. Transl. Oncol. 10.1016/j.tranon.2020.100851.
ChIP-Seq	H3K27me3, H3K4me3	Acta Neuropathol.	2019	Andre Silveira <i>et al.</i> H3.3 K27M depletion increases differentiation and extends latency of diffuse intrinsic pontine glioma growth in vivo. Acta Neuropathol. 137: 1021
ChIP-Seq	RNA Pol II	Agronomy	2019	Sonja Klemme <i>et al.</i> Selection of Salicylic Acid Tolerant Epilines in Brassica napus. Agronomy. 9: 92
ChIP-Seq	H3K4me1, H3K-4me3, H3K27ac, H3K27me3	BMC Biol.	2019	Lingzhao Fang <i>et al.</i> Functional annotation of the cattle genome through systematic discovery and characterization of chromatin states and butyrate-induced variations. BMC Biol. 17: 68
ATAC-Seq		Cancer Cell	2019	Ana P. Gomes <i>et al.</i> Dynamic incorporation of histone H3 variants into chromatin is essential for acquisition of aggressive traits and metastatic colonization. Cancer Cell. 36: 402
ChIP-Seq	H3K27me3	Cancer Cell	2019	Zulekha A. Qadeer <i>et al.</i> ATRX in-frame fusion neuroblastoma is sensitive to EZH2 inhibition via modulation of neuronal gene signatures. Cancer Cell. 36: 512
ChIP-Seq	H3K27me3, H3K4me1, H3K-4me3, H3K9me3, H3K36me3, EZH2, SMARCB1, SUZ12, SMARCA4, REST	Cancer Cell	2019	Serap Erkek <i>et al.</i> Comprehensive Analysis of Chromatin States in Atypical Teratoid/Rhabdoid Tumors Identified Diverging Roles for SWI/SNF and Polycomb in Gene Regulation. Cancer Cell. 35: 95
ChIP-Seq	H3K27ac	Cancer Res.	2019	Joesph Castillo <i>et al.</i> CBP/p300 drives the differentiation of Regulatory T cells through transcriptional and non-transcriptional mechanisms. Cancer Res. 79: 3916
ChIP-Seq	Estrogen Receptor	Cell	2019	"Jane Guan <i>et al.</i> Therapeutic ligands antagonize estrogen receptor function by impairing its mobility. Cell. 178: 15"
ChIP-Seq	H4 panAc	Cell Metab.	2019	Ramon C. Sun <i>et al.</i> Nuclear glycogenolysis modulates histone acetylation in human non-small cell lung cancers. Cell Metab. 30: 1

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K4me1, H3K-4me3, H3K27ac	Cell Metab.	2019	"Liming Du <i>et al.</i> IGF-2 Preprograms Maturing Macrophages to Acquire Oxidative Phosphorylation-Dependent Antiinflammatory Properties. <i>Cell Metab.</i> 29: 1"
Mod Spec		Cell Rep.	2019	"Banushree Kumar <i>et al.</i> Quantitative Multiplexed ChIP Reveals Global Alterations that Shape Promoter Bivalency in Ground State Embryonic Stem Cells. <i>Cell Reports.</i> 28: 3274–3284."
ChIP-Seq, ChIP-qPCR	Androgen Receptor, Estrogen Receptor	Cell Rep.	2019	"Karyn Schmidt <i>et al.</i> The lncRNA SLNCR Recruits the Androgen Receptor to EGR1-Bound Genes in Melanoma and Inhibits Expression of Tumor Suppressor p21. <i>Cell Reports.</i> 27: 2493"
ChIP-Seq, RNA-Seq Data Analysis	ASH2L, H3K4me3	Cell Rep.	2019	"Liang Li <i>et al.</i> The COMPASS Family Protein ASH2L Mediates Corticogenesis via Transcriptional Regulation of Wnt Signaling. <i>Cell Reports.</i> 28: 698"
ChIP-Seq, ChIP-qPCR	CBP, p300, H3K27ac	Cell Rep.	2019	Denise de Almeida Nagata <i>et al.</i> Regulation of Tumor-Associated Myeloid Cell Activity by CBP/EP300 Bromodomain Modulation of H3K27 Acetylation. <i>Cell Reports.</i> 27:269
ChIP-Seq	RNA Pol II	Cell Stem Cell	2019	Kimberley N. Babos <i>et al.</i> Mitigating antagonism between transcription and proliferation allows near-deterministic cellular reprogramming. <i>Cell Stem Cell.</i> 25: 486
ChIP-qPCR	BRD4, RelA	Clin Epigenetics	2019	"Laura Tsujikawa <i>et al.</i> Apabetalone (RVX-208) reduces vascular inflammation in vitro and in CVD patients by a BET-dependent epigenetic mechanism. <i>Clin Epigenet.</i> 11: 102"
RNA-Seq		EMBO J.	2019	"Kathryn A. Jacobs <i>et al.</i> Paracaspase MALT1 regulates glioma cell survival by controlling endo-lysosome homeostasis. <i>EMBO J.</i> doi: 10.15252/embj.2019102030"
MeDIP-Seq	5-Methylcytosine	Epigenomes	2019	Liliana Ferreira <i>et al.</i> Uncovering Differentially Methylated Regions (DMRs) in a Salt-Tolerant Rice Variety under Stress: One Step towards New Regulatory Regions for Enhanced Salt Tolerance. <i>Epigenomes.</i> 3:4
ATAC-Seq		Haematologica	2019	James Ropa <i>et al.</i> SETDB1 mediated histone H3 lysine 9 methylation suppresses MLL-fusion target expression and leukemic transformation. <i>Haematologica.</i> doi: 10.3324/haematol.2019.223883
ChIP-Seq, Spike-in	H3K27me3	Haematologica	2019	"Shao Xie <i>et al.</i> EZH2 inhibitors abrogate upregulation of trimethylation of H3K27 by CDK9 inhibitors and potentiate its activity against diffuse large B-cell lymphoma. <i>Haematologica.</i> doi:10.3324/haematol.2019.222935"
ChIP-qPCR	PPAR γ	Int J Obes (Lond).	2019	"Laura Butruille <i>et al.</i> Maternal high-fat diet during suckling programs visceral adiposity and epigenetic regulation of adipose tissue stearoyl-CoA desaturase-1 in offspring. <i>Int J Obes (London)</i> doi: 10.1038/s41366-018-0310-z"

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K27me3, H3K27ac, H3K4me3	Invest Ophthalmol Vis Sci.	2019	"Christina F. Tingle <i>et al.</i> Paradoxical Changes Underscore Epigenetic Reprogramming During Adult Zebrafish Extraocular Muscle Regeneration. <i>Invest Ophthalmol Vis Sci.</i> 2019;60:4991–4999"
ChIP-Seq, RNA-Seq	Glucocorticoid Receptor	J Am Heart Assoc.	2019	Elena Severinova <i>et al.</i> Glucocorticoid Receptor Binding and Transcriptome Signature in Cardiomyocytes. <i>J Am Heart Assoc.</i> 8: e011484.
ChIP-Seq	H4K20me1	J Autoimmun.	2019	Keqi Fan <i>et al.</i> CRL4DCAF2 is required for mature T-cell expansion via Aurora B-regulated proteasome activity. <i>J of Autoimmun.</i> 96:74
ChIP-Seq	SMC1a, H3K4me1, H3K4me3, H3K27ac	J Biol Chem.	2019	Sylvia C. Hewitt <i>et al.</i> A distal super enhancer mediates estrogen-dependent mouse uterine-specific gene transcription of Igf1 (Insulin-like growth factor 1). <i>J Biol Chem.</i> 294: 9746
ChIP-Seq, ChIP-qPCR	Progesterone Receptor	J Clin Endocrinol Metab.	2019	Ru-pin Alicia Chi <i>et al.</i> Human endometrial transcriptome and progesterone receptor cistrome reveal important pathways and epithelial regulators. <i>J Clin Endocrinol Metab.</i> doi: 10.1210/clinem/dgz117
ChIP-Seq	BRG1	J Clin Invest.	2019	Yufeng Ding <i>et al.</i> Chromatin remodeling ATPase BRG1 and PTEN are synthetic lethal in prostate cancer. <i>J Clin Invest.</i> 129: 759
ChIP-Seq	H3K9me3	J Exp Med.	2019	"Jun Hyung Sin <i>et al.</i> The epigenetic regulator ATF7ip inhibits IL2 expression, regulating Th17 responses. <i>J. Exp Med.</i> 216: 2024"
ChIP-Seq	Androgen Receptor	J Mol Med (Berl).	2019	Kinza Younas <i>et al.</i> Delayed endometrial decidualisation in polycystic ovary syndrome; the role of AR-MAGEA11. <i>J Mol Med</i> 2019 doi: 10.1007/s00109-019-01809-6
ChIP-Seq	RNA Pol II, H3K27ac	J Neurooncol.	2019	Isabel Tegeder <i>et al.</i> Functional relevance of genes predicted to be affected by epigenetic alterations in atypical teratoid/rhabdoid tumors. <i>J. Neurooncol.</i> 141: 43
ChIP-Seq	EGR1, H3K27ac	J Steroid Biochem Mol Biol.	2019	"Maria Szwarc <i>et al.</i> Early growth response 1 transcriptionally primes the human endometrial stromal cell for decidualization. <i>J Steroid Biochem.</i> 189: 283"
ChIP-Seq, RNA-Seq	H3K4me3, H3K27me3	JCI Insight	2019	Taiyi Huo <i>et al.</i> Induction of a cell-restricted Gc in dedifferentiating β cells contributes to stress-induced β -cell dysfunction. <i>J Clin Invest.</i> 5: 128351
ChIP-Seq, ChIP-qPCR	Androgen Receptor	Mol Cell Endocrinol.	2019	Claire Nash <i>et al.</i> Genome-wide analysis of AR binding and comparison with transcript expression in primary human fetal prostate fibroblasts and cancer associated fibroblasts. <i>Mol Cell Endocrinol.</i> 471: 1
ChIP-Seq	YAP1, NFIA	Nat Commun.	2019	Kristian W. Pajtler <i>et al.</i> YAP1 subgroup supratentorial ependymoma requires TEAD and nuclear factor I-mediated transcriptional programmes for tumorigenesis. <i>Nat Commun.</i> 10: 3914
ChIP-Seq	BRG1, H3K9Ac	Nat Commun.	2019	Min Liu <i>et al.</i> BRG1 attenuates colonic inflammation and tumorigenesis through autophagy-dependent oxidative stress sequestration. <i>Nat Commun.</i> 10: 4614
ChIP-Seq	BRD4	Nat Commun.	2019	"Nicolas Mercado <i>et al.</i> IRF2 is a master regulator of human keratinocyte stem cell fate. <i>Nat Commun.</i> 10: 4676"

Technique	Target	Journal	Year	Reference
ChIP-Seq	RUNX1	Nat Commun.	2019	Barbara Nicol <i>et al.</i> RUNX1 maintains the identity of the fetal ovary through an interplay with FOXL2. Nat. Commun. 10: 5116
ChIP-Seq	H3K27ac	Nat Commun.	2019	Jayaram Vijayakrishnan <i>et al.</i> Identification of four novel associations for B-cell acute lymphoblastic leukaemia risk. Nat Commun. 10: 5348
ChIP-Seq	H3K27ac	Nat Commun.	2019	"Liqing Tian <i>et al.</i> Long-read sequencing unveils IGH-DUX4 translocation into the silenced IGH allele in B-cell acute lymphoblastic leukemia. Nat. Commun. 10: 2789"
ChIP-Seq	SALL3, DNMT3B	Nat Commun.	2019	"Takuya Kuroda <i>et al.</i> SALL3 expression balance underlies lineage biases in human induced pluripotent stem cell differentiation. Nat Commun. 10: 2175"
ChIP-qPCR	SRF, CEBP	Nat Commun.	2019	"S. Sarcar <i>et al.</i> Next-generation muscle-directed gene therapy by in silico vector design. Nat Commun. 10: 492"
ChIP-Seq	H3K27ac, H3K4me3, CTCF, NR4A3	Nat Commun.	2019	Florian Haller <i>et al.</i> Enhancer hijacking activates oncogenic transcription factor NR4A3 in acinic cell carcinomas of the salivary glands. Nat Commun. 10: 336
ChIP-Seq	GFI1, LSD1	Nat Commun.	2019	Catherine Lee <i>et al.</i> Lsd1 as a therapeutic target in Gfi1-activated medulloblastoma. Nat Commun. 10: 332
Mod Spec		Nat Genet.	2019	"Sara Martire <i>et al.</i> Phosphorylation of histone H3.3 at serine 31 promotes p300 activity and enhancer acetylation. Nat. Genet. 51: 941"
ChIP-qPCR	NRF2	Nat Immunol.	2019	Anne-Valerie Burgener <i>et al.</i> SDHA gain-of-function engages inflammatory mitochondrial retrograde signaling via KEAP1–Nrf2. Nat Immunol. 20: 1311
ChIP-qPCR	ChREBP	Nat Metab.	2019	"Pauline Morigny <i>et al.</i> Interaction between hormone-sensitive lipase and ChREBP in fat cells controls insulin sensitivity. Nat. Metab. 1: 133"
ATAC-Seq, ChIP-Seq	H3K4me3, H3K27ac, H3K27me3	Nature	2019	Nina Oberbeck <i>et al.</i> The RIPK4–IRF6 signalling axis safeguards epidermal differentiation and barrier function. Nature. 574: 249
ChIP-Seq	BRD4, H3K9me3	Neuron	2019	Jackie L. Norrie <i>et al.</i> Nucleome dynamics during retinal development. Neuron. 104:1
RIME	PPARb/d	Nucleic Acids Res.	2019	Nathalie Legrand <i>et al.</i> PPARb/d recruits NCOR and regulates transcription reinitiation of ANGPTL4. Nucleic Acids Res. 47: 9573
Next-Gen Bisulfite-Seq		Plant Biotechnol J.	2019	"Aurine Verkest <i>et al.</i> Impact of differential DNA methylation on transgene expression in cotton (<i>Gossypium hirsutum</i> L.) events generated by targeted sequence insertion. Plant Biotechnol. J. 17: 1236"
ChIP-Seq	Progesterone Receptor, H3K27ac	Sci Rep.	2019	Diem T. Dinh <i>et al.</i> Tissue-specific progesterone receptor-chromatin binding and the regulation of progesterone-dependent gene expression. Sci Rep. 19: 9
ChIP-Seq	LSD1, H3K4me2, H3K27ac	Sci Signal.	2019	Arnaud Augert <i>et al.</i> Targeting NOTCH activation in small cell lung cancer through LSD1 inhibition. Sci.Signal. 12: eaau2922

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K27me3	Acta Neuropathol Commun.	2018	David Castel <i>et al.</i> Transcriptomic and epigenetic profiling of 'diffuse midline gliomas, H3 K27M-mutant' discriminate two subgroups based on the type of histone H3 mutated and not supratentorial or infratentorial location. Acta Neuropathol Commun. 6: 17
Mod Spec		Acta Neuropathol.	2018	Leah Katz <i>et al.</i> Loss of histone H3K27me3 identifies a subset of meningiomas with increased risk of recurrence Acta Neuropathol. 135:955
ChIP-Seq, Next-Gen Bisulfite-Seq	H3K4me3, RNA Pol II	Agronomy	2018	"Martin Schmidt <i>et al.</i> Methylome and epialleles in rice epilines selected for energy use efficiency. Agronomy. 8: 163"
ChIP-Seq	GATA2	Am J Hum Genet.	2018	Katelyn M. Mika <i>et al.</i> An ancient fecundability-associated polymorphism creates a GATA2 binding site in a distal enhancer of HLA-F. Am J Hum Genet. 103: 509
ChIP-qPCR	Glucocorticoid Receptor, TIF1	Arthritis Rheumatol.	2018	Yanhua Hu <i>et al.</i> Development of a molecular signature to monitor pharmacodynamic response mediated by in vivo administration of glucocorticoids. Arthritis Rheumatol. 70: 1331
ChIP-Seq	BRD4	Atherosclerosis	2018	Dean Gilham <i>et al.</i> Apabetalone downregulates factors and pathways associated with vascular calcification Atherosclerosis 280:75
ChIP-Seq	RNA Pol II, H2A.Z, CDK9, ANP32e	Biochim Biophys Acta Gene Regul Mech.	2018	Hyewon Shin <i>et al.</i> Transcriptional regulation mediated by H2A.Z via ANP32e-dependent inhibition of protein phosphatase 2A. Biochim Biophys Acta. 1861: 481
ChIP-Seq	BRD4	Blood	2018	Rebecca Kohnken <i>et al.</i> Diminished microRNA-29b level is associated with BRD4 mediated activation of oncogenes in cutaneous T-cell lymphoma. Blood. 131: 771
ChIP-Seq	BRD4, RNA Pol II	Cancer Cell	2018	Elizabeth Stewart <i>et al.</i> Identification of therapeutic targets in rhabdomyosarcoma through integrated genomic, epigenomic, and proteomic analyses. Cancer Cell. 34: 1
ChIP-Seq	HDAC2, H3K27ac, H3K27me3	Cancer Immunol Immunother.	2018	David Briere <i>et al.</i> The class I/IV HDAC inhibitor mocetinostat increases tumor antigen presentation, decreases immune suppressive cell types and augments checkpoint inhibitor therapy. Cancer Immunol Immunother. 67: 381
ChIP-Seq	MTA1	Cancer Med.	2018	Nasir A. Butt <i>et al.</i> Targeting MTA1/HIF-1 α signaling by pterostilbene in combination with histone deacetylase inhibitor attenuates prostate cancer progression. Cancer Med. 6: 2673
RNA-Seq		Cell Commun Signal.	2018	Keri Callegari <i>et al.</i> Pharmacological inhibition of LSD1 activity blocks REST-dependent medulloblastoma cell migration. Cell Commun Signal. 16: 60
ChIP-Seq	H3K27ac	Cell	2018	Xun Huang <i>et al.</i> Targeting epigenetic crosstalk as a therapeutic strategy for EZH2-aberrant solid tumors. Cell. 175: 1
ChIP-Seq	H3K36me2	Cell Rep.	2018	"Jingjing Chen <i>et al.</i> Methyltransferase Nsd2 Ensures Germinal Center Selection by Promoting Adhesive Interactions between B Cells and Follicular Dendritic Cells Cell Reports. 25: 3393"

Technique	Target	Journal	Year	Reference
ChIP-Seq	RNA Pol II, BRD4, H3K9me3	Cell Rep.	2018	Lu Wang <i>et al.</i> Retinal Cell Type DNA Methylation and Histone Modifications Predict Reprogramming Efficiency and Retinogenesis in 3D Organoid Cultures. Cell Reports. 22: 2601
ATAC-Seq		Cell Rep.	2018	Xiaolong Zhang <i>et al.</i> OX40 Costimulation inhibits Foxp3 expression and Treg induction via BATF3-dependent and independent mechanisms. Cell Reports. 24: 607
RIME	TRPS1	Cell Rep.	2018	Robert M. Witwicki <i>et al.</i> TRPS1 Is a lineage-specific transcriptional dependency in breast cancer. Cell Reports. 25: 1255
ChIP-Seq, RIME	p300, CBP, CDK9, BRD4, H3K27ac, H3K4me1, H3K4me3, H3K18Ac	Cell Rep.	2018	Ryan Raisner <i>et al.</i> Enhancer activity requires CBP/P300 bromodomain-dependent histone H3K27 acetylation. Cell Reports. 24: 1722
ChIP-Seq	ZFP24	Cell Rep.	2018	Benayahu Elbaz <i>et al.</i> Phosphorylation state of ZFP24 controls oligodendrocyte differentiation. Cell Reports. 23: 2254
ChIP-Seq, Histone PTM Quantitation	H3K9me3, H3K27me3	Cell Rep.	2018	Jessica Camacho <i>et al.</i> The memory of environmental chemical exposure in <i>C. elegans</i> is dependent on the jumonji demethylases <i>jmjd-2</i> and <i>jmjd-3/utx-1</i> . Cell Reports. 23: 2392
ChIP-Seq	H3K27ac	Clin Cancer Res.	2018	Shori Saito <i>et al.</i> Eradication of central nervous system leukemia of T-cell origin with a brain-permeable LSD1 inhibitor Clin Cancer Res. 25:1601
ChIP-Seq	H3K27me3	G3 (Bethesda)	2018	James Ferguson <i>et al.</i> PRC2 is dispensable in vivo for b-catenin-mediated repression of chondrogenesis in the mouse embryonic cranial mesenchyme. G3. 8: 491
ChIP-qPCR	p65, RelA	Hum Mol Genet.	2018	Jeffrey R. Gehlhausen <i>et al.</i> A proteasome-resistant fragment of NIK mediates oncogenic NF-κB signaling in schwannomas. Hum Mol Gen. 28:572
ChIP-Seq	FOXL2	Hum Mol Genet.	2018	"Barbara Nicol <i>et al.</i> Genome-wide identification of FOXL2 binding and characterization of FOXL2 feminizing action in the fetal gonads. Hum Mol Genet. Advance online publication doi: 10.1093/hmg/ddy312"
ChIP-Seq	NRF2	J Biol Chem.	2018	"Junsheng Fu <i>et al.</i> Hyperactivity of the transcription factor Nrf2 causes metabolic reprogramming in mouse esophagus. J. Biol. Chem. 294: 327."
ChIP-Seq, ChIP-qPCR	TRRAP	J Cell Biol.	2018	Zhao Wang <i>et al.</i> TRRAP is a central regulator of human multiciliated cell formation. J Cell Biol. 217: 1941
ChIP-Seq, RNA-Seq	H3K27ac, MED1, RNA Pol II	J Clin Invest.	2018	Maria Florencia Martinez <i>et al.</i> Super-Enhancers maintain renin-expressing cell identity and memory to preserve multi-systemhomeostasis. J. Clin Invest. Advance online publication doi: 10.1172/JCI121361
ChIP-Seq	EZH2	J Exp Med.	2018	Xingli Zhang <i>et al.</i> Macrophage/microglial Ezh2 facilitates autoimmune inflammation through inhibition of Socs3. J Exp Med. 215: 1365
ChIP-Seq	ELL2	J Immunol.	2018	Ashley M. Nelson <i>et al.</i> RNA splicing in the transition from B cells to antibody-secreting cells: The influences of ELL2, small nuclear RNA, and endoplasmic reticulum stress. J Immunol. doi: 10.4049/jimmunol.1800557

Technique	Target	Journal	Year	Reference
ChIP-Seq	SIRT1	J Neuroimmune Pharmacol.	2018	Nikki Bortell <i>et al.</i> Sirtuin 1-chromatin-binding dynamics points to a common mechanism regulating inflammatory targets in SIV infection and in the aging brain. J Neuroimmune Pharmacol. 13: 163
ChIP-Seq	BRD4, H3K27ac	JCI Insight	2018	Anisley Valenciaga <i>et al.</i> Transcriptional targeting of oncogene addiction in medullary thyroid cancer. JCI Insight. 3: e122225
ChIP-Seq	FOXG1	Mol Neurobiol.	2018	Stefan Weise <i>et al.</i> FOXG1 Regulates PRKAR2B Transcriptionally and Posttranscriptionally via miR200 in the Adult Hippocampus. Mol. Neurobiol. S12035-018-1444.
ChIP-Seq	MTA1	Mol Oncol.	2018	Avinash Kumar <i>et al.</i> MTA drives malignant progression and bone metastasis in prostate cancer. Mol. Oncol. 12:1596
ChIP-Seq	NRF2	Mol Pharmacol.	2018	Rance Nault <i>et al.</i> Comparison of hepatic NRF2 and AHR binding in 2,3,7,8- tetrachlorodibenzo-p-dioxin (TCDD) treated mice demonstrates NRF2- independent PKM2 induction. Mol Pharmacol. 94: 876
ChIP-Seq	SOX17	Nat Commun.	2018	Xiaoqiu Wang <i>et al.</i> SOX17 regulates uterine epithelial–stromal cross- talk acting via a distal enhancer upstream of Ihh. Nat Commun. 9: 4421
ChIP-Seq, RNA-Seq	NRF2, RNA Pol II	Nat Commun.	2018	David Olgarnier <i>et al.</i> Nrf2 negatively regulates STING indicating a link between antiviral sensing and metabolic reprogramming. Nat Commun. 9: 3506
ChIP-Seq	TET1, DNMT1	Nat Genet.	2018	Nipun Verma <i>et al.</i> TET proteins safeguard bivalent promoters from de novo methylation in human embryonic stem cells. Nat Genet. 50: 83
RNA-Seq		Pain	2018	Pradipta Ray <i>et al.</i> Comparative transcriptome profiling of the human and mouse dorsal root ganglia: an RNA-Seq based resource for pain and sensory neuroscience research Pain. 159:1325
ATAC-Seq, ChIP-Seq	H3K36me3, H3K27ac	PLoS One	2018	"Ryan K. Shultzaberger <i>et al.</i> Agnostic detection of genomic alterations by holistic DNA structural interrogation. PLoS ONE 13 (11):e02080554"
ChIP-Seq	BHLHE40	PLoS One	2018	Kelly A. Hamilton <i>et al.</i> Mice lacking the transcriptional regulator Bhlhe40 have enhanced neuronal excitability and impaired synaptic plasticity in the hippocampus. PLoS One. 13: e0196223
ChIP-Seq	RNA Pol II	PLoS One	2018	Liana Basova <i>et al.</i> Dopamine and its receptors play a role in the modulation of CCR5 expression in innate immune cells following exposure to Methamphetamine: Implications to HIV infection. PLoS One. 13: e0199861
ChIP-Seq	H3K4me3, H3K9me3, H3K27me3	Proc Biol Sci.	2018	Theresa K. Kelly <i>et al.</i> Epigenetic regulation of transcriptional plasticity associated with developmental song learning. Proc Biol Sci. 285: 20180160
ChIP-Seq	CREB, SWc	Sci Rep.	2018	Jason L. Larabee <i>et al.</i> Unique, Intersecting, and Overlapping Roles of C/EBP β and CREB in Cells of the Innate Immune System Sci Rep. 8:16931
MeDIP-Seq	5-Methylcytosine	Sci Rep.	2018	Jose P. silva <i>et al.</i> Analysis of diet-induced differential methylation, expression, and interactions of lncRNA and protein- coding genes in mouse liver. Sci Rep. 8: 11537

Technique	Target	Journal	Year	Reference
ChIP-Seq	STAT3, STAT5, ROCK2	Sci Rep.	2018	Wei Chen <i>et al.</i> ROCK2, but not ROCK1 interacts with phosphorylated STAT3 and co-occupies TH17/TFH gene promoters in TH17-activated human T cells. <i>Sci Rep.</i> 8: 16636
ChIP-Seq	HA-tagged JP2NT, TBP, MEF2C	Science	2018	Ang Guo <i>et al.</i> E-C coupling structural protein junctophilin-2 encodes a stress-adaptive transcription regulator. <i>Science.</i> 362:6421.
ChIP-Seq	G34R	Acta Neuropathol Commun.	2017	Farhana Haque <i>et al.</i> Evaluation of a novel antibody to define histone 3.3 G34R mutant brain tumours. <i>Acta Neuropathol Commun.</i> 5: 45
ChIP-Seq	Progesterone Receptor	Biol Reprod.	2017	Margeaux Wetendorf <i>et al.</i> Decreased epithelial progesterone receptor A at the window of receptivity is required for preparation of the endometrium for embryo attachment. <i>Biol Reprod.</i> 96: 313.
ATAC-Seq, ChIP-Seq	H3K4me3, H3K9me3	Cancer Cell	2017	Gulfem Dilek Guler <i>et al.</i> Repression of stress-induced LINE-1 expression protects cancer cell subpopulations from lethal drug exposure. <i>Cancer Cell.</i> 32: 1
ChIP-Seq, ChIP-qPCR	EZH2, H3K27me3, H3K27ac	Cancer Cell	2017	Eric E. Gardner <i>et al.</i> Chemosensitive relapse in small cell lung cancer proceeds through an EZH2-SLFN11 axis. <i>Cancer Cell.</i> 31: 286
RIME	Estrogen Receptor, Progesterone Receptor	Cancer Res.	2017	"Jessica Finlay-Schultz <i>et al.</i> Breast cancer suppression by progesterone receptors is mediated by their modulation of estrogen receptors and RNA polymerase III. <i>Cancer Res.</i> 77: 4934"
ChIP-Seq, Spike-in	p300, H3K27ac, Androgen Receptor	Cancer Res.	2017	Lingyan Jin <i>et al.</i> Therapeutic targeting of the CBP/p300 bromodomain blocks the growth of castration-resistant prostate cancer. <i>Cancer Res.</i> 77: 5564
ChIP-Seq	SLY1	Cell Death Differ.	2017	Charlotte Moretti <i>et al.</i> SLY regulates genes involved in chromatin remodeling and interacts with TBL1XR1 during sperm differentiation. <i>Cell Death Differ.</i> 24: 1029
ChIP-qPCR	H3K9Ac, RNA Pol II, ChREBP, PPAR α	Cell Rep.	2017	Alison Iroz <i>et al.</i> A specific ChREBP and PPAR α cross-talk is required for the glucose-mediated FGF21 response. <i>Cell Reports.</i> 21: 403
ChIP-Seq	TRIM28	Cell Rep.	2017	Per Ludvik Brattas <i>et al.</i> TRIM28 controls a gene regulatory network based on endogenous retroviruses in human neural progenitor cells. <i>Cell Reports</i> 18: 1
MeDIP-Seq	5-Methylcytosine	Cereb Cortex	2017	Daniela Grassi <i>et al.</i> Neuronal activity, TGF β -Signaling and unpredictable chronic stress modulate transcription of Gadd45 family members and DNA methylation in the hippocampus. <i>Cereb Cortex.</i> 27: 4166
ChIP-Seq	LMX1b	Development	2017	Endika Haro <i>et al.</i> Lmx1b-targeted cis-regulatory modules involved in limb dorsalization. <i>Dev.</i> 444: 2009
ChIP-Seq, ChIP-qPCR	Estrogen Receptor	Endocrinology	2017	Sylvia C. Hewitt <i>et al.</i> Role of ER α in mediating female uterine transcriptional Responses to IGF1. <i>Endocrinology.</i> 158: 2427
ChIP-Seq	Estrogen Receptor	Endocrinology	2017	Laurel A. Coons <i>et al.</i> DNA sequence constraints define functionally active steroid nuclear receptor binding sites in chromatin. <i>Endocrinology.</i> 158: 3212
ChIP-Seq	H3K27ac, H3K4me1	Epigenomics	2017	John P. Thomson <i>et al.</i> Defining baseline epigenetic landscapes in the rat liver. <i>Epigenomics.</i> 9: 1503

Technique	Target	Journal	Year	Reference
MeDIP-Seq		Genes Brain Behav.	2017	M. Fonteneau <i>et al.</i> Inhibition of DNA methyltransferases regulates cocaine self-administration by rats: a genome-wide DNA methylation study. <i>Genes Brain Behav.</i> 16:313
ChIP-qPCR	IRF5, NFkB (p65)	J Biol Chem.	2017	Leah Cushing <i>et al.</i> IRAK4 kinase activity controls Toll-like receptor induced inflammation through the transcription factor IRF5 in primary human monocytes. <i>J Biol Chem.</i> 292: 18689
ChIP-Seq	RNA Pol II	J Biol Chem.	2017	Leah A. Gates <i>et al.</i> Acetylation on histone H3 lysine 9 mediates a switch from transcription initiation to elongation. <i>J Biol Chem.</i> 292: 14456
ChIP-Seq	H3K36me3	J Clin Invest.	2017	Huairui Yuan <i>et al.</i> Histone methyltransferase SETD2 modulates alternative splicing to inhibit intestinal tumorigenesis. <i>J Clin Invest.</i> 127: 3375
ChIP-Seq	H3K36me2	J Clin Invest.	2017	Ni Li <i>et al.</i> AKT-mediated stabilization of histone methyltransferase WHSC1 promotes prostate cancer metastasis. <i>J Clin Invest.</i> 127: 1284
ChIP-Seq	SALL4	J Hematol Oncol.	2017	Lina Yang <i>et al.</i> The stem cell factor SALL4 is an essential transcriptional regulator in mixed lineage leukemia-rearranged leukemogenesis. <i>J Hematol Oncol.</i> 10: 159
Next-Gen Bisulfite-Seq, ChIP-Seq	KLF4	Leukemia	2017	"Y Shen <i>et al.</i> Inactivation of KLF4 promotes T-cell acute lymphoblastic leukemia and activates the MAP2K7 pathway. <i>Leukemia.</i> 31: 1314"
ChIP-qPCR	EZH2	Mech Ageing Dev.	2017	Khyobeni Mozhui <i>et al.</i> Conserved effect of aging on DNA methylation and association with EZH2 polycomb protein in mice and humans. <i>Mech Ageing Dev.</i> 162: 27
ChIP-Seq	H3K79me2	Mol Cancer Ther.	2017	Carly T. Campbell <i>et al.</i> Mechanisms of Pinometostat (EPZ-5676) treatment-emergent resistance in MLL-rearranged leukemia. <i>Mol Cancer Ther.</i> 16: 1669
ChIP-Seq	EZH2, H3K27me3, H3K4me3	Mol Cancer Ther.	2017	Dorothy Brach <i>et al.</i> EZH2 inhibition by tazemetostat results in altered dependency on B-cell activation signaling in DLBCL. <i>Mol Cancer Ther.</i> 16:2586
ChIP-Seq	Androgen Receptor	Mol Cell Endocrinol.	2017	Claire Nash <i>et al.</i> Genome-wide analysis of AR binding and comparison with transcript expression in primary human fetal prostate fibroblasts and cancer associated fibroblasts. <i>Mol Cell Endocrinol.</i> Advance online publication doi: 10.1016/j.mce.2017.05.006
ChIP-Seq	BACH2	Nat Commun.	2017	Nicolas Hipp <i>et al.</i> IL-2 imprints human naive B cell fate towards plasma cell through ERK/ELK1-mediated BACH2 repression. <i>Nat Commun.</i> 8: 1443
MeDIP-Seq	5-Methylcytosine	Nature	2017	Rama S. Akondy <i>et al.</i> Origin and differentiation of human memory CD8 T cells after vaccination. <i>Nature.</i> 552: 362
ChIP-Seq	10 different Histone mods, CTCF, RNA Pol II, BRD4	Nature	2017	Elizabeth Stewart <i>et al.</i> Orthotopic patient-derived xenografts of paediatric solid tumours. <i>Nature.</i> 549: 96
ChIP-Seq	BRD4, H3K9me3, RNA Pol II	Neuron	2017	Issam Aldiri <i>et al.</i> The dynamic epigenetic landscape of the retina during development, reprogramming and tumorigenesis. <i>Neuron.</i> 94: 550
ChIP-Seq	SRC-2	PLoS Genet.	2017	Shruthy Suresh <i>et al.</i> SRC-2-mediated coactivation of anti-tumorigenic target genes suppresses MYC-induced liver cancer. <i>PLoS Genet.</i> 13: e1006650

Technique	Target	Journal	Year	Reference
ChIP-Seq	EZH2	Proc Natl Acad Sci USA	2017	Yongfeng Liu <i>et al.</i> Epithelial EZH2 serves as an epigenetic determinant in experimental colitis by inhibiting TNF α -mediated inflammation and apoptosis. Proc Natl Acad Sci. 114: E3796
ChIP-Seq	ETV5	Proc Natl Acad Sci USA	2017	Zhen Zhang <i>et al.</i> Transcription factor Etv5 is essential for the maintenance of alveolar type II cells. Proc Natl Acad Sci. 114: E3903
ChIP-Seq Data Analysis	H3K27me3	Sci Rep.	2017	Brid O'Leary <i>et al.</i> Long non-coding RNA PARTICLE bridges histone and DNA methylation. Sci Rep. 7: 1790
ChIP-Seq	CBX3/HP1	Sci Rep.	2017	Michael Sun <i>et al.</i> Cbx3/HP1 γ deficiency confers enhanced tumor-killing capacity on CD8 $^+$ T cells. Sci Rep. 7: 42888
ChIP-Seq	H3K27ac	Sci Rep.	2017	Yao Shen <i>et al.</i> Epigenetic and genetic dissections of UV-induced global gene dysregulation in skin cells through multi-omics analyses. Sci Rep. 7: 42646
ChIP-qPCR	SIRT1	Sci Rep.	2017	Jung-Yoon Yoo <i>et al.</i> KRAS activation and over-expression of SIRT1/BCL6 contributes to the pathogenesis of endometriosis and progesterone resistance. Sci Rep. 7: 6765
Low Cell ChIP-Seq	H3K4me3, H3K27ac, H3K- 27me3	Biol Reprod.	2016	Kazadi Mutoji <i>et al.</i> TSPAN8 expression distinguishes spermatogonial stem cells in the prepubertal mouse testis. Biol Reprod. 95: 17.
ChIP-Seq	OTX2, MITF, BRD4, H3K27ac	Cancer Cell	2016	Pascal D. Johann <i>et al.</i> Atypical teratoid/rhabdoid tumors are comprised of three epigenetic subgroups with distinct enhancer landscapes. Cancer Cell. 29: 379
ChIP-Seq	HDAC1, H3K27ac	Cancer Discov.	2016	Anjali Mishra <i>et al.</i> Mechanism, consequences, and therapeutic targeting of abnormal IL15 signaling in cutaneous T-cell lymphoma. Cancer Discov. 6: 986
ChIP-Seq	H3K27me3	Cancer Res.	2016	John P. Thomson <i>et al.</i> Loss of Tet1 associated 5-hydroxymethylcytosine is concomitant with aberrant promoter hypermethylation in liver cancer. Cancer Res. 76: 3097
ChIP-Seq	GATA2, Progesterone Receptor	Cell Rep.	2016	Cory A. Rubel <i>et al.</i> A Gata2-dependent transcription network regulates uterine progesterone responsiveness and endometrial function. Cell Reports. 17: 1414
ChIP-qPCR	Androgen Receptor	Cell Rep.	2016	Karyn Schmidt <i>et al.</i> The lncRNA SLNCR1 mediates melanoma invasion through a conserved SRA1-like region. Cell Rep. 15: 2025
ChIP-Seq, Spike-in	KDM5C	Cell Rep.	2016	Shigeki Iwase <i>et al.</i> A mouse model of X-linked intellectual disability associated with impaired removal of histone methylation. Cell Reports. 14: 1
ChIP Libraries, ChIP-Seq, ChIP-Seq Data Analysis	H3K27me3	Dev Dyn.	2016	Oyvind Dahle <i>et al.</i> Inhibiting smad2/3 signaling in pluripotent mouse embryonic stem cells enhances endoderm formation by increasing transcriptional priming of lineage-specifying target genes. Dev Dyn. 245: 807
ChIP-Seq	PLZF, SALL4	Development	2016	Dawn L. Lovelace <i>et al.</i> The regulatory repertoire of PLZF and SALL4 in undifferentiated spermatogonia. Development. 143:1893
ChIP-Seq, ChIP-qPCR	WIZ	eLife	2016	Luke Isbel <i>et al.</i> Wiz binds active promoters and CTCF-binding sites and is required for normal behaviour in the mouse. Elife. 5: e15082

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K27ac, RNA Pol II, JDP2	Genome Res.	2016	"Petros Kolovos <i>et al.</i> Binding of nuclear factor κ B to noncanonical consensus sites reveals its multimodal role during the early inflammatory response. <i>Genome Res.</i> 26:1478"
ChIP-Seq	V5-tagged DUX4	Hum Mol Genet.	2016	Jocelyn O. Eidahl <i>et al.</i> Mouse Dux is myotoxic and shares partial functional homology with its human paralog DUX4. <i>Hum Mol Genet.</i> 25: 4577
ChIP-Seq	LL-37 (antimicrobial peptide)	J Cancer	2016	Mindy Munoz <i>et al.</i> Antimicrobial peptide LL-37 participates in the transcriptional regulation of melanoma cells. <i>J Cancer.</i> 26: 2341
ChIP-Seq, Spike-in	BRD4	J Clin Invest.	2016	Mark L. McClelland <i>et al.</i> CCAT1 is an enhancer-templated RNA that predicts BET sensitivity in colorectal cancer. <i>J Clin Invest.</i> 126: 639
ChIP-Seq	Myc-tagged PET-1	J Neurosci.	2016	Steven C. Wyler <i>et al.</i> Pet-1 switches transcriptional targets postnatally to regulate maturation of serotonin neuron excitability. <i>J Neurosci.</i> 36:1758
ChIP-Seq	Androgen Receptor, Estrogen Receptor	Mol Cancer Res.	2016	Nicholas D'Amato <i>et al.</i> Cooperative Dynamics of AR and ER Activity in Breast Cancer. <i>Mol Cancer Res.</i> 14:1054
ChIP-Seq	Androgen Receptor	Mol Cell	2016	Boyu Zhang <i>et al.</i> Non-Cell-Autonomous Regulation of Prostate Epithelial Homeostasis by Androgen Receptor. <i>Mol Cell.</i> 63: 976
ChIP-Seq	H3K27me3	Mol Cell Biol.	2016	Kyung Hyun Yoo <i>et al.</i> Histone demethylase KDM6A controls the mammary luminal lineage through enzyme-independent mechanisms. <i>Mol Cell Biol.</i> 36: 2108
ChIP-Seq, ChIP-qPCR	p53	Mol Oncol.	2016	Cheryl Chan <i>et al.</i> Global re-wiring of p53 transcription regulation by the hepatitis B virus X protein. <i>Mol Oncol.</i> 10: 1183
hMeDIP-Seq	5-Hydroxymethylcytosine	Nat Commun.	2016	Carolina M. Greco <i>et al.</i> DNA hydroxymethylation controls cardiomyocyte gene expression in development and hypertrophy. <i>Nat Commun.</i> 7: 12418
Low Cell ChIP-Seq	H3K27me3, CTCF	Nat Genet.	2016	Joachim Weischenfeldt <i>et al.</i> Pan-cancer analysis of somatic copy-number alterations implicates IRS4 and IGF2 in enhancer hijacking. <i>Nat Genet.</i> 49: 65
ChIP-Seq	H3K27ac	Nat Genet.	2016	Pratiti Bandopadhyay <i>et al.</i> MYB-QKI rearrangements in angiocentric glioma drive tumorigenicity through a tripartite mechanism. <i>Nat Genet.</i> 48: 273
ChIP-Seq	BRD4, H3K27ac, H3K4me1, H3K27me3, HLX, LHX2, LMX1A	Nature	2016	Charles Y. Lin <i>et al.</i> Active medulloblastoma enhancers reveal subgroup-specific cellular origins. <i>Nature.</i> 530: 57
ChIP-qPCR	H3K9me3, H3 pan-acetyl	Nucleic Acids Res.	2016	Xiaoyu Chen <i>et al.</i> Probing the impact of chromatin conformation on genome editing tools. <i>Nucleic Acids Res.</i> 44: 6482
ChIP-Seq, ChIP-qPCR	H3K36me3	Oncogene	2016	Thai H. Ho <i>et al.</i> High-resolution profiling of histone h3 lysine 36 trimethylation in metastatic renal cell carcinoma. <i>Oncogene.</i> 35: 1565
ChIP-Seq	MTA1	Oncotarget	2016	Swati Dhar <i>et al.</i> Dietary pterostilbene is a novel MTA1-targeted chemopreventive and therapeutic agent in prostate cancer. <i>Oncotarget.</i> 7: 18469

Technique	Target	Journal	Year	Reference
ChIP-Seq	PLZF	PLoS Genet.	2016	Ramakrishna Kommagani <i>et al.</i> The Promyelocytic leukemia zinc finger transcription factor Is critical for human endometrial stromal cell decidualization. PLoS Genet. 12: e1005937
ChIP-Seq, Spike-in	H3K27me3, H3K-4me3, H3K9me3	PLoS One	2016	Brian Egan <i>et al.</i> An alternative approach to ChIP-Seq normalization enables detection of genome- wide changes in histone H3 lysine 27 trimethylation upon EZH2 inhibition. PLoS One. 11: e0166438
ChIP-Seq	HDAC1, HDAC2, GATA2	PLoS One	2016	"Jeffrey R. Shearstone <i>et al.</i> Chemical Inhibition of histone deacetylases 1 and 2 Induces fetal hemoglobin through activation of GATA2. PLoS One. 11: e0153767"
ChIP-Seq	NR4A1	PLoS One	2016	Ryan P. Duren <i>et al.</i> Genome wide mapping of NR4A binding reveals cooperativity with ETS factors to promote epigenetic activation of distal enhancers in acute myeloid leukemia cells. PLoS One. 11: e0150450
ChIP-Seq, ChIP-qPCR	H3K9Ac	Stem Cells Transl Med.	2016	Dalia Ali <i>et al.</i> Epigenetic library screen identifies abexinostat as novel regulator of adipocytic and osteoblastic differentiation of human skeletal (mesenchymal) stem cells. Stem Cells Transl Med. 5: 1036
ChIP-Seq	FLAG-tagged TWIST2	Am J Hum Genet.	2015	Shannon Marchegiani <i>et al.</i> Recurrent mutations in the basic domain of TWIST2 cause ablepharon macrostomia and barber-say syndromes. Am J Hum Genet. 97: 99
ChIP-Seq	H3K4me1	BMC Biol.	2015	Sara K. Harten <i>et al.</i> The recently identified modifier of murine metastable epialleles, Rearranged L-Myc Fusion, is involved in maintaining epigenetic marks at CpG island shores and enhancers. BMC Biol. 13: 21
ChIP-Seq	FOXM1	Breast Cancer Res Treat.	2015	Christina Yau <i>et al.</i> FOXM1 cistrome predicts breast cancer metastatic outcome better than FOXM1 expression levels or tumor proliferation index. Breast Cancer Res Treat. 154: 23
ChIP-Seq	SIX1	Cancer Cell	2015	Jenny Wegert <i>et al.</i> Mutations in the SIX1/2 pathway and the DROSHA/DGCR8 miRNA microprocessor complex underlie high-risk blastemal type Wilms tumors. Cancer Cell. 27: 298
ChIP-Seq, ChIP-qPCR	b-Arrestin-1	Cancer Res.	2015	Smitha Pillai <i>et al.</i> β -arrestin-1 mediates nicotine-induced metastasis through E2F1 target genes that modulate epithelial-mesenchymal transition. Cancer Res. 75: 1009
ChIP-Seq	Progesterone Receptor	Cell Rep.	2015	Vincent J. Lynch <i>et al.</i> Ancient transposable elements transformed the uterine regulatory landscape and transcriptome during the evolution of mammalian pregnancy. Cell Rep. 10: 551
ChIP-Seq	H3K9Ac, RNA Pol II, TFIIB	Circ Heart Fail.	2015	Danish Sayed <i>et al.</i> Acute targeting of general transcription factor IIB restricts cardiac hypertrophy via selective inhibition of gene transcription. Circ Heart Fail. 8: 138
ChIP-Seq	HA-tagged SHOX2	Development	2015	Wenduo Ye <i>et al.</i> A common Shox-2-Nkx2-5 antagonistic mechanism primes the pacemaker cell fate in the pulmonary vein myocardium and sinoatrial node. Development. 142: 2521
ChIP-qPCR	EZH2, SUZ12, H3K27me3, RNA Pol II	EMBO J.	2015	Stephen G. Dann <i>et al.</i> Reciprocal regulation of amino acid import and epigenetic state through Lat1 and EZH2. EMBO J. 34: 1773

Technique	Target	Journal	Year	Reference
ChIP-Seq	Progesterone Receptor, FOSL2	Endocrinology	2015	Erik C. Mazur <i>et al.</i> Progesterone receptor transcriptome and cistrome in decidualized human endometrial stromal cells. <i>Endocrinology</i> . 156: 2239
ChIP-Seq	H3K4me3, H3K9me3	Epigenetics	2015	Nioka C. Chisholm <i>et al.</i> Histone methylation patterns in astrocytes are influenced by age following ischemia. <i>Epigenetics</i> . 10: 142
ChIP-Seq	LHX6	Hum Mol Genet.	2015	"Jeffrey M. Cesario <i>et al.</i> Lhx6 and Lhx8 promote palate development through negative regulation of a cell cycle inhibitor gene, p57Kip2. <i>Hum Mol Genet.</i> 24: 5024"
ChIP-Seq	H3K27me3, H3K4me1, H3K4me3, H3K27ac, p300, RAR α	Immunity	2015	Chrysothemis C. Brown <i>et al.</i> Retinoic acid is essential for Th1 cell lineage stability and prevents transition to a Th17 cell program. <i>Immunity</i> . 42: 1
Next-Gen Bisulfite-Seq		J Clin Invest.	2015	Coralie Hoareau-Aveilla <i>et al.</i> Reversal of microRNA-150 silencing disadvantages crizotinib-resistant NPM-ALK(+) cell growth. <i>J Clin Invest.</i> 125: 3505
ChIP-Seq	H3K4me3	J Natl Cancer Inst.	2015	Cristian Taccioli <i>et al.</i> Repression of esophageal neoplasia and inflammatory signaling by anti-miR-31 delivery in vivo. <i>J Natl Cancer Inst.</i> 107: djv220
ChIP-Seq	BRD4	Mol Cancer Ther.	2015	Ryan Lenhart <i>et al.</i> Sensitivity of small cell lung cancer to BET inhibition is mediated by regulation of ASCL1 gene expression. <i>Mol Cancer Ther.</i> 14: 2167
ChIP-Seq	BAF180/PBRM1	Mol Cell	2015	Bokai Zhu <i>et al.</i> Coactivator-dependent oscillation of chromatin accessibility dictates circadian gene amplitude via REV-ERB loading. <i>Mol Cell.</i> 60: 769
ChIP-Seq	ZNF384	Mol Endocrinol.	2015	Paul Childress <i>et al.</i> Genome-wide mapping and interrogation of the Nmp4 anti-anabolic bone axis. <i>Mol Endocrinol.</i> 29: 1269
ChIP-Seq	FOXO1, RNA Pol II	Mol Endocrinol.	2015	Yasmin M. Vasquez <i>et al.</i> FOXO1 is required for binding of PR on IRF4, novel transcriptional regulator of endometrial stromal decidualization. <i>Mol Endocrinol.</i> 29: 421
ChIP-Seq	EGR2	Nat Commun.	2015	Tomohisa Okamura <i>et al.</i> TGF- β 3-expressing CD4+CD25-LAG3+ regulatory T cells control humoral immune responses. <i>Nat Commun.</i> 6: 6329
ChIP-Seq	SOX9	Nucleic Acids Res.	2015	Zhongcheng Shi <i>et al.</i> Context-specific role of SOX9 in NF- κ B mediated gene regulation in colorectal cancer cells. <i>Nucleic Acids Res.</i> 43: 6257
ChIP-Seq	H3K27me3	Nucleic Acids Res.	2015	Kyung Hyun Yoo <i>et al.</i> Loss of EZH2 results in precocious mammary gland development and activation of STAT5-dependent genes. <i>Nucleic Acids Res.</i> 43: 8774
ChIP-Seq	CCND1 (Cyclin D1)	Oncotarget	2015	Mathew C. Casimiro <i>et al.</i> Kinase-independent role of cyclin D1 in chromosomal instability and mammary tumorigenesis. <i>Oncotarget.</i> 6: 8524
ChIP-Seq Data Analysis	KDM3A	PLoS Biol.	2015	Mo-bin Cheng <i>et al.</i> Specific phosphorylation of histone demethylase KDM3A determines target gene expression in response to heat shock. <i>PLoS Biol.</i> 12: e1002026
ChIP-Seq	TRIM33	PLoS Genet.	2015	Luke Isbel <i>et al.</i> Trim33 binds and silences a class of young endogenous retroviruses in the mouse testis; a novel component of the arms race between retrotransposons and the host genome. <i>PLoS Genet.</i> 11: e1005693

Technique	Target	Journal	Year	Reference
ChIP-Seq	EGR1	PLoS One	2015	Anthony A. Portale <i>et al.</i> Characterization of FGF23-dependent Egr-1 cistrome in the mouse renal proximal tubule. <i>PLoS One</i> . 10: e0142924
ChIP-Seq	RAR α , RAR β , RXR α	Am J Physiol Gastrointest Liver Physiol.	2014	Yuqi He <i>et al.</i> Biological functional annotation of retinoic acid alpha and beta in mouse liver based on genome-wide binding. <i>Am J Physiol Gastrointest Liver Physiol</i> . 307: G205
ChIP-Seq	H3K27me3	Bioinformatics	2014	Yanxiao Zhang <i>et al.</i> PePr: A peak-calling prioritization pipeline to identify consistent or differential peaks from replicated ChIP-Seq data. <i>Bioinformatics</i> . 30: 2568
ChIP-Seq	PPAR γ	Biol Reprod.	2014	Kelsey E. Brooks <i>et al.</i> Peroxisome proliferator activator receptor gamma (PPAR γ) regulates conceptus elongation in sheep. <i>Biol Reprod</i> . 92: 42.
Next-Gen Bisulfite-Seq		Cell Rep.	2014	Deepak Kumar <i>et al.</i> Fibroblast growth factor maintains chondrogenic potential of limb bud mesenchymal cells by modulating DNMT3A recruitment. <i>Cell Rep</i> . 8: 1419
ChIP-Seq	SRC-2	Cell Rep.	2014	Erin Stashi <i>et al.</i> SRC-2 Is an essential coactivator for orchestrating metabolism and circadian rhythm. <i>Cell Rep</i> . 6: 633
ChIP-Seq, ChIP-qPCR	ZBTB20	Cereb Cortex	2014	Jakob V. Nielsen <i>et al.</i> Zbtb20 Defines a hippocampal neuronal identity through direct repression of genes that control projection neuron development in the isocortex. <i>Cereb Cortex</i> . 24: 1216
ChIP-Seq	H3K9me3, H3K36me3, H3K27me3	Genome Res.	2014	Goran Kungulovski <i>et al.</i> Application of histone modification-specific interaction domains as an alternative to antibodies. <i>Genome Res</i> . 24: 1842
ChIP-Seq	H3K36me3	Genome Res.	2014	Jeremy M. Simon <i>et al.</i> Variation in chromatin accessibility in human kidney cancer links H3K36 methyltransferase loss with widespread RNA processing defects. <i>Genome Res</i> . 24: 241
ChIP-Seq	H3K27ac	J Biol Chem.	2014	Andre Landin Malt <i>et al.</i> Identification of a face enhancer reveals direct regulation of LIM homeobox 8 (Lhx8) by wingless-int (WNT)/ β -catenin Signaling. <i>J Biol Chem</i> . 289: 30289
ChIP-Seq, ChIP-qPCR	SRC-3	J Clin Invest.	2014	Jun Qin <i>et al.</i> Androgen deprivation–induced NCoA2 promotes metastatic and castration-resistant prostate cancer. <i>J Clin Invest</i> . 124: 5013
ChIP-Seq, ChIP-qPCR	ZBTB20	J Med Genet.	2014	Malene B. Rasmussen <i>et al.</i> Neurodevelopmental disorders associated with dosage imbalance of ZBTB20 correlate with the morbidity spectrum of ZBTB20 candidate target genes. <i>J Med Genet</i> . 51: 605
ChIP-qPCR	EZH2, SUZ12, H3K27me3	Mol Cancer Ther.	2014	Sarah K. Knutson <i>et al.</i> Selective inhibition of EZH2 by EPZ-6438 leads to potent antitumor activity in EZH2 mutant Non-Hodgkin lymphoma. <i>Mol Cancer Ther</i> . 13: 842
ChIP-Seq	Fish H3K4me3	Mol Ecol Resour.	2014	Claudius F. Kratochwil <i>et al.</i> Mapping active promoters by ChIP-seq profiling of H3K4me3 in cichlid fish – a first step to uncover cis-regulatory elements in ecological model teleosts. <i>Mol Ecol Resour</i> . 15: 761
ChIP-Seq	SRC-1	Mol Endocrinol.	2014	Mounia Tannour-Louet <i>et al.</i> Hepatic SRC-1 activity orchestrates transcriptional circuitries of amino acid pathways with potential relevance for human metabolic pathogenesis. <i>Mol Endocrinol</i> . 28: 1707

Technique	Target	Journal	Year	Reference
ChIP-Seq	Estrogen Receptor	Mol Endocrinol.	2014	Sylvia C. Hewitt <i>et al.</i> Novel DNA motif binding activity observed In vivo with an estrogen receptor a mutant mouse. Mol Endocrinol. 26: 899
ChIP-Seq	H3K4me3, H3K-27me3	Nat Commun.	2014	Brian C. Belyea <i>et al.</i> Identification of renin progenitors in the mouse bone marrow that give rise to B-cell leukaemia. Nat Commun. 5: 3273
Next-Gen Bisulfite-Seq		Nat Neurosci.	2014	Brian G. Dias <i>et al.</i> Parental olfactory experience influences behavior and neural structure in subsequent generations. Nat Neurosci. 17: 89
ChIP-Seq	H3K9ac, H3K27Ac	Nature	2014	Paul A. Northcott <i>et al.</i> Enhancer hijacking activates GFI1 family oncogenes in medulloblastoma. Nature. 511: 428
ChIP-Seq	H3K4me3, H3K-27me3	Nature	2014	Volker Hovestadt <i>et al.</i> Decoding the regulatory landscape of medulloblastoma using DNA methylation sequencing. Nature. 510: 537
ChIP-Seq, ChIP-qPCR	SRF	Physiol Genomics	2014	Sharolyn V. Kawakami-Schulz. Serum response factor: positive and negative regulation of an epithelial gene expression network in the destrin mutant cornea. Physiol Genomics. 46: 277
ChIP-Seq, LightSwitch	PXR, p300, H3K-4me1, H3K27ac	PLoS Genet.	2014	Robin P. Smith <i>et al.</i> Genome-wide discovery of drug-dependent human liver regulatory elements. PLOS Genet. 10: e1004648
ChIP-Seq	GATA2, FOXA1, p300, CBP, SRC-1, SRC-2, SRC-3, Androgen Receptor	Proc Natl Acad Sci USA	2014	Bin He <i>et al.</i> GATA2 facilitates steroid receptor coactivator recruitment to the androgen receptor complex. Proc Natl Acad Sci. 111: 18261
ChIP-Seq	Vitamin D Receptor	BMC Med.	2013	Adam E. Handel <i>et al.</i> Vitamin D receptor ChIP-seq in primary CD4+ cells: relationship to serum 25-hydroxyvitamin D levels and autoimmune disease. BMC Med. 11: 163
ChIP-Seq	H3K27me3	Cancer Cell	2013	Sebastian Bender <i>et al.</i> Reduced H3K27me3 and DNA hypomethylation are major drivers of gene expression in K27M mutant pediatric high-grade gliomas. Cancer Cell. 24: 660
ChIP-Seq	H3K36me3	Cancer Discov.	2013	Lynn Bjerke <i>et al.</i> Histone H3.3 Mutations Drive Pediatric Glioblastoma through Upregulation of MYCN. Cancer Discov. 3: 512
ChIP-Seq	CCND1 (Cyclin D1)	Cancer Res.	2013	Xiaoming Ju <i>et al.</i> Identification of a cyclin D1 network in prostate cancer that antagonizes epithelial-mesenchymal restraint. Cancer Res. 74: 508
ChIP-Seq	PPARα	Chem Biol Interact.	2013	Patrick D. McMullen <i>et al.</i> A map of the PPARα transcription regulatory network for primary human hepatocytes. Chem Biol Interact. 209: 14
Next-Gen Bisulfite-Seq		Clin Cancer Res.	2013	David S. Shames <i>et al.</i> Loss of NAPRT1 expression by tumor-specific promoter methylation provides a novel predictive biomarker for NAMPT inhibitors. Clin Cancer Res. 19: 6912
ChIP-Seq	COUP-TFII (NR2F2)	Dev Cell	2013	San-pin Wu <i>et al.</i> Atrial identity is determined by a COUP-TFII regulatory network. Dev Cell. 25: 417
ChIP-Seq	FOXA2	EMBO J.	2013	Justyna Filant <i>et al.</i> Integrated chromatin immunoprecipitation sequencing and microarray analysis identifies FOXA2 target genes in the glands of the mouse uterus. FASEB J. 28: 230

Technique	Target	Journal	Year	Reference
ChIP-Seq	RNA Pol II	J Allergy Clin Immunol.	2013	Purvi Mehrotra <i>et al.</i> PARP-14 and its enzyme activity regulates Th2 differentiation and allergic airway disease. J Allergy Clin Immunol. 131: 521
ChIP-qPCR	NFkB (p52 and p65)	J Biol Chem.	2013	Sarah L. Doyle <i>et al.</i> Nuclear factor kB2 p52 protein has a role in antiviral immunity through IkB kinase epsilon-dependent induction of Sp1 protein and interleukin 15. J Biol Chem. 288: 25066
ChIP-chip, ChIP-qPCR	C/EBPβ, RNA Pol II	J Biol Chem.	2013	Hana Vakili <i>et al.</i> CCAAT-enhancer-binding protein β (C/EBPβ) and downstream human placental growth hormone genes are targets for dysregulation in pregnancies complicated by maternal obesity. J Biol Chem. 288: 22849
ChIP-Seq	KDM2B	J Clin Invest.	2013	Alexandros Tzatsos <i>et al.</i> KDM2B promotes pancreatic cancer via Polycomb-dependent and -independent transcriptional programs. J Clin Invest. 123: 727
ChIP-Seq	ASXL1	J Exp Med.	2013	Omar Abdel-Wahab <i>et al.</i> Deletion of Asxl1 results in myelodysplasia and severe developmental defects in vivo. J Exp Med. 210: 2641
ChIP-Seq	NR2F2	Mol Endocrinol.	2013	Xilong Li <i>et al.</i> COUP-TFII regulates human endometrial stromal genes involved in inflammation. Mol Endocrinol. 27: 2041
ChIP-Seq	Progesterone Receptor	Mol Endocrinol.	2013	Ashlee R. Lain <i>et al.</i> Research Resource: Progesterone receptor targetome underlying mammary gland branching morphogenesis. Mol Endocrine. 27: 1743
ChIP-qPCR	H3K9me3, H3	Nat Immunol.	2013	Patrick M. Gubser <i>et al.</i> Rapid effector function of memory CD8+ T cells requires an immediate-early glycolytic switch. Nat Immunol. 14: 1064
ChIP-Seq	GFI1	Nat Immunol.	2013	Chauncey J. Spooner <i>et al.</i> Specification of type 2 innate lymphocytes by the transcriptional determinant Gfi1. Nat Immunol. 14: 1229
ChIP-Seq	p53	Physiol Genomics	2013	Yuwen Li <i>et al.</i> Genome-wide analysis of the p53 gene regulatory network in the developing mouse kidney. Physiol Genomics. 45: 948
ChIP-Seq	MYRF	PLoS Biol.	2013	Helena Bujalka <i>et al.</i> MYRF is a membrane-associated transcription factor that autoproteolytically cleaves to directly activate myelin genes. PLoS Biol. 11: e1001625
ChIP-Seq	p300	PLoS Genet.	2013	Aaron M. Wenger <i>et al.</i> The enhancer landscape during early neocortical development reveals patterns of dense regulation and co-option. PLoS Genet. 9: e1003728
ChIP-Seq	RNA Pol II	PLoS One	2013	Jonathan P. Riley <i>et al.</i> PARP-14 binds specific DNA sequences to promote Th2 cell gene expression. PLoS One. 8: e83127
ChIP-Seq, ChIP-qPCR	RXRα, RNA Pol II	PLoS One	2013	Astrid Kosters <i>et al.</i> Sexually dimorphic genome-wide binding of retinoid X receptor alpha (RXRα) determines male-female differences in the expression of hepatic lipid processing genes in mice. PLoS One. 8: e71538
ChIP-Seq	H3K4me3	PLoS One	2013	Alex Gutteridge <i>et al.</i> Novel pancreatic endocrine maturation pathways identified by genomic profiling and causal reasoning. PLoS One. 8: e56024.
ChIP-Seq	Farnesoid X Receptor	Am J Physiol Gastrointest Liver Physiol.	2012	Julia Yue Cui <i>et al.</i> Bile acids via FXR initiate the expression of major transporters involved in the enterohepatic circulation of bile acids in newborn mice. Am J Physiol Gastrointest Liver Physiol. 302: G979

Technique	Target	Journal	Year	Reference
ChIP-Seq	Androgen Receptor	BMC Genomics	2012	Zhou Zhu <i>et al.</i> Dose-dependent effects of small-molecule antagonists on the genomic landscape of androgen receptor binding. <i>BMC Genomics</i> . 13: 355
ChIP-Seq	RNA Pol II	Cardiovasc Res.	2012	Mingyue Han <i>et al.</i> GATA4 expression is primarily regulated via a miR-26b-dependent post-transcriptional mechanism during cardiac hypertrophy. <i>Cardiovasc Res</i> . 93: 645
ChIP-qPCR	PDX1, TCF3	Chem Biol.	2012	Alice Kiselyuk <i>et al.</i> HNF4a antagonists discovered by a high-throughput screen for modulators of the human insulin promoter. <i>Chem Biol</i> . 19:806
ChIP-Seq	RNA Pol II	J Biol Chem.	2012	Danish Sayed <i>et al.</i> Transcriptional regulation patterns revealed by high-resolution chromatin immunoprecipitation during cardiac hypertrophy. <i>J Biol Chem</i> . 288: 2546
ChIP-Seq	NKX3.1	J Clin Invest.	2012	Philip D. Anderson <i>et al.</i> Nkx3.1 and Myc crossregulate shared target genes in mouse and human prostate tumorigenesis. <i>J Clin Invest</i> . 122: 1907
ChIP-Seq	FLAG-tagged CCND1 (Cyclin D1)	J Clin Invest.	2012	Mathew C. Casimiro <i>et al.</i> ChIP sequencing of cyclin D1 reveals a transcriptional role in chromosomal instability in mice. <i>J Clin Invest</i> . 122: 833
ChIP-Seq	LSD1, H3K4me2	Mol Cell Biol.	2012	Venugopalan D. Nair <i>et al.</i> Involvement of histone demethylase LSD1 in short-time-scale gene expression changes during cell cycle progression in embryonic stem cells. <i>Mol Cell Biol</i> . 32: 4861
ChIP-Seq	Progesterone Receptor	Mol Endocrinol.	2012	Cory A. Rubel <i>et al.</i> Genome-Wide Profiling of Progesterone Receptor Binding in the Mouse Uterus. <i>Mol Endocrinol</i> . 26: 1428
ChIP-Seq	Estrogen Receptor, RNA Pol II	Mol Endocrinol.	2012	Sylvia C. Hewitt <i>et al.</i> Whole-Genome Estrogen Receptor Binding in Mouse Uterine Tissue Revealed by ChIP-Seq. <i>Mol Endocrinol</i> . 26: 887
ChIP-Seq Data Analysis	pTyr37 H2B	Nat Struct Mol Biol.	2012	Kiran Mahajan <i>et al.</i> H2B Tyr37 phosphorylation suppresses expression of replication-dependent core histone genes. <i>Nat Struct Mol Biol</i> . 19: 930
ChIP-Seq	EZH2	Nature	2012	Michael T. McCabe <i>et al.</i> EZH2 inhibition as a therapeutic strategy for lymphoma with EZH2-activating mutations. <i>Nature</i> . 492: 108
ChIP-Seq	Glucocorticoid Receptor	PLoS One	2012	Marie-José C. van Lierop <i>et al.</i> Org 214007-0: A novel non-steroidal selective glucocorticoid receptor modulator with full anti-inflammatory properties and improved therapeutic index. <i>PLoS One</i> . 7: e48385
ChIP-chip	B-MYB	PLoS One	2012	Ming Zhan <i>et al.</i> The B-MYB transcriptional network guides cell cycle progression and fate decisions to sustain self-renewal and the identity of pluripotent stem cells. <i>PLoS One</i> . 7: e42350
ChIP-chip	RORa	PLoS One	2012	Yongjun Wang <i>et al.</i> Regulation of p53 Stability and Apoptosis by a ROR Agonist. <i>PLoS One</i> . 7: e34921
ChIP-Seq	Progesterone Receptor	PLoS One	2012	Ping Yin <i>et al.</i> Genome-wide progesterone receptor binding: cell type-specific and shared mechanisms in T47D breast cancer cells and primary leiomyoma cells. <i>PLoS One</i> . 7: e29021
ChIP-Seq	FLAG-tagged BAP1, OGT, HCF1	Science	2012	Anwasha Dey <i>et al.</i> Loss of the tumor suppressor BAP1 causes myeloid transformation. <i>Science</i> . 337: 1541

Technique	Target	Journal	Year	Reference
ChIP-Seq	RBPJ	Stem Cells	2012	Yaochen Li <i>et al.</i> Genome-wide Analysis of N1ICD/RBPJ Targets in vivo Reveals Direct Transcriptional Regulation of Wnt, SHH, and Hippo Pathway Effectors by Notch1. <i>Stem Cells</i> . 30: 741
ChIP-chip	IRF4	Allergy	2011	Barrenas S Bruhn <i>et al.</i> Increased expression of IRF4 and ETS1 in CD4+ cells from patients with intermittent allergic rhinitis. <i>Allergy</i> . 67: 33
ChIP-qPCR	Progesterone Receptor	Biochem Pharmacol.	2011	Matthew R Yudt <i>et al.</i> Discovery of a novel mechanism of steroid receptor antagonism: WAY-255348 modulates progesterone receptor cellular localization and promoter interactions. <i>Biochem Pharmacol</i> . 82: 1709
MeDIP-chip	5-Methylcytosine	BMC Biol.	2011	Genevieve Lavoie <i>et al.</i> PKC isoforms interact with and phosphorylate DNMT1. <i>BMC Biol</i> . 9: 31
ChIP-qPCR	IRF4, cMyc	Br J Haematol.	2011	Antonia Lopez-Girona <i>et al.</i> (2011) Lenalidomide downregulates the cell survival factor, interferon regulatory factor-4, providing a potential mechanistic link for predicting response. <i>Br J Haematol</i> . 154: 325
ChIP-qPCR	CTCF	DNA Cell Biol.	2011	"Yan Jin <i>et al.</i> Enhancer-Blocking Activity Is Associated with Hypersensitive Site V Sequences in the Human Growth Hormone Locus Control Region. <i>DNA Cell Biol</i> . 30: 995"
ChIP-chip	Progesterone Receptor	EMBO J.	2011	Heather L. Franco <i>et al.</i> Epithelial progesterone receptor exhibits pleiotropic roles in uterine development and function. <i>FASEB J</i> . 26: 1218
ChIP-qPCR	PPARg	Invest Ophthalmol Vis Sci.	2011	Gerard A Rodrigues <i>et al.</i> Differential Effects of PPARg Ligands on Oxidative Stress-Induced Death of Retinal Pigmented Epithelial Cells. <i>Invest Ophthalmol Vis Sci</i> . 52: 890
ChIP-Seq	NELF-b (COBRA1)	J Biol Chem.	2011	"Jianlong Sun <i>et al.</i> Genetic and Genomic Analyses of RNA Polymerase II-pausing Factor in Regulation of Mammalian Transcription and Cell Growth. <i>J Biol Chem</i> . 286: 36248"
ChIP-chip	RNA Pol II, H3K9Ac	J Biol Chem.	2011	Hong Hao <i>et al.</i> The Transcription Factor NRL Controls Photoreceptor-specific Expression of Myocyte Enhancer Factor Mef2c from an Alternative Promoter. <i>J Biol Chem</i> . 286: 34893
ChIP-chip	LXR, RXR	J Biol Chem.	2011	"Qi Shen <i>et al.</i> Liver X Receptor-Retinoid X Receptor (LXR-RXR) Heterodimer Cistrome Reveals Coordination of LXR and AP1 Signaling in Keratinocytes. <i>J Biol Chem</i> . 286: 14554"
ChIP-chip	C/EBPb	J Mol Endocrinol.	2011	Aristides Lytras <i>et al.</i> Identification of functional CCAAT/ enhancer-binding protein and Ets protein binding sites in the human chorionic somatomammotropin enhancer sequences. <i>J Mol Endocrinol</i> . 47: 179
ChIP-Seq	MSGN1	Nat Commun.	2011	Ravindra B Chalamalasetty <i>et al.</i> The Wnt3a/ β -catenin target gene Mesogenin1 controls the segmentation clock by activating a Notch signalling program. <i>Nat Commun</i> . 2: 390
ChIP-Seq	FOXP3	Nucleic Acids Res.	2011	Fabian Birzele <i>et al.</i> Next-generation insights into regulatory T cells: expression profiling and FoxP3 occupancy in Human. <i>Nucleic Acids Res</i> . 39: 7946
ChIP-chip	DAF-12	PLoS Genet.	2011	Daniel Hochbaum <i>et al.</i> DAF-12 Regulates a Connected Network of Genes to Ensure Robust Developmental Decisions. <i>PLoS Genet</i> . 7: e1002179

Technique	Target	Journal	Year	Reference
Next-Gen Bisulfite-Seq		PLoS One	2011	Michela Deleidi <i>et al.</i> Oct4-Induced Reprogramming Is Required for Adult Brain Neural Stem Cell Differentiation into Midbrain Dopaminergic Neurons. PLoS ONE. 6: e19926
ChIP-qPCR	BRD4	Proc Natl Acad Sci USA	2011	Jennifer A. Mertz <i>et al.</i> Targeting MYC dependence in cancer by inhibiting BET bromodomains. Proc Natl Acad Sci. 108:16669
ChIP-chip	p73	Proc Natl Acad Sci USA	2011	Jennifer M Rosenbluth <i>et al.</i> Differential regulation of the p73 cistrome by mammalian target of rapamycin reveals transcriptional programs of mesenchymal differentiation and tumorigenesis. Proc Natl Acad Sci. 108: 2076
ChIP-qPCR	Aryl Hydrocarbon Receptor	Toxicol Appl Pharmacol.	2011	K Nadira De Abrew <i>et al.</i> Regulation of Bach2 by the aryl hydrocarbon receptor as a mechanism for suppression of B-cell differentiation by 2,3,7,8-tetrachlorodibenzo-p-dioxin. Toxicol Appl Pharmacol. 252: 150
ChIP-qPCR	NFkB (p50 and p65)	Cardiovasc Res.	2010	Alina G Sofronescu <i>et al.</i> FGF-16 is a target for adrenergic stimulation through NF-kB activation in postnatal cardiac cells and adult mouse heart. Cardiovasc Res. 87: 102
ChIP-Seq	Vitamin D Receptor	Genome Res.	2010	Sreeram V Ramagopalan <i>et al.</i> A ChIP-seq defined genome-wide map of vitamin D receptor binding: Associations with disease and evolution. Genome Res. 20: 1352
ChIP-qPCR	MAF1, RPC39, mTOR, BRF1, Raptor	J Biol Chem.	2010	Boris Shor <i>et al.</i> Requirement of the mTOR Kinase for the Regulation of Maf1 Phosphorylation and Control of RNA Polymerase III-dependent Transcription in Cancer Cells. J Biol Chem. 286: 15380
ChIP-qPCR	p73	Mol Cancer	2010	Kathryn G Eby <i>et al.</i> ISG20L1 is a p53 family target gene that modulates genotoxic stress-induced autophagy. Mol Cancer. 9: 95
ChIP-Seq	SRC-3, RNA Pol II	Mol Endocrinol.	2010	Rainer B Lanz <i>et al.</i> Global Characterization of Transcriptional Impact of the SRC-3 Coregulator. Mol Endocrinol. 24: 859
ChIP-chip	Androgen Receptor	Mol Endocrinol.	2010	Anastasia Wyce <i>et al.</i> The Androgen Receptor Modulates Expression of Genes with Critical Roles in Muscle Development and Function. Mol Endocrinol. 24: 1665
ChIP-chip	RNA Pol II	Mol Vis.	2010	Padmaja Tummala <i>et al.</i> Temporal ChIP-on-Chip of RNA-Polymerase-II to detect novel gene activation events during photoreceptor maturation. Mol Vis. 16: 252
ChIP-qPCR	Aryl Hydrocarbon Receptor, cMaf	Nat Immunol.	2010	Lionel Apetoh <i>et al.</i> The aryl hydrocarbon receptor interacts with c-Maf to promote the differentiation of type 1 regulatory T cells induced by IL-27. Nat Immunol. 11: 854
ChIP-qPCR	Myc-tagged PET-1	Nat Neurosci.	2010	Chen Liu <i>et al.</i> Pet-1 is required across different stages of life to regulate serotonergic function. Nat Neurosci. 13: 1190
ChIP-Seq	Pregnane X Receptor	Nucleic Acids Res.	2010	Julia Yue Cui <i>et al.</i> ChIPing the cistrome of PXR in mouse liver. Nucleic Acids Res. 38: 7943
ChIP-chip	O-GlcNAc, RNA Pol II	Proc Natl Acad Sci USA	2010	Dona C Love <i>et al.</i> Dynamic O-GlcNAc cycling at promoters of Caenorhabditis elegans genes regulating longevity, stress, and immunity. Proc Natl Acad Sci. 107: 7413
ChIP-Seq	Pregnane X Receptor	Toxicol Sci.	2010	Julia Yue Cui <i>et al.</i> Genetic and Epigenetic Regulation and Expression Signatures of Glutathione S-Transferases in Developing Mouse Liver. Toxicol Sci. 116: 32

Technique	Target	Journal	Year	Reference
ChIP-chip	Aryl Hydrocarbon Receptor	Toxicol Sci.	2010	K Nadira De Abrewet <i>et al.</i> An Integrated Genomic Analysis of Aryl Hydrocarbon Receptor-Mediated Inhibition of B-Cell Differentiation. <i>Toxicol Sci.</i> 118:454
ChIP-chip	STAT4	J Immunol.	2009	"Seth R Good <i>et al.</i> Temporal Induction Pattern of STAT4 Target Genes Defines Potential for Th1 Lineage-Specific Programming. <i>J Immunol.</i> 183: 3839"
ChIP-Seq	p73	Mol Cell Biol.	2008	Jennifer M Rosenbluth <i>et al.</i> A Gene Signature-Based Approach Identifies mTOR as a Regulator of p73. <i>Mol Cell Biol.</i> 28: 5951
ChIP-qPCR	Aryl Hydrocarbon Receptor	Nature	2008	Francisco J Quintana <i>et al.</i> Control of Treg and TH17 cell differentiation by the aryl hydrocarbon receptor. <i>Nature.</i> 453: 65