

## DiaMag anti-mouse IgG-coated magnetic beads

Cat. No. **C03010022-220** Format: 220  $\mu$ l

Cat. No. **C03010022-660** Format: 660  $\mu$ l

Cat. No. **C03010022-150** Format: 1.5 ml

### Description

These magnetic beads coated with anti-mouse IgG antibody have been extensively validated in chromatin immunoprecipitation assay (ChIP) and methylated DNA immunoprecipitation assays (MeDIP). The beads bind mouse IgG1, IgG2a, IgG2b. The beads should be washed before use.

### Format

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

### Storage

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.