



Goat anti-Mouse IgG (H&L) conjugated to Surelight™ 488

Product Number	D44-112	<u>Spectral Characteristics</u>	
Lot Number	SL488044-18-001	Visible absorption maxima	496
Concentration	4.40 mg/mL	Emission maximum	525
Amount	0.5 mg		
F:P	8.91		
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.

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.

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



COLUMBIA BIOSCIENCES

4985 Winchester Blvd. Frederick, MD 21703

310.732.5336 www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com