



Rabbit anti-HA (Hemagglutinin) IgG conjugated to R-Phycoerythrin

Product Number D5-1830
Amount 100 µg total protein
Store at 4°C

Form/ Storage

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

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.

Stability

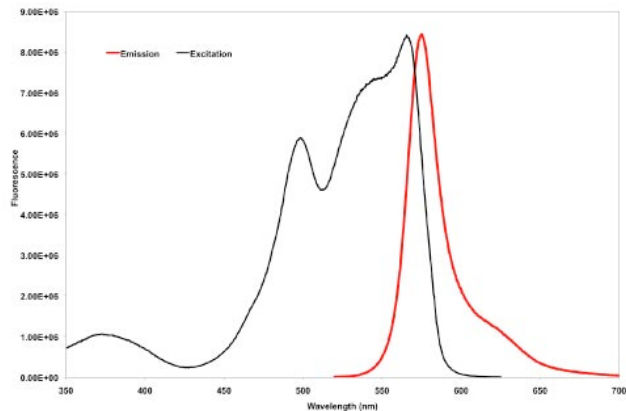
Lyophilized material is stable for one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 3 months.

Note

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

Antigen Info

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



Fluorescence excitation and emission spectra of R-phycoerythrin in 100 mM sodium phosphate (pH 7.2) + 1 mM EDTA and 1 mM sodium azide. Emission scan was taken with excitation at 498 nm. Excitation scan was taken with emission at 575 nm. Scans were normalized to equalize peak heights.

Spectral Characteristics

Visible absorption maxima 565>540>498
Emission maximum 578 nm

Concentration

After reconstitution to 1.0 ml

0.1 mg/mL

Fluor : Protein ~2:1

