



# HTRF® G9a HISTONE H3K27 DI-METHYLATION ASSAY (me0 → me2)

## TECHNICAL NOTE

**ABSTRACT** G9a Histone H3K27 di-methylation assay measures the dimethylation of a biotinylated histone H3(21-44) peptide at lysine 27.

The HTRF G9a Histone H3K27 dimethylation assay uses a H3(21-44) lysine 27 un-methylated biotinylated peptide (substrate), a Eu3+-cryptate labeled anti-H3K27 me2 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 dimethylated H3(21-44) peptide. The assays within this technical note were performed in a 384-well plate in a 20 µL final volume.

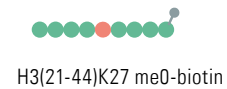
Enzyme	G9a
Substrate	H3(21-44)K27 me0-biotin ATKAARKSAPATGGVKK- PHRYRPG-GK(Biotin)
Detection Antibody	Anti-H3K27 me2-Eu(K)

## G9a HISTONE H3K27 DI-METHYLATION ASSAY AND REAGENTS

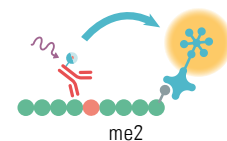
H3K27 me2-Eu(K) Ab.	Cisbio Bioassays	# 61KC2KAE
Streptavidin XL-665	Cisbio Bioassays	# 610SAXLA
Detection buffer	Cisbio Bioassays	# 62SDBRDD
G9a	BPS Bioscience	# 51000
Histone H3(21-44) lysine 27 un-methylated biotinylated peptide	AnaSpec	# 64440
S-(5'-Adenosyl)-L-methionine chloride (SAM)	Sigma	# A7007
Sinefungin	Sigma	# S8559
Enzymatic buffer	50 mM Tris-HCl, pH 8.8, 10 mM NaCl, 4 mM DTT, 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  $\mu$ L of inhibitor (2.5X) or enzymatic buffer
  - 2  $\mu$ L of G9a enzyme (5X)
  - Incubate for 5 min at room temperature
  - 4  $\mu$ L of H3(21-44)K27 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-H3K27 me2-Eu(K) 2X according to the product datasheet recommended final concentration and SA-XL665 at 50 nM in detection buffer. Final concentration of 25 nM for SA-XL665 corresponds to 0.25X the final concentration of peptide substrate.
- Add 10  $\mu$ L 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.

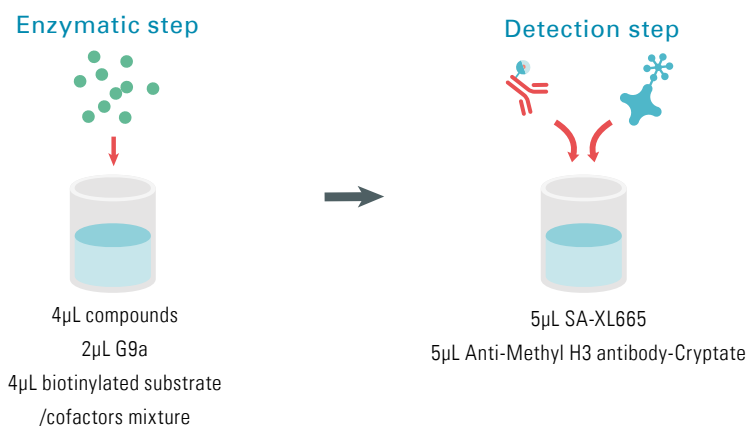
$$\text{HTRF Ratio} = (665\text{nm}/620\text{nm}) \times 10^4$$

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

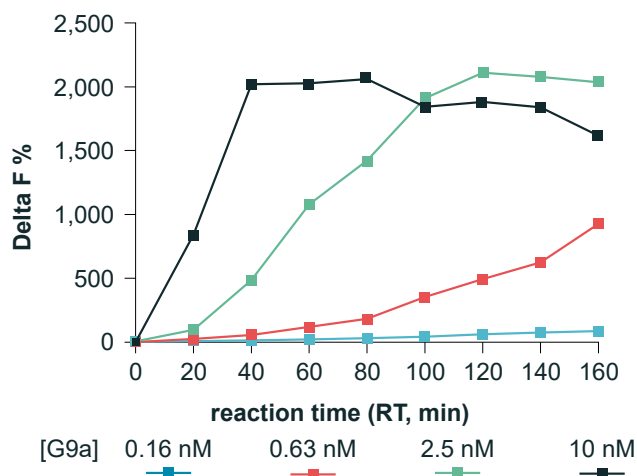
$$\text{Delta F\%} = (\text{Delta Ratio}/\text{Ratio Negative}) \times 100$$

### DISTRIBUTION: ENZYME INHIBITION STUDY

	ENZYMATIC STEP				DETECTION STEP	
	ENZYMATIC BUFFER	INHIBITOR	G9A	COFACTOR/SUBSTRATE MIXTURE	CRYPTATE-Ab	SA-XL 665
<b>SAMPLE</b>	-	4 $\mu$ L	2 $\mu$ L	4 $\mu$ L	5 $\mu$ L	5 $\mu$ L
<b>POSITIVE CONTROL</b>	4 $\mu$ L	-	2 $\mu$ L	4 $\mu$ L	5 $\mu$ L	5 $\mu$ L
<b>NEGATIVE CONTROL</b>	6 $\mu$ L	-	-	4 $\mu$ L	5 $\mu$ L	5 $\mu$ L

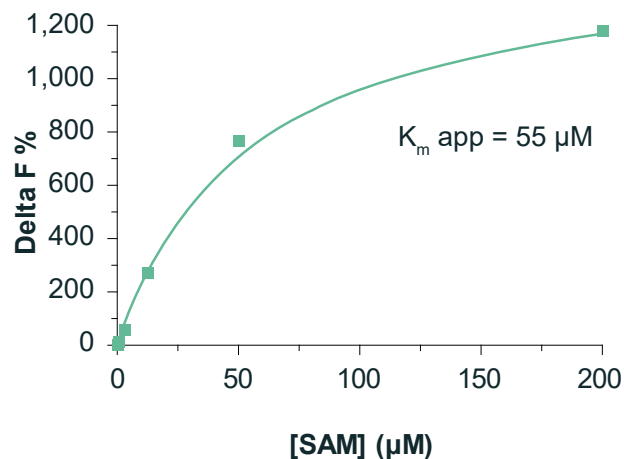


## 1. TIME COURSE AND ENZYME TITRATION



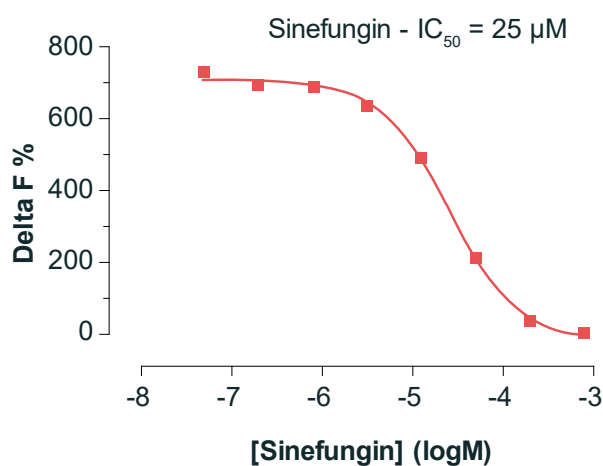
This step allows the optimal enzyme concentration and enzyme reaction time to be determined. Human recombinant G9a was serially diluted to the concentrations indicated in the figure (0.16, 0.63, 2.5, 10 nM) and the assay was carried out with 200 nM biotinylated H3(21-44)-me0 peptide substrate and 200  $\mu$ M SAM. Enzyme kinetics depends on the G9a specific activity and substrate concentrations. The enzymatic reaction was carried out at RT and then stopped by adding H3K27me2-K Ab and SA-XL665 (detection reagents) after each time point (20-160 min). For further experiments, a reaction time of 90min at RT and 2.5 nM enzyme were selected.

## 2. SAM TITRATION



This step enables the determination of  $K_m$  for SAM. The  $K_m$  value was determined with 2.5 nM G9a and 200 nM biotinylated H3 (21-44)me0 substrate in the enzymatic step. We recommend testing SAM concentrations ranging from 400  $\mu$ M to 3.1 nM (serial dilutions). The enzyme reaction was stopped at the optimal incubation period (RT, 90 min) by adding the detection reagents. The 55  $\mu$ M  $K_m$  value for SAM was determined from this experiment using a Michaelis-Menten plot.

## 3. ENZYME INHIBITION



G9a H3K27 dimethylation inhibitor assay was validated by measuring the activity of sinefungin inhibitor. This assay was performed using 100  $\mu$ M SAM and 2.5 nM G9a. Serial dilutions of sinefungin ranged from 13 nM to 1,000  $\mu$ M and were pre-incubated for 5 min with G9a. Enzymatic reaction was initiated by the addition of 200 nM biotinylated H3 (21-44) me0 peptide substrate plus 100  $\mu$ M SAM. The enzyme reaction was stopped with the detection reagents after 90 min incubation at RT.  $IC_{50}$  value calculated from the inhibition curve was 25  $\mu$ M.

For more information, please visit us at [www.htrf.com/epigenetic-toolbox-reagents](http://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 Japan +81-(0)43-306-8712

Visit [www.cisbio.com](http://www.cisbio.com) to find a list of our regional distributors



cisbio