



Product Data Sheet

Product Name: QXL[®] 490 acid, SE

Size: 5 mg

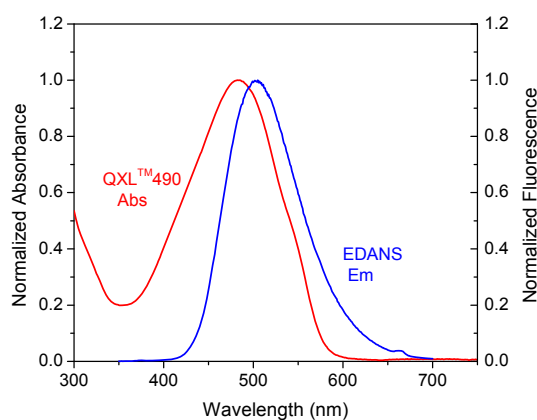
Catalog Number: AS-81826-5

Molecular Weight: 474.49

Solvent: DMF or DMSO

Product Description: QXL[®] 490 dyes are optimized quenchers for EDANS, AMCA and most coumarin fluorophores. Their absorption spectra perfectly overlap with the fluorescence spectra of EDANS and AMCA. This characteristic makes QXL[®] 490 dyes excellent quenchers for EDANS, AMCA and coumarin fluorophores. QXL[®] 490 acid, SE readily reacts with amino-containing compounds. Product purity is over 95%.

Spectra: Maximum Ex/Em wavelength is 495 / none nm.



Application: Fluorescence Resonance Energy Transfer (FRET)

Stock Solution Preparation: Prepare quencher stock solution in high quality anhydrous DMSO. Completely dissolve all quencher content by vortexing. Quencher solution should be prepared fresh immediately before an experiment. Extensive storage may reduce activity. Any solution containing the quencher should be protected from light.

Shelf Life and Storage: Shelf life is two years if stored at -20C, desiccated and protected from light.

This product is for *in vitro* research use only