



Rabbit anti-V5 IgG conjugated to Dylight® 488

Product Number D8-1834
Lot Number XXXX
Approximately 5-7
Amount 100 µg
Concentration 1.00 mg/mL
Dylight 488:α-V5 4.12
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 as a preservative.

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.

