



Anti-6xHistidine IgG conjugated to Dylight® 488

Product Number	D8-1711
Lot	XXXX
Amount	100 µg antibody
Concentration	1.0 mg/mL
Dylight 488/Antibody ratio	6.19
Store at	2-8°C

Form/Shipping & Storage

Supplied as a 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, 0.10M NaCl, pH7.4, 2 mM NaN₃

Stability

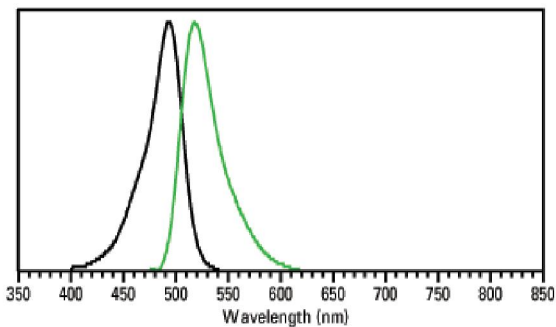
Product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

Note

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

Spectral Characteristic

Visible absorption maxima	488 nm
Emission maximum	518 nm



References

Puertollano R, Alonso MA. MAL, an integral element of the apical sorting machinery, is an itinerant protein that cycles between the trans-Golgi network and the plasma membrane. *Mol Biol Cell*. 1999 Oct;10(10):3435-3447.

Roy SF, Laporte SA, Escher E, Leduc R, Guillemette G. Epitope tagging and immunoreactivity of the human angiotensin II type 1 receptor. *Can J Physiol Pharmacol*. 1997 Jun;75(6):690-695.

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