



## Goat Anti-Human IgG conjugated to Dylight® 488

Product Number D8-110-1MG  
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% NaN<sub>3</sub>

### Stability

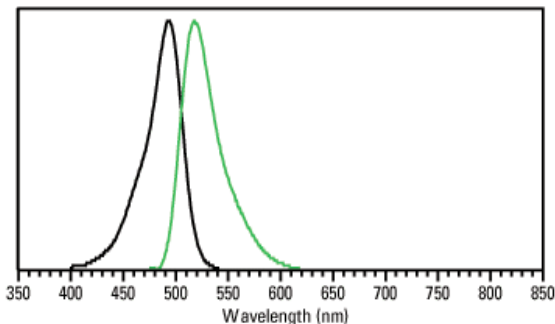
Product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

### Note

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

### Spectral Characteristic

Visible absorption maxima 488 nm  
Emission maximum 518 nm



**COLUMBIA BIOSCIENCES**