

SGC3

SGC3 is a coumarin styryl dye that binds to membranes in a manner similar to the FM $^{\odot}$ dyes. It becomes highly fluorescent upon insertion into lipid bilayers.



Call us: 800-304-5357

Product attributes

| Probe cellular localization | Membrane/vesicular |
|-----------------------------|-----------------------|
| For live or fixed cells | For live/intact cells |
| Assay type/options | Real-time imaging |
| Excitation/Emission | 484/563 nm |

Product Description

SGC3 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.

- λ_{Ex}/λ_{Em}: 484/563 nm (in MeOH)
- Dark Orange solid soluble in water
- MW: 625

FM is a registered trademark of Thermo Fisher Scientific.

References

- 1. Biophysical J 97, 101 (2009).
- 2. Biochemistry (2011), doi: 10.1021/bi2006288

This datasheet was generated on November 26, 2025 at 07:14:01 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/sgc3/