

Control Magnetic Beads

Cat. # UM400M

Background: Control Magnetic beads is generated by blocking an activated matrix (identical to that used for TUBEs conjugation) with ethanolamine.

Product Information

Physical State: Control magnetic beads (1 mL) is supplied as a slurry in PBS, pH 7.4.

Storage: +4°C. Storage below this temperature may result in freezing and/or loss of function.

Applications: For use in "pre-clearing" cell lysates, as well as a negative control for the pulldown of polyubiquitylated proteins in conjunction with magnetic beads-conjugated TUBEs (UM401M and UM402M).

General recommendations: use 100µl of control magnetic beads slurry with 1ml of cell lysate (3-5mg/ml total protein) by incubation for 1hr (4C). Individual conditions may have to be optimized by the end user.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2011 LifeSensors, Inc. All Rights Reserved