



Europium Chelate W-1024 conjugated to Anti-6xHistidine IgG (clone AD1.1.10)

Product Number D17-1711
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

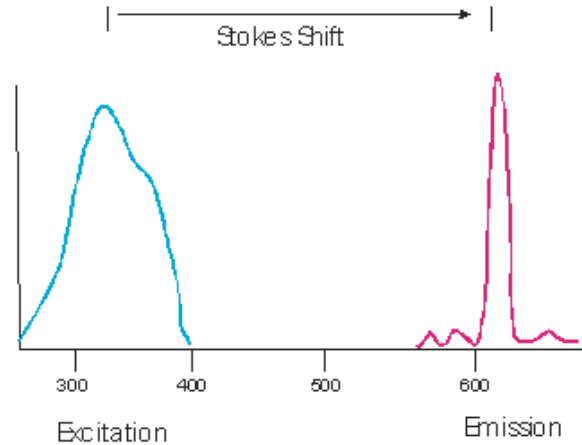
The product is in 50 mM Tris (pH 7.8) + 0.9% NaCl + 0.05% NaN₃.

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



COLUMBIA BIOSCIENCES