

Rat IgG

Cat. No. C15420001

Type: Polyclonal

Source: Rat

Size: 25 µg/25 µl

Concentration: 1 µg/µl

Specificity: NA

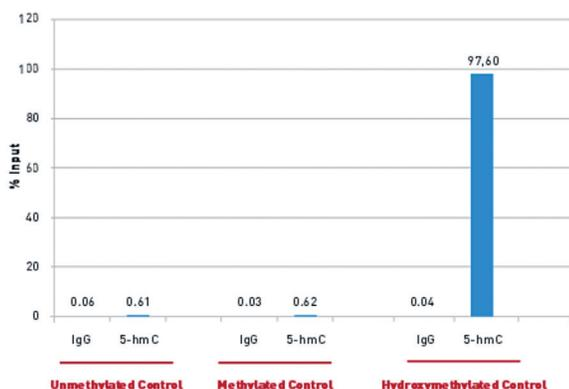
Purity: Purified by protein A chromatography. In 10 mM phosphate, 150 mM NaCl, pH 7.8; 0.08% sodium azide. Contains sucrose for stabilization.

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

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

Product description: The negative control IgG from rat has been extensively validated in methylated DNA immunoprecipitation (MeDIP). It contains a spectrum of the IgG subclasses present in serum of healthy rats. This IgG preparation is intended for use as a negative control in MeDIP, ChIP, IF and other experiments performed with specific antibodies made in rat. The negative control IgG from rat should be used in parallel with the specific antibody at the same concentration. It is also included in our rat hMeDIP kit.

Results



hMeDIP with the Diagenode rat IgG negative control antibody

hMeDIP assays were performed using the Diagenode rat monoclonal antibody against 5-hmC (cat. No. C15220001) and the "hMeDIP" kit (rat, cat. No. C02010030) on 1 µg of sheared DNA from mouse ES cells. The DNA was spiked with 0.025 ng of each of the DNA standards from the 5-hmC, 5-mC & cytosine DNA standard pack (cat. No. C02040011), prior to the IP. Rat IgG (cat. No. C15420001) was used as a negative IP control. 2.5 µg of antibody per ChIP experiment was used for both antibodies. Quantitative PCR was performed with the primers specific for the unmethylated, methylated and hydroxymethylated controls present in the kit. 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 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: June 7, 2016