AM2-10

AM2-10 is a fixable analog of SynaptoGreen™ C2.



Product attributes

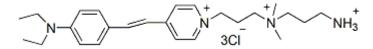
Probe cellular localization	Membrane/vesicular
For live or fixed cells	For live/intact cells
Assay type/options	Real-time imaging
Fixation options	Fix after staining (formaldehyde), Permeabilize after staining
Colors	Green
Evoltation/Emission	490/E00 pm (in mombrance)

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

Product Description

AM2-10 is a fixable analog of $\underline{\text{SynaptoGreen}^{\text{TM}} \text{C2}}$.

- Dark red solid soluble in water
- C₂₅H₃₇Cl₃N₄
- MW: 503.5



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) Background Reducers	70021 70050	View full list of SynaptoRed™ Dyes	
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 TM dyes	
Nerve Terminal Staining Kits			
Nerve Terminal Staining Kit I	70030	Includes SynaptoGreen TM 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 2, 2025 at 05:08:38 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/am2-10/