

Mouse anti-HA.11 (haemagglutinin) IgG conjugated to Alkaline Phosphatase

Product Number AP-1722 Lot Number XXXX Amount 100 µg Store at 4°C

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at 4° C in the dark. Conjugates are sensitive to freeze-thaw cycles; after reconstitution, store at $2-8^{\circ}$ C in the dark, do not freeze.

Handling

Prior to use, reconstitute to 1 mL with ddH_2O providing a concentration of 100 μ g/mL, vortex gently and allow material to reconstitute for 20 minutes. Avoid exposure to heat.

Buffer

Upon rehydration with 1 mL of ddH_2O , the product is in 25mM Tris (pH 7.4), 150mM sodium chloride, 5 mM $MgCl_2$, 0.12 mM $ZnCl_2$, and 2mM sodium azide as a preservative.

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.

Antibody/Antigen Information

Mouse IgG₁ clone 16B12 versus sequence CYPYDVPDYASL. Recognizes HA epitopes located in the middle of protein sequences as well as at the N- or C-terminus.

Note

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