



Goat anti-Human IgG F(ab')₂ fragment, conjugated to SureLight™ 488

Product Number D44-110-Fab
Lot# SL488061-19-0001
Amount 1 mg
Clone N/A
Store at 4°C

Form/Shipping&Storage

Supplied as a stock solution. Upon receipt, store at 4°C in the dark. Do not freeze.

Handling

Prior to use, dilute to the appropriate concentration in protein-containing buffered solution pH 7-8 (e.g. PBS + 0.5% BSA). Avoid exposure to heat and light.

Buffer

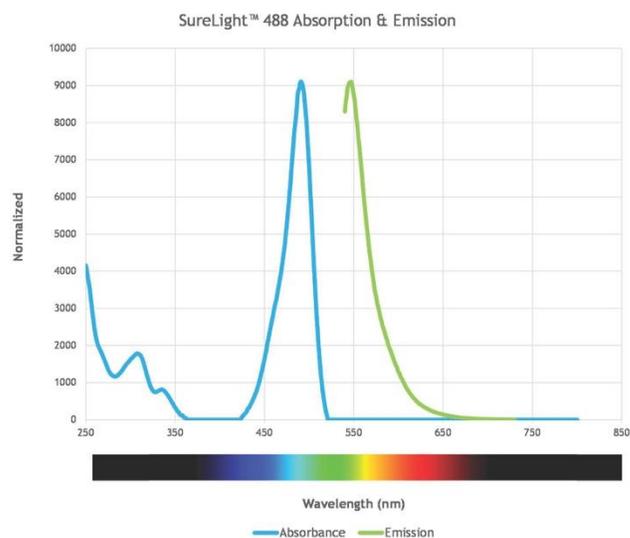
The product is in 100 mM sodium phosphate (pH 7.4), 100 mM sucrose, 150 mM sodium chloride, 0.1% BSA as a stabilizer, and 2 mM sodium azide as a preservative. The concentration is 1 mg/mL total protein.

Stability

The material is stable for up to one year if stored at 2-8°C in the dark. 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	494 ± 3 nm
Emission maximum	518 ± 4 nm

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com