

# Histone H3 phospho Ser10 ELISA

Catalog No. 53111

# Histone H3 phospho Ser28 ELISA

Catalog No. 53100

(version A)

## **Active Motif North America**

1914 Palomar Oaks Way, Suite 150

Carlsbad, California 92008, USA

Toll free: 877 222 9543

Telephone: 760 431 1263

Fax: 760 431 1351

## **Active Motif Europe**

104 Avenue Franklin Roosevelt

B-1330 Rixensart, Belgium

UK Free Phone: 0800 169 31 47

France Free Phone: 0800 90 99 79

Germany Free Phone: 0800 181 99 10

Telephone: +32 (0)2 653 0001

Fax: +32 (0)2 653 0050

## **Active Motif Japan**

Azuma Bldg, 7th Floor

2-21 Ageba-Cho, Shinjuku-Ku

Tokyo, 162-0824, Japan

Telephone: +81 3 5225 3638

Fax: +81 3 5261 8733

Copyright 2010 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.

© 2010 Active Motif, Inc., 1914 Palomar Oaks Way, Suite 150; Carlsbad, CA 92008. All rights reserved.

All trademarks, trade names, service marks or logos referenced herein belong to their respective companies.

| <b>TABLE OF CONTENTS</b>                                    | <b>Page</b> |
|---|-------------|
| <b>Overview</b> .....                                       | 1           |
| <b>Introduction</b> .....                                   | 2           |
| <b>Kit Performance and Benefits</b> .....                   | 3           |
| <b>Kit Components and Storage</b>                           |             |
| Additional Materials Required .....                         | 4           |
| <b>Protocols</b>  |             |
| Buffer Preparation and Recommendations .....                | 5           |
| Quick Chart for Preparing Buffers .....                     | 6           |
| ELISA Protocol .....  | 7           |
| <b>References</b> .....                                     | 9           |
| <b>Appendix</b>   |             |
| Section A. Preparation of Acid Extract/Crude Histones ..... | 10          |
| Section B. Troubleshooting Guide .....                      | 12          |
| Section C. Related Products .....                           | 13          |
| <b>Technical Services</b> .....                             | 16          |

## Overview

---

The addition or removal of modifications such as phospho-, methyl- and acetyl- functional groups to histones can have a profound effect on nuclear signaling as these dynamic modifications are critical in the regulation of transcription, chromosome packaging, DNA damage repair and functional genomics. Screening extracts for specific histone modifications is a simple way to assess cell health and the effect of treatment compounds on cell division.

Phosphorylation of serine 10 and serine 28 in the tail of histone H3 occur early in mitosis and serve as excellent mitotic markers. Histone H3 is phosphorylated on serine 10 during late S phase or G2 phase, while the phosphorylation of serine 28 occurs during prophase. In contrast to serine 10, serine 28 phosphorylation has never been observed in interphase.

The Histone H3 phospho Ser10 ELISA is a simple solution for screening purified core histones or extracts for levels of phosphorylated serine 10 on histone H3 in human and other (wide range predicted) systems. These kits are sandwich ELISAs that utilize a C-terminal Histone H3 antibody to capture histone H3 from your sample and a mouse monoclonal antibody phosphorylated at serine 10 on histone H3 for specific detection. A secondary antibody conjugated to horseradish peroxidase (HRP) and developing solutions provide a sensitive colorimetric readout that is easily quantified by spectrophotometry on a standard plate reader. The assay is performed in a convenient 96-stripwell plate, enabling you to simultaneously screen from 1 to 96 samples in a single experiment. It works with acid extracts from tissue or cell samples and is able to detect phospho-histone H3 (Ser10) in as little as 156 nanograms of acid extract. For added convenience, negative and positive control acid extract is also included.

The Histone H3 phospho Ser28 ELISA is a simple solution for screening purified core histones or extracts for levels of phosphorylated serine 28 on histone H3 in human, mouse, hamster and bovine systems. These kits are sandwich ELISAs that utilize a C-terminal Histone H3 antibody to capture histone H3 from your sample and a rat monoclonal antibody phosphorylated at serine 28 on histone H3 for specific detection. A secondary antibody conjugated to horseradish peroxidase (HRP) and developing solutions provide a sensitive colorimetric readout that is easily quantified by spectrophotometry on a standard plate reader. The assay is performed in a convenient 96-stripwell plate, enabling you to simultaneously screen from 1 to 96 samples in a single experiment. It works with acid extracts from tissue or cell samples and is able to detect phospho-histone H3 (Ser28) in as little as 156 nanograms of acid extract. For added convenience, negative and positive control acid extract is also included.

| product                        | format      | catalog no. |
|--------------------------------|-------------|-------------|
| Histone H3 phospho Ser10 ELISA | 1 x 96 rxns | 53111       |
| Histone H3 phospho Ser28 ELISA | 1 x 96 rxns | 53100       |

## Introduction

---

### Histone H3 phospho Serine 10 and Serine 28

The basic structural unit of chromatin is the nucleosome, which consists of 146 base pairs (bp) of DNA wrapped around a histone octamer. The histone octamer consists of two copies each of the core histone H2A-H2B dimers and a tetramer of H3-H4. A linker histone, histone H1, binds chromatin outside the nucleosome unit to regulate chromatin structure.

Histone modifications such as phosphorylation, acetylation, and methylation at specific amino acid residues on the histone tails that extend beyond the core nucleosome have been found to influence and regulate transcription, chromosome packaging, and DNA damage repair. Many of these specific histone modifications are conserved throughout eukaryotes. While the biological significance of some histone modifications remains to be understood, some have been demonstrated to correlate very closely with specific cellular states like mitosis.

Cell division is a complex, tightly regulated process that is marked by mitosis. Two significant mitotic events include microtubule formation and chromosome condensation. Phosphorylation of serine 10 and serine 28 in the tail of histone H3 occur early in mitosis when chromosomes begin to condense and during premature chromosome condensation induced in S-phase cells. Histone H3 is phosphorylated on serine 10 during late S phase or G2 phase, while the phosphorylation of serine 28 occurs during prophase<sup>1,2</sup>. In contrast to serine 10, serine 28 phosphorylation has never been observed in interphase<sup>1</sup>. Both phosphorylations occur during mitotic chromatin condensation before nuclear division occurs, which make phosphorylation of histone H3 at Ser10 or Ser28 important markers for cells undergoing mitosis.

Phosphorylated serine 28 on histone H3 has recently been shown to associate with destabilized nucleosomes in transcribed chromatin, making this an interesting indicator of both mitotic activity and transcriptional activation<sup>3</sup>.

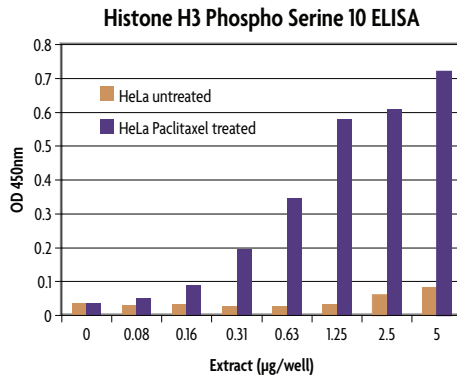
## Kit Performance and Benefits

**Detection limit:** < 5 µg acid extract/well.

**Range of detection:** Histone phospho Ser10 and Ser28 ELISAs provide quantitative results from 156 ng to 5 µg of histones isolated by acid extraction.

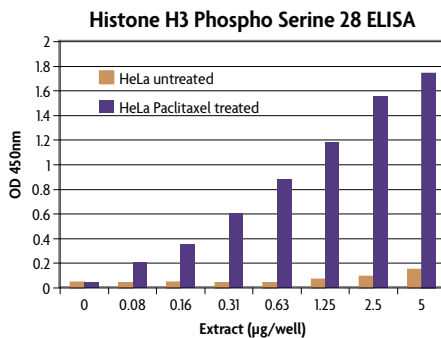
**Cross-reactivity:** Histone H3 phospho Ser10 Kits work with human and a wider range of species reactivity is predicted due to the high degree of sequence homology of histone H3. Histone H3 phospho Ser28 Kits work with human, mouse, hamster and bovine systems.

**Assay time:** 3.5 hours.



### Histone H3 phospho Ser 10 levels in paclitaxel and untreated HeLa acid extracts.

78 ng to 5 µg of Paclitaxel treated (Catalog No. 36201) and untreated HeLa acid extract (Catalog No. 36200), prepared as stated in Appendix A, were assayed in the Histone H3 phospho Ser10 ELISA. Data shown are the results from wells assayed in duplicate. These results are provided for demonstration only.



### Histone H3 phospho Ser 28 levels in paclitaxel and untreated HeLa acid extracts.

78 ng to 5 µg of Paclitaxel treated (Catalog No. 36201) and untreated HeLa acid extract (Catalog No. 36200), prepared as stated in Appendix A, were assayed in the Histone H3 phospho Ser28 ELISA. Data shown are the results from wells assayed in duplicate. These results are provided for demonstration only.

## Kit Components and Storage

---

Histone H3 phospho Ser10 Kits and Histone H3 phospho Ser28 Kits are for research use only. Not for use in diagnostic procedures. All components are guaranteed stable for 6 months from date of receipt when stored properly.

| Reagents  | Quantity                 | Storage    |
|---|--------------------------|------------|
| Histone H3 phospho Ser10 Mouse mAb, or<br>Anti-phospho Histone H3 Ser28 Rat mAb | 10 $\mu$ l<br>11 $\mu$ l | 4°C<br>4°C |
| HRP-conjugated anti-mouse IgG, or<br>HRP-conjugated anti-rat IgG                | 10 $\mu$ l<br>10 $\mu$ l | 4°C<br>4°C |
| Histone H3 Capture Plate  | 1                        | 4°C        |
| Plate sealer  | 1                        | 4°C        |
| Assay Dilution Buffer AM2   | 15 ml                    | 4°C        |
| 20X Wash Buffer   | 25 ml                    | 4°C        |
| Developing Solution   | 10 ml                    | 4°C        |
| Stop Solution   | 10 ml                    | 4°C        |
| Positive control HeLa acid extract<br>(Paclitaxel treated)                      | 25 $\mu$ g/50 $\mu$ l    | -80°C      |
| Negative control HeLa acid extract  | 25 $\mu$ g/50 $\mu$ l    | -80°C      |

### Additional materials required

- Multi-channel pipettor
- Multi-channel pipettor reservoirs
- Rocking platform
- Acid extracted samples or purified core histones
- Microplate spectrophotometer capable of reading at 450 nm (655 nm as optional reference wavelength)

## Protocols

---

### Buffer Preparation and Recommendations

#### Assay Dilution Buffer AM2

Assay Dilution Buffer AM2 is provided as a 1X solution and is ready for use once thawed.

#### 20X Wash Buffer

Prepare the amount of 1X Wash Buffer required for the assay as follows: For every 100 ml of 1X Wash Buffer required, dilute 5 ml 20X Wash Buffer with 95 ml distilled water (see Quick Chart for preparing buffers in this section). Mix gently to avoid foaming. The 1X Wash Buffer may be stored at 4°C for one week. The Tween 20 contained in the 20X Wash Buffer may form clumps, therefore homogenize the buffer by incubating at 50°C for 2 minutes and mixing prior to use.

#### Preparation of antibodies (see Quick Chart for preparing buffers in this section)

Dilute the Histone H3 phospho Ser 10 mouse monoclonal antibody 1:2000 with Assay Dilution Buffer AM2. Use 50 µl per well. Dilute the anti-mouse HRP-conjugated secondary 1:5000 with Assay Dilution Buffer AM2. Use 50 µl per well.

Dilute the Histone H3 phospho Ser28 rat monoclonal antibody 1:500 with Assay Dilution Buffer AM2. Use 50 µl per well. Dilute the anti-rat HRP-conjugated secondary 1:1000 with Assay Dilution Buffer AM2. Use 50 µl per well.

#### Developing Solution

100 µl are needed per well. The Developing Solution should be warmed to room temperature before use. The Developing Solution is light sensitive, therefore, we recommend avoiding direct exposure to intense light during storage. The Developing Solution may develop a yellow hue over time. This does not affect product performance. A blue color present in the Developing Solution indicates that it has been contaminated and must be discarded. Prior to use, place the Developing Solution at room temperature for at least 1 hour. Transfer the amount of Developing Solution required for the assay into a secondary container before aliquoting into the wells (see Quick Chart for preparing buffers in this section). After use, discard remaining Developing Solution.

#### Stop Solution

100 µl are needed per well. Prior to use, transfer the amount of Stop Solution required for the assay into a secondary container (see Quick Chart for preparing buffers in this section). After use, discard remaining Stop Solution.

**WARNING:** The Stop Solution is corrosive. Wear personal protective equipment when handling, *i.e.* safety glasses, gloves and labcoat.



## Diluting HeLa acid extracts

The HeLa acid extracts (paclitaxel treated and untreated) are both provided as controls for comparing levels of phosphorylated Serine 10 or Serine 28 on Histone H3. Sufficient extract is supplied for 1 control curve of each. The paclitaxel treated HeLa acid extract is optimized to give a strong signal when used at 156 ng to 5 µg/well. The untreated HeLa acid extract is provided as a negative control. Avoid multiple freeze/thaw cycles of the extract.

## Preparation of sample extracts

Histones isolated by acid extraction of tissue or cell samples or purified core histones can be used in the assay. Histone phospho Ser10 and Ser28 ELISAs provide quantitative results from 156 ng to 5 µg. It is recommended initially to use a range of sample concentrations to determine the amount of sample necessary to fall within the linear range of the assay. Once the appropriate amount of sample has been determined, perform the rest of the assays within the linear range. To prepare histones isolated by acid extraction, follow the protocol in Appendix Section A.

## Quick Chart for Preparing Buffers

| Reagents to prepare | Components                    | For 1 well      | For 1 strip (8 wells) | For 6 strips (48 wells) | For 12 strips (96 wells) |
|---------------------|-------------------------------|-----------------|-----------------------|-------------------------|--------------------------|
| Primary Antibody    | Histone H3 phospho Ser10 Ab   | 0.026 µl        | 0.23 µl               | 1.35 µl                 | 2.7 µl                   |
|                     | Assay Dilution Buffer         | 52 µl           | 450 µl                | 2.7 ml                  | 5.4 ml                   |
|                     | <b>TOTAL REQUIRED</b>         | <b>52 µl</b>    | <b>450 µl</b>         | <b>2.7 ml</b>           | <b>5.4 ml</b>            |
| OR                  |                               |                 |                       |                         |                          |
|                     | Histone H3 phospho Ser28 Ab   | 0.1 µl          | 0.9 µl                | 5.4 µl                  | 10.8 µl                  |
|                     | Assay Dilution Buffer         | 51.9 µl         | 450 µl                | 2.7 ml                  | 5.4 ml                   |
|                     | <b>TOTAL REQUIRED</b>         | <b>52 µl</b>    | <b>450 µl</b>         | <b>2.7 ml</b>           | <b>5.4 ml</b>            |
| Secondary Antibody  | HRP-conjugated anti-mouse IgG | 0.01 µl         | 0.09 µl               | 0.54 µl                 | 1.1 µl                   |
|                     | Assay Dilution Buffer         | 52 µl           | 450 µl                | 2.7 ml                  | 5.4 ml                   |
|                     | <b>TOTAL REQUIRED</b>         | <b>52 µl</b>    | <b>450 µl</b>         | <b>2.7 ml</b>           | <b>5.4 ml</b>            |
| OR                  |                               |                 |                       |                         |                          |
|                     | HRP-conjugated anti-rat IgG   | 0.05 µl         | 0.45 µl               | 2.7 µl                  | 5.4 µl                   |
|                     | Assay Dilution Buffer         | 52 µl           | 450 µl                | 2.7 ml                  | 5.4 ml                   |
|                     | <b>TOTAL REQUIRED</b>         | <b>52 µl</b>    | <b>450 µl</b>         | <b>2.7 ml</b>           | <b>5.4 ml</b>            |
| 1X Wash Buffer      | Distilled water               | 1.9 ml          | 17.1 ml               | 95 ml                   | 190 ml                   |
|                     | 20X Wash Buffer               | 100 µl          | 0.9 ml                | 5 ml                    | 10 ml                    |
|                     | <b>TOTAL REQUIRED</b>         | <b>2 ml</b>     | <b>18 ml</b>          | <b>100 ml</b>           | <b>200 ml</b>            |
| Developing Solution | <b>TOTAL REQUIRED</b>         | <b>112.5 µl</b> | <b>900 µl</b>         | <b>5.4 ml</b>           | <b>10.8 ml</b>           |
| Stop Solution       | <b>TOTAL REQUIRED</b>         | <b>112.5 µl</b> | <b>900 µl</b>         | <b>5.4 ml</b>           | <b>10.8 ml</b>           |

## ELISA Protocol

Determine the appropriate number of microwell strips required for testing samples, controls and blanks in duplicate. If less than 8 wells in a strip need to be used, cover the unused wells with a portion of the plate sealer while you perform the assay. The content of these wells is stable at room temperature if kept dry and, therefore, can be used later for a separate assay. Store the unused strips in the aluminum pouch at 4°C. Use the strip holder for the assay.

Prepare the 1X Wash Buffer as described above in the section Buffer Preparation and Recommendations. Multi-channel pipettor reservoirs may be used for dispensing the Wash Buffer, Assay Dilution Buffer, Developing Solution and Stop Solution into the wells being used.

### Control Curve Preparation (Optional)

If a control curve is not desired, skip this portion and proceed with Step 1 below.

|   | Untreated |          | Paclitaxel treated |          |   |   | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
|---|-----------|----------|--------------------|----------|---|---|---|---|---|---|---|----|----|----|
|   | 1         | 2        | 3                  | 4        |   |   |   |   |   |   |   |    |    |    |
| A | 5 µg      | 5 µg     | 5 µg               | 5 µg     | – | – | – | – | – | – | – | –  | –  | –  |
| B | 2.5 µg    | 2.5 µg   | 2.5 µg             | 2.5 µg   | – | – | – | – | – | – | – | –  | –  | –  |
| C | 1.25 µg   | 1.25 µg  | 1.25 µg            | 1.25 µg  | – | – | – | – | – | – | – | –  | –  | –  |
| D | 0.625 µg  | 0.625 µg | 0.625 µg           | 0.625 µg | – | – | – | – | – | – | – | –  | –  | –  |
| E | 312 ng    | 312 ng   | 312 ng             | 312 ng   | – | – | – | – | – | – | – | –  | –  | –  |
| F | 156 ng    | 156 ng   | 156 ng             | 156 ng   | – | – | – | – | – | – | – | –  | –  | –  |
| G | 78 ng     | 78 ng    | 78 ng              | 78 ng    | – | – | – | – | – | – | – | –  | –  | –  |
| H | 0 µg      | 0 µg     | 0 µg               | 0 µg     | – | – | – | – | – | – | – | –  | –  | –  |

1. Dilute the untreated and paclitaxel treated HeLa acid extracts to a starting concentration of 0.1 µg/µl. The control extracts provided with the kit contain 25 µg/50 µl (0.5 µg/µl) per vial. Using the reported concentration, add 200 µl Assay Dilution Buffer AM2 to the vial and pipette up and down to mix thoroughly. The result will be 25 µg HeLa acid extract in a total volume of 250 µl = 0.1 µg/µl.
2. Add 100 µl of the diluted untreated HeLa extracts to wells A1 and A2.
3. Add 100 µl of the diluted paclitaxel treated HeLa extracts to wells A3 and A4.
4. Add 50 µl of Assay Dilution Buffer to wells B1 through H4.
5. Perform a serial two-fold dilution of the extracts by transferring 50 µl of the extracts in row A to the wells in row B.

- Mix the contents of row B by pipetting up and down 3-5 times. Do not change pipette tips between well transfers.
- Transfer 50  $\mu$ l of the contents of row B to row C and mix, as previous.
- Continue this process until row G is reached.
- When row G is reached, discard 50  $\mu$ l of the well contents so that the final volume is 50  $\mu$ l.
- Row H will serve as the blank wells and receives no acid extracts.

## Step 1: Binding of H3 to the Capture Plate

- In duplicate, prepare the amount of desired sample. It is recommended to try a range of sample concentrations in order to determine the amount of sample necessary for optimal detection. Add desired amount of sample in 50  $\mu$ l volume to plate.
- Incubate plate containing the optional control curve and samples for 1 hour at room temperature with mild agitation (100 rpm on a rocking platform).
- After the incubation, wash the wells 3 times with 200  $\mu$ l of Wash Buffer.

## Step 2: Binding of Primary Antibody

- Histone H3 phospho Ser10:** Dilute the Histone H3 phospho Ser10 antibody 1:2000 in Assay Dilution Buffer AM2 and mix thoroughly.  
**Histone H3 phospho Ser28:** Dilute the anti-phospho-Histone H3 Ser28 antibody 1:500 in Assay Dilution Buffer AM2 and mix thoroughly.
- Add 50  $\mu$ l of diluted primary antibody to each well.
- Incubate at room temperature for 1 hour with agitation.
- After the incubation, wash the wells 3 times with 200  $\mu$ l of Wash Buffer.

## Step 3: Binding of Secondary Antibody

- Histone H3 phospho Ser10:** Dilute the HRP-conj. anti-mouse IgG 1:5000 in Assay Dilution Buffer AM2 and mix thoroughly.  
**Histone H3 phospho Ser28:** Dilute the HRP-conj. anti-rat IgG 1:1000 in Assay Dilution Buffer AM2 and mix thoroughly.
- Add 50  $\mu$ l of the diluted secondary antibody solution to each well.
- Incubate at room temperature for 1 hour with agitation.
- After the incubation, wash the wells 3 times with 200  $\mu$ l of wash buffer.

## Step 4: Colorimetric Reaction

12. Remove as much of the final wash as possible by blotting the plate on paper towels.
13. Add 100  $\mu$ l of room temperature Developing Solution to each well.
14. Incubate under low light conditions for several minutes while monitoring color development. Please read the Certificate of Analysis supplied with this kit for optimal development time associated with this lot number. Monitor the blue color development until the samples turn medium to dark blue. For the optional control curve, monitor the wells containing 5  $\mu$ g paclitaxel treated HeLa acid extract until they turn medium to dark blue.
15. Add 100  $\mu$ l of Stop Solution to all the wells. In the presence of the acid, the blue color turns yellow.
16. Read absorbance on a spectrophotometer within 5 minutes at 450nm with an optional reference wavelength of 655 nm. Blank the plate reader according to the manufacturer's instructions using the blank wells.

Reading the reference wavelength is optional. Most microtiter plate readers are equipped to perform dual wavelength analysis and with the appropriate software, will automatically subtract the reference wavelength absorbance from the test wavelength absorbance. If your plate reader does not have this capability, you may read the plate twice, once at 450 nm and once at 655 nm then manually subtract the 655 nm OD from the 450 nm OD values.

## References

---

1. Goto, H., *et al.* (1999) *J. Biol. Chem.* 274(36): 25543-25549
2. Hooser, A. V. *et al.* (1998) *J. of Cell Science* 111: 3497-3506
3. Sun, J-M., *et al* (2007) *Nucleic Acids Res.* 35(19): 6640-6647

## Appendix

---

### Section A. Preparation of Acid Extract/Crude histone proteins

This procedure can be used for a 150 mm plate that is 70% confluent. The yield is approximately 0.15 mg of nuclear proteins from  $9 \times 10^6$  cells.

1. Grow HeLa cells to 70% confluency in DMEM with 10% FBS.
2. Treat cells as desired.
3. Wash cells with 1X PBS and aspirate.
4. Add 3-5 ml of PBS per 150 mm plate.
5. Scrape cells from the plate and transfer to a 50 ml conical tube.
6. Pellet the cells by centrifugation in a pre-cooled 4°C rotor at 200 x g for 5-10 minutes.
7. Aspirate as much of the PBS as possible without disturbing the cell pellet.
8. Resuspend the cell pellet in 5 volumes of Lysis Buffer (see below).
9. Incubate on ice for 30 minutes and occasionally invert the tube to mix.
10. Centrifuge the lysate at 11,000 x g for 10 minutes at 4°C.
11. Collect the supernatant fraction containing acid soluble proteins, and discard the acid-insoluble pellet.
12. Immediately neutralize the acid extracted proteins by adding 2/5 the total volume of Neutralization Buffer (see below).
13. Quantify the protein concentration of your acid extraction using either gel electrophoresis or a Bradford Assay.

Gel electrophoresis is a more sensitive technique to determine histone concentration as histones are most effectively stained by Coomassie dye in a gel matrix. To determine the protein concentration run a BSA or histone standard curve on the gel.

A Bradford Assay can be used to determine total protein concentration, not just the concentration of crude histone proteins. A total protein determination, however, is sufficient for use in the Histone Phospho Serine ELISA Kits. The quantity of acid extract tested in the Histone Phospho Serine ELISA Kits are based on total protein determination values.

14. Immediately aliquot the extract in small volumes to avoid multiple freeze/thaws.
15. Store the protein at -80°C for long-term stability.

**Lysis Buffer:**

0.4 M HCl

**Neutralization Buffer:**

1 M Sodium Phosphate, dibasic, pH 12.5. Use 5M NaOH to adjust pH.

2.5 mM DTT

10 mM PMSF

**Note:** DTT and PMSF must be added immediately prior to use.

## Section B: Troubleshooting Guide

| Problem/question   | Possible cause                             | Recommendation  |
|--|--|---|
| No signal or weak signal                                 | Omission of key reagent                    | Check that all reagents have been added in all wells in the correct order   |
|  | Substrate or conjugate is no longer active | Test conjugate and substrate for activity by mixing a small aliquot of HRP and Developing Solution together   |
|  | Enzyme inhibitor present                   | Sodium azide will inhibit the peroxidase reaction. Follow our recommendations to prepare buffers  |
|  | Plate reader settings not optimal          | Verify the wavelength and filter settings in the plate reader   |
|  | Incorrect assay temperature                | Bring Developing Solution and Stop Solution to room temperature before using  |
|  | Inadequate volume of Developing Solution   | Check to make sure that correct volume is delivered by pipette  |
| High background in all wells                             | Developing time too long                   | Stop enzymatic reaction as soon as the positive wells turn medium-dark blue   |
|  | Concentration of antibodies is too high    | Increase antibody dilutions   |
|  | Inadequate washing                         | Ensure all wells are filled with Wash Buffer and follow washing recommendations   |
| Uneven color development                                 | Incomplete washing of wells                | Ensure all wells are filled with Wash Buffer and follow washing recommendations   |
|  | Well cross-contamination                   | Follow washing recommendations  |
| High background in sample wells                          | Too much sample per well                   | Decrease amount of sample per well. For acid extracts, dilute down to 1-2 µg/well   |
|  | Concentration of antibodies is too high    | Perform antibody titration to determine optimal working concentration. For Histone H3 phospho Ser10, start using 1:5000 for primary antibody and 1:5000 for the secondary antibody. For Histone H3 phospho Ser28, start using 1:1000 for primary antibody and 1:2000 for the secondary antibody. The sensitivity of the assay will be decreased |
| No signal or weak signal in sample wells                 | Not enough sample per well                 | For extracts, make sure you are using an acid extract by following the protocol recommended in Appendix Section A. Increase amount of acid extract to 10 µg/well  |
| No signal or weak signal in positive control curve wells | Incomplete mixing of control extracts      | Upon addition of Assay Dilution Buffer AM2 to the positive and negative control HeLa acid extract vials, pipette up and down to thoroughly mix the extract. Follow the protocol to generate an optional control curve   |

## Section C. Related Products

| Histone ELISAs                   | Format      | Catalog No. |
|----------------------------------|-------------|-------------|
| Histone H3 dimethyl Lys9 ELISA   | 1 x 96 rxns | 53108       |
| Histone H3 trimethyl Lys9 ELISA  | 1 x 96 rxns | 53109       |
| Histone H3 trimethyl Lys27 ELISA | 1 x 96 rxns | 53106       |
| Total Histone H3 ELISA           | 1 x 96 rxns | 53110       |

| Recombinant Methylated Histones         | Format | Catalog No. |
|---|--------|-------------|
| Recombinant Histone H2A                 | 50 µg  | 31251       |
| Recombinant Histone H3 (C110A)          | 50 µg  | 31207       |
| Recombinant Histone H3 monomethyl Lys4  | 50 µg  | 31208       |
| Recombinant Histone H3 dimethyl Lys4    | 50 µg  | 31209       |
| Recombinant Histone H3 trimethyl Lys4   | 50 µg  | 31210       |
| Recombinant Histone H3 monomethyl Lys9  | 50 µg  | 31211       |
| Recombinant Histone H3 dimethyl Lys9    | 50 µg  | 31212       |
| Recombinant Histone H3 trimethyl Lys9   | 50 µg  | 31213       |
| Recombinant Histone H3 monomethyl Lys27 | 50 µg  | 31214       |
| Recombinant Histone H3 dimethyl Lys27   | 50 µg  | 31215       |
| Recombinant Histone H3 trimethyl Lys27  | 50 µg  | 31216       |
| Recombinant Histone H3 monomethyl Lys36 | 50 µg  | 31217       |
| Recombinant Histone H3 dimethyl Lys36   | 50 µg  | 31218       |
| Recombinant Histone H3 trimethyl Lys36  | 50 µg  | 31219       |
| Recombinant Histone H3 monomethyl Lys79 | 50 µg  | 31220       |
| Recombinant Histone H3 dimethyl Lys79   | 50 µg  | 31221       |
| Recombinant Histone H3 trimethyl Lys79  | 50 µg  | 31222       |
| Recombinant Histone H4                  | 50 µg  | 31223       |
| Recombinant Histone H4 monomethyl Lys20 | 50 µg  | 31224       |
| Recombinant Histone H4 dimethyl Lys20   | 50 µg  | 31225       |
| Recombinant Histone H4 trimethyl Lys20  | 50 µg  | 31226       |

| Control Acid Extracts                       | Format | Catalog No. |
|---|--------|-------------|
| HeLa acid extract                           | 100 µg | 36200       |
| HeLa acid extract (Paclitaxel treated)      | 100 µg | 36201       |
| HeLa acid extract (Sodium Butyrate treated) | 100 µg | 36202       |
| HeLa acid extract (Etoposide treated)       | 100 µg | 36203       |
| HeLa acid extract (Anacardic acid treated)  | 100 µg | 36204       |

| DNA Methylation             | Format  | Catalog No. |
|-----------------------------|---------|-------------|
| MethylDetector™             | 50 rxns | 55001       |
| MethylCollector™            | 25 rxns | 55002       |
| MethylCollector™ Ultra      | 30 rxns | 55005       |
| UnMethylCollector™          | 30 rxns | 55004       |
| Fully Methylated Jurkat DNA | 10 µg   | 55003       |

| Histone Purification          | Format  | Catalog No. |
|-------------------------------|---------|-------------|
| Histone Purification Kit      | 10 rxns | 40025       |
| Histone Purification Mini Kit | 20 rxns | 40026       |



| <b>Chromatin Assembly</b>                                 | <b>Format</b> | <b>Catalog No.</b> |
|---|---------------|--------------------|
| Chromatin Assembly Kit                                    | 10 rxns       | 53500              |
| HeLa Core Histones  | 36 µg         | 53501              |
| <b>Histone Acetyltransferase and Deacetylase Activity</b> | <b>Format</b> | <b>Catalog No.</b> |
| HAT Assay Kit (Fluorescent)                               | 1 x 96 rxns   | 56100              |
| Recombinant p300 protein, catalytic domain                | 5 µg          | 31205              |
| HDAC Assay Kit (Fluorescent)                              | 1 x 96 rxns   | 56200              |
| HDAC Assay Kit (Colorimetric)                             | 1 x 96 rxns   | 56210              |
| <b>Co-Immunoprecipitation</b>                             | <b>Format</b> | <b>Catalog No.</b> |
| Nuclear Complex Co-IP Kit                                 | 50 rxns       | 54001              |
| Universal Magnetic Co-IP Kit                              | 25 rxns       | 54002              |
| <b>SUMOylation</b>  | <b>Format</b> | <b>Catalog No.</b> |
| SUMOLink™ SUMO-1 Kit                                      | 20 rxns       | 40120              |
| SUMOLink™ SUMO-2/3 Kit                                    | 20 rxns       | 40220              |
| <b>ChIP-IT™ Kits</b>                                      | <b>Format</b> | <b>Catalog No.</b> |
| ChIP-IT™ Express  | 25 rxns       | 53008              |
| ChIP-IT™ Express Enzymatic                                | 25 rxns       | 53009              |
| ChIP-IT™ Express HT                                       | 96 rxns       | 53018              |
| ChIP-IT™ Protein G Magnetic Beads                         | 25 rxns       | 53014              |
| Re-ChIP-IT™   | 25 rxns       | 53016              |
| ChIP-IT™  | 25 rxns       | 53001              |
| ChIP-IT™ w/o controls                                     | 25 rxns       | 53004              |
| ChIP-IT™ Shearing Kit                                     | 10 rxns       | 53002              |
| ChIP-IT™ Enzymatic  | 25 rxns       | 53006              |
| ChIP-IT™ Enzymatic w/o controls                           | 25 rxns       | 53007              |
| Enzymatic Shearing Kit                                    | 10 rxns       | 53005              |
| Salmon Sperm DNA/Protein G agarose                        | 25 rxns       | 53003              |
| ChIP-IT™ Control Kit – Human                              | 5 rxns        | 53010              |
| ChIP-IT™ Control Kit – Mouse                              | 5 rxns        | 53011              |
| ChIP-IT™ Control Kit – Rat                                | 5 rxns        | 53012              |
| Ready-to-ChIP HeLa Chromatin                              | 10 rxns       | 53015              |
| Ready-to-ChIP Hep G2 Chromatin                            | 10 rxns       | 53019              |
| Ready-to-ChIP K-562 Chromatin                             | 10 rxns       | 53020              |
| Ready-to-ChIP NIH/3T3 Chromatin                           | 10 rxns       | 53021              |

Active Motif also offers a growing list of application validated antibodies, including antibodies for histones and histone modifications, transcription factor antibodies, DNA methylation-related antibodies and ChIP validated antibodies. For a complete list go to [www.activemotif.com/abs](http://www.activemotif.com/abs)

| <b>Transcription Factor ELISAs</b> | <b>Format</b>      | <b>Catalog No.</b> |
|------------------------------------|--------------------|--------------------|
| TransAM™ AP-1 Family               | 2 x 96-well plates | 44296              |
| TransAM™ AP-1 c-Jun                | 1 x 96-well plate  | 46096              |
| TransAM™ GR                        | 1 x 96-well plate  | 45496              |
| TransAM™ HIF-1                     | 1 x 96-well plate  | 47096              |
| TransAM™ IRF-3 (Human)             | 1 x 96-well plate  | 48396              |
| TransAM™ IRF-7                     | 1 x 96-well plate  | 50196              |
| TransAM™ NFATc1                    | 1 x 96-well plate  | 40296              |
| TransAM™ NFκB Family               | 2 x 96-well plates | 43296              |
| TransAM™ NFκB p50                  | 1 x 96-well plate  | 41096              |
| TransAM™ NFκB p52                  | 1 x 96-well plate  | 48196              |
| TransAM™ NFκB p65                  | 1 x 96-well plate  | 40096              |
| TransAM™ p53                       | 1 x 96-well plate  | 41196              |
| TransAM™ STAT Family               | 2 x 96-well plates | 42296              |

For a complete list of the over 40 TransAM Kits available, please visit [www.activemotif.com/transam](http://www.activemotif.com/transam)

## Technical Services

---

If you need assistance at any time, please call Active Motif Technical Service at one of the numbers listed below.

### Active Motif North America

1914 Palomar Oaks Way, Suite 150

Carlsbad, CA 92008

USA

Toll Free: 877 222 9543

Telephone: 760 431 1263

Fax: 760 431 1351

E-mail: [tech\\_service@activemotif.com](mailto:tech_service@activemotif.com)

### Active Motif Europe

104 Avenue Franklin Roosevelt

B-1330 Rixensart, Belgium

UK Free Phone: 0800 169 31 47

France Free Phone: 0800 90 99 79

Germany Free Phone: 0800 181 99 10

Telephone: +32 (0)2 653 0001

Fax: +32 (0)2 653 0050

E-mail: [eurotech@activemotif.com](mailto:eurotech@activemotif.com)

### Active Motif Japan

Azuma Bldg, 7th Floor

2-21 Ageba-Cho, Shinjuku-Ku

Tokyo, 162-0824, Japan

Telephone: +81 3 5225 3638

Fax: +81 3 5261 8733

E-mail: [japantech@activemotif.com](mailto:japantech@activemotif.com)

Visit Active Motif on the worldwide web at <http://www.activemotif.com>

At this site:

- Read about who we are, where we are, and what we do
- Review data supporting our products and the latest updates
- Enter your name into our mailing list to receive our catalog, MotifVations newsletter and notification of our upcoming products
- Share your ideas and results with us
- View our job opportunities

**Don't forget to bookmark our site for easy reference!**