



Streptavidin conjugated to Alkaline Phosphatase

Product Number AP-2212
Lot Number XXXX
Amount 100 µg
Store at 4°C

Form/Shipping

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°C in the dark, do not freeze.

Suggested Dilution Range

1:5000-1:50,000 for ELISA and Western blotting

Handling

Prior to use, reconstitute to 1 mL with ddH₂O providing a concentration of 100 µg/mL, vortex gently and allow material to reconstitute for 20 minutes. Avoid exposure to heat.

Buffer

Upon rehydration with 1 mL of ddH₂O, the product is in 25mM Tris (pH 7.4), 150mM sodium chloride, 5 mM MgCl₂, 0.12 mM ZnCl₂, 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. For extended storage, add an equal volume of glycerol (ACS grade) for a final concentration of 50%, and store at -20 degrees Celsius. Prepare a working dilution daily.

Country of Origin

USA.

Antibody/Antigen Information

Biotin-Streptavidin

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