

## SynaptoGreen™ C1

SynaptoGreen<sup>TM</sup> C1 has the shortest (one-carbon) lipophilic tail and is the most water soluble of the nerve terminal dye series. Thus it is expected to have the slowest on-rate and the fastest off-rate among the SynaptoGreen<sup>TM</sup> dyes.



## 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> C1 has the shortest (one-carbon) lipophilic tail and is the most water soluble of the nerve terminal dye series. Thus it is expected to have the slowest on-rate and the fastest off-rate among the SynaptoGreen<sup>TM</sup> dyes.

- Red solid soluble in water
- C24H37Br2N3
- MW: 528

## **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) 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	<u>70030</u>	Includes SynaptoGreen™ 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)	70031-1	Includes AM1-43 and SCAS
Nerve Terminal Staining Kit III	70032	Includes SynaptoGreen <sup>TM</sup> C4 and Sulforhodamine 101
Nerve Terminal Staining Kit V	<u>70034</u>	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.

This datasheet was generated on November 23, 2025 at 10:50:59 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/synaptogreentm-c1/">https://biotium.com/product/synaptogreentm-c1/</a>