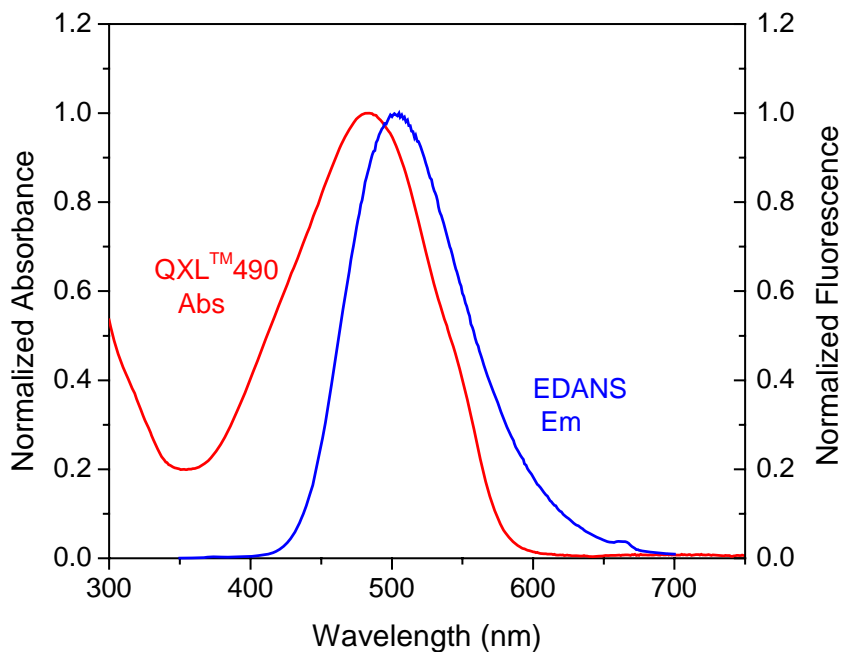




Product Data Sheet

Product Name:	QXL™ 490 hydrazide
Size:	5 mg
Catalog Number:	81846
Price (US\$):	\$225
Molecular Weight:	391.45
Solvent:	DMSO or DMF
Product Description:	QXL™ 490 dyes are the 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 hydrazide readily reacts with carbonyl-containing compounds. Product purity is over 95%.
Spectra:	Maximum Ex/Em wavelength is 485 / none nm.



Stock solution preparation: Make quencher stock solution in high quality anhydrous DMSO. Completely dissolve all quencher content by vortexing. Quencher solution should be prepared fresh right before 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.

Datasheet download: [Product data sheet](#)

MSDS:

This product is for *in vitro* research use only.