



Rabbit Anti-V5 IgG conjugated to Dylight® 650

Product Number	D10-1834
Lot#	XXXX
Approximately	5-7
Conc	1.0 mg/mL
Dye: Protein Ratio	2.25
Amount	100 µg total protein

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, 0.1M NaCl, pH7.4, 2 mM 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	650 nm
Emission maximum	672 nm