



# Build a better epigenetic assay.

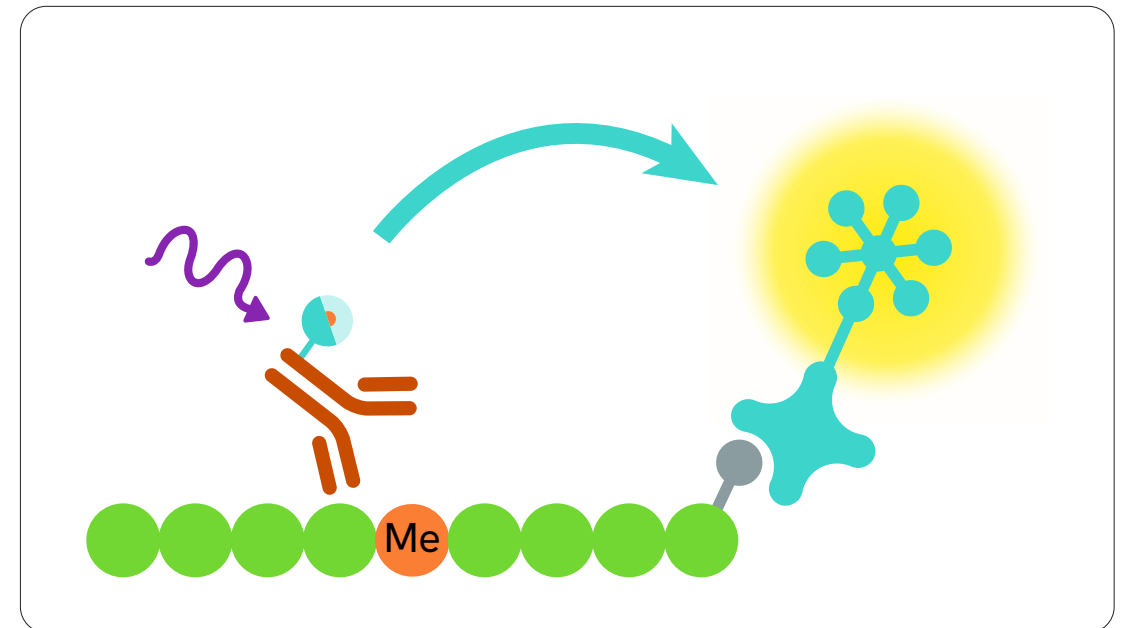
## EPIgeneous™ Toolbox reagents

Build your own robust HTRF® assay using Revvity'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 Eu3+-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 and acetylation level	K4	K9	K27	K36
Me0 and Ac0	▪	▪		
Me1	▪	▪	▪	▪
Me2	▪	▪	▪	▪
Me3	▪		▪	

## Easy to implement validated assays

Revvity 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 and H3K9Me0), HDAC6 (H3K9Me0), HDAC8 (H3K9Me0), SIRT1 (H3K9Me0), SIRT2 (H3K9Me0), SIRT3 (H3K9Me0) and SIRT6 (H3K9Me0).

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

## Ordering information

### Anti-histone H3 antibodies labeled with cryptate

Methylation and 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

