

DiaMag protein A-coated magnetic beads

Cat. No. **C03010020-220** Format: 220 μ l

Cat. No. **C03010020-660** Format: 660 μ l

Cat. No. **C03010020-150** Format: 1.5 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.