



R-Phycoerythrin

Product Number D5-000
Concentration 21.26 mg/mL
Store at 2-8°C

Form/Shipping & Storage

Supplied as an ammonium sulfate precipitate. Upon receipt, store at 2-8°C in the dark - do not freeze. Phycobiliproteins are sensitive to freeze-thaw cycles.

Handling

Avoid exposure to heat and light.

Buffer

60% ammonium sulfate in 100mM sodium phosphate + 100mM sodium chloride buffer pH 7.0 with 2mM-sodium azide.

Note

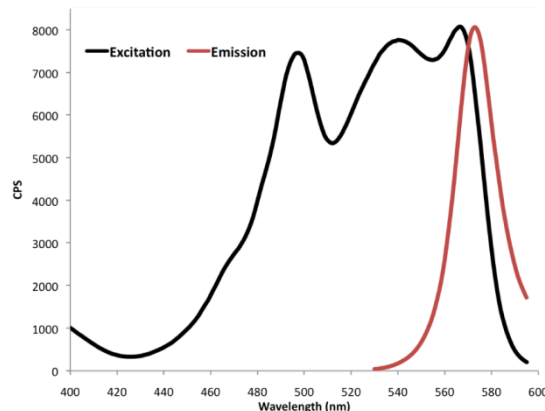
For research use only, not for diagnostic or therapeutic use.

Method

The material concentration was measured in 100 mM phosphate buffer pH 7.0 with 2mM sodium azide. The reported result is the average over all dilutions and all readings.

Specifications: _____ Actual Result:

A_{566} / A_{280} :	> 5.00	5.27
A_{620} / A_{566} :	< 0.010	0.002
A_{566} / A_{498} :	< 1.50	1.35
Ex at 498 nm	peak 571-575	573.0
HPLC	>95%	98.38%



Fluorescence excitation and emission spectra of R-phycoerythrin in 100 mM sodium phosphate (pH 7.2) + 1 mM EDTA and 1 mM sodium azide. Emission scan was taken with excitation at 498 nm. Excitation scan was taken with emission at 575 nm.

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com