

4-Sulforhodamine 101 Cadaverine, Trifluoroacetate Salt

4-Sulforhodamine 101 cadaverine can be coupled to activated carboxy groups or sulfonyl chlorides.



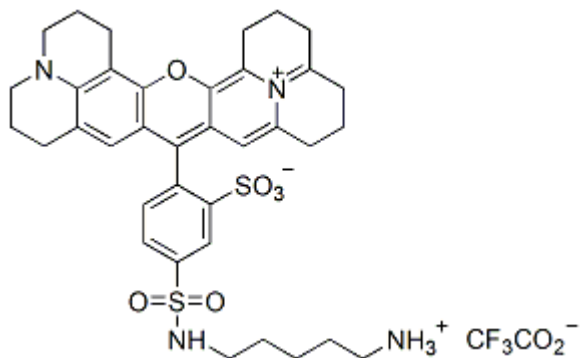
Product attributes

Excitation/Emission	583/603 nm

Product Description

4-Sulforhodamine 101 cadaverine can be coupled to activated carboxy groups or sulfonyl chlorides.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 583/603 nm
- Dark purple solid soluble in DMSO or DMF
- Store at 4°C and protect from light
- $C_{38}H_{43}F_3N_4O_8S_2$
- MW: 804.90



This datasheet was generated on December 7, 2025 at 01:51:30 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/4-sulforhodamine-101-cadaverine-trifluoroacetate-salt/>