AM4-66

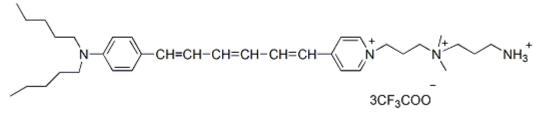
A fixable red fluorescent nerve terminal probe that is spectrally identical to SynaptoRed™ C2 (also called FM®4-64).



Product Description

AM4-66 is a fixable red fluorescent nerve terminal probe. The dye is spectrally identical to SynaptoRedTM C2 (also called FM®4-64), but it has an additional amine group at the hydrophilic end of the molecule, making the dye fixable with formaldehyde or gluteraldehyde. The long emission wavelength makes the dye suitable to be used with GFP in two-color imaging.

- Red fluorescent nerve terminal probe that is fixable with formaldehyde or gluteraldehyde
- $\lambda_{Ex}/\lambda_{Em}$ (membranes) = ~510/750 nm*
- Dark purple solid soluble in water
- MW: 872.85



^{*} 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 AM4-66.

Product attributes

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

	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
	Colors	Far-red
	Excitation/Emission	510/750 nm (see product description)

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	<u>70029</u>	Sulfonated cyclodextrin that aids in removal of free dye during washes		
SCAS	<u>70037</u>	Quenches extracellular fluorescences with fewer wash steps than ADVASEP-7		
Sulforhodamine 101	80101	Red fluorescent dye that quenches extracellular fluorescence of SynaptoGreen TM 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)	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.

Features

This datasheet was generated on November 1, 2025 at 11:20:28 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/am4-66/