

Product Information

HRP Goat Anti-Rabbit IgG (H+L)

Cat. No.	Unit Size
L153A	50 µg
L153B	250 µg

Storage upon receipt:

- 2-8 °C
- Protect from light

Physical Form: Lyophilized powder

Reconstitution and Storage: Store lyophilized powder at 2-8°C. When ready to use, rehydrate with dH₂O (50 µL for 50 µg antibody or 250 µL for 250 µg antibody) to make 1 mg/mL solution and centrifuge if not clear. Product is stable for about 6 months at 2-8°C as an undiluted liquid. Prepare working dilution fresh each day. For extended storage after rehydration, add an equal volume of glycerol for a final concentration of 50%, and store at -20 °C as a liquid.

Warning: Confirm that glycerol is ACS grade or better before adding to this product. Lower grades of glycerol may seriously inhibit peroxidase enzyme activity.

Antibody Concentration: 1 mg/mL.

Buffer: 10 mM Sodium Phosphate, 250 mM NaCl, pH 7.5, 15 mg/mL BSA.

Peroxidase: Purified from horseradish root.

Purity: The antibody was purified from antisera by immunoaffinity chromatography using antigens coupled to agarose beads.

Antibody Specificity: Based on immunoelectrophoresis and/or ELISA, the antibody reacts with whole molecule rabbit IgG. It also reacts with the light chains of other rabbit immunoglobulins. No antibody was detected against non-immunoglobulin serum proteins. The antibody may cross-react with immunoglobulins from other species.

Suggested Dilution Range:

1:500 – 1:5,000 for immunohisto/cytochemistry

1:5,000 – 1:100,000 for ELISA and Western blotting with chromogenic substrates

1:10,000 – 1:200,000 for Western blotting with ECL substrates

Country of Origin: USA

Note: For in vitro research use only, not for diagnostic or therapeutic use.