



Anti-DYKDDDDK conjugated to Dylight® 650

Product Number	D10-1718
Lot Number	XXXX
Approximately	5-7
Amount	100 µg total protein
Concentration	1.0 mg/ml
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, 0.1M NaCl, pH7.4, 2mM 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	655 nm
Emission maximum	672 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