AM3-25

AM3-25 is a lipophilic styryl dye that has an amine group at the hydrophilic head. The amine group reacts with formaldehyde or glutaldehyde to make the dye fixable.



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	480/600 nm (in membranes)

Call us: 800-304-5357 Email: btinfo@biotium.com

glutaldehyde to make the dye fixable.

AM3-25 is a lipophilic styryl dye that has an amine group at the hydrophilic head. The amine group reacts with formaldehyde or glutaldehyde to make the dye fixable.

- Ex/Em (MeOH) = 510/625 nm; Ex/Em (liposomes) = 480/600 nm
- Dark red soluble in MeOH or DMSO
- Store at 4°C and protect from light
- C57H105N4Cl3N4

Product Description

• MW: 952.83

Product

Features

Nerve Terminal Dyes and Kits

Catalog No

Product	Catalog No.	Features	
SynaptoGreen™ Dyes (Ex/Em ~480/660 nm in membranes)	70020 70053	<u>View full list of</u> <u>SynaptoGreen™ Dyes</u>	
SynaptoRed Dyes™ (Ex/Em ~510/750 nm in membranes)	70021 70050	View full list of SynaptoRed™ Dyes	
Background Reducers			
ADVASEP-7	<u>70029</u>	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)	<u>70031</u>	Includes AM1-43 and ADVASEP-7	
Nerve Terminal Staining Kit II (B)	<u>70031-1</u>	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 November 15, 2025 at 03:05:22 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/am3-25/