

EPIgeneous™ Binding Domain Assays



ROBUST ASSAY TOOLS TO DECIPHER KEY EPIGENETIC INTERACTIONS

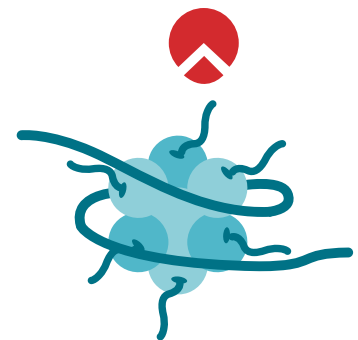
Natural relationships can reveal more than meets the eye.

Cisbio EPIgeneous™ assays help you study epigenetic targets with a complete solution that detects and screens the vast variety of epigenetic binding domain interactions with histones.

Validated on critical bromo-, chromo-, tudor- and MBT domains, these assays provide high-level, reliable output for high-throughput biochemical screening and downstream investigation, with minimal compound interference.

Leverage a leading reference in screening technology for reader protein investigation

- Validated for 28 key reader proteins
- Cisbio Discovery Assay Kit to assess your own GST-tagged binding protein
- DMSO tolerance compatible with compound screening conditions
- Low consumption of binding domain protein



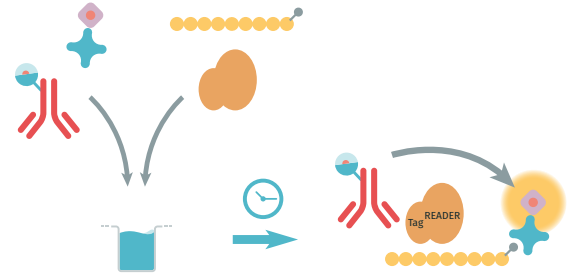
EPIgeneous™ Binding Domain Assays enable straightforward detection of reader protein activities interacting with specific chromatin sites modified by epigenetic-related enzymes.

EPIGENETICS

HOW IT WORKS HIGHLY SENSITIVE, ROBUST ASSAY KITS FOR STUDYING EPIGENETICS READERS

EPIgeneous™ Binding Domain assay kits provide fast and easy assessment of reader domain binding. Binding to the biotinylated peptide substrate, the GST-tagged reader domain protein is detected by a conjugate comprised of a donor Cryptate-labeled antibody raised against the donor GST tag, and a streptavidin-acceptor, to detect biotin-peptide. Reader domain interaction with the substrate brings the donor and acceptor dyes into close proximity, enabling FRET to occur upon light excitation. The specific signal at 665 nm is inhibited when a specific compound prevents the reader domain protein from binding to its substrate.

Performed in a single assay plate without transfer, the assay procedure is straightforward and does not require any washing or separation steps.

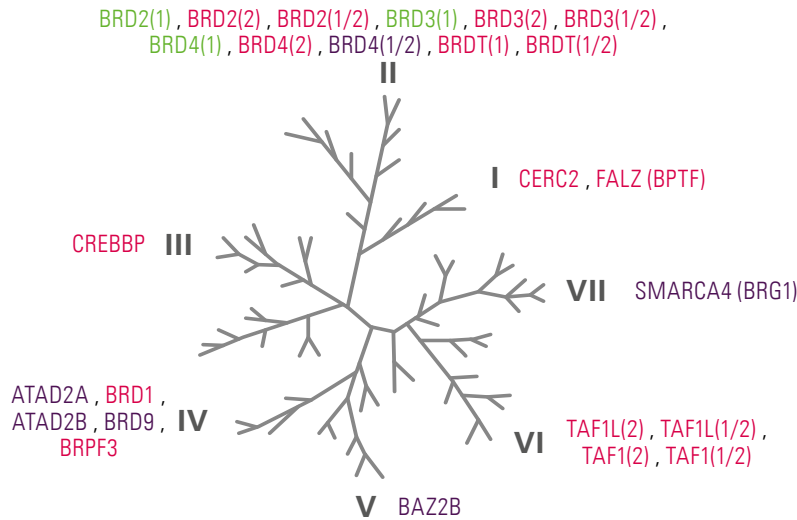


Simple procedure saves time without compromising robustness.

28 BINDING DOMAINS ALREADY VALIDATED WITH THREE EPIGENEOUS™ BINDING DOMAIN KITS

Cisbio offers three distinct kits that have been validated successfully with 28 binding domains. We continually add to this list. For other binding domains, pre-screen with the Binding Domain Discovery Kit to determine which kit to use. Please contact us for the latest details on validated domains.

Bromodomain



... and more:

MBT Domain: L3MBTL1

Chromodomain: CBX1

Tudordomain: UHRF1

■ Binding Domain Kit A ■ Binding Domain Kit B
■ Binding Domain Kit C

ORDERING INFORMATION

PRODUCT	PART#	PRODUCT	PART#
EPIgeneous™ Binding Discovery Kit – 500 tests	62BDDPEG	EPIgeneous™ Binding Domain Kit B – 10,000 tests	62BDBPEH
EPIgeneous™ Binding Domain Kit A – 500 tests	62BDAPEG	EPIgeneous™ Binding Domain Kit C – 500 tests	62BDCPEG
EPIgeneous™ Binding Domain Kit A – 10,000 tests	62BDAPEH	EPIgeneous™ Binding Domain Kit C – 10,000 tests	62BDCPEH
EPIgeneous™ Binding Domain Kit B – 500 tests	62BDBPEG		

EXPERT SERVICES TO SATISFY YOUR EXACT NEEDS



In addition to providing ready-to-use kits, Cisbio's team of experienced experts works hand-in-hand with its clients to create custom solutions. If your requirements involve binding domain proteins other than those already validated, Cisbio is the ideal trusted partner to identify the exact kit you need and save you time.

Keep your epigenetic research moving forward with faster, more robust, physiologically relevant results. Call us or visit www.htrf.com/epigenetic-screening

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 to find a list of our regional distributors



cisbio