

## H4K20me2 monoclonal antibody - Classic

Cat. No. **C15200205**

Type: Monoclonal

Source: Mouse

Lot #: 41582

Size: 100 µl

Concentration: 1 µg/µl

**Specificity:** Human

**Purity:** Protein G purified

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

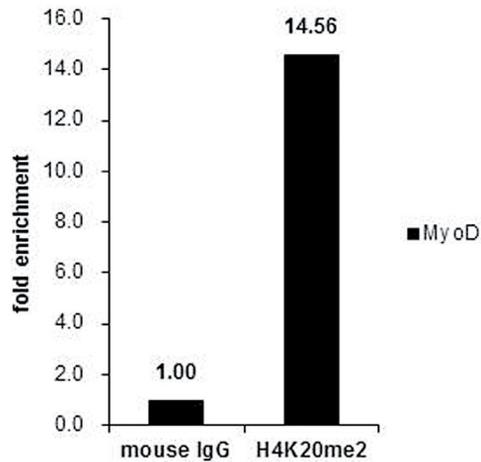
**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

### Applications

	Suggested dilution*	Results
ChIP assay	5 µg/ChIP	Figure 1
ICC/IF	1:100-1:1,000	Figure 3
WB	1:1,000-1:10,000	Figure 2

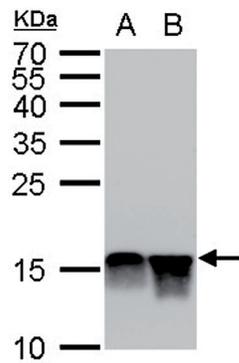
\*Optimal dilutions/concentrations should be determined by the researcher.

### Results



**Figure 1. H4K20me2 ChIP results**

ChIP was performed with HeLa chromatin extract and 5 µg of either control mouse IgG (Cat. No. C15200001) or H4K20me2 antibody. The precipitated DNA was detected by PCR with primer set targeting to MyoD.



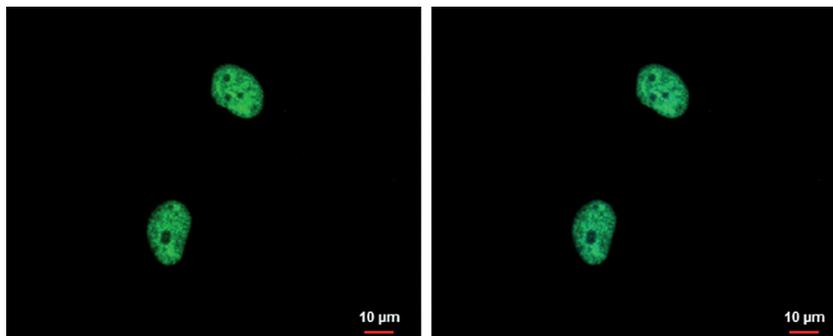
**Figure 2. H4K20me2 Western Blot**

H4K20me2 antibody detects Histone H4 protein dimethylated at Lys20 by western blot analysis.

A. 30 µg HeLa whole cell extract

B. 30 µg HeLa nuclear extract

H4K20me2 antibody dilution: 1:5,000



**Figure 3. H4K20me2 IF results**

H4K20me2 antibody detects Histone H4 dimethylated at Lys20 at nucleus by immunofluorescent analysis.

Sample: HeLa cells were fixed in 4% paraformaldehyde at RT for 15 min.

Green: Histone H4 (dimethylated at Lys20) stained by H4K20me2 antibody diluted 1:500.