



**Goat Anti-GST conjugated to Dylight® 488**

Product Number D8-1310  
Lot# XXXX  
Approximately 5-7  
Conc 1.0 mg/mL  
F/P 5.08  
Amount 100 µg total protein  
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 NaN<sub>3</sub>

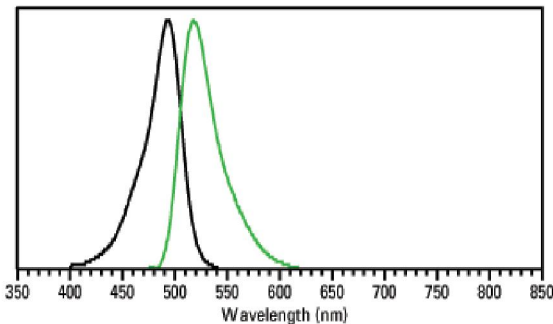
Stability

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.

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