

## SynaptoGreen™ C5 (Equivalent to FM®1-84)

SynaptoGreen™ C5 is similar to SynaptoGreen™ C4 ([70020](#)), except that the lipophilic tail is one carbon longer.



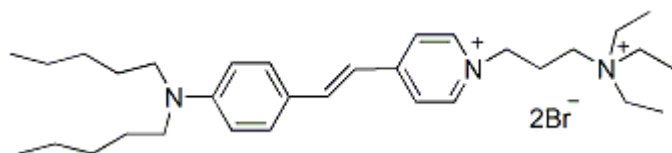
### Product attributes

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™ C5 (equivalent to FM®1-84) is similar to SynaptoGreen™ C4 ([70020](#)), except that the lipophilic tail is one carbon longer. Thus SynaptoGreen™ C5 is less water soluble than SynaptoGreen™ C4 and is predicted to have a faster staining rate but slower de-staining rate.

- Ex/Em: 510/625 nm (in MeOH)
- Red solid soluble in water
- C<sub>32</sub>H<sub>53</sub>Br<sub>2</sub>N<sub>3</sub>
- MW: 640



## Nerve Terminal Dyes and Kits

Product	Catalog No.	Features
<b>SynaptoGreen™ Dyes</b> (Ex/Em ~480/660 nm in membranes)	70020... 70053	<a href="#">View full list of SynaptoGreen™ Dyes</a>
<b>SynaptoRed Dyes™</b> (Ex/Em ~510/750 nm in membranes)	70021... 70050	<a href="#">View full list of SynaptoRed™ Dyes</a>
<b>Background Reducers</b>		
<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
<b>Nerve Terminal Staining Kits</b>		
<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.