

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

SynaptoGreen<sup>TM</sup> 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.

# Biotlum

## Product attributes

Call us: 800-304-5357

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<sup>TM</sup> 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
- C58H105Br2N3
- MW: 1004

$$[CH_3(CH_2)_{17}]_2N$$
 —  $CH=CH$   $N$  —  $(CH_2)_3$  —  $N$   $(CH_2CH_3)_3$ 

### **Nerve Terminal Dyes and Kits**

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) Background Reducers	70021 70050	View full list of SynaptoRed™ Dyes
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

#### References

1. Mol Pain (1997) doi: 10.1186/1744-8069-3-1

This datasheet was generated on November 23, 2025 at 10:49:53 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/synaptogreen-c18-also-known-as-fm3-25/">https://biotium.com/product/synaptogreen-c18-also-known-as-fm3-25/</a>

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.