SGC₅

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.



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

Call us: 800-304-5357 Email: btinfo@biotium.com

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.

- λ_{Ex}/λ_{Em}: 484/563 nm (in MeOH)
- Dark Orange solid soluble in water
- MW: 682

FM is a registered trademark of Thermo Fisher Scientific.

References

- 1. Biophysical J, 97, 101 (2009), <u>DOI: 10.1016/j.bpj.2009.04.028</u> 2. Biochemistry, 50, 35 (2011), <u>DOI: 10.1021/bi2006288</u>

This datasheet was generated on November 2, 2025 at 12:56:45 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/sgc5/