



Goat anti-Rabbit IgG conjugated to Alkaline Phosphatase

Product Number AP-114
Lot Number AP004-21-003
Amount 1.0 mg
Store at 4°C

Form/Shipping & Storage

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

Handling

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.

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.

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com