



### Mouse anti-C-Myc IgG conjugated to Dylight® 550

Product Number	D9-1714
Lot#	XXXX
Approximately	5-7
Amount	100 µg
Concentration	0.514 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. 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 rehydration with 1.0 ml distilled deionized water; 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	655
Emission maximum	672
F/P Ratio	4.66

4985 Winchester Blvd. Frederick, MD 21703  
301.732.5415

[www.columbiabiosciences.com](http://www.columbiabiosciences.com)

For technical inquiries: [info@columbiabiosciences.com](mailto:info@columbiabiosciences.com)

For sales inquiries: [sales@columbiabiosciences.com](mailto:sales@columbiabiosciences.com)