

## Rabbit IgG antibody

**Cat. No.** C15410206

**Type:** Polyclonal

**Source:** Rabbit

**Lot:** RIG002

**Size:** 250 µg

**Concentration:** 1 µg/µl

**Specificity:** NA

**Purity:** Purified by protein A chromatography. In 2 mM phosphate, 30 mM NaCl, pH 7.8; 0.02% sodium azide. Contains sucrose for stabilization.

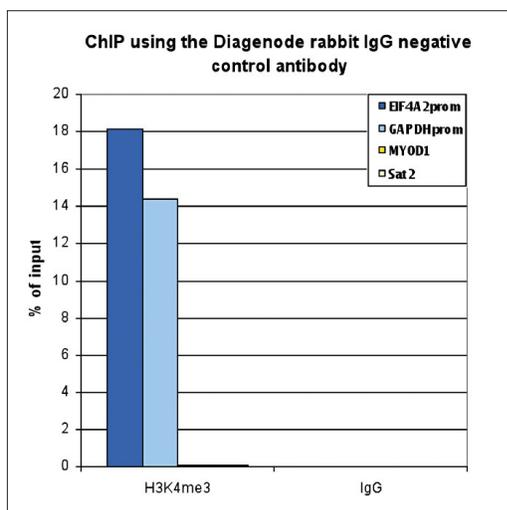
**Storage:** Store at 4°C/-20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

**Storage buffer:** PBS containing 0.05% azide.

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

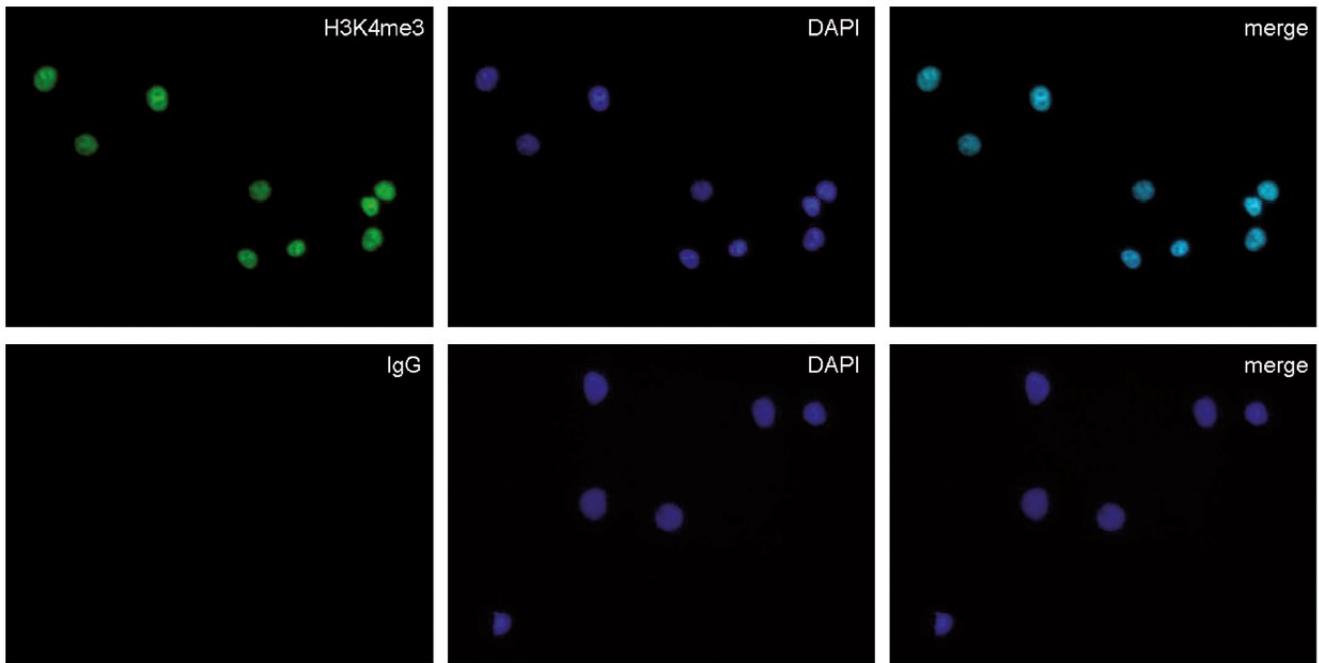
**Description:** The negative control IgG from rabbit has been extensively validated in chromatin immunoprecipitation (ChIP). It contains a spectrum of the IgG subclasses present in serum of healthy rabbits. This IgG preparation is intended for use as a negative control in ChIP, MeDIP, IF and other experiments performed with specific antibodies made in rabbit. The negative control IgG from rabbit should be used in parallel with the specific antibody at the same concentration. It is also included in many of our ChIP and MeDIP kits.

## Results



**Figure 1. ChIP with the Diagenode rabbit IgG negative control antibody**

ChIP assays were performed using the Diagenode rabbit polyclonal antibody against H3K4me3 (Cat. No. C15410003) and the "iDeal ChIP-seq" kit (Cat. No. C01010051) on sheared chromatin from 1 million HeLa cells. Rabbit IgG (cat. No. C15410206) was used as a negative IP control. One µg of antibody per ChIP experiment was used for both antibodies. 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)



**Figure 2. Immunofluorescence with the Diagenode rabbit IgG negative control antibody**

HeLa cells were stained with the Diagenode rabbit polyclonal antibody against H3K4me3 (Cat. No. C15410003) (top) and with DAPI. Rabbit IgG (Cat. No. C15410206) was used as a negative control (bottom). Cells were fixed with 4% formaldehyde for 10' and blocked with PBS/TX-100 containing 5% normal goat serum and 1% BSA. The cells were immunofluorescently labeled with the H3K4me3 or rabbit IgG negative control antibody (left) diluted 1:200 in blocking solution followed by an anti-rabbit antibody conjugated to Alexa488. The middle panel shows staining of the nuclei with DAPI. A merge of the two stainings is shown on the right.