



Glowing products for science

SynaptoGreen™ C2 (Equivalent to FM®2-10)

SynaptoGreen™ C2 is more water soluble than SynaptoGreen™ C4 and thus has a faster de-staining rate than the latter.



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

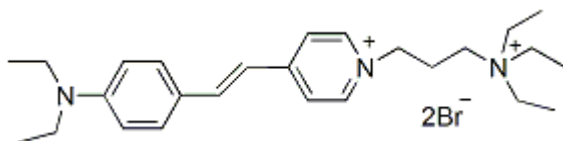
Product attributes

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™ C2 is more water soluble than SynaptoGreen™ C4 and thus has a faster de-staining rate than the latter.

- Red solid soluble in water
- C₂₆H₄₁Br₂N₃
- MW: 556



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
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

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. Proc
R Soc Lond B
255, 61 (1994).

This datasheet was generated on December 21, 2025 at 03:38:05 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/synaptogreentm-c2-also-known-as-fm2-10-a-trademark-of-molecular-probes-inc/>