520 Mercury Drive, Sunnyvale, CA 94085 Tel: 408-733-1055; Fax: 408-733-1304

www.aatbio.com; E-Mail: info@aatbio.com

Product Information Sheet

Ordering Information

Product Number: 610

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

Unit Size: 1g

Storage Conditions: <-15 °C and kept from light and moisture

Expiration Date: 24 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight: 266.32

Chemical Structure:

Appearance: Off-white solid Soluble in: DMSO or DMF

Excitation Wavelength: 335 nm Emission Wavelength: 493 nm

Application Notes

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.