

Rabbit IgG

Cat. No. C15410206

Lot:	/
Size:	250 µg
Type:	Polyclonal, ChIP-grade, CUT&Tag-grade
Isotype:	NA
Source:	Rabbit
Concentration:	1 µg/µl

Specificity:	NA
Purity:	Purified by protein A chromatography.
Storage buffer:	2 mM phosphate, 30 mM NaCl, pH 7.8; 0.04% sodium azide. Contains sucrose for stabilization.

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 rabbit-derived negative control IgG has been extensively validated in chromatin immunoprecipitation (ChIP). It contains a spectrum of the IgG subclasses present in the serum of healthy rabbits. This IgG preparation is intended for use as a negative control in ChIP, CUT&Tag, MeDIP, IF and other experiments performed with specific antibodies made in rabbit. The rabbit-derived negative control IgG should be used in parallel with the specific antibody at the same concentration. It is also included in the Antibody package (anti-rabbit), Cat. No. C01070022.

Results

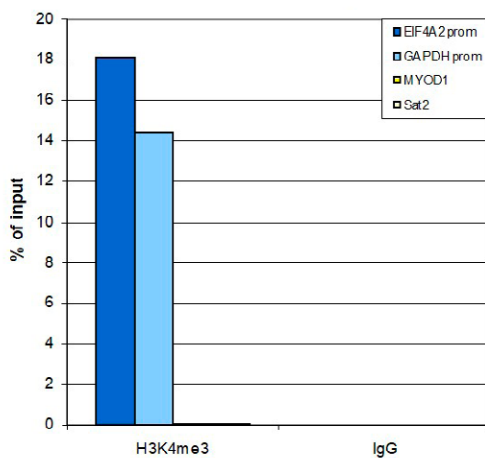


Figure 1: ChIP with the rabbit IgG negative control antibody

ChIP assays were performed using the 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 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).

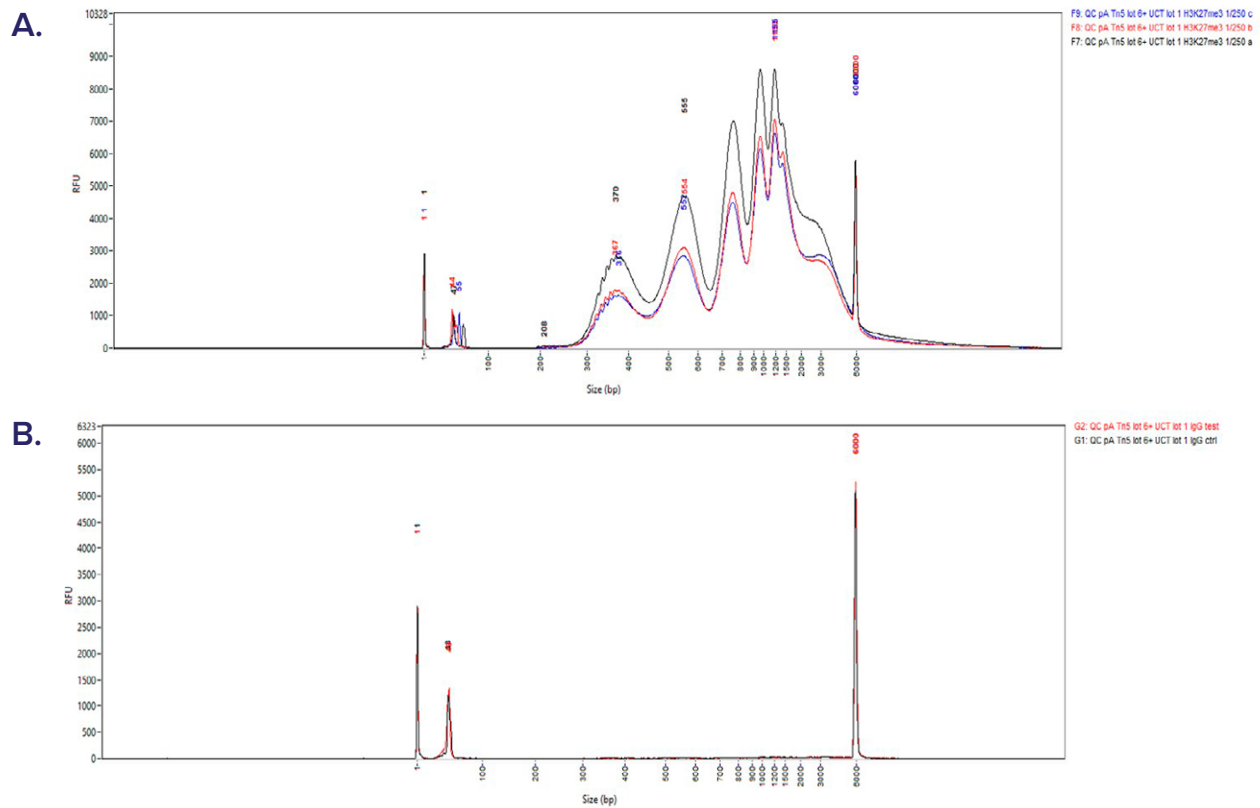


Figure 2: CUT&Tag results obtained with the rabbit IgG negative control antibody

CUT&Tag was performed on 50,000 K562 cells using 0.2 µg of the rabbit polyclonal antibody against H3K27me3 (cat. No. C15410195) and the Universal CUT&Tag kit (cat. No. C01070024). Rabbit IgG (cat. No. C15410206) was used as a negative control. The resulting libraries were analysed on a Fragment Analyzer. Figure 2 shows the results with the H3K27me3 antibody (3 replicates, top) and the IgG negative control (2 replicates, bottom).

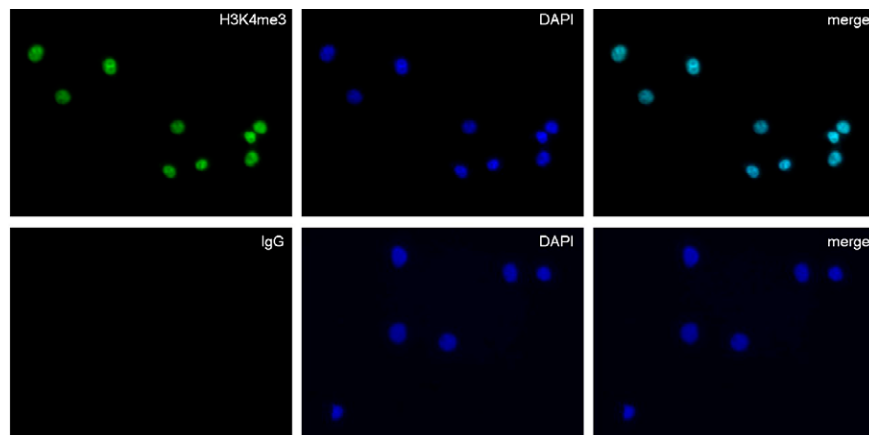


Figure 3: Immunofluorescence with the rabbit IgG negative control antibody

HeLa cells were stained with the 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 minutes 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.