

Protein G Agarose Beads

Catalog No: 37499

Format: 1.2 ml

Quality Control: Protein G Agarose (part number 104889) is used for the immunoprecipitation of IgG from mammalian samples (including human, mouse, rabbit, goat, sheep, horse, cow, rat, mouse and guinea pig). It does not bind IgM, IgA, IgE or serum albumin. Protein G agarose can be used in chromatin immunoprecipitation (ChIP) or Co-immunoprecipitation (Co-IP) reactions.

Protein G agarose beads are supplied as a 50% slurry in PBS, pH 7.0 and 20% ethanol. The beads have a binding capacity of 10 µg IgG/µl bead. We recommend using 30 µl Protein G agarose beads per reaction.

Storage and Guarantee: Protein G agarose beads are shipped at room temperature and should be stored at 4°C upon arrival. 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.