



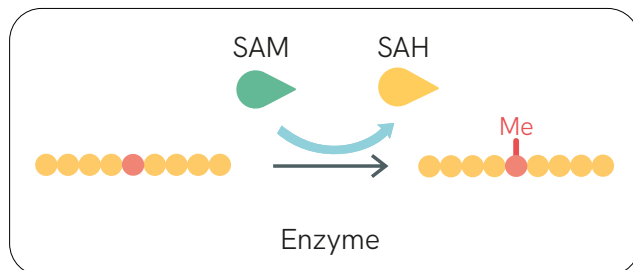
# Adapt to anything with a truly universal assay.

## EPIgeneous™ Methyltransferase Assay

Imagine being able to use the same biochemical assay for every methyltransferase, reducing assay development time and costs. That's the simplicity and flexibility you get with the new EPIgeneous Methyltransferase Assay from Revvity.

Built around our proprietary HTRF® (Homogeneous Time-Resolved Fluorescence) technology, this fast, simple, mix-and-read assay offers:

- A universal solution for all methyltransferases and substrates (peptides, nucleosomes, full histones, DNA, RNA, p53 and small molecules)
- Exceptional sensitivity without radioactivity
- Easy miniaturization for high throughput screening and profiling
- Eliminate false positives with direct detection of SAH (s-adenosylhomocysteine) from SAM (s-adenosylmethionine) conversion
- Enzymatic flexibility with a wide range of SAM concentrations (0.4  $\mu$ M – 200  $\mu$ M)
- Comprehensive ready-to-use kit

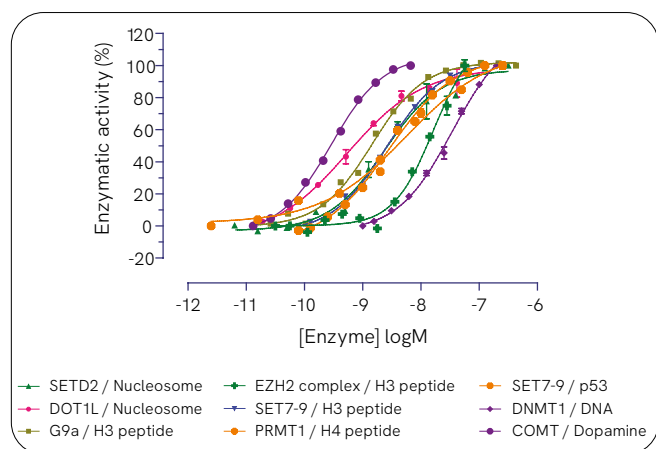


## Assay principal

Methyltransferase activity is assessed through SAM to SAH conversion. Direct measurement of SAH is done with an anti SAH antibody.

## Universally excellent results

No need for site-specific antibodies to detect substrate modifications, this truly universal assay delivers exceptional results for all enzymes within the methyltransferase families that produce s-adenosylhomocysteine (SAH).



EPIgenetic Methyltransferase Assay has been validated with a large set of enzymes and different types of substrates. Enzymatic reactions were carried out at RT or 30 °C for 1 to 2h with 0.5 to 10 $\mu$ M of SAM and related substrates according to the enzymes tested. Reactions were stopped with the addition of detection reagents in 384 low volume plate (20 $\mu$ L) and then read on a BMG LABTECH PHERAstar FS (flash lamp) after 1h incubation.

## Ordering information

Enzymatic assay conditions compatible with 0.4 $\mu$ M-200 $\mu$ M SAM at a 10% conversion of SAM

Product	Tests	Part #
EPIgenetic Methyltransferase Kit	1,000 tests	62SAHPEB
(for assays with < 20 $\mu$ M SAM)	10,000 tests	62SAHPEH

## If assay requires 20 $\mu$ M - 200 $\mu$ M of SAM also order

Ultrapure SAM Cofactor (2.4 $\mu$ mol)	1,000 tests	62SAHZLC
Ultrapure SAM Cofactor (22 $\mu$ mol)	10,000 tests	62SAHZLD
SAH-d2 Conjugate	1,000 tests	62SAHDLA
	10,000 tests	62SAHDLB

## Additional reagents available

SAH Calibrator		62SAHCLB
EPIgenetic Detection Buffer One (24mL)		62SAHRDD
EPIgenetic Detection Buffer Two (130mL)		62SAHFDD

## Interaction is everything

From the chemistry of our assays to the collaborative nature of our relationship with customers, Revvity is founded on the belief that "interaction is everything". Talk to us. And discover a partner dedicated to providing exactly what you need:

- Convenient ready-to-use kits
- Custom assay capabilities
- Expert technical support
- Complete confidentiality
- Trusted and proven HTRF technology

revvity