Nerve Terminal Staining Kits

Nerve Terminal Staining Kits contain Biotium's nerve terminal dyes and accessories for washing away or quenching extracellular dye fluorescence.



Product attributes

Probe cellular localization	Membrane/vesicular
For live or fixed cells	For live/intact cells
Assay type/options	Real-time imaging
Colors	Green, Far-red

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

Product Description

Nerve Terminal Staining Kits contain Biotium's nerve terminal dyes and background reducing agents for washing away or quenching extracellular dye fluorescence. Kits are available with SynaptoGreen™ C4 (equivalent to FM®1-43), the fixable green nerve terminal dye AM1-43, or SynaptoRed™ C2 (equivalent to FM®4-64).

The kits are available with a choice of background reducing agents. ADVASEP-7 is a sulfonated cyclodextrin derivative that has been reported to reduce background fluorescence when using SynaptoGreen™ C4 to stain brain slices. SCAS (4-Sulfonato calix[8]arene, sodium salt) is a quencher developed by Biotium to reduce background fluorescence. Unlike ADVASEP-7, SCAS dramatically lowers background fluorescence as soon as it is added to the preparation without the need for repeated wash steps. Sulforhodamine 101 has been used as a quencher dye to reduce SynaptoGreen™ C4 fluorescence.

See the Neuronal Probes Flyer for detailed information.

Nerve Terminal Dyes and Kits

Product	Catalog No.	Features	
SynaptoGreen™ Dyes (Ex/Em ~480/660 nm in membranes)	70020 70053	<u>View full list of</u> <u>SynaptoGreen™ Dyes</u>	
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	<u>80101</u>	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)	<u>70031</u>	Includes AM1-43 and ADVASEP-7	
Nerve Terminal Staining Kit II (B)	<u>70031-1</u>	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.

This datasheet was generated on November 11, 2025 at 10:18:32 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/nerve-terminal-staining-kits/