

EDANS C5 maleimideCatalog number: 619
Unit size: 5 mg**Product Details**

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

Chemical Properties

Appearance	Solid
Molecular Weight	417.44
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. EDANS C5 maleimide is a thiol-reactive EDANS derivative. It is water-soluble, and has the same spectral properties with EDANS.