

## AM1-43 (Equivalent to FM®1-43FX)

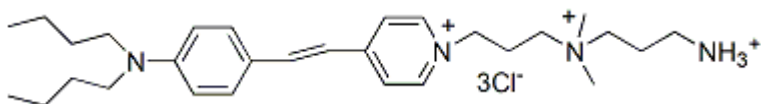
AM1-43 (equivalent to FM®1-43FX) is a fixable, activity-dependent fluorescent nerve terminal probe. AM1-43 functions similarly to SynaptoGreen C4 (catalog # 70020), but with the additional property of being fixable.



### Product Description

AM1-43 (equivalent to FM®1-43FX) is a fixable, activity-dependent fluorescent nerve terminal probe. AM1-43 functions similarly to SynaptoGreen™ C4 ([70020](#)), but with the additional property of being fixable. Thus, AM1-43 is a useful tool for synapse studies where subsequent fluorescent immunocytochemistry is desired. AM1-43 and FM®1-43FX have identical excitation and emission wavelengths. Also see AM1-44 (catalog no. [70038](#)), which is similar to AM1-43, but has longer lipophilic tails, resulting in better fixability.

- Ex/Em: 510/625 nm (in MeOH); 480/598 nm (in membranes)
- Red solid soluble in water
- C<sub>29</sub>H<sub>49</sub>Cl<sub>3</sub>N<sub>4</sub>
- MW: 559.5



### 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	Green
Excitation/Emission	480/598 nm (in membranes)

# 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.

## References

1. PLoS One, 5(8), e12386.
2. Anat Cell Biol 44(4), 274 (2011).

This datasheet was generated on June 15, 2026 at 03:09:58 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/am1-43/>