

2-Sulforhodamine 101 Cadaverine, Trifluoroacetate Salt

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



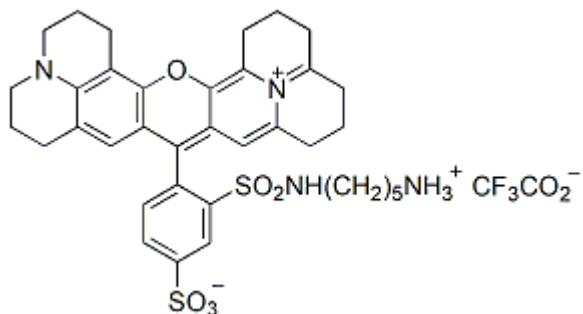
Product attributes

Excitation/Emission	583/603 nm
---------------------	------------

Product Description

2-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 January 5, 2026 at 06:48:54 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/2-sulforhodamine-101-cadaverine-trifluoroacetate-salt/>