



Glowing products for science

## SynaptoGreen™ C4 (Equivalent to FM®1-43)

A widely used green fluorescent dye for following synaptic activity by staining synaptic vesicles at the synapse or neuromuscular junctions. When used in combination with the red fluorescent dye SynaptoRed™ (), synapses or neuromuscular junctions can be imaged independently in two colors.

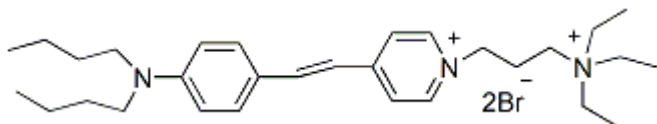


## Product Description

A widely used green fluorescent dye for following synaptic activity by staining synaptic vesicles at the synapse or neuromuscular junctions. When used in combination with the red fluorescent dye SynaptoRed™ ([70021](#))&comma; synapses or neuromuscular junctions can be imaged independently in two colors. Research results from the Tsien and Kay labs showed that brain slices can be stained with SynaptoGreen™ when SR101 (Sulforhodamine 101&comma; [80101](#)) or ADVASEP-7 ([70029](#)) is used to reduce the background staining. Biotium offers several [Nerve Terminal Staining Kits](#) for staining brain slices. Also see AM1-43 ([70024](#))&comma; which is similar to SynaptoGreen™&comma; but with a formaldehyde-fixable amine group.

SynaptoGreen™ also can be used to stain bacteria (gram-positive or gram-negative) and yeast&comma; although in yeast it rapidly internalizes to vacuolar membranes. See our [Cellular Stains Table](#) for more information on how our dyes stain various organisms.

- $\lambda_{\text{Ex}}/\lambda_{\text{Em}}$ : 510/625 nm (in MeOH); 480/598 nm (in membranes)
- Red solid soluble in water
- $\text{C}_{30}\text{H}_{49}\text{Br}_2\text{N}_3$
- MW: 612



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

### Product attributes

|                             |                           |
|-----------------------------|---------------------------|
| CAS number                  | 149838-22-2               |
| 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) |

# Nerve Terminal Dyes and Kits

| Product  | Catalog No.             | Features   |
|--|-------------------------|--|
| SynaptoGreen™ Dyes<br>(Ex/Em ~480/660 nm in membranes) | 70020... 70053          | <a href="#">View full list of SynaptoGreen™ Dyes</a>                               |
| SynaptoRed Dyes™<br>(Ex/Em ~510/750 nm in membranes)   | 70021... 70050          | <a href="#">View full list of SynaptoRed™ Dyes</a>                                 |
| Background Reducers                                    |                         |  |
| <a href="#">ADVASEP-7</a>                              | <a href="#">70029</a>   | Sulfonated cyclodextrin that aids in removal of free dye during washes             |
| <a href="#">SCAS</a>                                   | <a href="#">70037</a>   | Quenches extracellular fluorescences with fewer wash steps than ADVASEP-7          |
| <a href="#">Sulforhodamine 101</a>                     | <a href="#">80101</a>   | Red fluorescent dye that quenches extracellular fluorescence of SynaptoGreen™ dyes |
| Nerve Terminal Staining Kits                           |                         |  |
| <a href="#">Nerve Terminal Staining Kit I</a>          | <a href="#">70030</a>   | Includes SynaptoGreen™ C4 and ADVASEP-7  |
| <a href="#">Nerve Terminal Staining Kit II (A)</a>     | <a href="#">70031</a>   | Includes AM1-43 and ADVASEP-7  |
| <a href="#">Nerve Terminal Staining Kit II (B)</a>     | <a href="#">70031-1</a> | Includes AM1-43 and SCAS   |
| <a href="#">Nerve Terminal Staining Kit III</a>        | <a href="#">70032</a>   | Includes SynaptoGreen™ C4 and Sulforhodamine 101                                   |
| <a href="#">Nerve Terminal Staining Kit V</a>          | <a href="#">70034</a>   | 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 January 16, 2026 at 03:27:08 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/synaptogreentm-c4-also-known-as-fm1-43-a-trademark-of-molecular-probes-inc/>