

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: SGC5**CATALOG #:** 70057**MOLECULAR INFORMATION:** C₃₃H₅₁Cl₂N₃O₂
MW: 593**PROPERTIES:**

Color & Form	Dark orange solid
Purity	≥ 95% by HPLC
Solubility	Soluble in DMSO or DMF
Absorption/Emission	492/576 nm (in MeOH)
Extinction Coefficient	≈ 58,000

STORAGE AND HANDLING: Store desiccated at ≤ 4 °C. Protect from light, especially in solution.**APPLICATION:** SGC5 is a styryl dye that binds to membranes in a manner similar to the FM dyes. It becomes highly fluorescent upon insertion into lipid bilayers.**References:**Wu, et al., Biophysical Characterization of Styryl Dye-Membrane Interactions,
Biophysical Journal, 97, 101-109

Biotium has the largest collection of high quality nerve terminal dyes available to you at a fraction of the prices of our competitor's products. A comprehensive list of nerve terminal dyes are shown in Table 1.

*FM is a trademark of Molecular Probes, Inc.

Table 1. Related Nerve Terminal Probes

Cat. #	Product Name	Equivalent FM™ dye	$\lambda_{\text{abs}}/\lambda_{\text{em}}$ (nm, in MeOH) ^a	$\lambda_{\text{abs}}/\lambda_{\text{em}}$ (nm, in membranes)	Fixability
70024	AM1-43	none	510/625	479/598	Yes
70036	AM2-10	none	502/625	d	Yes
70042, 70043	SynaptoGreen C1	none	510/625	d	No
70044, 70045	SynaptoGreen C2	FM2-10	505/620	d	No
70023, 70026	SynaptoGreen C3	none	510/625	d	No
70020, 70022	SynaptoGreen C4	FM1-43	510/625	479/598	No
70046, 70047	SynaptoGreen C5	FM1-84	510/625	d	No
70040, 70041	SynaptoRed C1	none	543/b	c	No
70021, 70027	SynaptoRed C2	FM4-64	543/b	c	No

^a The spectra of styryl dyes are known to undergo a large blue shift (30 to 40 nm) when going from polar media (MeOH) to nonpolar media (membranes). ^b AM4-64, SynaptoRed C1 and SynaptoRed C2 have similar absorption spectra and their emission spectra are too weak to measure in MeOH. ^c Data have not been determined yet. However, excitation/emission wavelength setting at 515/640 nm has been used for detection of yeast vacuole membrane staining with SynaptoRed C2 (FM4-64) [J. Cell. Biol. 128, 779(1995)], and this setting should also be applicable to SynaptoRed C1. ^d Data are not available, but expected to be similar to that for AM1-43 or SynaptoGreen C4.

TOXICITY: Unknown.

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--