

## AM4-65

AM4-65 is a fixable red fluorescent nerve terminal probe. The dye is spectrally identical to SynaptoRed™ C2 (equivalent to FM®4-64), but it has an additional amine group at the hydrophilic end of the molecule, making the dye fixable with formaldehyde or glutaraldehyde.



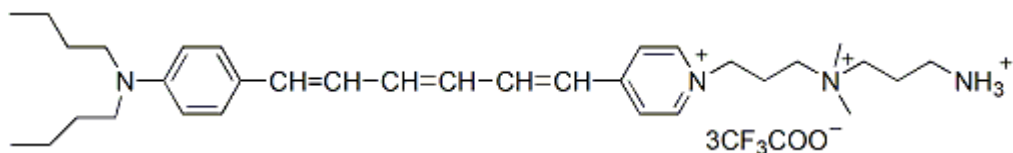
### Product attributes

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

## Product Description

AM4-65 is a fixable red fluorescent nerve terminal probe. The dye is spectrally identical to SynaptoRed™ C2 (equivalent to FM®4-64), but it has an additional amine group at the hydrophilic end of the molecule, making the dye fixable with formaldehyde or glutaraldehyde. The long emission wavelength makes the dye suitable to be used with GFP in two-color imaging.

- $\lambda_{Ex}/\lambda_{Em}$  (membranes) = ~510/750 nm\*
- Dark purple solid soluble in water
- MW: 844.85



\* 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 AM4-65.

# 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.

This datasheet was generated on January 9, 2026 at 03:31:02 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/am4-65/>