



## SureLight™ PerCP-695

Product Number	D27-000
Lot Number	Cy5.5PerCP000-20-017
Concentration	10.55 mg/ml
Expiration Date	8/12/2021
Store at	2-8° C

### Product Specifications

Emission Ratio 695/676	0.3 - 0.5
Absorbance Ratio 483/676	2.0 - 2.6
Concentration	≥10 mg/mL

### Lot Specific Values

Emission Ratio 695/676	0.31
Absorbance Ratio 483/676	2.53
Concentration	10.55 mg/mL

### Form/Shipping & Storage

Supplied as liquid. Upon receipt, store at 2-8° C. DO NOT FREEZE. PerCP is sensitive to freeze-thaw cycle.

### Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Light Sensitive. Please use an AU/mg of 11.6 at 483nm to determine PerCP concentration.

### Buffer

10mM sodium phosphate (pH 7.0) and 15mM sodium azide

### Note

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

### Method

The PerCP concentration was measured in 100 mM Sodium Phosphate, 2 mM Sodium Azide (pH7.4)

### 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)



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)