

Streptavidin conjugated to Dylight® 650

Product Number	D10-2212
Lot#	XXXX
Approximately	5-7
Amount	1.0 mg
Concentration	1.0 mg/mL
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.

<u>Buffer</u>

The product is in 100 mM sodium phosphate (pH 7.4), 100 mM sodium chloride, 2mM sodium azide as a preservative.

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

Spectral Characteristics

Visible absorption maxima	655
Emission maximum	672
F/P Ratio	2.84