## **HM1-43**

Product

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



## **Product attributes**

Probe cellular localization	Membrane/vesicular
For live or fixed cells	For live/intact cells
Assay type/options	Real-time imaging
Fixation options	Fix after staining (formaldehyde), Permeabilize after staining
Excitation/Emission	480/598 nm (in membranes)

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

## **Product Description**

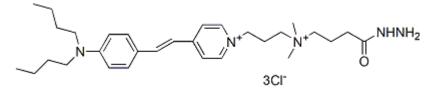
HM1-43 is a fixable, activity-dependent fluorescent nerve terminal probe.

Catalog No.

• Functions similarly to SynaptoGreen™ (70020), but with the additional property of being fixable

**Features** 

- Useful for synapse studies where subsequent fluorescent immunochemistry is desired
- HM1-43 and FM®1-43 have identical excitation and emission wavelengths
- λEx/λEm (in membranes) = 480/598 nm



## **Nerve Terminal Dyes and Kits**

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 <sup>TM</sup> 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 <sup>TM</sup> 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	

This datasheet was generated on November 12, 2025 at 12:32:35 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/hm1-43/

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.