



Goat Anti-Mouse IgG, Fcy Fragment Specific conjugated to HRP

Product Number HRP-112-Fc
Lot # XXX
Amount 1.0 mg
Store at 2-8°C

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at 2-8°C in the dark, do not freeze. Proteins are sensitive to freeze-thaw cycles.

Handling

Reconstitute to 1.1 ml with distilled deionized water vortex gently and allow vial to sit on ice for 30 minutes. We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light.

Buffer

Upon rehydration with 1.1 ml distilled deionized water; The product is in 10 mM sodium phosphate (pH 7.2), 150 mM sodium chloride, 1.0% BSA, 0.1% Kathon CG

Stability

Lyophilized material is stable for up to one year. After product has been reconstituted, product should be stored at 2-8°C in the dark. For long term storage after reconstitution, dilute with 50% glycerol and store at -20 °C as a liquid. Prepare working dilution prior to use and then discard..

Note

For research use only, not for diagnostic or therapeutic use.