

Product Information

SGC5

Catalog Number: 70057

Unit Size: 1 mg

Storage and Handling

Store desiccated at 4°C. Protect from light, especially in solution. Product is stable for at least 12 months from date of receipt when stored as recommended.

Molecular Information: C₃₃H₅₁Br₂N₃O₂

Molecular Weight: 682

Color and Form: Dark orange solid

Solubility: Soluble in DMSO or DMF

Absorption/Emission: 484/563 nm (in MeOH)

Extinction Coefficient: 58,000

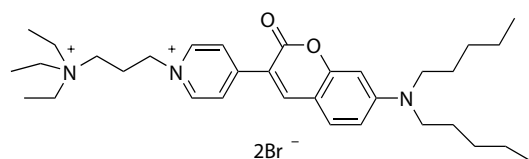


Figure 1. Structure of SGC5.

Product Description

SGC5 is a coumarin styryl dye that binds to membranes in a manner similar to the FM® dyes. It becomes highly fluorescent upon insertion into lipid bilayers.

References

1. Biophysical J 97, 101 (2009).

Related Products

Cat. No.	Product
70056	SGC3
70024	AM1-43
70038	AM1-44
70036	AM2-10
70051	AM3-25
70025	AM4-64
70039	AM4-65
70050	AM4-66
70053	HM1-43
70042	SynaptoGreen™ C1
70044	SynaptoGreen™ C2 (equivalent to FM®2-10)
70023	SynaptoGreen™ C3
70020	SynaptoGreen™ C4 (equivalent to FM®1-43)
70046	SynaptoGreen™ C5 (equivalent to FM®1-84)
70048	SynaptoGreen™ C18 (equivalent to FM®3-25)
70040	SynaptoRed™ C1
70021	SynaptoRed™ C2 (equivalent to FM®4-64)
70019	SynaptoRed™ C2M (equivalent to FM®5-95)

Please visit our website at www.biotium.com for information on our membrane potential dyes, nerve terminal dyes, and other reagents for neuroscience research.

FM is a registered trademark of Thermo Fisher Scientific. Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.