



Sulfo-SMCC-Activated SureLight® R-Phycoerythrin (R-PE)

Product Number D5-010
Activation Sulfo-SMCC
Amount: 1.0 mg
Concentration 10 mg/mL upon reconstitution

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt please, store at -20° C. After resuspension store a 4°C.

Handling

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

Buffer

100mM Sodium Phosphate (pH 7.0) + 5mM EDTA + 50 mM sucrose + 0.05% sodium azide after resuspension with 0.5 mL of deionized water

Stability

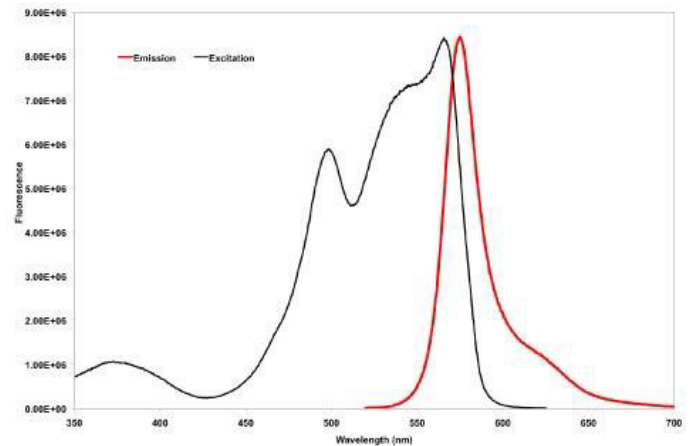
Product should be stored at 2-8°C in the dark. Product is stable for up to 1 year. Please use within 1 month after resuspension

Note

Succinimidyl 4-[N-maleimidomethyl]cyclohexane-1-carboxylate (SMCC) activated R-PE for research use only, not for diagnostic or therapeutic use.

General Information on RPE:

- Approximate molecular weight 250,000 kDa
- Subunit structure (αβ)6γ
- Absorbance/excitation maximum 565 > 540 >498 nm
- Emission maximum 572 nm



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. Scans were normalized to equalize peak heights.

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com



Sulfo-SMCC-Activated R-Phycoerythrin (R-PE)

Product Number D5-010
Activation Sulfo-SMCC
Amount: 5.0 mg
Concentration 10 mg/mL upon reconstitution

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt please, store at -20° C. After resuspension store at 4° C.

Handling

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

Buffer

100mM Sodium Phosphate (pH 7.0) + 5mM EDTA + 50 mM sucrose + 0.05% sodium azide after resuspension with 0.5 mL of deionized water

Stability

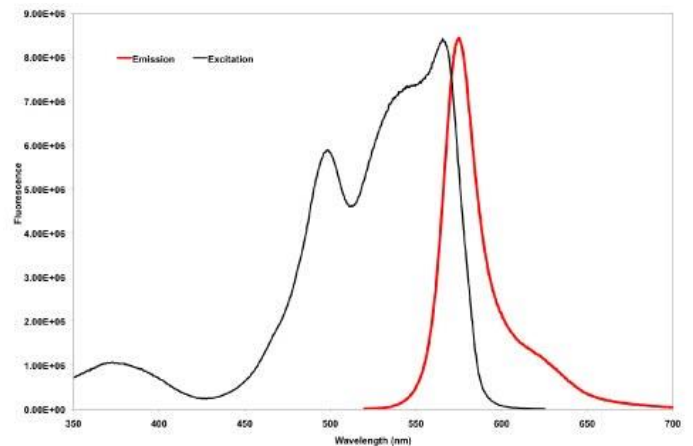
Product should be stored at 2-8°C in the dark. Product is stable for up to 1 year. Please use within 1 month after resuspension

Note

Succinimidyl 4-[N-maleimidomethyl]cyclohexane-1-carboxylate (SMCC) activated R-PE for research use only, not for diagnostic or therapeutic use.

General Information on RPE:

- Approximate molecular weight 250,000 kDa
- Subunit structure $(\alpha\beta)_6$
- Absorbance/excitation maximum 565 > 540 > 498 nm
- Emission maximum 572 nm



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. Scans were normalized to equalize peak heights.

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com



Sulfo-SMCC-Activated R-Phycoerythrin (R-PE)

Product Number D5-010
Activation Sulfo-SMCC
Amount: 10 mg
Concentration 10 mg/mL upon reconstitution

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt please, store at -20° C. After resuspension store a 4° C.

Handling

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

Buffer

100mM Sodium Phosphate (pH 7.0) + 5mM EDTA + 50 mM sucrose + 0.05% sodium azide after resuspension with 0.1 mL of deionized water

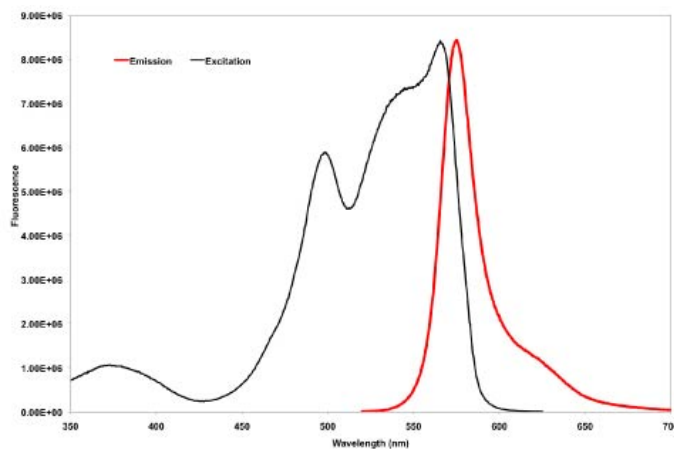
Stability

Product should be stored at 2-8°C in the dark. Product is stable for up to 1 year. Please use within 1 month after resuspension

Note

Succinimidyl 4-[N-maleimidomethyl]cyclohexane-1-carboxylate (SMCC) activated R-PE for research use only, not for diagnostic or therapeutic use.

Specifications:	Actual Result:
A ₅₆₆ / A ₂₈₀ :	≥ 5.00 5.41
A ₆₂₀ / A ₅₆₆ :	≤ 0.010 -0.00015
A ₅₆₆ / A ₄₉₆ :	≤ 1.50 1.45
Ex at 498 nm	peak 571-575 574.0
HPLC	≥95% 98.7%



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. Scans were normalized to equalize peak heights.

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com