

## 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.



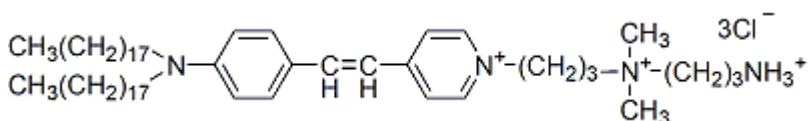
### Product attributes

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)

### Product Description

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
- C<sub>57</sub>H<sub>105</sub>N<sub>4</sub>Cl<sub>3</sub>N<sub>4</sub>
- MW: 952.83



### 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.