



Rabbit Anti-T7 conjugated to Dylight® 488

Product Number D8-1832
Lot# XXXX
Approximately 5-7
Dye:Ab ratio 3.49
Conc 1.0 mg/ml
Amount 100 µg total protein

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, 2 mM NaN₃

Stability

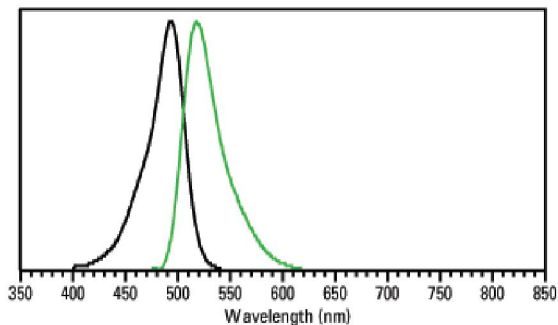
Product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

Note

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

Spectral Characteristic

Visible absorption maxima 488 nm
Emission maximum 518 nm



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