



Mouse anti-6xHistidine IgG conjugated to Alkaline Phosphatase

Product Number AP-1711
Lot Number AP010-21-006
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, do not freeze.

Handling

Avoid exposure to heat.

Buffer

After reconstitution with 1 mL ddH₂O, the product is in 25mM Tris (pH 7.4), 150mM sodium chloride, 5 mM MgCl₂, 0.12 mM ZnCl₂, and 2mM sodium azide.

Stability

Product should be stored at 4°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 AD1.1.10.

Note

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

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com