



SureLight® PE-777 (Phycoerythrin Cyanine 7.0)

Product Number	D16-000
Lot Number	CY7RPE000-21-020
Concentration:	10.75 mg/ml
Mfg. Date	10/05/2021
Exp. Date	10/05/2023
Store at	2-8° C

Minimum Specifications

Emission Ratio 776/572	0.5
Concentration	10 mg/mL

Lot Specific Values

Emission Ratio 776/572	0.725
Concentration	10.75 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. 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

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com