

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 attributes

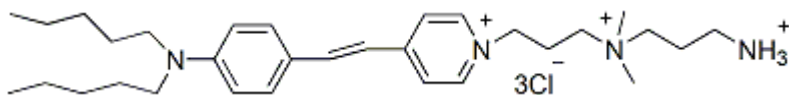
Probe cellular localization	Membrane/vesicular
For live or fixed cells	For live/intact cells
Assay type/options	Real-time imaging
Fixation options	Fix after staining (formaldehyde), Permeabilize after staining
Colors	Green
Excitation/Emission	480/598 nm (in membranes)

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.

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

- Soluble in MeOH, DMSO and water
- $C_{31}H_{52}Cl_3N_4$
- MW: 587.5



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 June 9, 2026 at 05:02:56 PM. Visit product page to check for updated information before use.

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