

SynaptoRed™ C1

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



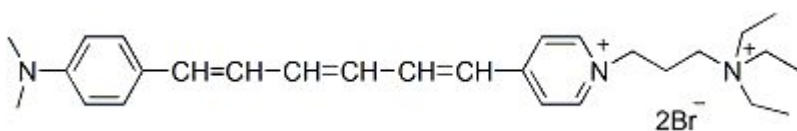
Product attributes

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)

Product Description

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

Product	Catalog No.	Features
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)	70031	Includes AM1-43 and ADVASEP-7
Nerve Terminal Staining Kit II (B)	70031-1	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

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.

This datasheet was generated on June 15, 2026 at 01:29:34 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/synaptoRedtm-c1/>

