



## Goat anti-Rabbit IgG conjugated to Dylight® 488

Product Number D8-114  
Lot Number DY488004-18-001  
Amount 1 mg  
Concentration 1.0 mg/mL  
F:P 5.42  
Store at 2-8°C

### Form/Shipping & Storage

Supplied as a liquid Upon receipt, store at 2-8°C in the dark, 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 100 mM sodium phosphate (pH 7.4), 100 mM sodium chloride and 2 mM sodium azide.

### Note

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

### Spectral Characteristics

Visible absorption maxima 488  
Emission maximum 518

### Note

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

