

Product Information

HRP Goat Anti-Human IgG (H+L)

Cat. No.	Unit Size
L154A	50 µL
L154B	250 µL

Storage upon receipt:

- -20 °C
- Protect from light
- Avoid freeze-thaw cycles

Antibody Concentration: 0.8 mg/mL.

Buffer: 10 mM Sodium Phosphate, 250 mM NaCl, 50% glycerol, 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 human IgG. It also reacts with the light chains of other human 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.