



CUT&Tag-IT[®] Core Assay Kit

Catalog No. 53176 (16 rxns)

(Version A1)

US Pat. No. 10,689,643, EP Pat. No. 2999784
and related patents and applications

Copyright 2024 Active Motif, Inc.

Information in this manual is subject to change without notice and does not constitute a commitment on the part of Active Motif, Inc. It is supplied on an “as is” basis without any warranty of any kind, either explicit or implied. Information may be changed or updated in this manual at any time. This documentation may not be copied, transferred, reproduced, disclosed, or duplicated, in whole or in part, without the prior written consent of Active Motif, Inc. This documentation is proprietary information and protected by the copyright laws of the United States and international treaties. The manufacturer of this documentation is Active Motif, Inc.

US Pat. No. 10,689,643, EP Pat. No. 2999784 and related patents and applications

© 2024 Active Motif, Inc., 1914 Palomar Oaks Way, Suite 100; Carlsbad, CA 92008. All rights reserved.

All trademarks, trade names, service marks or logos referenced herein belong to their respective companies. Oligonucleotide sequences © 2019 Illumina, Inc. All rights reserved.

The guinea pig anti-rabbit secondary antibody is provided under license from Rockland.

For Research Use Only. Not for use in diagnostic procedures.

Permitted Use; Resale Prohibited

In the absence of an express written agreement to the contrary, all products are sold and services deliverables are provided by Active Motif for (a) internal in vitro research purposes only and may not be used for services or any other commercial purpose (b) the exclusive use of the original purchaser, and are not to be resold. You agree not to reverse engineer or otherwise attempt to discover the structure or composition of products or services deliverables unless we otherwise agree in writing.

Contents

Overview	1
Kit Components and Storage	2
CUT&Tag-IT® Core Assay Kit Protocol	4
Prepare Fresh Cells (30 minutes)	4
Bind Cells to Concanavalin A Beads (30 minutes)	4
Bind Primary Antibody (2 hours to overnight)	5
Bind Secondary Antibody (60 minutes)	5
Bind CUT&Tag-IT® Assembled pA-Tn5 Transposomes (60 minutes)	6
Tagmentation (60 minutes)	6
DNA Extraction (60 minutes)	7
References	8
Technical Services	9

Overview

Cleavage Under Targets and Tagmentation (CUT&Tag) is a method to investigate genomic localization of histone modifications and some transcription factors that reveals interactions between proteins and DNA or identifies DNA binding sites for proteins of interest¹.

Unlike MNase-Seq or ATAC-Seq methods that target open chromatin and are therefore dependent on chromatin accessibility, CUT&Tag utilizes an antibody-based enzyme tethering strategy to target specific histone modifications or proteins to reveal chromatin-binding information that is specific to those sites or proteins of interest.

CUT&Tag is based on the same principles as ChIP-Seq, but with several changes to the protocol that are advantageous in certain situations. Instead of the sonication of fixed chromatin and immunoprecipitation steps performed in ChIP-Seq protocols, in CUT&Tag, fresh (not frozen) unfixed cells are bound to concanavalin A beads and the antibody incubation is performed with cells in their native state. Directly following antibody binding, the chromatin is sheared and NGS libraries are prepared in a single step by tagmentation using the protein A-Tn5 (pA-Tn5) transposome enzyme that has been pre-loaded with sequencing adapters.

The CUT&Tag-IT[®] Core Assay Kit is optimized for 50,000 to 500,000 cells per reaction, although 200,000 cells are recommended. This Kit provides optimized reagents and protocol to perform 16 CUT&Tag reactions and yields purified DNA that can be amplified and indexed for NGS sequencing. This kit does not include indexing primers or DNA polymerase for library amplification.

CUT&Tag-IT[®] Core Assay Kit Advantages

- Compatible with as few as 5,000 cells
- Low background signal enables lower sequencing depth
- Use your own indexing primers and DNA polymerase for library creation
- No artifacts caused by formaldehyde crosslinking

Product	Format	Catalog No.
CUT&Tag-IT [®] Core Assay Kit	16 rxns	53176



Kit Components and Storage

The kit contains sufficient reagents for 16 CUT&Tag Assay reactions. The reagents in this kit have multiple storage temperatures. Please store components according to the storage conditions below. All reagents are guaranteed stable for 6 months from date of receipt when stored properly.

Reagents	Quantity	Storage
5% Digitonin	610 μ L	-20°C
Concanavalin A Beads	320 μ L	4°C
CUT&Tag-IT™ Assembled pA-Tn5 Transposomes	16 μ L	-20°C
Tagmentation Buffer	2 mL	4°C or -20°C
1X Binding Buffer	55 mL	4°C
1X Wash Buffer	40 mL	4°C
Dig-Wash Buffer	50 mL	4°C
Antibody Buffer	800 μ L	4°C
Dig-300 Buffer	50 mL	4°C
Guinea Pig Anti-Rabbit Antibody	16 μ L	-20°C
Rabbit Anti-Mouse Antibody	16 μ L	-20°C
Protease Inhibitor Cocktail	1.42 mL	-20°C
0.5M EDTA	68 μ L	RT
10% SDS	20 μ L	RT
10 μ g/ μ L Proteinase K	18 μ L	-20°C
DNA Purification Columns	16 columns	RT
DNA Purification Binding Buffer	10 mL	RT
DNA Purification Wash Buffer	12 mL	RT
DNA Purification Elution Buffer	560 μ L	RT
3M Sodium Acetate	128 μ L	RT

Additional Materials Required

- Molecular grade nuclease free water
- 100% Ethanol
- 80% Ethanol
- Magnetic Rack for 1.5/2 mL tubes
- Magnetic Rack for 200 μ L 8-well PCR strip tubes
- Nutator or orbital rotating mixer
- Vortexer
- Thermal cycler
- NGS indexing primers
- DNA Polymerase for library creation
- 1.5 mL Low-bind Microcentrifuge tubes
- 2 mL Low-bind Microcentrifuge tubes
- Microcentrifuge for 8-well PCR strip tubes with caps, or for 0.2 and 1.5 mL tubes
- Multichannel pipette (20 - 200 μ L)
- Filter tips for pipettes

CUT&Tag-IT[®] Core Assay Kit Protocol

This Kit is optimized for 50,000 to 500,000 cells, although 200,000 cells are recommended.

Note: This protocol is optimized for non-adherent cells. If adherent cells are to be used, avoid using Trypsin or cell detachment solutions that digest cell membrane proteins. Cell membrane proteins are necessary to bind the cells to the Concanavalin A Beads in Section B of the assay. To lift adherent cells from growth plates, use an enzyme-free dissociation method such as scraping with a rubber policeman.

Prepare Fresh Cells (30 minutes)

Note: Before using 1X Wash Buffer in Step 3 below, you must add Protease Inhibitor Cocktail to the 1X Wash Buffer. Add 10 μ L Protease Inhibitor Cocktail for every 1 mL of 1X Wash Buffer. 2.5 mL total are used per cell sample. Keep on ice. Wash Buffer can be stored at 4°C for up to 1 week once Protease Inhibitor Cocktail has been added.

1. Harvest fresh culture(s) at room temperature and count cells. Place the desired amount of cells into a 1.5 mL microcentrifuge tube.
2. Centrifuge for three minutes at 600 x *g* at room temperature, then remove supernatant.
3. Resuspend in 1 mL 1X Wash Buffer at room temperature, then centrifuge for another three minutes at 600 x *g* at room temperature, and remove supernatant.
4. Resuspend in 1.5 mL 1X Wash Buffer and transfer to a 2 mL microcentrifuge tube. Place tube with resuspended pellet on ice while the beads are prepared in the next steps.

Bind Cells to Concanavalin A Beads (30 minutes)

Note: This Kit is optimized for 50,000 to 500,000 cells. Best results have been obtained with 200,000 cells. 500,000 cells can become viscous in the steps below (**Bind Cells to Concanavalin A Beads, Steps 5-9**) and DNA purification.

5. Gently resuspend and withdraw enough of the Concanavalin A Bead slurry such that there will be 20 μ L per each sample. Use 20 μ L for up to 500,000 cells. The following steps are written for 1 sample of a 500,000 cells reaction. Do not adjust the volumes for lower cell numbers.
6. Into a 2 mL tube, transfer 20 μ L Concanavalin A bead slurry into 1.6 mL 1X Binding Buffer and mix by pipetting. Place the tube on a magnetic stand to clear (30 seconds to 2 minutes).
7. Withdraw the liquid completely and remove the tube from the magnetic stand. Add 1.5 mL 1X Binding Buffer, mix by pipetting, and collect any liquid from the sides or cap with a quick pulse spin in a microcentrifuge.

8. Place tube on magnetic stand to clear (30 seconds to 2 minutes), and remove liquid by pipette. Resuspend in 20 μL 1X Binding Buffer (20 μL per 500,000 cells sample) and keep the tube containing prepared Concanavalin A Beads at room temperature until the fresh cells are ready.
9. Slowly add the bead slurry from Step 8 and invert tube to mix. Place the tube containing the cells and Beads on an end-over-end rotator for 10 minutes.

Bind Primary Antibody (2 hours to overnight)

Note: Before using Antibody Buffer in Step 11 below, you must add Protease Inhibitor Cocktail and Digitonin. For every 1 mL of buffer to be used, add 10 μL of Protease Inhibitor Cocktail and 10 μL of 5% Digitonin. Keep on ice. Antibody Buffer can be kept at 4°C for up to 2 days once Protease Inhibitor Cocktail and Digitonin have been added. For each reaction, you will need 50 μL supplemented with 0.5 μL Protease Inhibitor Cocktail and 0.5 μL of 5% Digitonin.

10. After a quick spin (<100 x *g*) to remove liquid from the cap, place the tubes on a magnetic stand to clear (30 seconds to 2 minutes), and remove the liquid via a pipette.
11. Resuspend cells in a volume of ice-cold Antibody Buffer such that each sample is in a 50 μL volume, not to exceed 500,000 cells per 50 μL . Gently vortex. Place the tube on ice.
12. Add 1 μL (or at least 1 μg) undiluted primary antibody to each sample, and gently vortex or pipette to mix.

Note: Use a rabbit or mouse primary antibody. We recommend using a 1:50 to 1:100 dilution or the antibody manufacturer's recommendation for the dilution used in applications such as immunofluorescence. Histone H3K27me3 ([Catalog No. 39157](#)) is a common mark that can be a positive control.

13. Incubate 2 hours to overnight at 4°C with orbital mixing, make certain liquid stays together in the bottom of the tube so that it all mixes together.

Bind Secondary Antibody (60 minutes)

Note: Before using Dig-Wash Buffer in Step 15 below, you must add Protease Inhibitor Cocktail and Digitonin. For every 1 mL of buffer, add 10 μL Protease Inhibitor Cocktail and 10 μL of 5% Digitonin. 3.1 mL total are used per sample. Keep on ice. Dig-Wash Buffer can be kept at 4°C for up to 2 days once Protease Inhibitor Cocktail and Digitonin have been added.

Note: Choose the right secondary antibody for the reactions! If a mouse primary antibody was used, then use the Rabbit Anti-Mouse Secondary Antibody. If a rabbit primary antibody was used, then use the Guinea Pig Anti-Rabbit Secondary Antibody.

14. After a quick spin (<100 x g), place each tube on a magnetic stand to clear (30 seconds to 2 minutes), and then remove the liquid via a pipette.
15. Dilute the secondary antibody 1:100 in Dig-Wash Buffer. 100 µL of the diluted secondary are needed per reaction. Add 100 µL of this diluted secondary antibody to each sample while gently vortexing to dislodge the beads from the sides of the tube.
16. Place the tubes on an orbital rotator at room temperature for 60 minutes.
17. After a quick spin, place the tubes on a magnetic stand to clear (30 seconds to 2 minutes), then remove the liquid via a pipette.
18. Add 1 mL Dig-Wash Buffer to each reaction. Gently vortex or pipette to dislodge any beads that have aggregated together.
19. Repeat Step 18 twice, for a total of three washes.

Bind CUT&Tag-IT® Assembled pA-Tn5 Transposomes (60 minutes)

Note: Before using Dig-300 Buffer in Step 20 below, you must add Protease Inhibitor Cocktail and Digitonin. For every 1 mL of buffer, add 10 µL Protease Inhibitor Cocktail and 2 µL of 5% Digitonin. 3.1 mL total are used per sample. Keep on ice. Dig-300 Buffer can be kept at 4°C for up to 2 days once Protease Inhibitor Cocktail and Digitonin have been added.

20. Dilute the CUT&Tag-IT Assembled pA-Tn5 Transposomes with Dig-300 Buffer to a final concentration of 1:100. 100 µL will be needed per reaction. For example, for one reaction, add 1 µL of the CUT&Tag-IT™ Assembled pA-Tn5 Transposomes to 100 µL of Dig-300 Buffer.
21. After a quick spin (<100 x g), place each tube containing the immunoprecipitated samples on a magnetic stand to clear (30 seconds to 2 minutes), and then remove the liquid via a pipette.
22. Add 100 µL of the diluted CUT&Tag-IT™ Assembled pA-Tn5 Transposomes from Step 20 while gently pipetting to make sure it mixes together.
23. Incubate the reactions at room temperature on an orbital rotator for 60 minutes.
24. After a quick spin (<100 x g), place the tubes on a magnetic stand to clear (30 seconds to 2 minutes) and remove the liquid with a pipette.
25. Add 1 mL Dig-300 Buffer. Gently vortex or pipette to dislodge the beads.
26. Repeat Step 25 twice for a total of three washes.

Tagmentation (60 minutes)

Note: Before using Tagmentation Buffer, you must add Protease Inhibitor Cocktail and Digitonin. For every 1 mL of Tagmentation Buffer to be used, add 10 µL Protease Inhibitor Cocktail and 2 µL of 5% Digitonin.

27. After a quick spin (<100 x g), place the tubes on a magnetic stand (30 seconds to 2 minutes) and remove the liquid with a pipette.
28. Add 125 µL Tagmentation Buffer while gently vortexing or pipetting to mix each tube.
29. Incubate tubes at 37°C for 60 minutes.

DNA Extraction (60 minutes)

- 30.** To stop tagmentation and solubilize DNA fragments, add to each sample:
 - 4.2 μL 0.5 M EDTA
 - 1.25 μL 10% SDS
 - 1.1 μL Proteinase K (10 mg/mL)
- 31.** Mix by vortexing at full speed briefly for approximately two seconds, and incubate for 60 minutes at 55°C to digest.

Note: It is typical for the beads to form a large clump during incubation with Proteinase K and SDS due to the viscoelasticity of DNA. However, for abundant genome-wide epitopes, large-scale fragmentation of the genome will normally result in reduced clumping and release of beads into suspension, turning the liquid brownish relative to negative controls.

- 32.** After a quick spin (<100 x *g*), place the tubes on a magnetic stand (30 seconds to 2 minutes) and transfer the liquid with a pipette into a new 1.5 mL microcentrifuge tube.

Note: If the sample is too viscous, preventing the supernatant from being transferred, add 625 μL DNA Purification Binding Buffer, pipette up and down, place again on the magnetic stand and transfer to a new tube. Proceed to **step 34**.

- 33.** Add 625 μL DNA Purification Binding Buffer to each sample and mix by pipetting up and down. If the color indicator turns violet or orange, add 8 μL of 3 M Sodium Acetate.
- 34.** For each sample, place a labeled DNA Purification Column into a collection tube.
- 35.** Transfer each sample to its corresponding column, close the cap, and centrifuge at 17,000 x *g* (~14,000 rpm) for 1 minute.
- 36.** Discard the flow-through and return the collection tube to the column.

Note: 100% Ethanol must be added to the DNA Purification Wash Buffer before the first use, to a final concentration of 80% (40 mL of 100% Ethanol to the DNA Purification Wash Buffer bottle).

- 37.** Add 750 μL of DNA Purification Wash Buffer (prepared above) to the column and cap the column. Centrifuge at 17,000 x *g* for 1 minute.
- 38.** Discard the flow-through and return the collection tube to the column. Centrifuge the empty tube at 17,000 x *g* for 2 minutes to remove any remaining DNA Purification Wash Buffer.
- 39.** Transfer each column to a new microcentrifuge tube and add 35 μL of DNA Purification Elution Buffer to the center of the column matrix, cap the column, and incubate at room temperature for 1 minute.

40. Centrifuge at 17,000 x *g* for 1 minute. Discard the column. The DNA purification is complete, and is ready for next-gen library creation compatible with dual-index primers (Nextera) for Illumina® sequencing. The DNA can be stored at -20°C or you may proceed to PCR Amplification. You will need indexing primers, and high-fidelity DNA polymerase. For an example protocol please refer to our complete CUT&Tag-IT Assay Kit manuals (Catalog Nos. [53160](#) or [53165](#)).

References

1. Kaya-Okur, H.S., *et al.* (2019) *Nature comm.* 10:1930 (1).

Technical Services

If you need assistance at any time, please call or send an e-mail to Active Motif Technical Service at one of the locations listed below.

North America	Toll free: 877 222 9543 Direct: 760 431 1263 Fax: 760 431 1351 Email: tech_service@activemotif.com
Europe	Direct: +32 (0)2 653 0001 Fax: +32 (0)2 653 0050 Email: eurotech@activemotif.com
Japan	Direct: +81 (0)3 5225 3638 Fax: +81 (0)3 5261 8733 Email: japantech@activemotif.com
China	Direct: (86)-21-20926090 Cell Phone: 18521362870 Email: techchina@activemotif.com