



## Goat anti-Human IgG conjugated to DyLight® 650

Product Number D10-110-1MG  
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 NaPO<sub>4</sub> (pH 7.4), 100 mM NaCl, 2 mM Sodium Azide

### Stability

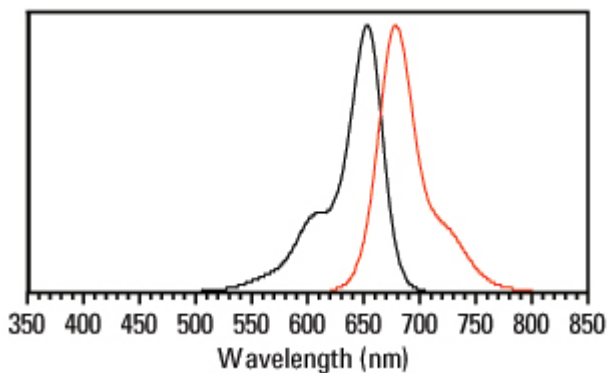
Product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

### Note

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

### Spectral Characteristic

Visible absorption maxima 652 nm  
Emission maximum 672 nm



**COLUMBIA BIOSCIENCES**