



Europium Chelate W-1024 conjugated to Goat anti-Rabbit IgG

Product Number D17-114-1MG
Lot# Eu1024004-19-002
Amount 1 mg
Concentration 1.0 mg/mL
Eu/Ab 7.8
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

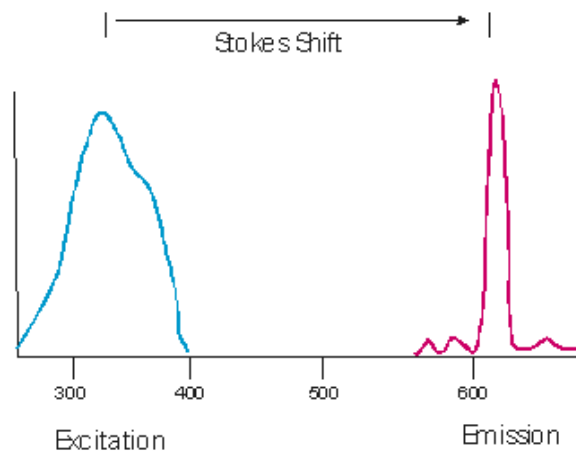
The product is in 50 mM Tris (pH 7.8) + 0.9% NaCl + 0.05% NaN₃.

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

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com