



Rabbit Anti-HA (haemagglutinin) conjugated to PerCP

Product Number D11-1830
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 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

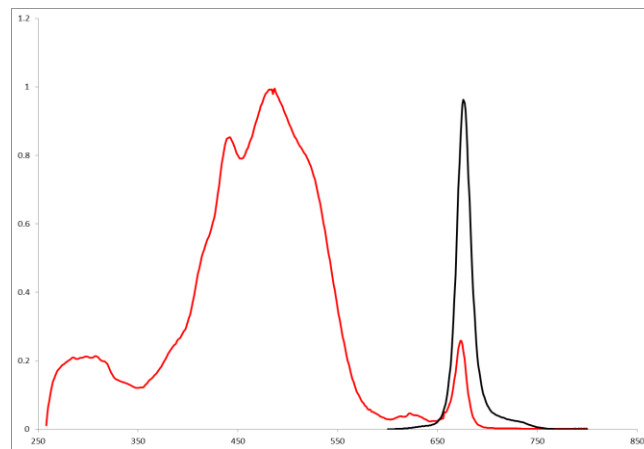
Lyophilized material is stable for one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 3 months.

Note

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

Antigen Info

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



Spectral Characteristics

Visible absorption maxima 483
Emission maximum 673 nm

Concentration 0.1 mg/mL
(after reconstitution)

References:

Nygard NR, Bono C, Brown LR, Gorka J, Giacometto KS, Schaiff WT, Graham MB, McCourt DW, Kabeer M, Braciale VL, et al. Antibody recognition of an immunogenic influenza hemagglutinin-human leukocyte antigen class II complex. J Exp Med. 1991 Jul 1;174(1):243-51.

Rogers BE, Chaudhuri TR, Reynolds PN, Della Manna D, Zinn KR. Non-invasive gamma camera imaging of gene transfer using an adenoviral vector encoding an epitope-tagged receptor as a reporter. Gene Ther. 2003 Jan;10(2):105-14. PubMed PMID: 12571639.



COLUMBIA BIOSCIENCES