



Mouse anti-6His IgG conjugated to Dylight® 550

Product Number	D9-1711
Lot#	XXXX
Approximately	5-7
Conc	1.00 mg/mL
Amount	100 µg total protein
Clone	AD1.1.10
Store at	2-8°C

Form/Shipping & Storage

Supplied liquid. Upon receipt, store at 2-8 °C. Fluorescent dyes are sensitive to light. Please store in the dark and avoid exposure to sunlight.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application.

Buffer

0.1M sodium phosphate, 0.1M NaCl, pH7.4, 2 mM NaN₃

Stability

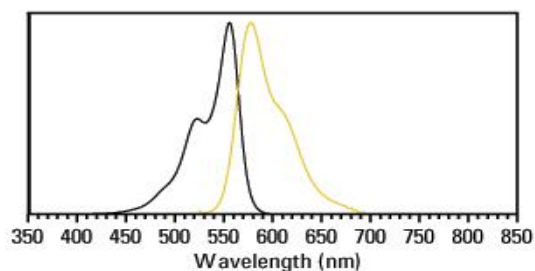
Product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

Note

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

Spectral Characteristic

Visible absorption maxima	562 nm
Emission maximum	576 nm
F/P Ratio	4.07



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