

## AM1-44

AM1-44 is a version of AM1-43 (catalog no. 70024) with improved fixability. The dye can also be used as a general probe for following endocytosis.

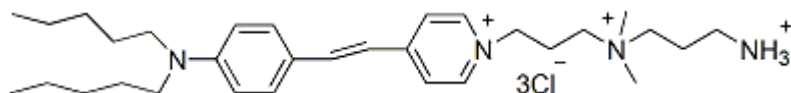


## Product Description

AM1-44 is a version of AM1-43 (catalog no. [70024](#)) with improved fixability. The dye can also be used as a general probe for following endocytosis. AM1-44 probe functions similarly as AM1-43, except that AM1-44 has lipophilic tails that are each one carbon longer than those of AM1-43, resulting in better fixability.

$\lambda_{Ex}$ : 480/598 nm (in MeOH); 480/598 nm (in membranes)  $\lambda_{Em}$ : 510/625 nm (in MeOH); 480/598 nm (in membranes)

- $C_{31}H_{52}Cl_3N_4$
- MW: 587.5



## 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 January 16, 2026 at 05:52:01 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/am1-44/>