

Anti-Mouse HRP

Goat Polyclonal Antibody

Catalog # **G32-62G-1000**

Lot # X574-6

Cited Applications

- Western blot (1:2000) - ELISA
(2% skim milk blocking solution is recommended)

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

Specificity

Recognizes mouse immunoglobulins

Cross Reactivity

- Western blot using mouse primary antibodies

Other immunoglobulins may also be detectable

Host

Goat

Immunogen

Affinity purified IgG from mouse

Formulation

PBS pH 7.2, 50% Glycerol

Stability

Store at 4°C for several months and at -20°C for longer periods; do not add NaN₃ as this will inhibit the HRP enzyme. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

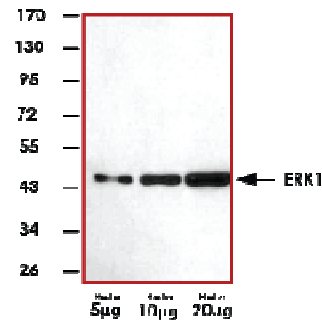
Scientific Background

Affinity purified mouse immunoglobulins were used to raise goat antibodies specific to the mouse IgG. The affinity purified antibodies are then conjugated to horse radish peroxidase (HRP) by reductive amination in a 2:1 molar ratio. The HRP conjugated secondary antibody can recognize the primary immunoglobulin and give a signal in the presence of a chemiluminescent substrate. The signal can be detected using an electronic imager or autoradiography film (1).

References

1. Nakane PK, Pierce GB: Enzyme-labeled antibodies: preparation and application for localization of antigens, *J Histochem Cytochem* 1966, 14:929

Sample Data



Representative western blot with a mouse monoclonal ERK1 antibody and Anti-Mouse HRP(1:2000) using 5 µg, 10 µg, and 20 µg HeLa cell lysate.

Anti-Mouse HRP

Goat Polyclonal Antibody

Catalog Number	G32-62G-1000
Quantity	1ml
Specific Lot Number	X574-6
Purification	Affinity purified
Format	1mg in 1 ml
Concentration	1.0 µg/µL
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.