



### Anti-GST conjugated to Dylight® 650

Product Number D10-1310  
Conc: 2.54 mg/mL  
Lot# XXXX  
Approximately 5-7  
Amount 100 µg total protein  
Store at 2-8°C

#### Immunogen

Glutathione-S-transferase from *S. japonicum*

#### 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

100 mM NaPO<sub>4</sub> (pH 7.4), 100 mM NaCl, 2mM Sodium Azide

#### Stability

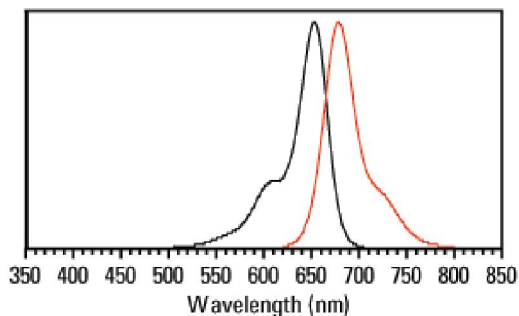
Product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

#### Note

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

#### Spectral Characteristic

Visible absorption maxima 652 nm  
Emission maximum 672 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