Glowing products for science

SynaptoRed™ C1

SynaptoRedTM C1 is similar to the widely used red fluorescent nerve terminal dye SynaptoRedTM C2 (also called FM®4-64) except that the lipophilic tail of SynaptoRedTM C1 is one carbon shorter than that of the SynaptoRedTM C2.



Product attributes

Probe cellular localization	Membrane/vesicular
For live or fixed cells	For live/intact cells
Assay type/options	Real-time imaging
Colors	Far-red
Excitation/Emission	510/750 nm (see product description)

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

SynaptoRedTM C1 is similar to the widely used red fluorescent nerve terminal dye <u>SynaptoRedTM C2</u> (also called FM®4-64) except that the lipophilic tail of SynaptoRedTM C1 is one carbon shorter than that of the SynaptoRedTM C2. Thus, SynaptoRedTM C1 is expected to have a faster off-rate compared to SynaptoRedTM C2.

- $\lambda_{Ex}/\lambda_{Em}$ (membranes) = ~510/750 nm*
- Dark purple solid soluble in water
- C₂₈H₄₁Br₂N₃
- MW: 580

^{*} Excitation/emission settings of 515/640 nm have been used for detection of yeast vacuole staining with SynaptoRed™ C2 (FM®4-64), and should be applicable to SynaptoRed™ C1.

Nerve Terminal Dyes and Kits

Catalog No.

Product

SynaptoGreen™ Dyes (Ex/Em ~480/660 nm in membranes)	70020 70053	View full list of SynaptoGreen™ Dyes			
SynaptoRed Dyes™ (Ex/Em ~510/750 nm in membranes)	70021 70050	View full list of SynaptoRed™ Dyes			
Background Reducers					
ADVASEP-7	70029	Sulfonated cyclodextrin that aids in removal of free dye during washes			
SCAS	70037	Quenches extracellular fluorescences with fewer wash steps than ADVASEP-7			
Sulforhodamine 101	80101	Red fluorescent dye that quenches extracellular fluorescence of SynaptoGreen™ dyes			
Nerve Terminal Staining Kits					
Nerve Terminal Staining Kit I	70030	Includes SynaptoGreen™ C4 and ADVASEP-7			
Nerve Terminal Staining Kit II (A)	<u>70031</u>	Includes AM1-43 and ADVASEP-7			
Nerve Terminal Staining Kit II (B)	<u>70031-1</u>	Includes AM1-43 and SCAS			
Nerve Terminal Staining Kit III	70032	Includes SynaptoGreen™ C4 and Sulforhodamine 101			
Nerve Terminal Staining Kit V	70034	Includes SynaptoRed™ C2 and ADVASEP-7			

Features

This datasheet was generated on November 15, 2025 at 03:38:29 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/synaptoredtm-c1/

m is the number of carbons in the lipophilic tail and n is the number of double bonds linking the two aromatic rings in the dye.

**The positively-charged end of SynaptoRed C2M is a trimethylammonium group.FM is a registered trademark of Thermo Fisher Scientific.