

Sulfo-SMCC Activated SureLight™ APC

Product Number D3-010 Lot # XXXX

Amount 1.00 mg total protein

Conc 10 mg/mL after resuspension

Store at -20°C

Form/Shipping & Storage

Supplied as freeze dried material. After reconstitution with 100 μL of ddH2O the concentration is 10mg/mL.

Handling

We recommend that the investigator determine the appropriate working concentration and molar excess needed for conjugation.

Buffer

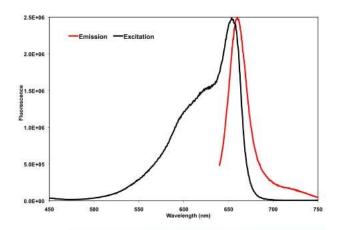
The product is in 100 mM Sodium Phosphate (pH 7.0) + 5 mM EDTA + 0.05% Sodium Azide +50mM Sucrose.

Stability

Upon receipt as freeze dried material, store at -20°C. After reconstitution store material at 4°C. Phycobiliproteins are sensitive to freeze-thaw cycles. Once reconstituted the material should be used within 1 month.

Note

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



Fluorescence excitation and emission spectra of APC in 100 mM sodium phosphate (pH 7.2) + 1 mM EDTA and 1 mM sodium azide. The emission scan was taken with excitation at 630 nm. The excitation scan was taken with an emission at 660 nm. The curves were normalized to equalize peak being the

Spectral Characteristics

Visible absorption maximum 652 Emission maximum 657

For technical inquiries: info@columbiabiosciences.com For sales inquiries: sales@columbiabiosciences.com