

## Fluorescein Hexylamine

Fluorescein hexylamine can be coupled to activated carboxy groups or sulfonyl chlorides. This product could be used as a direct replacement to fluorescein cadaverine (cat#: 92000).



### Product attributes

Excitation/Emission	493/517 nm (pH 9)
---------------------	-------------------

### Product Description

Fluorescein hexylamine can be coupled to activated carboxy groups or sulfonyl chlorides. This product could be used as a direct replacement to fluorescein cadaverine (cat#: 92000)

- $\lambda_{Ex}/\lambda_{Em}$  (pH 9) = 493/517 nm
- Yellow solid soluble in DMSO, DMF, or pH>6.5 buffers
- Store at 4 °C and protect from light
- $C_{29}H_{27}F_3N_2O_8$
- MW: 588.54

