

Anti-MOUSE, Biotin Conjugated

Goat Polyclonal Antibody

Catalog # M32-62BG

Lot # B3216-2

Cited Applications

WB, ELISA, IHC

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the IgG (H&L)

Cross Reactivity

Mouse

Host/Isotype/Clone#

Goat, IgG

Immunogen

The antibody was produced against mouse IgG whole molecule in goat.

Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH7.2 + 0.01% (w/v) sodium azide.

Stability

1yr at -20°C from date of shipment

Scientific Background

Anti-Mouse IgG (H&L) biotin conjugated antibody generated in goat detects specifically mouse IgG (H&L). This secondary conjugated antibody anti-mouse is ideal for investigators who routinely perform titration assays, western-blot, immunoprecipitation and more generally immunoassays.

Anti-MOUSE, Biotin Conjugated

Goat Polyclonal Antibody

Catalog #	M32-62BG
Lot #	B3216-2
Purification	Affinity chromatography
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.