

## Product Information Sheet

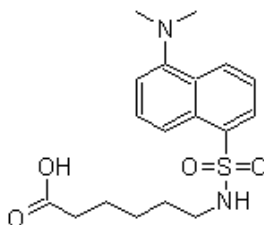
### Ordering Information

Product Number: 812  
Product Name: Dansyl-X, acid  
Unit Size: 5 g  
Storage Conditions: <-15 °C and kept from light and moisture  
Expiration Date: 12 months upon receiving

### Chemical and Spectral Properties

Molecular Weight: 364.46

Chemical Structure:



Soluble in: DMSO  
Excitation Wavelength: 333 nm  
Emission Wavelength: 518 nm

### Application Notes

Compared to Dansyl chloride, Dansyl-X acid gives better yield for labeling peptides and nucleotides. 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.