

Anti-Mouse HRP

Goat Polyclonal Antibody

Catalog # **G32-62G**

Lot # E4368-6

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.

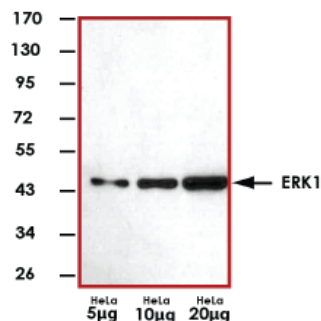
Scientific Background

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 #	G32-62G
Lot #	E4368-6
Purification	Affinity purified
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.

MATERIAL SAFETY DATA SHEET

Article 1 - Product Identification and Use

Product Name: Anti-Mouse HRP

Catalog # G32-62G-1000

This product is sold only for research use by qualified laboratory personnel, and is not to be used as a drug, medical device, food additive, cosmetic, nor household chemical. It is not to be used in diagnostic, therapeutic, consumer, agricultural, nor pesticidal applications.

Manufacturer's Name: SignalChem Pharmaceuticals Inc.
Street Address: 550-5600 Parkwood Way
City, Prov. Postal Code: Richmond, BC, V6V 2M2
Fax: 604-232-4601
EMERGENCY PHONE: 604-232-4600

Article 2 - Hazardous Ingredients

NOT AVAILABLE. We are not aware of any hazards associated with this product or its ingredients, but the chemical, physical, and toxicological properties of this product have not been investigated thoroughly. Observe normal laboratory precautions.

Article 3 - Physical Data

This product consists of purified antibody in PBS buffer shipped on ice packs. The physical properties of this product have not been investigated thoroughly.

Article 4 - Fire and Explosion Hazard

NOT APPLICABLE

Article 5 - Reactivity Data

NOT APPLICABLE

Article 6 - Toxicologically Data

May be harmful by inhalation, ingestion, or skin absorption. The toxicological properties of this product have not been investigated thoroughly. Exercise due caution.

Article 7 - Preventative Measures

Wear chemical safety goggles and compatible chemical-resistant gloves. Avoid inhalation, contact with eyes, skin or clothing.

*****MULTIPLE COMPONENT SPILL OR LEAK PROCEDURES*****

- Wear protective equipment.
- Absorb on sand or vermiculite and place in closed containers for disposal.
- Observe all federal, state and local environmental regulations.

Article 8 - First Aid Measures

- If swallowed, wash out mouth with water, provided person is conscious. Call a physician.
- In case of skin contact, flush with copious amounts of water for at least 15 minutes. Remove contaminated clothing and shoes. If a rash or other irritation develops, call a physician.
- If inhaled, remove to fresh air. If breathing becomes difficult, call a physician.
- In case of eye contact, flush with copious amounts of water for at least 15 minutes while separating the eyelids with fingers. Call a physician.

Article 9 - Preparation

Prepared By: Jun Yan

Phone #: 1-866-954-6273

The above information is believed to be correct but does not purport to be all-inclusive and shall be used only as a guide. SignalChem shall not be held liable for any damage resulting from handling or from contact with the above product. See the Technical Specification, Packing Slip, Invoice, and Product Catalogue for additional terms and conditions of sale.

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.