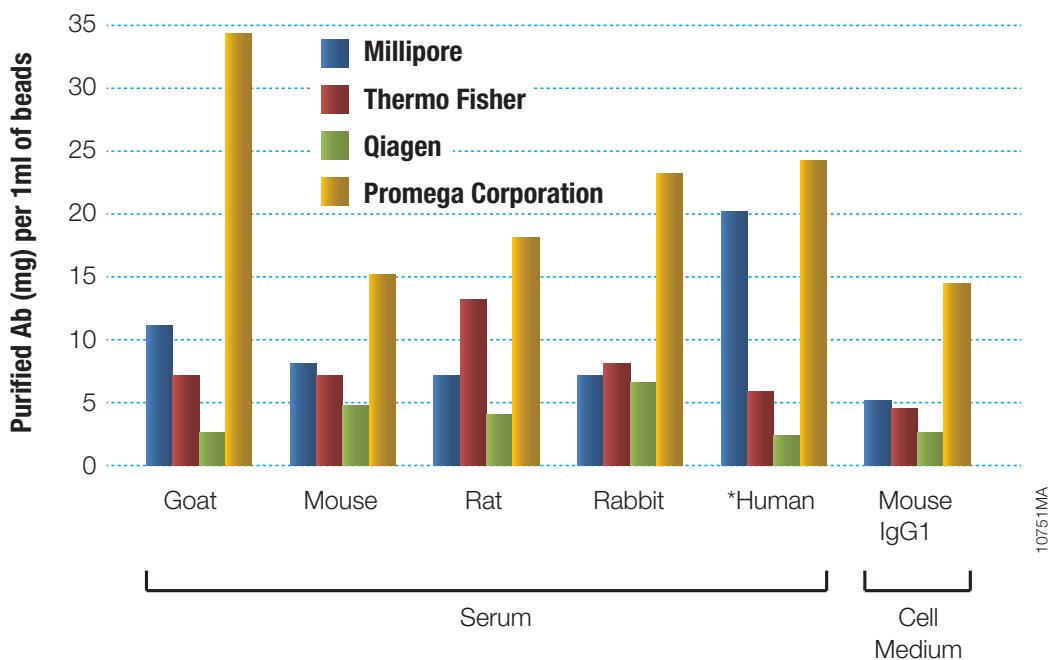


Magne™ Protein G and Protein A Beads

How effective is your antibody purification strategy?

Competitor Comparison for Magnetic Antibody Purification Resins



Comparison of antibody yield using the Magne™ Protein G Beads versus competitors' products from diverse sample types. *Delipidated human gamma globulin serum fraction was used as the source material.

High Capacity and Specificity: Magne™ Protein G and Magne™ Protein A Beads allow exceptional antibody yields (capacities in excess of 25mg per 1ml of beads depending on species and isotype) from diverse sample types such as serum, ascites and cell medium. The binding selectivity for immunoglobulins prevents copurification of albumin and other protein contaminants.

Optimized Performance: Robust magnetic response makes the Magne™ Protein G and Magne™ Protein A Beads ideal for high-throughput, automated applications. Validated protocols are available for microscale (20µl) to medium-scale (50ml) sample volumes.

Efficient Recovery: The magnetic method minimizes antibody losses encountered during the column chromatography, dialysis and centrifugation concentration steps found in traditional antibody purification protocols.

Cost Effectiveness: Magne™ Protein G and Magne™ Protein A Beads enable high-yield, high-purity antibody purification without using expensive liquid chromatography equipment.

Experience the magnetic antibody purification approach today!

Ordering Information

Product	Size	Cat.#
Magne™ Protein G Beads, 20% Slurry	1ml	G7471
	5ml (5 × 1ml)	G7472
	50ml	G7473
Magne™ Protein A Beads, 20% Slurry	1ml	G8781
	5ml (5 × 1ml)	G8782
	50ml	G8783

Patent Pending.
Magne is a trademark of Promega Corporation.

