

AM4-64

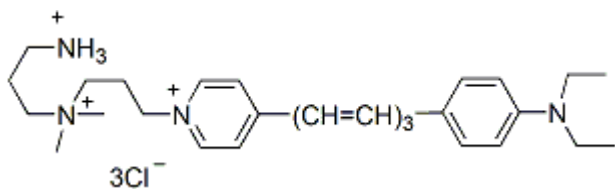
AM4-64 is a fixable, activity-dependent fluorescent nerve terminal probe. AM4-64 functions similarly as SynaptoRed™ C2 (#70021) (equivalent to FM®4-64), but with the additional property of being fixable.



Product Description

AM4-64 is a fixable, activity-dependent fluorescent nerve terminal probe. AM4-64 functions similarly as SynaptoRed™ C2 ([70021](#)) (equivalent to FM®4-64), but with the additional property of being fixable. Thus, AM4-64 is a useful tool for synapse studies where subsequent fluorescent immunocytochemistry is desired. AM4-64 and FM®4-64 have identical excitation and emission wavelengths.

- $\lambda_{Ex}/\lambda_{Em}$ (membranes) = ~510/750 nm*
- Dark purple solid soluble in water
- MW: 555.5



* 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 AM4-64.

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)

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 January 14, 2026 at 04:05:40 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/am4-64/>