

## H4R3me1 polyclonal antibody - Classic

**Cat. No.** C15410299

**Type:** Polyclonal

**Source:** Rabbit

**Lot #:** 001

**Size:** 50 µg

**Concentration:** 0.43 µg/µl

**Specificity:** Human, mouse, rat

**Purity:** Affinity purified

**Storage:** Store at -20°C; for long storage, store at -80°C.  
Avoid multiple freeze-thaw cycles.

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

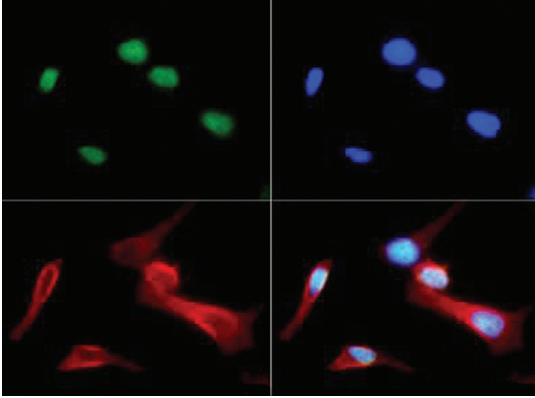
### Applications

	Suggested dilution	Results
ChIP	2-5 µg/million cells	
IF	1:50	Figure 1
Western blot	1:1,000	Figure 2
Immunochemistry	1:50	
Dot Blot	1:1,000	

### Target description

Histone H4 is an 11.4 kD nuclear protein that is a component of an octamer containing pairs of each of four core histones (H2A, H2B, H3, H4). The core histones create nucleosome structure of chromosomal fiber in eukaryotes and are dynamic in gene regulation. Histone H4 shows transcriptional regulation that is both cell cycle-dependent and -independent. Histone H4 is modified by phosphorylation, acetylation, ribosylation, and methylation. Histone H4 has been shown to interact with other histone acetyltransferases, histone deacetylases, and PARP.

## Results



**Figure 1. Immunofluorescence**

Immunofluorescence of H4R3me1 antibody. Tissue: Nonmitotic, prophase, and telophase HeLa cells. Fixation: 0.5% PFA. Primary antibody used at a 1:50 dilution for 1 h at RT. Secondary antibody: FITC secondary antibody at 1:10,000 for 45 min at RT. Localization: Histone H4R3me1 is nuclear and chromosomal. Staining: Histone H4R3me1 is expressed in green, nuclei are counterstained with DAPI (blue).



**Figure 2. Western Blot**

Western Blot of H4R3me1 antibody. 30 µg of NIH-3T3 histone extracts. Primary antibody used at a 1:1,000 dilution overnight at 4°C. Secondary antibody: IRDye800™ rabbit secondary antibody at 1:10,000 for 45 min at RT. Predicted/Observed size: ~13 kDa. Other band(s): None.

**Diagenode sa. BELGIUM | EUROPE**

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

**Diagenode Inc. 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:** September 15, 2014