



Mouse anti-HA.11 (haemagglutinin) IgG conjugated to Biotin

Product Number B1-1722
Amount 100 µg
Store at 4°C

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at 4°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

Prior to use, reconstitute to 1 ml with 50% glycerol, vortex gently and allow material to reconstitute for 20 minutes. Avoid exposure to heat.

Buffer

Upon rehydration with 1 ml of 50% glycerol, the product is in 25mM Tris (pH 7.2), 150mM sodium chloride, 50 mM sucrose and 2mM sodium azide as a preservative.

Stability

Lyophilized material is stable for up to one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 6 months. If further dilution of the conjugate is required, use diluted material within one week.

Antibody/Antigen Information

Mouse IgG₁ clone 16B12 versus sequence CYPYDVPDYASL. Recognizes HA epitopes located in the middle of protein sequences as well as at the N- or C-terminus.

Note

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

References:

Nygaard 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.

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415
www.columbiabiosciences.com
For technical inquiries: info@columbiabiosciences.com
For sales inquiries: sales@columbiabiosciences.com