

## Dansyl chloride [5-Dimethylaminonaphthalene-1-sulfonyl chloride] \*CAS 605-65-2\*

Catalog Number: 811

Unit Size: 100 mg

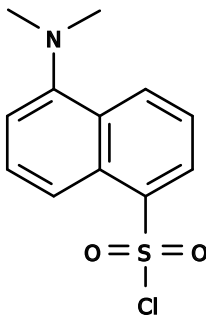
### Product Details

---

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

### Chemical Properties

---

Appearance	Solid yellow
Molecular Weight	269.75
Soluble In	DMF
Chemical Structure	

### Spectral Properties

---

Excitation Wavelength	335 nm
Emission Wavelength	518 nm

### Applications

---

Dansyl chloride is non-fluorescent until it reacts with amines. The resulting Dansyl amides have environmentally sensitive fluorescence quantum yields and emission maxima along with large Stokes shifts. This environment-sensitive fluorescence property has made Dansyl chloride an important tool for biophysical studies. It is 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.