



Rabbit Anti-Haemagglutinin (HA) IgG conjugated to Dylight® 550

Product Number	D9-1830
Lot#	XXXX
Approximately	5-7
Amount	100 µg
Concentration	1.0 mg/mL
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, 2mM sodium azide as a preservative.

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

Spectral Characteristics

Visible absorption maxima	557
Emission maximum	576
F/P Ratio	5.70

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

www.columbiabiosciences.com
For technical inquiries: info@columbiabiosciences.com
For sales inquiries: sales@columbiabiosciences.com