

## 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.



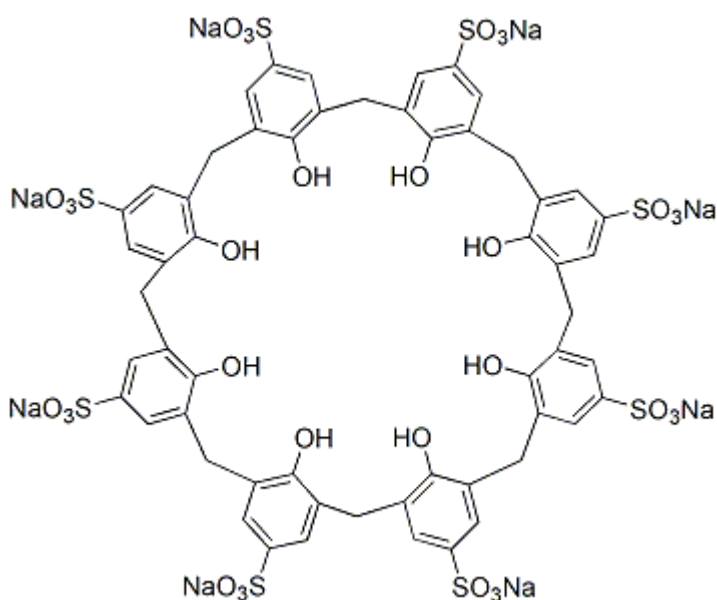
### Product attributes

| 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
- $C_{56}H_{40}Na_8O_{32}S_8$
- MW 1665.34



# Nerve Terminal Dyes and Kits

| Product  | Catalog No.             | Features   |
|--|-------------------------|--|
| SynaptoGreen™ Dyes<br>(Ex/Em ~480/660 nm in membranes) | 70020... 70053          | <a href="#">View full list of SynaptoGreen™ Dyes</a>                               |
| SynaptoRed Dyes™<br>(Ex/Em ~510/750 nm in membranes)   | 70021... 70050          | <a href="#">View full list of SynaptoRed™ Dyes</a>                                 |
| Background Reducers                                    |                         |  |
| <a href="#">ADVASEP-7</a>                              | <a href="#">70029</a>   | Sulfonated cyclodextrin that aids in removal of free dye during washes             |
| <a href="#">SCAS</a>                                   | <a href="#">70037</a>   | Quenches extracellular fluorescences with fewer wash steps than ADVASEP-7          |
| <a href="#">Sulforhodamine 101</a>                     | <a href="#">80101</a>   | Red fluorescent dye that quenches extracellular fluorescence of SynaptoGreen™ dyes |
| Nerve Terminal Staining Kits                           |                         |  |
| <a href="#">Nerve Terminal Staining Kit I</a>          | <a href="#">70030</a>   | Includes SynaptoGreen™ C4 and ADVASEP-7  |
| <a href="#">Nerve Terminal Staining Kit II (A)</a>     | <a href="#">70031</a>   | Includes AM1-43 and ADVASEP-7  |
| <a href="#">Nerve Terminal Staining Kit II (B)</a>     | <a href="#">70031-1</a> | Includes AM1-43 and SCAS   |
| <a href="#">Nerve Terminal Staining Kit III</a>        | <a href="#">70032</a>   | Includes SynaptoGreen™ C4 and Sulforhodamine 101                                   |
| <a href="#">Nerve Terminal Staining Kit V</a>          | <a href="#">70034</a>   | 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) [DOI: 10.1074/jbc.M706499200](#)
2. Dev. Cell, 44(1), 56(2018), [DOI: 10.1016/j.devcel.2017.12.014](#)
3. J. Cell Sci, 127(1), 250(2014), [DOI: 10.1242/jcs.140996](#)
4. J. Cell Sci, 134(2), (2021), [DOI: 10.1242/jcs.25773](#)

This datasheet was generated on January 25, 2026 at 01:38:07 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/scas4-sulfonato-calix8arene-sodium-salt/>