

SGC3

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.



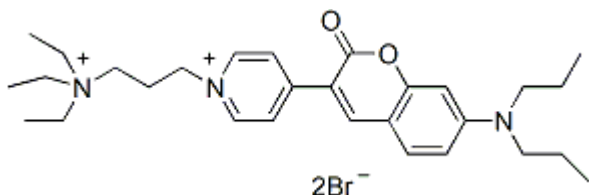
Product attributes

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.

- $\lambda_{Ex}/\lambda_{Em}$: 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 June 22, 2026 at 08:15:23 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/sgc3/>