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

SynaptoGreen  $^{\text{TM}}$  C2 is more water soluble than SynaptoGreen  $^{\text{TM}}$  C4 and thus has a faster de-staining rate than the latter.



#### **Product attributes**

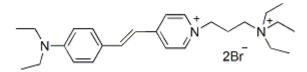
Probe cellular localization	Membrane/vesicular
For live or fixed cells	For live/intact cells
Assay type/options	Real-time imaging
Colors	Green
Excitation/Emission	480/598 nm (in membranes)

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

## **Product Description**

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

- Red solid soluble in water
- C<sub>26</sub>H<sub>41</sub>Br<sub>2</sub>N<sub>3</sub>
- 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	
<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 <sup>TM</sup> 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. Proc

R Soc Lond B

255, 61 (1994).

This datasheet was generated on November 2, 2025 at 05:08:42 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/synaptogreentm-c2-also-known-as-fm2-10-a-trademark-of-molecular-probes-inc/">https://biotium.com/product/synaptogreentm-c2-also-known-as-fm2-10-a-trademark-of-molecular-probes-inc/</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.