

## EDANS acid [5-((2-Aminoethyl)amino)naphthalene-1-sulfonic acid] \*CAS 50402-56-7\*

Catalog Number: 610

Unit Size: 1 g

### Product Details

---

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

### Chemical Properties

---

Appearance	Solid off-white
Molecular Weight	266.32
Soluble In	DMSO
Chemical Structure	

### Spectral Properties

---

Excitation Wavelength	336 nm
Emission Wavelength	455 nm

### Applications

---

EDANS is one of the most popular donors for developing FRET-based nucleic acid probes (such as Molecular Beacons) and protease substrates. EDANS is often paired with DABCYL or DABSYL in FRET-based probes. Its fluorescence is environment-sensitive.