



## Mouse anti-DYKDDDDK IgG conjugated to Alkaline Phosphatase

Product Number AP-1718  
Lot Number AP013-19-002  
Amount 100 µg  
Store at 4°C

### Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at 4°C in the dark. Phycobiliproteins are sensitive to freeze-thaw cycles; after reconstitution, store at 2-8°C in the dark, do not freeze.

### Handling

Prior to use, reconstitute to 1 ml with ddH<sub>2</sub>O, vortex gently and allow material to reconstitute for 20 minutes. Avoid exposure to heat.

### Buffer

Upon rehydration with 1 ml of ddH<sub>2</sub>O, the product is in 25mM Tris (pH 7.2), 150mM sodium chloride, 5 mM MgCl<sub>2</sub>, 0.12 mM ZnCl<sub>2</sub> 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 IgG<sub>1</sub> clone M2.

### Note

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