

Anti-GOAT, HRP Conjugated

Chicken Polyclonal Antibody

Catalog # G34-62DC

Lot # B3216-1

Cited Applications

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

Specificity

Recognizes IgG (H&L)

Cross Reactivity

Human

Host/Isotype/Clone#

Chicken, IgG

Immunogen

The antibody was produced against goat IgG whole molecule

Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH7.2 + 0.01% (w/v) Gentamicin Sulfate. Do NOT add Sodium Azide!

Stability

1yr at -20°C from date of shipment

Anti-GOAT, HRP conjugated

Chicken Polyclonal Antibody

Catalog #	G34-62DC
Lot #	B3216-1
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.