## 2-Sulforhodamine 101 Cadaverine, Trifluoroacetate Salt

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

## nictum 6

## **Product attributes**

Excitation/Emission 5

583/603 nm

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

## **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<sub>38</sub>H<sub>43</sub>F<sub>3</sub>N<sub>4</sub>O<sub>8</sub>S<sub>2</sub>
- MW: 804.90

This datasheet was generated on November 4, 2025 at 10:32:58 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/2-sulforhodamine-101-cadaverine-trifluoroacetate-salt/">https://biotium.com/product/2-sulforhodamine-101-cadaverine-trifluoroacetate-salt/</a>