

Mouse IgG IP Control Beads

Cat. # CIG02-Beads
Lot:

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 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 covalently 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

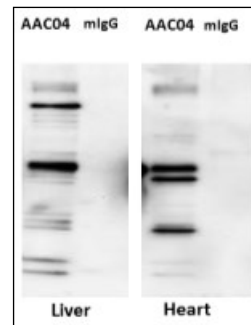
Shipped at ambient temperature. The lyophilized IP control beads can be stored desiccated at 4°C for 6 months. For reconstitution, the product tube should be briefly centrifuged to collect the lyophilized beads at the bottom of the tube. Reconstitute each tube in 500 µl of 50% glycerol and store in -20°C. 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 at -20°C.

Applications

Immuno-precipitation (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).