



Rabbit Anti-T7 IgG conjugated to PerCP

Product Number D11-1832
Lot Number PerCP015-12-001
Amount 100 µg total protein

Form/ Storage

Conjugate supplied as a lyophilized powder. Upon receipt, store at 2-8°C in the dark.

Handling

Avoid exposure to heat and light. Prior to use reconstitute to 1 ml with distilled deionized water, vortex and allow it to sit on ice for 20 minutes.

Buffer

Upon reconstitution, the product is in 100 mM sodium phosphate (pH 7.4), 50 mM sucrose, 100 mM sodium chloride, 0.1% BSA as a stabilizer, and 2 mM sodium azide as a preservative. The concentration of the conjugate is 100 µg/ml.

Stability

Lyophilized material is stable for one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 3 months.

Note

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

Spectral Characteristics

Visible absorption maxima 483 nm
Emission maximum 673 nm

Concentration (after reconstitution)

0.1 mg/mL
anti-T7 0.443 µM
PerCP 1.16 µM

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com
For sales inquiries: sales@columbiabiosciences.com