



SureLight® Allophycocyanin (APC)

Product Number D3-000
Lot Number APC000-21-070
Concentration 20.36 mg/ml
Exp. Date 09/21/2023
Store at 2-8° C

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. Avoid exposure to heat and light.

Buffer

50mM Sodium Phosphate (pH 7.0) +
50% Ammonium Sulfate

Stability

Product should be stored at 2-8°C in the dark.
Product is stable for up to 4 years depending on treatment by end-user.

Note: For research use only; not for diagnostic or therapeutic use.

General Information on APC:

- Approximate molecular weight 105,000 kDa
- Subunit structure $(\alpha\beta)_3$
- Absorbance/excitation maximum 652 nm $\pm$ 2nm
- Emission maximum 660 nm $\pm$ 2nm

Minimum ratio values:

$A_{650}/A_{280} \geq 4.6$ 5.54
 $A_{650}/A_{620} \geq 1.50$ 1.56

Lot Specific values:

Absorbance/Excitation maximum 653.5 nm
Emission maximum 659.5 nm

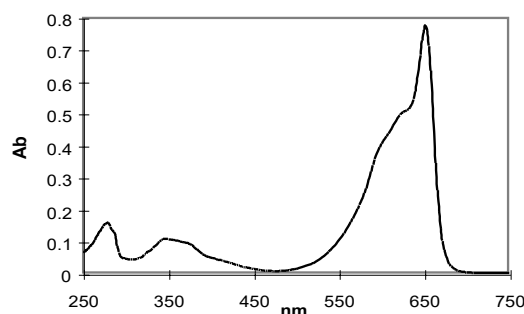


Figure 1. Absorbance spectrum of 0.1 mg/mL in 100 mM sodium phosphate (pH 7.0) and 0.05% sodium azide.

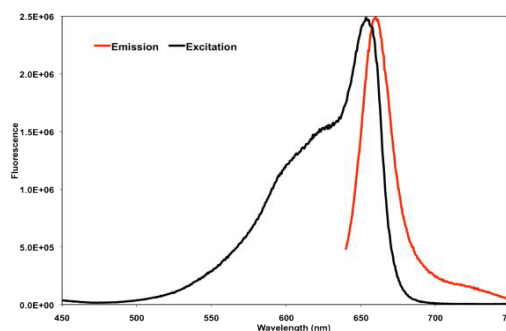


Figure 2. Normalized fluorescence emission and excitation spectra of sample at 0.167 µg/mL in 100 mM sodium phosphate (pH 7.2) and 0.05% sodium azide.

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com