Glowing products for science

SynaptoGreen™ C3

SynaptoGreenTM C3 is similar to SynaptoGreenTM C4 (70020), except that the lipophilic tail is one carbon shorter and also the hydrophilic end is a trimethylammonium group.



Product attributes

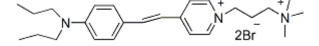
Probe cellular localization	Membrane/vesicular
For live or fixed cells	For live/intact cells
Assay type/options	Real-time imaging
Colors	Green
Excitation/Emission	480/598 nm (in membranes)

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

Product Description

SynaptoGreenTM C3 is similar to SynaptoGreenTM C4 (70020), except that the lipophilic tail is one carbon shorter and also the hydrophilic end is a trimethylammonium group. Thus SynaptoGreenTM C3 is slightly more water soluble than SynaptoGreenTM C4. Please also see the fixable nerve terminal dye AM1-43 (70024).

- λ_{Ex}/λ_{Em}: 510/625 nm (in MeOH); 480/598 nm (in membranes)
- Red solid soluble in water
- C₂₅H₃₉Br₂N₃
- MW: 542



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	View full list of SynaptoRed™ Dyes	
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 TM 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	<u>70034</u>	Includes SynaptoRed™ C2 and ADVASEP-7	

References

- 1. J Neurosci 12, 363 (1992).
- 2. Science 255, 200 (1992).
- 3. Neuron 24, 803 (1999).
- 4. Neuron 24, 809 (1999).

This datasheet was generated on November 2, 2025 at 05:08:15 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/synaptogreentm-c3/

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.