

Product Information Sheet

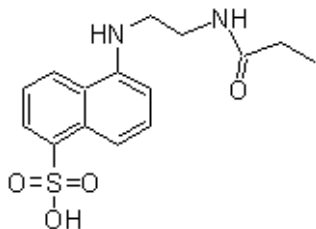
Ordering Information

Product Number: 618
Product Name: EDANS iodoacetamide
Unit Size: 25 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 434.25

Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 335 nm
Emission Wavelength: 493 nm

Application Notes

EDANS iodoacetamide [5-((((2-Iodoacetyl)amino)ethyl)amino)naphthalene-1-sulfonic acid] is a thiol-reactive EDANS derivative. It is water-soluble, and has the same spectral properties with EDANS.