

OmicsLink™ Peroxidase Conjugated Anti-Mouse IgG Antibody
(Goat)

Catalog number

CGAB-GOA-0025 (25 µg)

Immunogen

Mouse IgG whole molecule

Host

Goat

Label

Horseradish Peroxidase (HRP)

Specificity

Minimum cross reactivity to bovine, chicken, goat, guinea pig, hamster, horse, human, rabbit, rat and sheep serum proteins.

Purity

This product was prepared from monospecific antiserum by immunoaffinity chromatography using mouse IgG coupled to agarose beads followed by solid phase adsorption(s) to remove any unwanted reactivities. Assay by immunoelectrophoresis resulted in a single precipitin arc against anti-Peroxidase, anti-goat Serum, mouse IgG and mouse serum. Minimum cross reactivity to bovine, chicken, goat, guinea pig, hamster, horse, human, rabbit, rat and sheep serum proteins.

Formulation

0.25 mg/ml in PBS, pH 7.2 with 50% (v/v) glycerol, 2.5 mg/ml Bovine Serum Albumin (BSA) IgG and Protease free. Preservative: 0.02% Thimerosal.

Application

Immunoblotting (western or dot blot), ELISA, immunoperoxidase electron microscopy, immunohistochemistry as well as other peroxidase-antibody based enzymatic assays.

Storage/Stability

Stable for 1 year at -20°C from shipment date. Briefly centrifuge the original vial prior to opening the cap to collect the product at the bottom of the tube.

Starting dilution recommendations

Working dilution recommended for ELISA is 1:12,000 – 60,000. Optimal working dilution for other application should be determined by the user.

GeneCopoeia, Inc.
9620 Medical Center Drive, #101
Rockville, Maryland 20850
Tel: 301-762-0888 Fax: 301-762-8333
Email: inquiry@genecopoeia.com
Web: www.genecopoeia.com