



Anti-6xHistidine conjugated to SureLight™ APC

Product Number D3-1711
Clone AD1.1.10
Amount 100 µg total protein
Store at 2-8°C

Form/Shipping & Storage

Supplied lyophilized. Upon receipt, store at -20°C. Reconstitute with 1 ml of ddH₂O and store at 2-8°C. Phycobiliproteins are sensitive to freeze-thaw cycles.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light.

Buffer

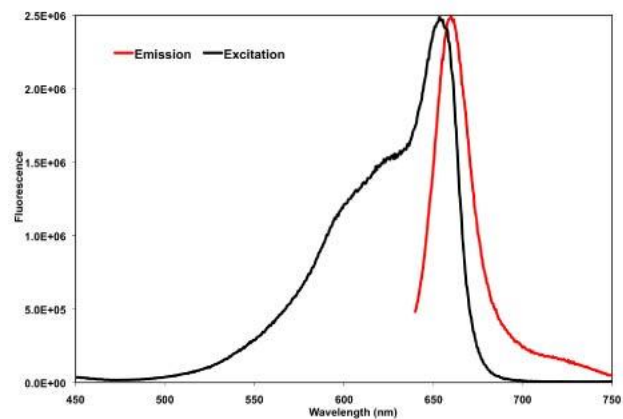
Upon reconstitution, the product is in 10 mM tris (pH 8.2) + 150 mM NaCl + 50 mM Sucrose + 1 µg/mL pentachlorophenol + 0.1% BSA.

Stability

Product should be stored at 2-8°C in the dark and be used within 6 months. If further dilution of the conjugate is required, use diluted material within one week.

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 heights.

Spectral Characteristics

Visible absorption maxima	652
Emission maximum	657

Concentration

Upon reconstitution with 1 ml ddH₂O:

[α-6xHistidine] = 0.518 µM
[SureLight™ APC] = 0.293 µM

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

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



Anti-6xHistidine conjugated to SureLight™ APC

Product Number D3-1711
Clone AD1.1.10
Amount 1.0 mg total protein
Store at 2-8°C

Form/Shipping & Storage

Supplied lyophilized. Upon receipt, store at -20°C. Reconstitute with 1 ml of ddH₂O and store at 2-8°C. Phycobiliproteins are sensitive to freeze-thaw cycles.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light.

Buffer

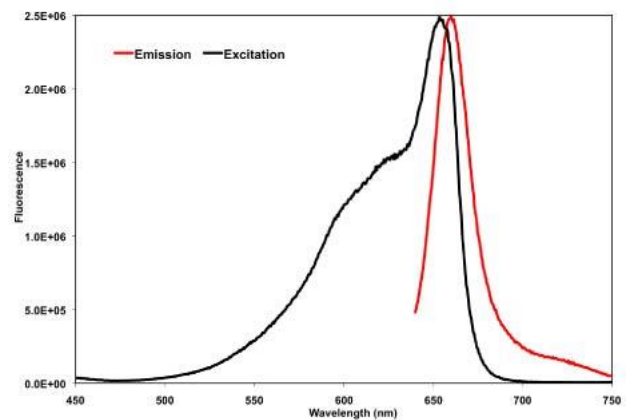
Upon reconstitution, the product is in 10 mM tris (pH 8.2) + 150 mM NaCl + 50 mM Sucrose + 1 µg/mL pentachlorophenol + 0.1% BSA.

Stability

Product should be stored at 2-8°C in the dark and be used within 6 months. If further dilution of the conjugate is required, use diluted material within one week.

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 heights.

Spectral Characteristics

Visible absorption maxima	652
Emission maximum	657

Concentration

Upon reconstitution with 1 ml ddH₂O:
[α-6xHistidine] = 5.78 µM
[SureLight™ APC] = 2.96 µM

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com