



## Europium Chelate W-1024 conjugated to Goat anti-Mouse IgG

Product Number D17-112-20, D17-112-50,  
D17-112-1MG  
Amount 20 µg, 50 µg, 1 mg  
Concentration ≈ 1mg/mL  
Store at 2-8°C  
EU/Ab 5-9 per Ab

### 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

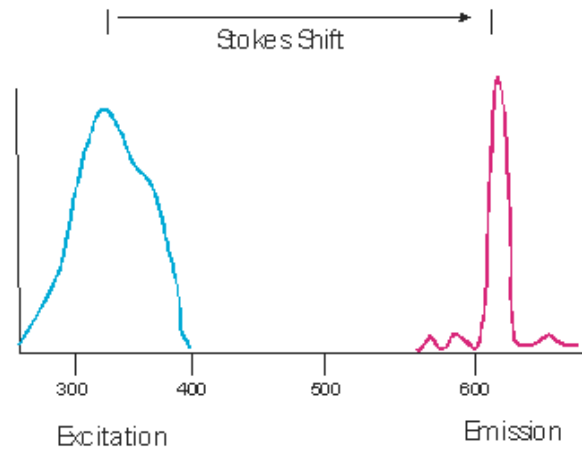
The product is in 50 mM Tris (pH 7.8) + 0.9% NaCl + 0.05% NaN<sub>3</sub>.

### Stability

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

### Note

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



### Spectral Characteristics

Visible absorption maxima	340
Emission maximum	615



**COLUMBIA BIOSCIENCES**