

# EPIgeneous™ biochemical and cellular assays



**DISCOVER AN EPIGENETICS ASSAY  
PORTFOLIO YOU'LL NEVER OUTGROW**



cisbio

## UNLOCKING THE POTENTIAL OF EPIGENETICS

Epigenetic mechanisms have been strongly linked to a broad array of illnesses ranging from cancer to mental disorders. The need to decipher key epigenetic interactions

and modifications has never been greater. Let Cisbio help you target more effectively and research more successfully with a broad portfolio of EPIgeneous assays and reagents offering uncompromising flexibility and sensitivity.

### HOW CAN WE HELP?

At Cisbio, our assay experience, expertise and product line spans every epigenetics-related application. Talk to us about your specific needs.

- Primary screening
- Secondary screening
- Target identification
- Target validation
- Inhibitor analysis
- SAR



### FEATURES AND BENEFITS

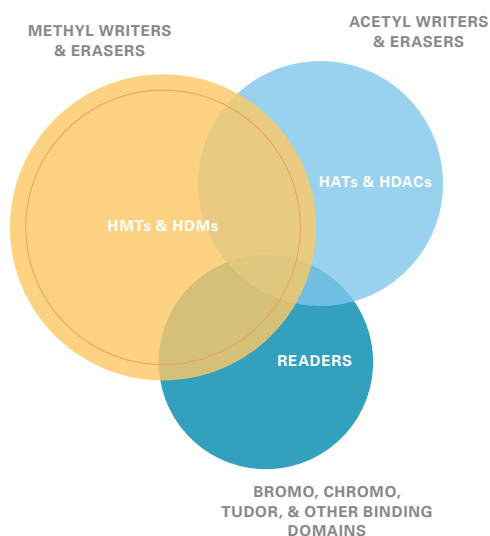
#### HTRF epigenetics assays

- Robust HTRF technology delivers reliable, reproducible results
- Multiple assay formats (universal, semi-universal, specific and cell-based) offer greater flexibility and applicability
- Assay-specific materials—including posters, technical notes, results, validated enzymes and binding domains—available online for added simplicity and confidence
- Compatible with existing HTRF plate readers for easy, economical adoption
- Homogeneous assays with no wash steps save time and money
- Easily miniaturized to take advantage of HTS cost efficiencies

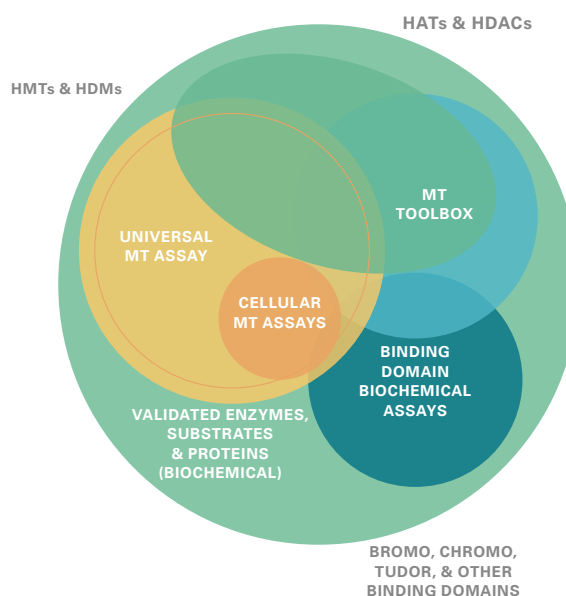
# INVESTIGATE MORE FREELY WITH A BROAD PORTFOLIO OF EPIGENETICS ASSAYS.

Whether your research concerns readers, writers, or erasers, you'll find the ideal solution with Cisbio's EPIgeneous product line. Our full portfolio features validated HTRF assay kits and toolbox reagents for reliable high throughput screening and investigation of epigenetic targets. We also provide custom labeling and assay development services that create new and innovative epigenetic screening approaches including cellular screening, full histone assays, protein:protein interaction assays and more.

## TARGETS



## SOLUTIONS



With assays for Histone Deacetylases (HDACs), Methyltransferases, Histone Demethylases, Histone Kinases, Bromodomains, Chromodomains, Tudor Domains and MBT Domains, Cisbio offers the ideal solution for anyone working with today's most popular targets in Epigenetics.



**ANSWERS ARE JUST A CLICK AWAY** For more information or help selecting the ideal solution for your specific needs, please visit [www.cisbio.com/epigenetic-screening](http://www.cisbio.com/epigenetic-screening)

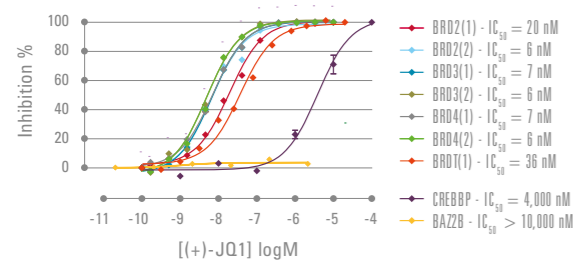
# AN EPIGENETICS ASSAY PORTFOLIO THAT EVOLVES WITH YOUR NEEDS.

For research to continually evolve and innovate, research tools must do the same. Which is why our EPIgeneous product line is constantly evolving and expanding, meeting the needs of today, and fueling the discoveries of tomorrow.

## BINDING DOMAIN ASSAYS

Cisbio simplifies the study of epigenetics readers with three assay kits already validated and optimized for a broad selection of 28 key binding domains. A fourth kit is also available to help select the best possible solution for non-validated reader domains.

PRODUCT LIST: Binding Domain Kit A Binding Domain Kit B Binding Domain Kit C Binding Domain Discovery Kit



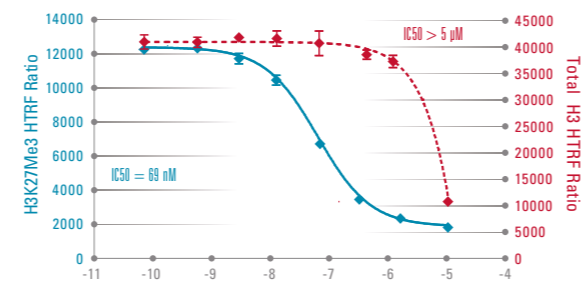
JQ1 profiling using EPIgeneous Binding Domain Kits (A & B) showing clear, reproducible results across multiple bromodomains.



## EPIGENETIC CELLULAR ASSAYS

Cisbio's specialized selection of cellular kits gives you a better, faster, more robust and reliable way to perform inhibitor studies.

PRODUCT LIST: H3K4Me2 H3K27Me3 H3K36Me2 Total H3 Phospho-Histone H3 (Ser10)

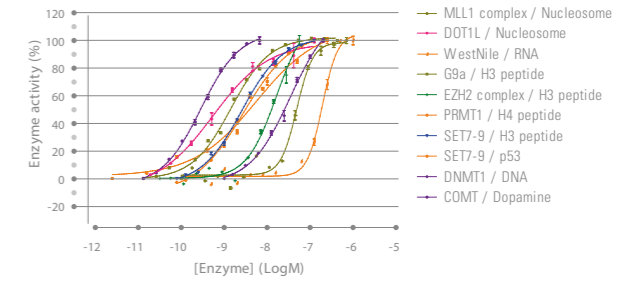


GSK126 compound titration using Cisbio's H3K27Me3 kit along with a normalization tool for precise total H3 measurement.

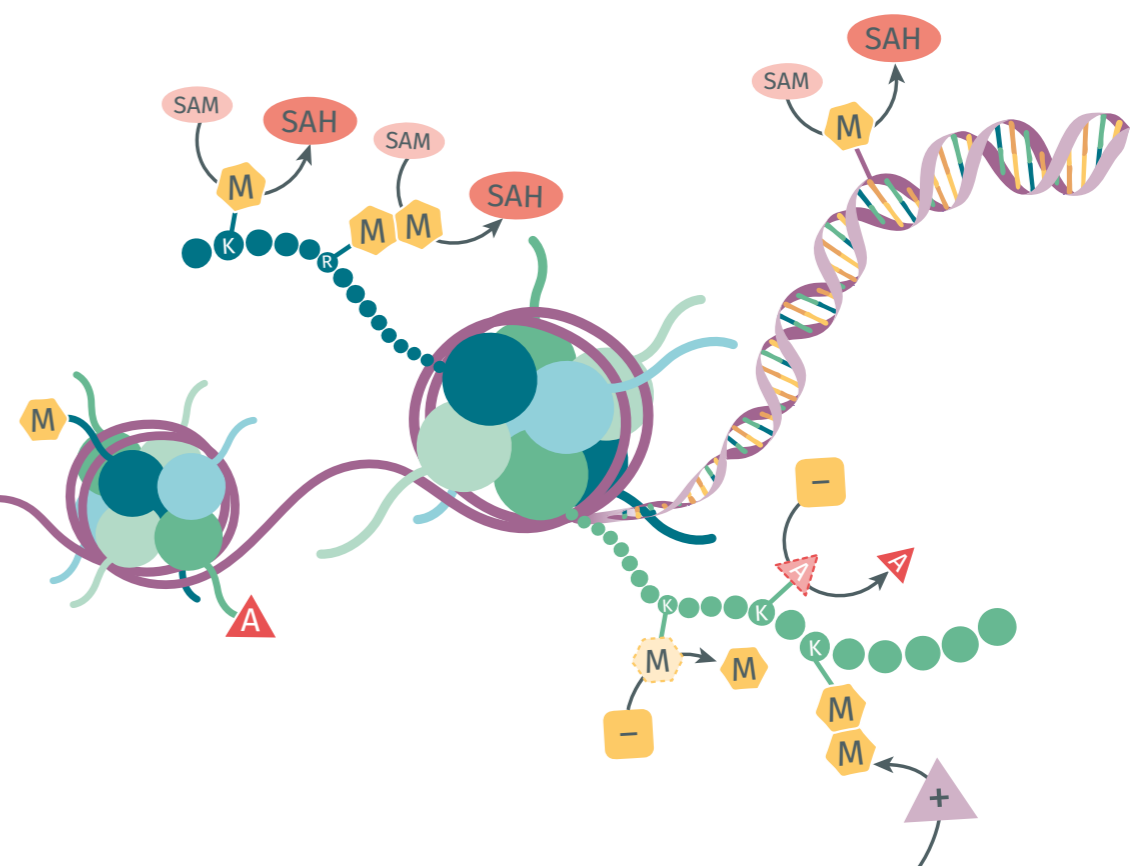
## UNIVERSAL METHYLTRANSFERASE ASSAY

This unique assay gives you the freedom to use all kinds of methyltransferases with any kind of substrate. Now you can work with the same single, reliable solution every time, simplifying assay development, avoiding conversion steps, and minimizing costs.

PRODUCT LIST: EPIgeneous Methyltransferase Assay Kit



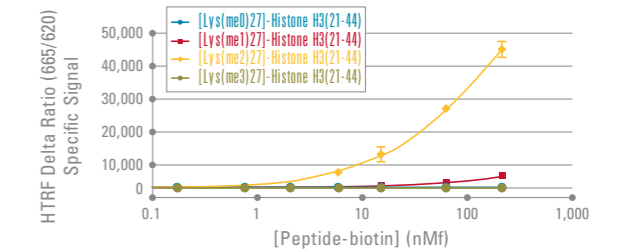
The universal EPIgeneous Methyltransferase Assay gives you the freedom and flexibility to work with different enzymes and different substrates.



## EPIGENETIC TOOLBOX REAGENTS

Customize and optimize your epigenetic screening by building your own robust HTRF assays using our toolbox reagents. Choose from an extensive line of labeled antibodies offering the highest levels of specificity and sensitivity, or select one of our enzyme-specific validated assays for fast, easy implementation.

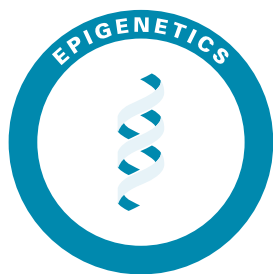
PRODUCT LIST: H3K4 H3K9 H3K4Me1,2,3 H3K9Me1,2 H3K27Me1,2,3 H3K36Me1,2



A strong signal and exceptional specificity of the H3K27Me2 antibody is shown in contrast to other levels of methylation on the same lysine.

*"There's nothing I like better than my daily interaction with customers, being able to provide them with the ideal assay or solving some technical challenge. Perhaps it's my lifelong scuba obsession, but I always dive full force into everything I do, always searching deeper, recognizing—and relishing—the fact that I'll never run out of new areas to explore. For me, that's the true beauty of what we do here at Cisbio."*

*- Arturo Gonzalez Moya,  
Custom Services Manager*



For more information or to discuss your specific labeling or custom epigenetics assay requirements, please contact your local Cisbio representative or visit [www.cisbio.com](http://www.cisbio.com)

**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  
[www.cisbio.com](http://www.cisbio.com)



cisbio