



Mouse Anti-DYKDDDDK IgG conjugated to HRP

Product Number	HRP-1718
Lot #	XXXX
Amount	100 µg
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 1.0 ml with distilled deionized water vortex gently and allow vial to sit on ice for 20 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 1.0 ml distilled deionized water; The product is in 100 mM NaPO₄ (pH 7.4) + 100 mM NaCl + 1.5% BSA, 0.01% Thimerosal + 50mM Sucrose

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 and be used within 6 months. If further dilution of the conjugate is required, use diluted material within one week.

Note

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

Concentration

After reconstitution - 0.1 mg/mL

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415
www.columbiabiosciences.com
For technical inquiries: info@columbiabiosciences.com
For sales inquiries: sales@columbiabiosciences.com