



Rabbit Anti-HA (haemagglutinin) conjugated to SureLight® APC

Product Number D3-1830
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 10 mM Tris (pH 8.2), 50 mM sucrose, 150 mM sodium chloride, 0.1% BSA as a stabilizer, and 1 µg/mL pentachlorophenol as a preservative. The concentration of the conjugate is 100 µg/ml.

Stability

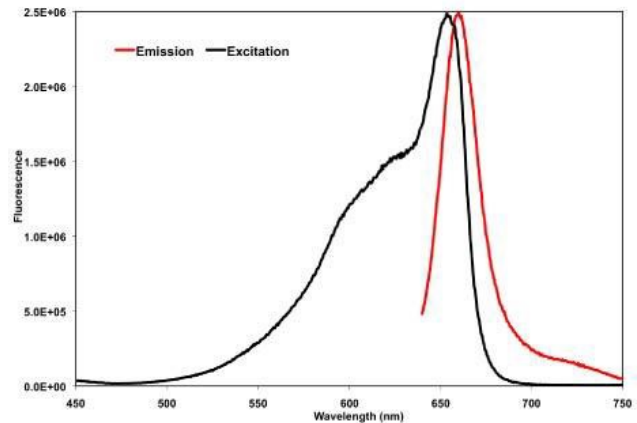
Lyophilized material is stable for one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 3 months.

Antigen Info

Antibody was raised against HA-tag with amino acid sequence CYPYDVPDYASL.

Note

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



Fluorescence excitation and emission spectra of APC in 100 mM sodium phosphate (pH 7.2) + 1 mM EDTA and 1 mM sodium azide. The emission scan was taken with excitation at 630 nm. The excitation scan was taken with an emission at 660 nm. The curves were normalized to equalize peak heights.

Spectral Characteristics

Visible absorption maxima 650
Emission maximum 656 nm

Concentration 0.1 mg/mL
(after reconstitution)

SureLight® APC 0.457 µM
Anti-HA 0.384 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