

## **SCAS**

SCAS (4-Sulfonato calix[8] arene, sodium salt) is a quencher developed by Biotium to reduce background fluorescence when using our fixable AM dyes, SynaptoGreen and SynaptoRed dyes. Unlike ADVASEP-7, SCAS dramatically lowers background fluorescence as soon as it is added to the preparation without the need for repeated wash steps.

Call us: 800-304-5357

#### **Product attributes**

Probe cellular localization Membrane/vesicular For live or fixed cells For live/intact cells

## **Product Description**

SCAS (4-Sulfonato calix[8]arene, sodium salt) is a quencher developed by Biotium to reduce background fluorescence when using our fixable AM dyes, SynaptoGreen and SynaptoRed dyes. Unlike ADVASEP-7, SCAS dramatically lowers background fluorescence as soon as it is added to the preparation without the need for repeated wash steps.

- Off-white solid soluble in water
- C56H40Na8O32S8
- MW 1665.34

# **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
SCAS	70037	Quenches extracellular fluorescences with fewer wash steps than ADVASEP-7
Sulforhodamine 101	80101	Red fluorescent dye that quenches extracellular fluorescence of SynaptoGreen <sup>TM</sup> dyes
Nerve Terminal Staining Kits		
Nerve Terminal Staining Kit I	<u>70030</u>	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.

#### References

- 1. J. Biol. Chem, 283(14), 9289(2008) <u>DOI: 10.1074/jbc.M706499200</u> 2. Dev. Cell, 44(1), 56(2018), <u>DOI: 10.1016/j.devcel.2017.12.014</u> 3. J. Cell Sci, 127(1), 250(2014), <u>DOI: 10.1242/jcs.140996</u> 4. J. Cell Sci, 134(2), (2021), <u>DOI: 10.1242/jcs.25773</u>

This datasheet was generated on November 24, 2025 at 04:08:38 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/scas4-sulfonato-calix8arene-sodium-salt/