Sulforhodamine G

Similar to other sulforhodamine dyes, the highly water-soluble sulforhodamine G can also be used as a polar tracer for the studies of cell morphology and neuronal cell-cell communication (1-3).



Product attributes

CAS number	5873-16-5
Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Red
Excitation/Emission	529/548 nm

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

Product Description

Similar to other sulforhodamine dyes, the highly water-soluble sulforhodamine G can also be used as a polar tracer for the studies of cell morphology and neuronal cell-cell communication (1-3).

- $\lambda_{Ex}/\lambda_{Em} = 529/548 \text{ nm}$
- Orange red solid soluble in water
- Store at room temperature
- C25H25N2NaO7S2
- MW: 530.6
- [5873-16-5]

References

- 1. Meth Enzymol 221, 234 (1993).
- 2. J Biol Chem 267, 18424 (1992).
- 3. Plant Cell Environ 17, 257 (1994).

This datasheet was generated on November 14, 2025 at 06:30:29 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/sulforhodamine-g/