

Fluorescein cadaverine

Fluorescein cadaverine can be coupled to activated carboxy groups or sulfonyl chlorides.



Call us: 800-304-5357 Product attributes

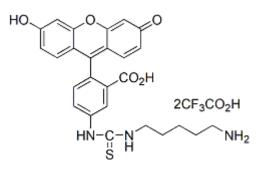
Excitation/Emission

493/517 nm (pH 9)

Product Description

Fluorescein cadaverine can be coupled to activated carboxy groups or sulfonyl chlorides.

- $\lambda Ex/\lambda Em(pH9) = 493/517 nm$
- Yellow solid soluble in DMSO, DMF, or pH>6.5 buffers
- Store at 4°C and protect from light
- C30H27F6N3O9S
- MW: 719.61



This datasheet was generated on November 28, 2025 at 07:17:54 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/fluorescein-cadaverine/