

Mouse IgG IP Control Beads

Cat. # CIG02-Beads
Lot: 045

Upon arrival, store at 4°C (desiccated)
See datasheet for storage after reconstitution

Form:	Lyophilized powder
Amount of material:	1 x 500 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 co-valently 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

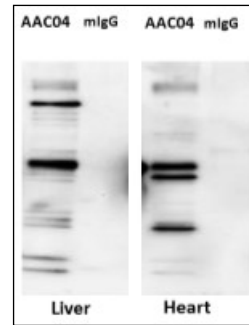
Shipped at ambient temperature. The lyophilized IP control beads can be stored desiccated at 4°C for 6 months. For reconstitution, the product tube should be briefly centrifuged to collect the lyophilized beads at the bottom of the tube. Reconstitute each tube in 330ul of Milli-Q water to achieve 50% 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 50-60 µl per IP. Sufficient for 8-10 IP reactions. See Figure 1 for representative data.

Fig. 1: Isolation and detection of acetylated proteins from mouse tissue



Mouse tissue extracts (liver and heart) were obtained with BlastR buffer. IP was performed using AAC04 beads (60ug) or CIG02-beads, mIgG control beads (60ug) in 1mg of tissue lysate. Enriched acetylated proteins were separated by SDS-PAGE and analyzed by western blot with AAC03-HRP (1:3000).