SynaptoGreen™ C5 (Equivalent to FM®1-84)

SynaptoGreen™ C5 is similar to SynaptoGreen™ C4 (70020), except that the lipophilic tail is one carbon longer.

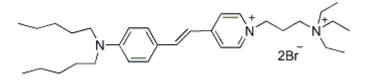
Product Description

SynaptoGreen™ C5 (equivalent to FM®1-84) is similar to SynaptoGreen™ C4 (70020), except that the lipophilic tail is one carbon longer. Thus SynaptoGreen™ C5 is less water soluble than SynaptoGreen™ C4 and is predicted to have a faster staining rate but slower de-staining rate.

Features

- Ex/Em: 510/625 nm (in MeOH)
- Red solid soluble in water
- C₃₂H₅₃Br₂N₃
- MW: 640

Product



Catalog No

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	<u>View full list of</u> <u>SynaptoRed™ Dyes</u>
Background Reducers		
ADVASEP-7	70029	Sulfonated cyclodextrin that aids in removal of free dye during washes
<u>SCAS</u>	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 TM 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 TM C2 and ADVASEP-7

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

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)

This datasheet was generated on November 8, 2025 at 08:07:14 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/synaptogreentm-c5-also-known-as-fm1-84-a-trademark-of-molecular-probes-inc/

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.