

# ADAPT TO ANYTHING WITH A TRULY UNIVERSAL ASSAY

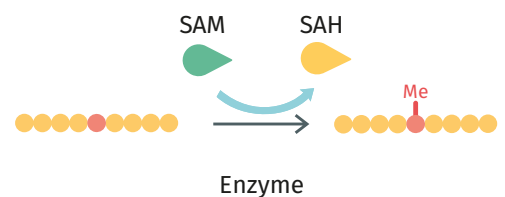
## EPIgeneous™ Methyltransferase Assay



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 Cisbio.

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

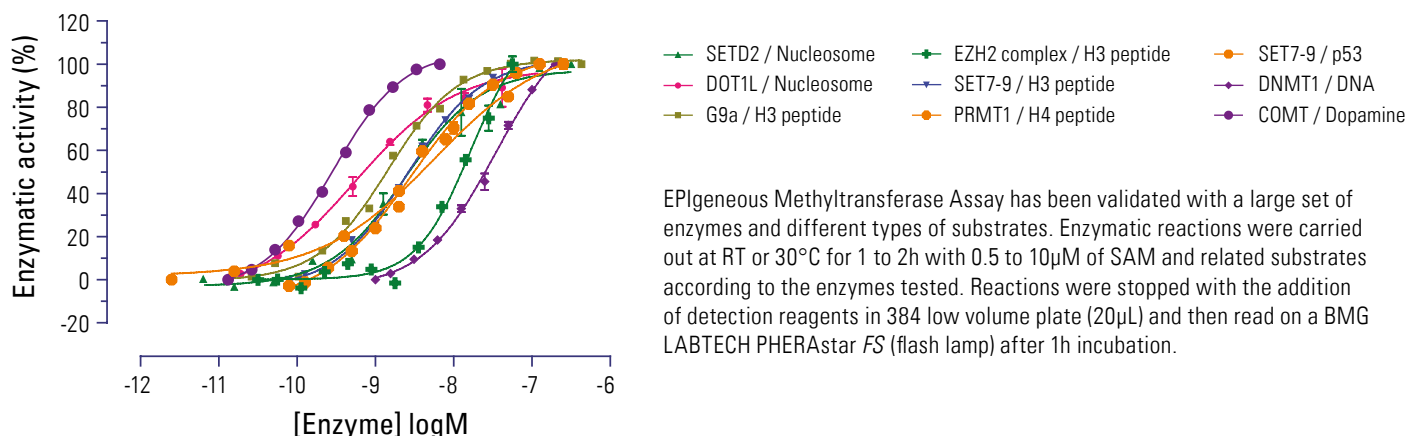


### ASSAY PRINCIPAL

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).



### ORDERING INFORMATION:

Enzymatic assay conditions compatible with 0.4µM-200µM SAM at a 10% conversion of SAM

#### PRODUCT

#### TESTS

#### PART#

EPIgeneous Methyltransferase Kit	1,000 tests	62SAHPEB
<i>(for assays with &lt; 20µM SAM)</i>	10,000 tests	62SAHPEH

If assay requires 20µM - 200µM of SAM also order:

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

Additional reagents available:

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

For more information, please visit us at [www.htrf.com/methyltransferaseassay](http://www.htrf.com/methyltransferaseassay)

## INTERACTION IS EVERYTHING

From the chemistry of our assays to the collaborative nature of our relationship with customers, Cisbio 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
- Complete confidentiality
- Custom assay capabilities
- Trusted and proven HTRF technology
- Expert technical support

#### 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