

EDANS acid [5-((2-Aminoethyl)amino)naphthalene-1-sulfonic acid]

Catalog Number: 611

Unit Size: 10 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

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.