



Mouse anti-HA (hemagglutinin) tag (YPYDVPDYA) IgG conjugated to Dylight® 650

Product Number	D10-1722
Lot Number	XXXX
Approximately	5-7
Amount	100 µg
Clone	16B12
Concentration	1.0 mg/mL
Dylight 650:α-HA	2.61
Store at	2-8°C

Form/Shipping & Storage

Supplied as a liquid. Upon receipt, store at 2-8°C in the dark, do not freeze.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light.

Buffer

The product is in 100 mM sodium phosphate (pH 7.4), 100 mM sodium chloride and 2 mM sodium azide.

Note

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

Spectral Characteristics

Visible absorption maxima	655
Emission maximum	672

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.