



Goat anti-Rabbit IgG conjugated to SureLight® P3

Product Number D7-114
Amount 1 mg
Store at 4°C

Form/Shipping & Storage

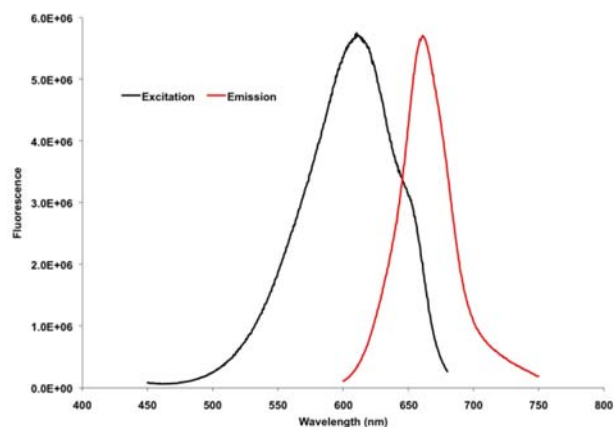
Supplied as a lyophilized powder. Upon receipt, store at 4°C in the dark. Phycobiliproteins are sensitive to freeze-thaw cycles; after reconstitution, store at 2-8°C in the dark, do not freeze.

Handling

Prior to use, reconstitute to 1 ml with distilled deionized water, vortex gently and allow vial to sit on ice for 20 minutes. Avoid exposure to heat and light.

Buffer

Upon rehydration with 1 ml distilled deionized water, the product is in 100 mM sodium phosphate (pH 7.4), 200 mM sucrose, 100 mM sodium chloride, 1 mg/ml BSA as a stabilizer, and 0.05% sodium azide as a preservative.



Stability

Lyophilized material is stable for up to one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 3 months. 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	614>650 (sh) nm
Emission maximum	662 nm



COLUMBIA BIOSCIENCES