



Goat anti-Mouse IgG conjugated to DyLight® 650

Product Number D10-112-1MG
Amount 1 mg total protein
Concentration 1mg/ml
Store at 2-8°C

Form/Shipping & Storage

Supplied as a liquid. Upon receipt, store at 2-8 °C. Fluorescent dyes are sensitive to light. Please store in the dark and avoid exposure to sunlight.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application.

Buffer

100 mM NaPO₄ (pH 7.4), 100 mM NaCl, 2 mM Sodium Azide

Stability

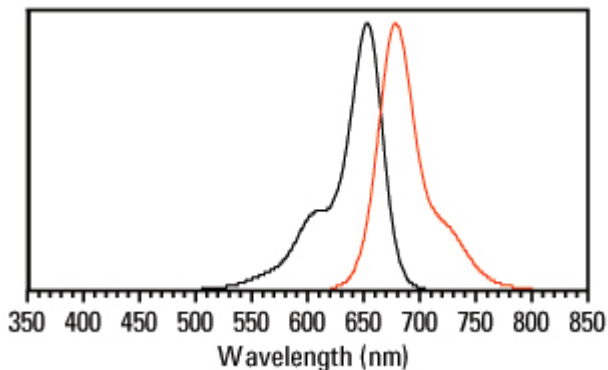
Product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

Note

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

Spectral Characteristic

Visible absorption maxima 652 nm
Emission maximum 672 nm



COLUMBIA BIOSCIENCES