

PRODUCT NAME	
DiaMag protein A-coated magnetic beads	
Cat. No. <b>kch-802-220</b>	Format: 220 µl
Cat. No. <b>kch-802-660</b>	Format: 660 µl
Cat. No. <b>kch-802-150</b>	Format: 1.5 ml
Cat. No. <b>kch-802-005</b>	Format: 5 ml

**Product description**

The protein A-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 performed. The beads are suitable for immunoprecipitation of rabbit polyclonal Abs, mouse IgG2a, IgG2b and IgA, guinea pig IgG, dog IgG, pig IgG. The beads should be washed before use.

**Format**

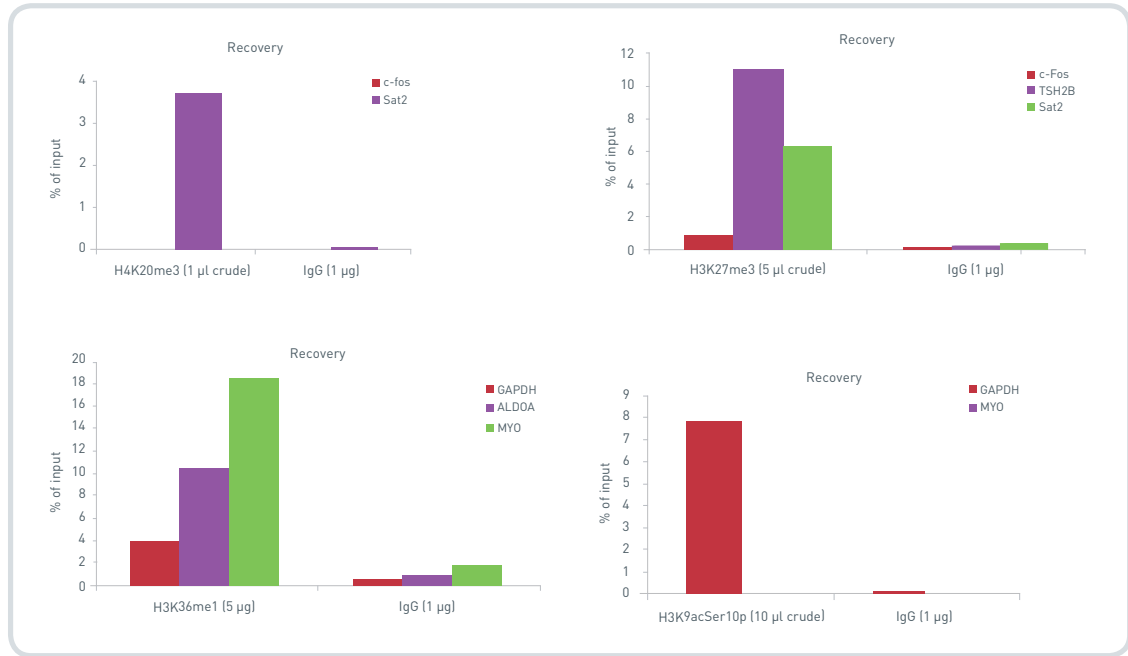
Supplied as a suspension in PBS (pH 7.4), with 0.1% Tween-20 and 0.02% sodium azide.

**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 A-coated magnetic beads.**

ChIP experiments were performed using chromatin of 10000 HeLA cells per immunoprecipitation reaction. Diagenode antibodies selected for the experiment were H4K20me3, H3K27me3, H3K36me3 and H3K9acSer10 rabbit polyclonal histone antibodies. Protein A-coated magnetic beads from Diagenode were selected as they are compatible with rabbit polyclonal IgG antibodies. Experiments were performed on the SX-8G IP-Star® Automated System. Different genomic control regions were selected according to the antibody tested. Data shown as % of input.