



Glowing products for science

## SynaptoRed™ C2M (Equivalent to FM®5-95)

SynaptoRed™ C2M is a red fluorescent probe used to follow synaptic vesicles.



Call us : [800-304-5357](tel:800-304-5357)

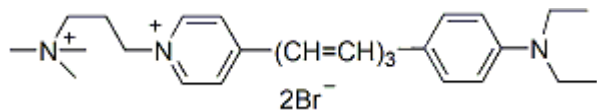
### Product attributes

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™ C2M (also known as FM®5-95), is a red fluorescent probe used to follow synaptic vesicles. SynaptoRed™ C2M is slightly more water soluble than SynaptoRed™ C2 (equivalent to FM®4-64).

- $\lambda_{Ex}/\lambda_{Em}$  (membranes) = ~510/750 nm\*
- Dark purple solid soluble in water
- MW: 565.43



\* 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™ C2M.

# Nerve Terminal Dyes and Kits

Product	Catalog No.	Features
SynaptoGreen™ Dyes (Ex/Em ~480/660 nm in membranes)	70020... 70053	<a href="#">View full list of SynaptoGreen™ Dyes</a>
SynaptoRed Dyes™ (Ex/Em ~510/750 nm in membranes)	70021... 70050	<a href="#">View full list of SynaptoRed™ Dyes</a>
Background Reducers		
<a href="#">ADVASEP-7</a>	<a href="#">70029</a>	Sulfonated cyclodextrin that aids in removal of free dye during washes
<a href="#">SCAS</a>	<a href="#">70037</a>	Quenches extracellular fluorescences with fewer wash steps than ADVASEP-7
<a href="#">Sulforhodamine 101</a>	<a href="#">80101</a>	Red fluorescent dye that quenches extracellular fluorescence of SynaptoGreen™ dyes
Nerve Terminal Staining Kits		
<a href="#">Nerve Terminal Staining Kit I</a>	<a href="#">70030</a>	Includes SynaptoGreen™ C4 and ADVASEP-7
<a href="#">Nerve Terminal Staining Kit II (A)</a>	<a href="#">70031</a>	Includes AM1-43 and ADVASEP-7
<a href="#">Nerve Terminal Staining Kit II (B)</a>	<a href="#">70031-1</a>	Includes AM1-43 and SCAS
<a href="#">Nerve Terminal Staining Kit III</a>	<a href="#">70032</a>	Includes SynaptoGreen™ C4 and Sulforhodamine 101
<a href="#">Nerve Terminal Staining Kit V</a>	<a href="#">70034</a>	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.

## References

1. Science 297, 1525 (2002).
2. Cell 107, 605 (2001).

This datasheet was generated on January 14, 2026 at 03:50:27 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/synaptoredtmc2m-also-known-as-fm5-95-a-trademark-of-molecular-probes-inc/>