

Mouse anti-6His IgG conjugated to Dylight® 650

Product Number	D10-1711
Clone	AD1.1.10
Lot#	XXXX
Approximately	5-7
Amount	100 µg
Concentration	1.00 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.

<u>Handling</u>

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	2.81