

## Mouse IgG

**Cat. No. C15400001**

Lot:	/	Specificity:	NA
Size:	15 µg / 100 µg	Purity:	Purified by protein A chromatography.
Type:	Polyclonal, <b>ChIP-grade, CUT&amp;Tag-grade</b>	Storage buffer:	5 mM phosphate, 75 mM NaCl, pH 7.8; 0.06% sodium azide. Contains sucrose for stabilization
Isotype:	NA		
Source:	Rabbit		
Concentration:	1 µg/µl		

**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.

**Description:** The negative control IgG derived from mouse has been extensively validated in chromatin immunoprecipitation (ChIP). It contains a range of the IgG subclasses present in the serum of healthy mice. This IgG preparation is intended for use as a negative control in ChIP, CUT&Tag, MeDIP, IF and other experiments performed with mouse-derived antibodies. The negative control IgG from mouse should be used in parallel with the specific antibody at the same concentration. It is also included in the Antibody package (anti-mouse), Cat. No. C01070023.

## Results

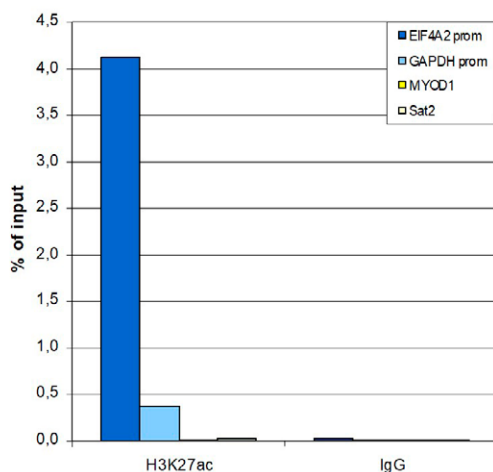


Figure 1: ChIP with the mouse IgG negative control antibody

ChIP assays were performed on sheared chromatin from 1 million HeLa cells using the mouse monoclonal antibody against H3K27ac (cat. No. C15200184) and the Hologic Diagenode ChIP kit. Mouse IgG (cat. No. C15400001) was used as a negative IP control. One microgram (µg) of each antibody was used per ChIP experiment. Quantitative PCR was performed with primers specific for the promoters of the active GAPDH and EIF4A2 genes, and for the inactive MYOD1 gene and the Sat2 satellite repeat. Figure 1 shows the recovery, expressed as a % of input (the relative amount of immunoprecipitated DNA compared to input DNA after qPCR analysis).

### 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: May, 2026

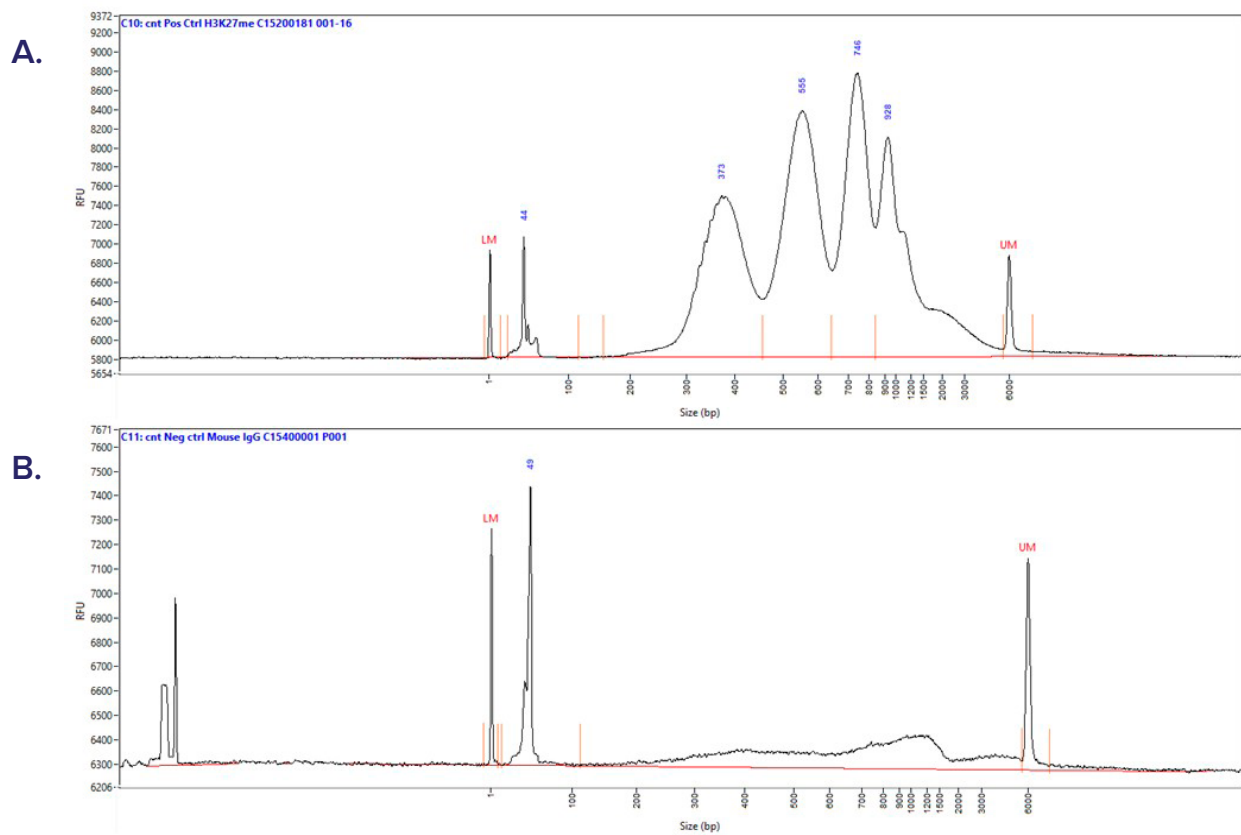


Figure 2: CUT&Tag results obtained with the mouse IgG negative control antibody

CUT&Tag was performed on 50,000 K562 cells using 0.5  $\mu$ g of the rabbit monoclonal antibody against H3K27me3 (cat. No. C15200181) and the Universal CUT&Tag kit (cat. No. C01070024). Mouse IgG (cat. No. C15400001) was used as a negative control. The resulting libraries were analyzed on a Fragment Analyzer. Figure 2 shows the results with the H3K27me3 antibody on top and the IgG negative control at the bottom.

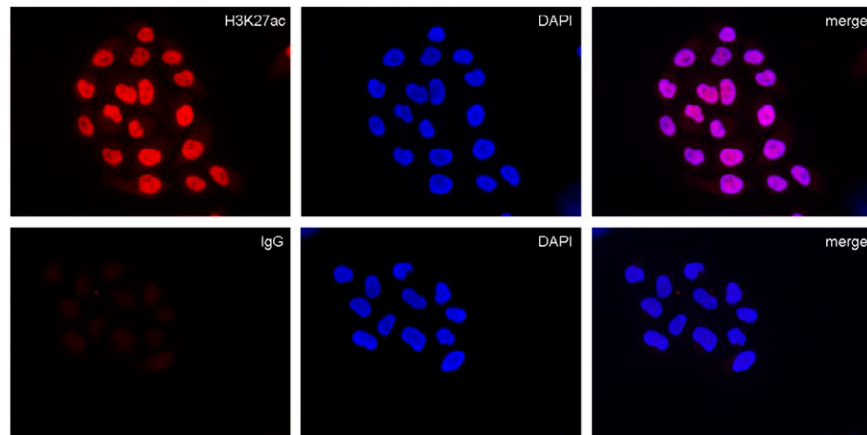


Figure 3: Immunofluorescence with the mouse IgG negative control antibody

HeLa cells were stained with the mouse monoclonal antibody against H3K27ac (cat. No. C15200184) (top) and with DAPI. Mouse IgG (cat. No. C15400001) was used as a negative control (bottom). Cells were fixed with 4% formaldehyde for 10 minutes and blocked with PBS/TX-100 containing 5% normal goat serum and 1% BSA. The cells were immunofluorescently labeled with either the H3K27ac antibody or the mouse IgG negative control antibody (left), each diluted 1:500 in blocking solution; followed by an anti-mouse antibody conjugated to Alexa594. The middle panel shows staining of the nuclei with DAPI. A merge of the two stainings is shown on the right.