

BUILD A BETTER EPIGENETIC ASSAY



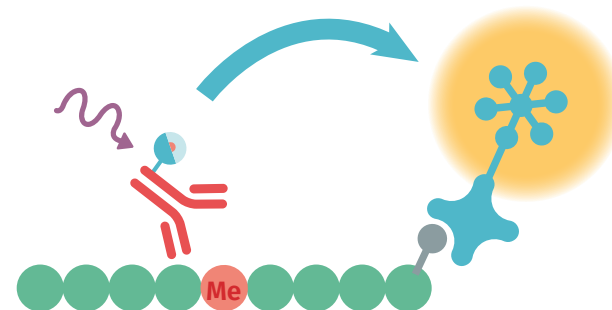
EPiGeneous™ Toolbox Reagents

Build your own robust HTRF® assay using PerkinElmer's anti-histone cryptate labeled antibodies. These antibodies provide assay specificity and flexibility to support your epigenetic screening.

Assay specificity is derived from the combination of the chosen Eu^{3+} cryptate labeled anti-histone antibody as well as the biotinylated peptide substrate.

Built around our proprietary HTRF (Homogeneous Time-Resolved Fluorescence) technology, this tool box offers:

- An extensive line of labeled antibodies
- The highest levels of specificity
- Validated assays for key histone methyltransferases (HMTs), demethylases (HDMs) and deacetylases (HDACs)
- Exceptional sensitivity without washes and radioactivity



Epigenetic Modification Profiling and Screening

HTRF signal is proportional to the concentration of methyl- or unacetyl specific peptide (e.g. unmethylated and unacetylated, mono-, di- or tri-methylated peptide).

An Array of Eu3+- Cryptate Labeled Anti-Histone H3 Antibodies

Each antibody is specific for a single H3 lysine residue and methylation or acetylation state. These antibodies used in conjunction with a biotinylated peptide substrate and XL-665 conjugated Streptavidin create custom, high specificity histone methyltransferases (HMTs), demethylases (HDMs) or deacetylases (HDACs) assay.

Methylation & Acetylation Level	K4	K9	K27	K36
Me0 & Ac0	.	.		
Me1
Me2
Me3	.		.	

Easy to Implement Validated Assays

PerkinElmer scientists have developed assays for a number of histone methyltransferases, demethylases and deacetylases. For histone methyltransferases and demethylases: EZH2 (H3K27Me3), SET7/9 (H3K4Me1), G9a (H3K9Me1), G9a (H3K9Me2), G9a (H3K27Me1), G9a (H3K27Me2), MLL1(H3K4Me2), MLL1 (HK4Me3), LSD1 (H3K4Me0), JMJD2A (H3K36Me2), JMJD2C (H3K9Me2), JHDM1A (H3K36Me1) and JMJD1A (H3K9Me0). For histone deacetylases: HDAC1 (H3K9Me0), HDAC2 (H3K4Me0 & H3K9Me0), HDAC6 (H3K9Me0), HDAC8 (H3K9Me0), SIRT1 (H3K9Me0), SIRT2 (H3K9Me0), SIRT3 (H3K9Me0) and SIRT6 (H3K9Me0).

Our list of validated assays is always expanding. Please visit perkinelmer.com for the most up-to-date list of assays.

Ordering Information

Anti-Histone H3 antibodies labeled with Cryptate

Methylation & Acetylation Level	K4	K9	K27	K36
Me0 / Ac0	61KA0KAE*	61KB0KAE		
	61KA0KAD**	61KB0KAD		
Me1	61KA1KAE*	61KB1KAE*	61KC1KAE*	61KD1KAE*
	61KA1KAD**	61KB1KAD**	61KC1KAD**	61KD1KAD**
Me2	61KA2KAE*	61KB2KAE*	61KC2KAE*	61KD2KAE*
	61KA2KAD**	61KB2KAD**	61KC2KAD**	61KD2KAD**
Me3	61KA3KAE*		61KC3KAE*	
	61KA3KAD**		61KCC3KAD**	

*500 tests, **10,000 tests based on 20µL assay volume

	5,000 tests	20,000 tests
Streptavidin-XL665	610SAXLA	610SAXLB