

## AM4-64

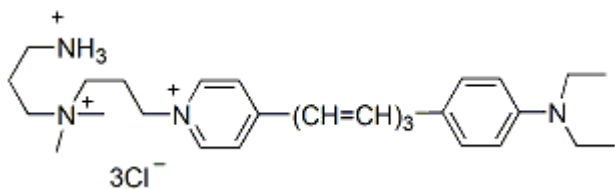
AM4-64 is a fixable, activity-dependent fluorescent nerve terminal probe. AM4-64 functions similarly as SynaptoRed™ C2 (#70021) (equivalent to FM®4-64), but with the additional property of being fixable.



## 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 immunochemistry 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

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

## Nerve Terminal Dyes and Kits

Product	Catalog No.	Features
<b>SynaptoGreen™ Dyes</b> (Ex/Em ~480/660 nm in membranes)	70020... 70053	<a href="#">View full list of SynaptoGreen™ Dyes</a>
<b>SynaptoRed Dyes™</b> (Ex/Em ~510/750 nm in membranes)	70021... 70050	<a href="#">View full list of SynaptoRed™ Dyes</a>
<b>Background Reducers</b>		
<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
<b>Nerve Terminal Staining Kits</b>		
<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 June 17, 2026 at 01:07:35 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/am4-64/>