

## DiaMag protein G-coated magnetic beads

Cat. No. **C03010021-220** Format: 220  $\mu$ l

Cat. No. **C03010021-660** Format: 660  $\mu$ l

Cat. No. **C03010021-150** Format: 1.5 ml

### Product description

The protein G-coated magnetic beads have been extensively validated in chromatin immunoprecipitation assay (ChIP). These beads are intended for isolation of immune complexes (chromatin and specific antibody) in ChIP experiments. The beads are suitable for immunoprecipitation of mouse IgG1, IgG2a, IgG2b and IgG3, rat IgG1, IgG2a, IgG2b and IgG3 rabbit and goat polyclonal Abs and human IgG1, IgG2, IgG3 and IgG4. The beads should be washed before use.

### Format

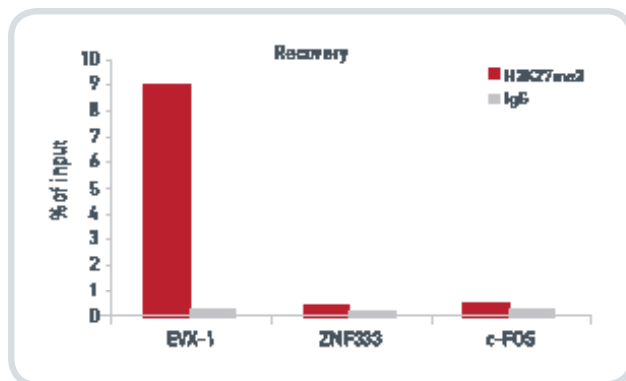
Supplied as a suspension in PBS (pH 7.4), with 0.1% Tween-20 and 0.02% sodium azide.

### Storage and stability

Store at 4°C. Do not freeze. Keep the beads in liquid suspension during storage as drying will result in reduced performance.

### Precautions:

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



**Figure 1: ChIP results obtained with Diagenode Auto ChIP kit (Cat. No: AB-Auto01-A100) including protein G-coated magnetic beads.**

ChIP experiments were performed using chromatin of 100.000 Ntera1 cells per immunoprecipitation reaction. H3K27me3 was used and the enrichment of a positive genomic control region (EVX-1) and two negative genomic control regions (ZNF333 and c-fos) were analyzed. Data are shown as % of input. Experiments were performed using protein G-coated magnetic beads from Diagenode as they are compatible with rabbit polyclonal IgG antibodies. The ChIP was run on Diagenode's SX-8G IP-Star® Automated System.