

**Rabbit Anti-S tag conjugated to PerCP**

Product Number D11-1831  
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, 150 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.

Antigen Info

Antibody was raised against S-tag with amino acid sequence KETAAAKFERQHMS.

**Spectral Characteristics**

Visible absorption maxima	483
Emission maximum	673 nm

**COLUMBIA BIOSCIENCES**