

HM1-43

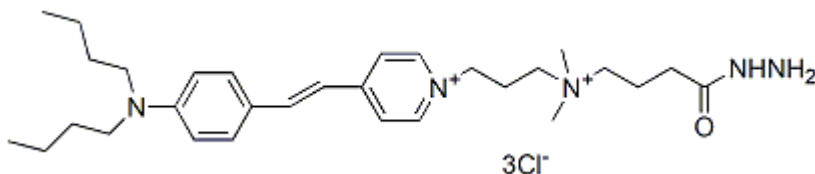
HM1-43 is a fixable, activity-dependent fluorescent nerve terminal probe.



Product Description

HM1-43 is a fixable, activity-dependent fluorescent nerve terminal probe.

- Functions similarly to SynaptoGreen™ (70020), but with the additional property of being fixable
- Useful for synapse studies where subsequent fluorescent immunochemistry is desired
- HM1-43 and FM®1-43 have identical excitation and emission wavelengths
- $\lambda_{\text{Ex}}/\lambda_{\text{Em}}$ (in membranes) = 480/598 nm



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.

This datasheet was generated on January 25, 2026 at 01:39:23 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/hm1-43/>