

## EDANS sodium salt [5-((2-Aminoethyl)aminonaphthalene-1-sulfonic acid, sodium salt] \*CAS 100900-07-0\*

Catalog Number: 615, 616

Unit Size: 1 g, 10 g

### Product Details

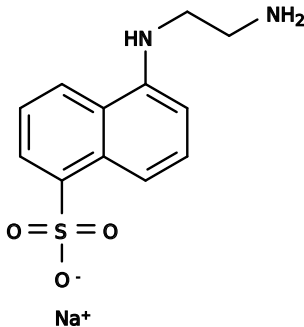
---

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

### Chemical Properties

---

Appearance	Solid light yellow
Molecular Weight	288.30
Soluble In	DMSO
Chemical Structure	



### Spectral Properties

---

Excitation Wavelength	336 nm
Emission Wavelength	455 nm

### Applications

---

EDANS sodium salt is much more water-soluble than EDANS acid. It 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.