MTS-Sulforhodamine 101 (Texas Red®)

MTS-sulforhodamine 101 is a fast-reacting and highly selective thiol-reactive rhodamine dye.



Product attributes

Excitation/Emission 583/603 nm

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

Product Description

MTS-sulforhodamine 101 is a fast-reacting and highly selective thiol-reactive rhodamine dye.

- λEx/λEm (MeOH) = 583/603 nm
- Dark Red solid soluble in DMF or DMSO
- Store at -20°C and protect from light
- C34H37N3O8S4
- MW: 743.94

This datasheet was generated on November 4, 2025 at 04:28:22 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mts-sulforhodamine-101-mts-texas-red/