



Rabbit anti-Goat IgG conjugated to Dylight® 488

Product Number D8-116-1MG
Conc 1 mg/mL

Amount 1mg total protein
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

0.1M sodium phosphate, 0.15M NaCl, pH 7.4, 1% BSA, 0.02% NaN₃

Stability

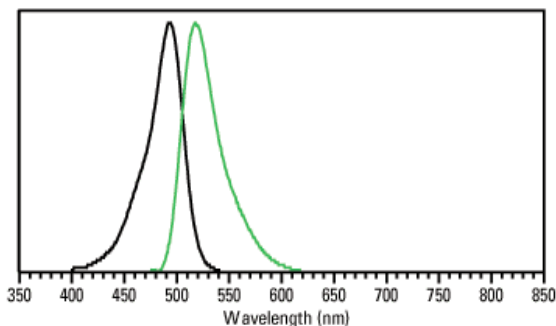
Product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

Note

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

Spectral Characteristic

Visible absorption maxima 488 nm
Emission maximum 518 nm



COLUMBIA BIOSCIENCES