

H3R8me2(sym) antibody

Cat. No. C15410287

Lot:	003
Size:	50 µg
Type:	Polyclonal, ChIP-grade
Isotype:	NA
Source:	Rabbit
Concentration:	0.84 µg/µl

Specificity:	Human, mouse, C. elegans, rat, chicken, Xenopus, Drosophila, plant
Purity:	Affinity purified polyclonal antibody
Storage buffer:	NA

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Description: Polyclonal antibody raised in rabbit against H3 (sym-dimethyl Arg8), using a KLH-conjugated synthetic peptide.

Applications

Applications	Suggested dilution	References
ChIP*	2 – 5 µg/million cells	Fig 1
Western blotting	2 µg/mL	Fig 2, 3

*Please note that the optimal antibody amount per IP should be determined by the end-user. We recommend testing 1–5 µg per ChIP.

Target description

Chromatin is the arrangement of DNA and proteins from which chromosomes are formed. Accordingly, chromatin is formed from nucleosomes, which are composed of a set of four histone proteins (H2A, H2B, H3, and H4) wrapped around DNA. Chromatin is a very dynamic structure in which numerous post-translational modifications work together to activate or repress the accessibility of DNA to be copied, transcribed, or repaired. These marks determine which DNA is open and commonly active (euchromatin) or tightly wound, preventing access and activation (heterochromatin). Common histone modifications include methylation of lysine and arginine, acetylation of lysine, phosphorylation of threonine and serine, and sumoylation, biotinylation, and ubiquitylation of lysine. Specifically, methylation of arginine 8 on histone H3 (H3 R8me2s) is associated with transcriptional repression and is modified by PRMT5 but not CARM1.

Diagenode SA BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue du Bois Saint-Jean, 3
4102 Seraing - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders.diagenode@hologic.com
support.diagenode@hologic.com

Diagenode LLC USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: April, 2026

Results

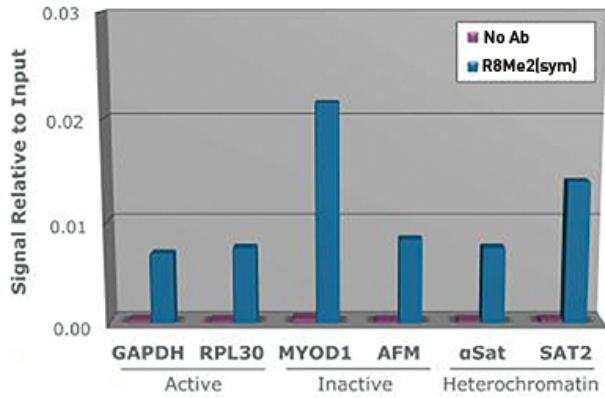


Figure 1. H3R8me2(sym) antibody ChIP results

Chromatin immunoprecipitation using the H3R8me2(sym) antibody. Chromatin from one million formaldehyde cross-linked HeLa cells was used with 2 µg of H3R8me2(sym) antibody to immunoprecipitate DNA from fixed HeLa cells, alongside a no antibody (No Ab) control. DNA was measured by PCR and normalized to total input.

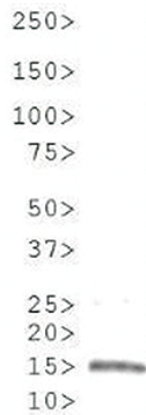


Figure 2. H3R8me2(sym) antibody Western blot results

Western blot of the H3R8me2(sym) antibody. 30 µg of *C. elegans* embryo lysate. Primary antibody: diluted 1:500 overnight at 4°C. Secondary antibody: IRDye800™ rabbit secondary antibody at 1:10,000 for 45 minutes at room temperature (RT). Predicted/observed size: ~15 kDa. Other bands: None.

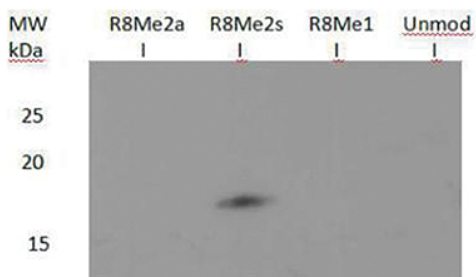


Figure 3. H3R8me2(sym) antibody Western blot results

Western blot of the H3R8me2(sym) antibody. Lane 1: R8Me2a. Lane 2: R8Me2s. Lane 3: R8Me1. Lane 4: Histone H3 R8. Load: 30 µg per lane. Primary antibody incubated at 2 µg/mL overnight at 4°C. Secondary antibody: IRDye800™ rabbit secondary antibody at 1:10,000 for 45 minutes at room temperature (RT). Predicted/observed size: ~15 kDa. Other bands: None.