



## Europium Chelate W-1024 conjugated to Goat anti-GST IgG (Glutathione-S-Transferase)

Product Number D17-1310-50  
Lot Number Eu1024011-14-001  
Amount 50 µg  
Concentration 1 mg/mL  
Eu1024/Ab 11.6  
Store at 2-8°C

### Form/Shipping & Storage

Supplied as liquid. Upon receipt, store at 2-8 °C.  
DO NOT FREEZE.

### Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light.

### Buffer

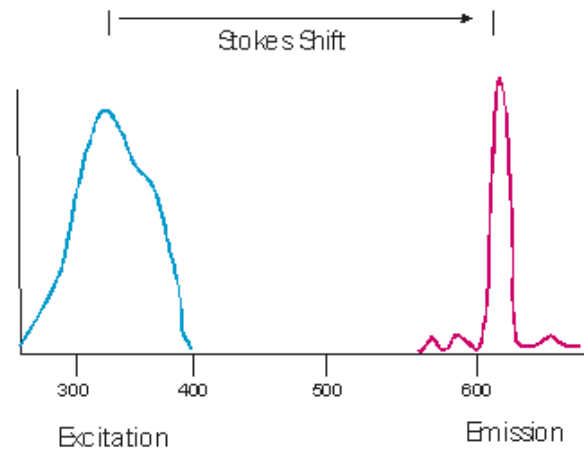
The product is in 50 mM Tris (pH 7.8) + 0.9% NaCl + 0.05% NaN<sub>3</sub>.

### Stability

Product should be stored at 2-8°C in the dark and used within 6 months.

### Note

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



### Spectral Characteristics

Visible absorption maxima	340
Emission maximum	615

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

[www.columbiabiosciences.com](http://www.columbiabiosciences.com)

For technical inquiries: [info@columbiabiosciences.com](mailto:info@columbiabiosciences.com)

For sales inquiries: [sales@columbiabiosciences.com](mailto:sales@columbiabiosciences.com)