



## SureLight® P3

Product Number D7-000  
Amount 5 mg  
Concentration >20 mg/mL  
Store at 4°C

### Form/Shipping & Storage

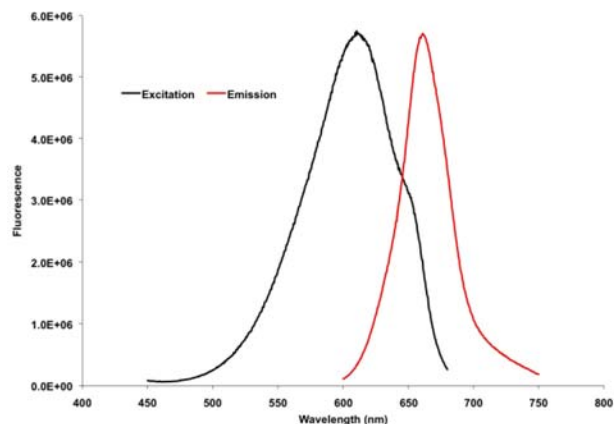
Supplied as ammonium sulfate precipitate. Upon receipt, store at 4°C in the dark. Phycobilisomes are sensitive to freeze-thaw cycles; do not freeze.

### Handling

Prior to use, dialyze into an appropriate physiological buffer. Avoid exposure to heat and light.

### Buffer

Ammonium sulfate (40%), 100 mM sodium phosphate (pH 7.4), and 2 mM sodium azide as a preservative.



### Stability

Material is stable for up to one year. Product should be stored at 2-8°C in the dark and be used within 12 months.

### Note

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

Approximate molecular weight 10,000 kDa  
AU/mg @ 615 nm 5.7

### Spectral Characteristics

Visible absorption maxima 614>650 (sh) nm  
Emission maximum 662 nm



**COLUMBIA BIOSCIENCES**