



HTRF® SET7/9 HISTONE H3K4 MONO-METHYLATION ASSAY (me0 → me1)

TECHNICAL NOTE

ABSTRACT SET7/9 Histone H3K4 mono-methylation assay measures the monomethylation of a biotinylated histone H3(1-21) peptide at lysine 4.

The HTRF SET7/9 Histone H3K4 monomethylation assay uses a H3(1-21) lysine 4 un-methylated biotinylated peptide (substrate), a Eu3+-cryptate labeled anti-H3K4 me1 detection antibody and XL665-conjugated Streptavidin (SA-XL665).

The assay is performed in a single well and run in two steps: the enzymatic step and the detection step. HTRF signal is proportional to the concentration of monomethylated H3(1-21) peptide. The assays within this technical note were performed in a 384-well plate in a 20 µL final volume.

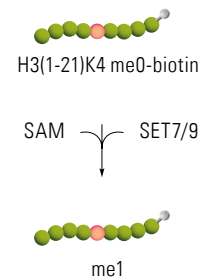
| | |
|--------------------|--|
| Enzyme | SET7/9 |
| Substrate | H3(1-21)K4 me0-biotin ARTKQTARKSTGGKAPRKQ- LA-GG-K(Biotin) |
| Detection Antibody | Anti-H3K4 me1-Eu(K) |

SET7/9 HISTONE H3K4 MONO-METHYLATION ASSAY AND REAGENTS

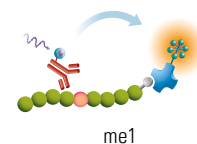
| | | |
|--|--|------------|
| H3K4 me1-Eu(K) Ab. | Cisbio Bioassays | # 61KA1KAE |
| Streptavidin XL-665 | Cisbio Bioassays | # 610SAXLA |
| Detection buffer | Cisbio Bioassays | # 62SDBRDD |
| SET7/9 | BPS Bioscience | # 51010 |
| Histone H3(1-21) lysine 4 un-methylated biotinylated peptide | AnaSpec | # 61702 |
| S-(5'-Adenosyl)-L-methionine chloride (SAM) | Sigma | # A7007 |
| Sinefungin | Sigma | # S8559 |
| Enzymatic buffer | 50 mM Tris-HCl, pH 8.8, 10 mM NaCl, 4 mM DTT, 4 mM MgCl ₂ , 0.01% Tween20 | |

Data shown on this application note has been obtained using Greiner # 784075, 384-well white microplates. For more information on the white plates we recommend, please visit <http://www.htrf.com/htrf-technology/microplate-recommendations>.

Enzymatic step



Detection step



ASSAY PROTOCOL

ENZYMATIC STEP

- Prepare working solutions of enzyme, peptide substrate, cofactors and inhibitor in enzymatic buffer just prior to use.
- Add to a 384-well small volume plate in the following order:
 - 4 µL of inhibitor (2.5X) or enzymatic buffer
 - 2 µL of SET7/9 enzyme (5X)
 - Incubate for 5 min at room temperature
 - 4 µL of H3(1-21)K4 me0-biotin peptide/ SAM pre-mixture (2.5X)
- Cover the plate with a plate sealer and incubate at room temperature.

DETECTION STEP

- Prepare detection mixture containing the anti-H3K4 me1-Eu(K) 2X according to the product datasheet recommended final concentration and SA-XL665 at 20 nM in detection buffer. Final concentration of 10 nM for SA-XL665 corresponds to 0.25X the final concentration of peptide substrate.
- Add 10 µL of detection mixture (2X) to the plate.
- Cover the plate with a plate sealer and incubate 1h at room temperature.
- Remove plate sealer and read fluorescence emission at 665nm and 620nm wavelengths on an HTRF compatible reader.

$$\text{HTRF Ratio} = (665\text{nm}/620\text{nm}) \times 10^4$$

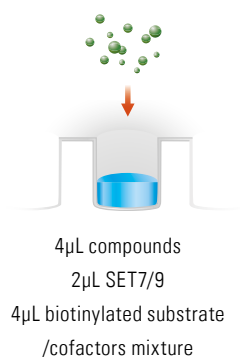
$$\text{Delta Ratio} = \text{Sample Ratio} - \text{Ratio negative}$$

$$\text{Delta F\%} = (\text{Delta Ratio}/\text{Ratio Negative}) \times 100$$

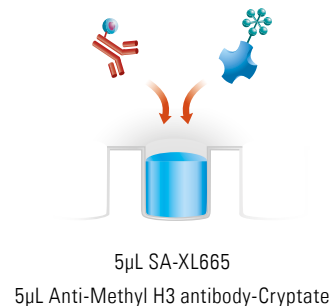
DISTRIBUTION: ENZYME INHIBITION STUDY

| | ENZYMATIC STEP | | | | DETECTION STEP | |
|-------------------------|------------------|-----------|--------|----------------------------|----------------|-----------|
| | ENZYMATIC BUFFER | INHIBITOR | SET7/9 | COFACTOR/SUBSTRATE MIXTURE | CRYPTATE-Ab | SA-XL 665 |
| SAMPLE | - | 4 µL | 2 µL | 4 µL | 5 µL | 5 µL |
| POSITIVE CONTROL | 4 µL | - | 2 µL | 4 µL | 5 µL | 5 µL |
| NEGATIVE CONTROL | 6 µL | - | - | 4 µL | 5 µL | 5 µL |

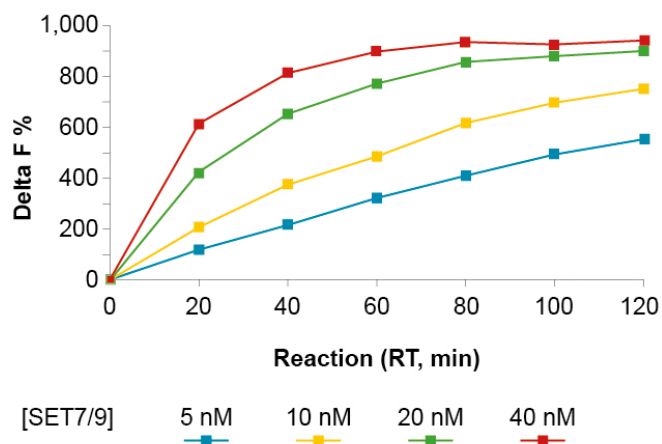
Enzymatic step



Detection step

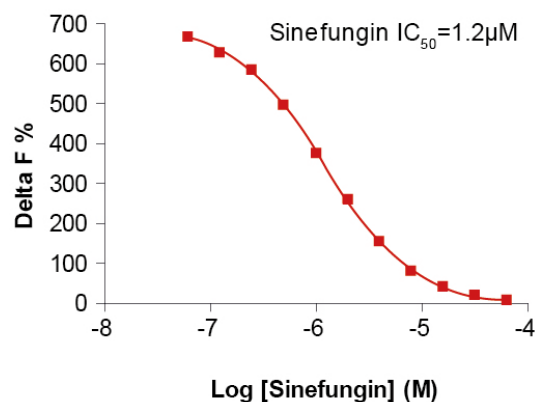


1. TIME COURSE AND ENZYME TITRATION



This step allows the optimal enzyme concentration and enzyme reaction time to be determined. Human recombinant SET7/9 was serially diluted to the concentrations indicated in the figure (5, 10, 20, 40 nM), and the assay was carried out with 40 nM biotinylated H3(1-21)me0 peptide substrate and 200 μ M SAM. Enzyme kinetics depends on the SET7/9 specific activity and substrate concentrations. The enzymatic reaction was carried out at RT and then stopped by adding H3K4me1-K Ab and SA-XL665 (detection reagents) after each time point (20, 40, 60, 80, 100, 120 min). A 60 min reaction time using 40 nM SET7/9 was selected for other experiments.

2. ENZYME INHIBITION



SET7/9 H3K4 monomethylation inhibitor assay was validated by measuring activity of the sinefungin inhibitor. This assay was performed using 200 nM SAM and 40 nM SET7/9. Serial dilutions of sinefungin were ranged from 6.1 nM to 100 μ M and pre-incubated for 5 min with SET7/9. Enzymatic reaction was initiated by the addition of 80 nM biotinylated H3(1-21) me0 peptide substrate plus 200 nM SAM. The enzyme reaction was stopped with the detection reagents after 60 min incubation at RT. IC_{50} value calculated from the inhibition curve was 1.2 μ M.

For more information, please visit us at www.htrf.com/epigenetic-toolbox-reagents

RELATED ARTICLES

EPIgeneous™ Methyltransferase assay: a new HTRF Universal, SAH detection assay to assess methyltransferase activity.

Roux T, Douayry N, Junique S, Sergeant L, Donsimoni G, Bourrier E, Trinquet E, LaRose R, Degorce F. - EpiCongress 2013, Boston, MA, USA.

High-Throughput, Homogeneous Histone Demethylase JARID1A, and JARID1C Enzymatic applications with HTRF Technology.

Adachi K, Tokuda C, Roux T, Trinquet E, Degorce F - Miptec 2013, Basel, Switzerland.

High-Throughput, Homogeneous Histone H3 Methyltransferase, (HMT) and Demethylase (HDM) Enzyme Assays using HTRF®, Technology: G9a H3K-27dimethylation assay example.

Roux T, Adachi K, Tokuda C, Verdi J, Junique S, Trinquet E, Gonzalez-Moya A, Degorce F - SLAS 2013, Orlando, USA.

High-Throughput, Homogeneous Histone H3 Methyltransferase (HMT) and Demethylase (HDM) Enzyme Assays using HTRF Technology.

Adachi K, Tokuda C, Chevallier F, Roux T, Gonzalez-Moya A, Degorce F. - Discovery on Target 2012, Boston, MA, USA.

Development of a panel of HTRF assay reagents for epigenetic targets.

Chevallier F, Jean A, Raynaldy D, Romier M, Servent F, Tokuda C, Adachi K. - Miptec 2011, Basel, Switzerland.

Development of G9a (Histone H3K9 methyltransferase) assay using HTRF technology.

Adachi K, Tokuda C, Chevallier F, Preaudat M. - SBS 2011, Orlando, USA.

FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567 China +86 10 8080 9288

Visit www.htrf.com to find a list of our regional distributors

