



Mouse anti-6His IgG conjugated to SureLight™ 488

Product Number	D44-1711
Lot Number	SL488010-19-001
Amount	100 µg
Concentration	1 mg/ml
F:P	4.24
Store at	2-8°C

Form/Shipping & Storage

Supplied as a liquid. Store material at 4-8C 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. The concentration of the conjugate is 1 mg/ml.

Stability

Product should be stored at 2-8C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one week.

Note

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

Spectral Characteristics

Visible absorption maxima	496nm
Emission maximum	518nm