



Glowing products for science

## SynaptoGreen™ C18 (Equivalent to FM®3-25)

SynaptoGreen™ C18 (equivalent to FM®3-25) is a membrane stain with a chemical structure similar to other nerve terminal dyes, except that it has the longest lipophilic tail (C18). Because of its large size, it is used as a control for ion channel permeability of nerve terminal dyes.



Call us : [800-304-5357](tel:800-304-5357)

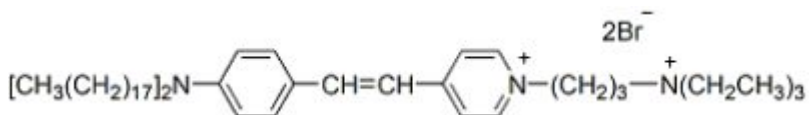
### Product attributes

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

## Product Description

SynaptoGreen™ C18 (equivalent to FM®3-25) is a membrane stain with a chemical structure similar to other nerve terminal dyes, except that it has the longest lipophilic tail (C18). Because of its large size, it is used as a control for ion channel permeability of nerve terminal dyes.

- Red solid soluble in DMSO or DMF
- Ex/Em: 510/625 nm (in MeOH); 480/598 nm (in membranes)
- Store desiccated at 4 °C and protect from light, especially in solution
- C<sub>58</sub>H<sub>105</sub>Br<sub>2</sub>N<sub>3</sub>
- MW: 1004



# Nerve Terminal Dyes and Kits

Product	Catalog No.	Features
SynaptoGreen™ Dyes (Ex/Em ~480/660 nm in membranes)	70020... 70053	<a href="#">View full list of SynaptoGreen™ Dyes</a>
SynaptoRed Dyes™ (Ex/Em ~510/750 nm in membranes)	70021... 70050	<a href="#">View full list of SynaptoRed™ Dyes</a>
Background Reducers		
<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
Nerve Terminal Staining Kits		
<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. Mol Pain (1997) doi: 10.1186/1744-8069-3-1

This datasheet was generated on January 14, 2026 at 10:33:36 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/synaptogreen-c18-also-known-as-fm3-25/>