

PRODUCT NAME

Pre-blocked Protein A/G beads

Cat. No.: **mcg-606-300**

Format: 300 µl

Product description

Pre-blocked Protein A/G beads are designed to capture immunoglobulins (Ig) for downstream immunoprecipitation of methylated DNA (MeDIP). The beads are supplied as a 50% suspension in 0.1 % BSA, containing azide.

Storage and stability

Store at 4°C. Do not freeze.

Applications

Methylated DNA Immunoprecipitation (MeDIP).

Quality control

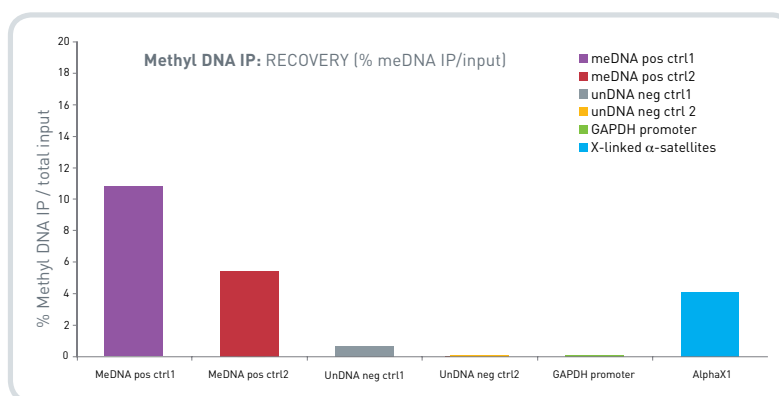


Figure 1: Methyl DNA IP results obtained with the Diagenode MeDIP Kit (Cat. No. mc-green-03), including Pre-blocked Protein A/G beads.

Methyl DNA IP assays were performed using DNA from NB4 cells, the Diagenode antibody directed against 5-methylcytosine and optimized PCR primer pairs for qPCR. The DNA was prepared with the GenDNA module. The IP on the human DNA sample, together with the internal kit controls. The internal positive and negative DNA controls included in the IP assay are methylated DNA (meDNA) and unmethylated DNA (unDNA). The DNA is then purified from the IP'd material and analysed by PCR using the primer pairs included in the kit (see below).

Each "primer pair" targets a specific DNA and expected results are as follows:

- Internal DNA controls

"meDNA pos ctrl1": meDNA control (positive signal is obtained for methylation).

"meDNA pos ctrl2": meDNA control (positive signal is obtained for methylation).

"unDNA neg ctrl1": unDNA control (no signal is obtained for 0% methylation).

"unDNA neg ctrl2": unDNA control (no signal is obtained for 0% methylation).

- Human DNA sample:

"GAPDH promoter": no signal is expected as this region is not methylated.

"X-linked alpha-satellites": a positive signal is expected as it is a methylated region.

The Methyl DNA IP controls reveal IP efficiency.