

ChIP-IT Protein G Magnetic Beads, 40 rxns

Catalog No: 53033

Format: 40 rxns

Quality Control: ChIP-IT Protein G Magnetic Beads are ideal for chromatin immunoprecipitation (ChIP) of protein/nucleic acid complexes and immunoprecipitation (or co-immunoprecipitation) of proteins and protein complexes.

ChIP-IT Protein G Magnetic Beads are validated using ChIP. An antibody directed against the protein of interest is used to capture formaldehyde cross-linked protein/DNA complexes from fragmented chromatin samples. Twenty-five microliters of the ChIP-IT Protein G Magnetic Beads are added to each capture reaction to bind the Fc-region of the antibody. A magnet is then used to collect the beads to the side of the reaction tube enabling easy removal of the unbound chromatin in the supernatant. The ChIP-enriched DNA is eluted and purified following reversal of the cross-links and Proteinase K treatment.

ChIP Protocol Reference: We recommend using a published and validated protocol such as the Myers Lab ChIP-Seq Protocol, <http://myers.hudsonalpha.org>

Contents:

ChIP-IT Protein G Magnetic Beads are 2.8 µm superparamagnetic beads supplied in PBS, pH 7.4, with 0.01% Tween-20 and 0.09% sodium azide. Each bead contains recombinant Protein G covalently coupled to the surface. Beads are provided at a concentration of 30 mg/ml.

Guidelines:

- Fully resuspend ChIP-IT Protein G Magnetic beads by pipetting before use. Beads settle quickly and should be resuspended more than once if working with a large number of samples.
- Use 25 µl resuspended beads per ChIP reaction.
- Use 25 µl resuspended beads per IP or Co-IP reaction.
- The binding capacity of the ChIP-IT Protein G Magnetic Bead is approximately 8 µg human IgG per mg beads. (Use up to 6 µg antibody per 25 µl beads (0.75 mg)).
- When working with antibodies with low binding affinity, increase the incubation time or pre-incubate sample and antibody prior to capture with the ChIP-IT Protein G Magnetic Beads.

Storage and Guarantee: Store at 4°C. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.