



## Europium Chelate W-1024 conjugated to Mouse Anti-c-myc IgG

Product Number D17-1714-20, D17-1714-50,  
D17-1714-1MG  
Amount 20  $\mu$ g, 50  $\mu$ g, 1 mg  
Concentration  $\approx$  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

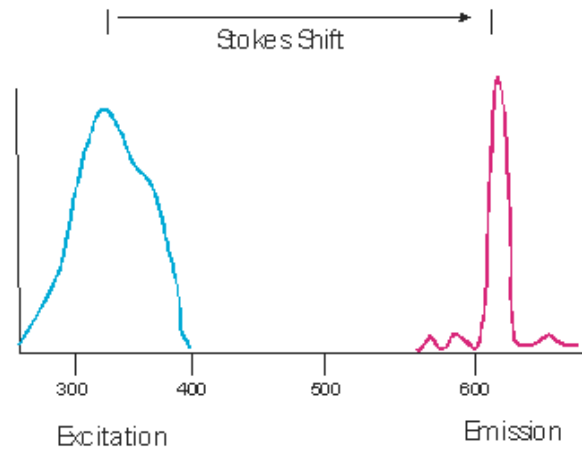
The product is in 50 mM Tris (pH 7.8) + 0.9% NaCl + 0.05% NaN<sub>3</sub>.

### 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**