

## Dansyl cadaverine [5-Dimethylaminonaphthalene-1-(N-(5-aminopentyl))sulfonamide]

Catalog Number: 810

Unit Size: 25 mg

### Product Details

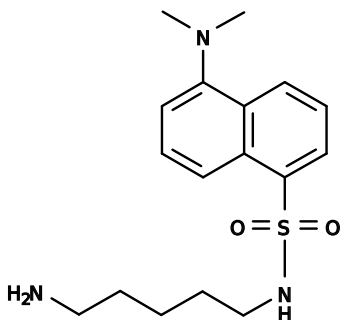
---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Solid yellow-green
Molecular Weight	335.46
Soluble In	DMSO
Chemical Structure	



### Spectral Properties

---

Excitation Wavelength	335 nm
Emission Wavelength	518 nm

### Applications

---

Dansyl cadaverine is an amine-containing Dansyl building block. Dansyl cadaverine is also the most popular fluorescent transglutaminase substrate to label proteins by transamidation. It has also been used to prepare other small fluorescent biomolecules via amidation or reductive amination. It should be noted that the Dansyl fluorescence is highly environment-sensitive. Dansyl dyes have environmentally sensitive fluorescence quantum yields and emission maxima along with large Stokes shifts. This environment-sensitive fluorescence property has made Dansyl dyes an important tool for biophysical studies. They are particularly useful for preparing fluorescent drug or ligand analogs that are expected to bind to hydrophobic sites in proteins or membranes or other biological receptors. Dansyl protein conjugates have fluorescence lifetimes of 10-20 nanoseconds.