



# HTRF® EZH2(Y641F) HISTONE H3K27 TRI-METHYLATION ASSAY (me0 → me3)

## TECHNICAL NOTE

**ABSTRACT** EZH2(Y641F) Histone H3K27 tri-methylation assay measures the trimethylation of a biotinylated histone H3(1-50) peptide at lysine 27.

The HTRF EZH2(Y641F) Histone H3K27 trimethylation assay uses a H3(1-50) lysine 27 un-methylated biotinylated peptide (substrate), a Eu<sup>3+</sup>-cryptate labeled anti-H3K27 me<sub>3</sub> 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 trimethylated H3(1-50) peptide. The assays within this technical note were performed in a 384-well plate in a 20 µL final volume.

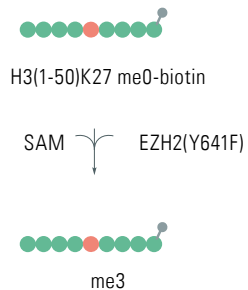
Enzyme	EZH2(Y641F)
Substrate	H3(1-50)K27 me <sub>0</sub> -biotin ARTKQTARKSTGG- KAPRKQLATKAARKSA- PATGGVKKPHRYRPGTVAL- REGG-K(Biotin)
Detection Antibody	Anti-H3K27 me <sub>3</sub> -Eu(K)

## EZH2(Y641F) HISTONE H3K27 TRI-METHYLATION ASSAY AND REAGENTS

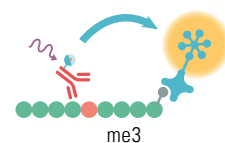
H3K27 me <sub>3</sub> -Eu(K) Ab.	Cisbio Bioassays	# 61KC3KAE
Streptavidin XL-665	Cisbio Bioassays	# 610SAXLA
Detection buffer	Cisbio Bioassays	# 62SDBRDD
EZH2(Y641F)	BPS Bioscience	# 51017
Histone H3(1-50) lysine 27 un-methylated biotinylated peptide	AnaSpec	# 65366
EZH2 complex	BPS Bioscience	# 51004
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 <sub>2</sub> , 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 EZH2(Y641F) enzyme (5X)
  - Incubate for 5 min at room temperature
  - 4  $\mu$ L of H3(1-50)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 me3-Eu(K) 2X according to the product datasheet recommended final concentration and SA-XL665 at 100 nM in detection buffer. Final concentration of 50 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.

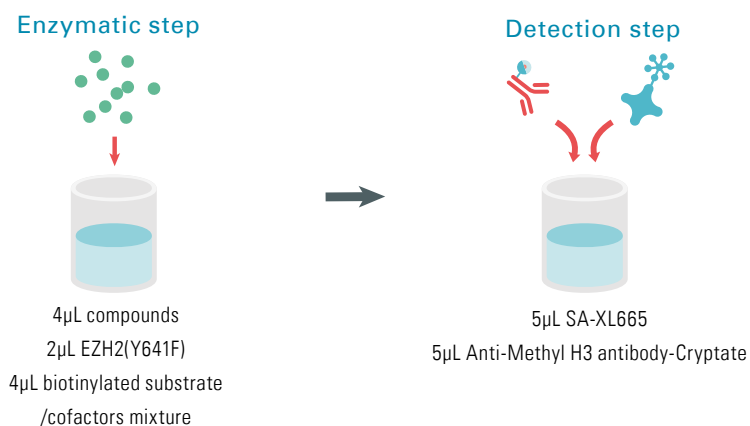
$$\text{HTRF Ratio} = (665\text{nm}/620\text{nm}) \times 10^4$$

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

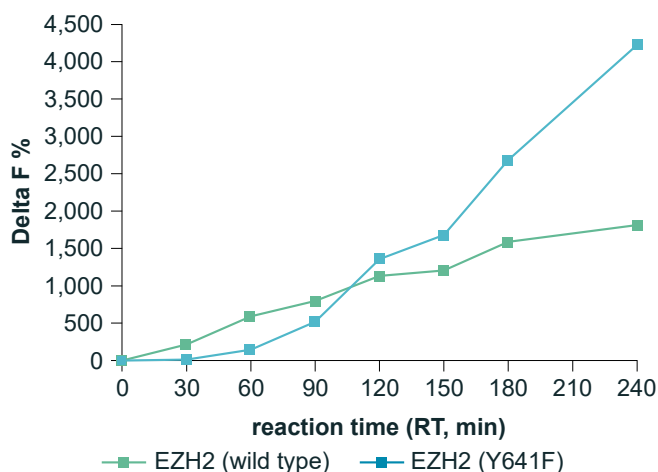
$$\text{Delta F\%} = (\text{Delta Ratio}/\text{Ratio Negative}) \times 100$$

### DISTRIBUTION: ENZYME INHIBITION STUDY

	ENZYMATIC STEP				DETECTION STEP	
	ENZYMATIC BUFFER	INHIBITOR	EZH2(Y641F)	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

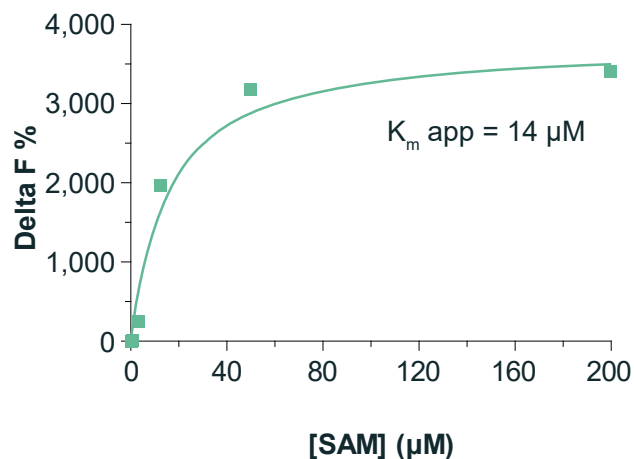


## 1. TIME COURSE



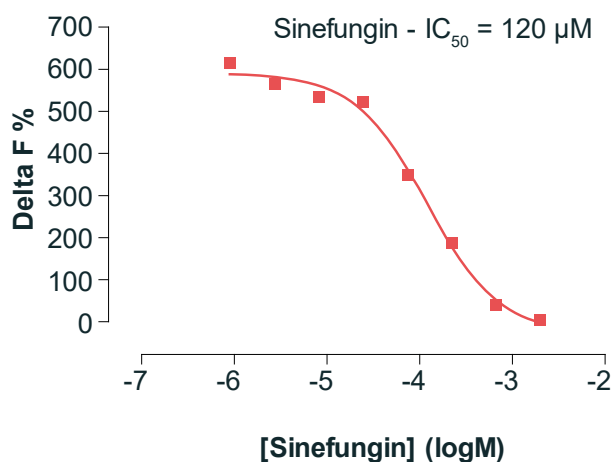
In this experiment, time course reaction of two types of human recombinant EZH2 complex (wild type and Y641 mutant) were compared at 30 °C. Both complex enzymes were added at 50 ng/well. This assay was carried out with 400 nM biotinylated H3(1-50)me0 peptide substrate and 200  $\mu$ M SAM and the reaction was then stopped by adding H3K27me3-K Ab and SA-XL665 (detection reagents) after each time point (30, 60, 90, 120, 150, 180, 210, 240 min). For further experiments, a reaction time of 180 min at RT and 40 ng/well enzyme complex were selected.

## 2. SAM TITRATION



This step enables the determination of  $K_m$  for SAM. The  $K_m$  value was determined with 40 ng/well EZH2(Y641) complex and 400 nM biotinylated H3(1-50)me0 substrate in the enzymatic step. We recommend testing SAM concentrations ranging from 200  $\mu$ M to 0.195  $\mu$ M (serial dilutions). The enzyme reaction was stopped at the optimal incubation period (RT, 180 min) by adding the detection reagents. The 14  $\mu$ M  $K_m$  value for SAM was determined from this experiment using a Michaelis-Menten plot.

## 3. ENZYME INHIBITION



EZH2 H3K27 trimethylation inhibitor assay was validated by measuring the activity of sinefungin inhibitor. This assay was performed using 15  $\mu$ M SAM and 40 ng/well EZH2(Y641) complex. Serial dilutions of sinefungin ranged from 1  $\mu$ M to 2 mM and were pre-incubated for 5 min with EZH2(Y641) complex. Enzymatic reaction was initiated by the addition of 400 nM biotinylated H3 (1-50) peptide substrate plus 15  $\mu$ M SAM. The enzyme reaction was stopped with the detection reagents after 180 min incubation at RT.  $IC_{50}$  value calculated from the inhibition curve was 120  $\mu$ 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