

Mouse IgG IP Control Beads

Cat. # CIG01-Beads

Lot:

Upon arrival, store at 4°C (desiccated)

See datasheet for storage after reconstitution

Form:	Lyophilized powder
Amount of material:	1 x 330 µl when reconstituted
Validated applications:	IP control for Ab affinity reagents
Species reactivity:	na
Host/Isotype:	Mouse/polyclonal
Clone:	na

Background Information

Many of Cytoskeleton Inc's Signal-Seeker™ affinity enrichment beads are based on mouse monoclonal antibody reagents covalently bound to beads. Mouse IgG IP Control Beads provide an ideal negative control and should be included in an IP experiment to control for non-specific binding in any antibody based affinity immunoprecipitation reaction, see **Figure 1**.

Material

Normal whole mouse IgG from non-immunized animals has been co-valently linked to agarose affinity beads. Antibody binding is in the range of 0.3-0.8 mg antibody per ml of bead slurry which is a similar range to Signal-Seeker™ affinity reagents.

Storage and Reconstitution

Shipped at ambient temperature. The lyophilized IP control beads can be stored desiccated at 4°C for 6 months. For reconstitution, the product tube should be briefly centrifuged to collect the lyophilized beads at the bottom of the tube. Reconstitute each tube in 330ul of Milli-Q water to achieve 50% slurry and store at 4°C. Alternatively, reconstitute in 330 µl of 50% glycerol and store in -20°C. In both cases, allow beads to rehydrate completely before use (15-20 minutes). Final buffer composition is 200 mM PIPES, 5% sucrose, and 1% dextran. When stored and reconstituted as described, the product is stable for at least 6 months in 4°C and 12 months in -20°C.

Applications

Immunoprecipitation (IP) Application

Use an equivalent volume of control bead slurry as that being used for an enrichment IP assay. This is generally in the region of 30-40 µl per IP. Sufficient for 8-10 IP reactions. See **Figure 1** for representative data.

Figure 1: Enrichment of SUMOylated proteins from cell lysates

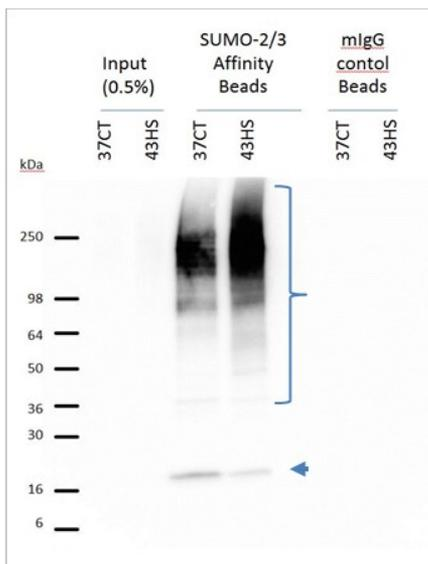


Figure Legend: A431 cell lysates were prepared from HS43: Heat Shock treated (43°C for 10min) and CT37: untreated cells. SUMO-2/3 conjugates were enriched from 1mg of lysate and immuno-blotted by anti-SUMO-2/3 antibody (Cytoskeleton Cat # ASM23) along with mouse IgG (mIgG) control beads (Cat # CIG01-Beads).

The level of total SUMO-2/3 conjugates in heat shock treated cell SUMO 2/3 Affinity Beads (Lane 43HS) is stronger than control cells (SUMO 2/3 Affinity Beads Lane 37CT). Total SUMOylated protein signal is delineated by the blue bracket, the blue arrowhead indicates the position of free SUMO 2/3. Lack of signal in the mouse IgG (mIgG) lanes demonstrates the specificity of the SUMO-2/3 bead reagent.