



## SureLight® PE-695 (Cyanine 5.5)

Product Number	D15-000
Lot Number	CY5.5RPE000-21-015
Concentration:	10.36 mg/mL
Store at	2-8° C

### Minimum Specifications

Emission Ratio 694/572	≥ 4.0
Concentration	10-11 mg/mL

### Lot Specific Values

Emission Ratio 694/572	9.95
Concentration	10.36 mg/mL

### Form/Shipping & Storage

Supplied as liquid. Upon receipt, store at 2-8°C. DO NOT FREEZE.

### Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. This product is light sensitive.

### Buffer

10mM Sodium Phosphate (pH 7.4) and 15mM sodium azide

### Stability

Product should be stored at 2-8°C in the dark

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

[www.columbiabiosciences.com](http://www.columbiabiosciences.com)

For technical inquiries: [info@columbiabiosciences.com](mailto:info@columbiabiosciences.com)

For sales inquiries: [sales@columbiabiosciences.com](mailto:sales@columbiabiosciences.com)