



## Anti-HA (hemagglutinin) IgG conjugated to SureLight®-Allophycocyanin

Product Number D3-1722-100, D3-1722-1MG  
Amount 100 µg, 1 mg  
Store at 2-8°C

### Form/Shipping & Storage

Supplied lyophilized. Upon receipt, store at -20 °C. Reconstitute with 1 ml of ddH<sub>2</sub>O and store at 2-8 °C. Phycobiliproteins are sensitive to freeze-thaw cycles.

### Handling

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

### Buffer

Upon reconstitution, the product is in 10 mM tris (pH 8.2), 150 mM NaCl, 1 µg/mL pentachlorophenol, 0.1% BSA and 50mM Sucrose.

### Stability

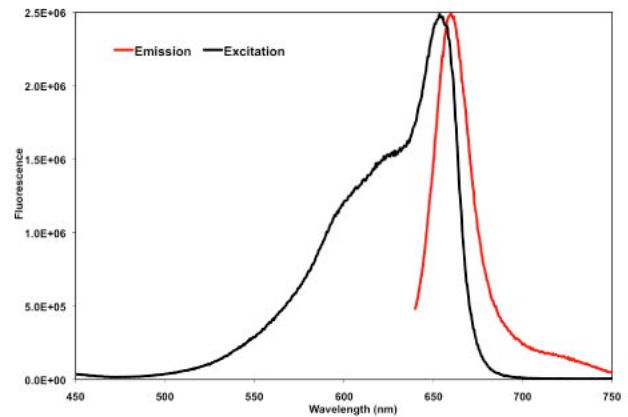
Product should be stored at 2-8°C in the dark and be used within 3 months. If further dilution of the conjugate is required, use diluted material within one week.

### Note

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

### Clone

Mouse monoclonal anti-HA (Hemagglutinin) IgG1, clone 16B12



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.

### Specificity

Monoclonal antibody HA.11 was raised against the twelve amino acid peptide CYPYDVPDYASL. The HA.11 antibody recognizes the influenza hemagglutinin epitope (YPYDVPDYA) which has been used extensively as a general epitope tag in expression vectors. The HA.11 antibody recognizes HA epitopes located in the middle of protein sequences as well as at the N- or C-terminus.

### Spectral Characteristics

Visible absorption maxima	652
Emission maximum	657

Fluor : Protein	~1:1
-----------------	------



**COLUMBIA BIOSCIENCES**