



Product Data Sheet

Product Name: QXL[®] 610 vinyl sulfone *Thiol-Reactive*

Size: 25 mg

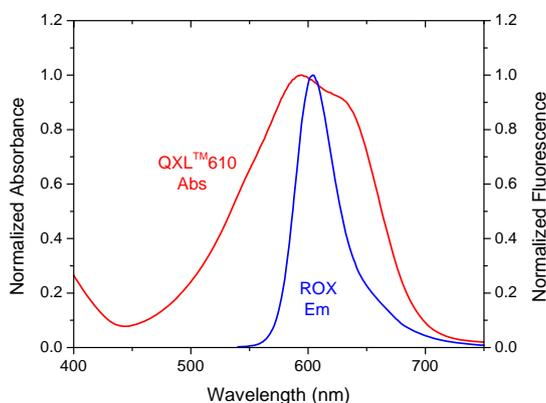
Catalog Number: 81817

Molecular Weight: 506.49

Solvent: Water or DMSO

Product Description: QXL[®] 610 dyes are optimized quenchers for ROX and Texas Red[®] fluorophores. Their absorption spectra perfectly overlap with the fluorescence spectra of ROX and Texas Red[®] dyes. This characteristic makes QXL[®] 610 dyes excellent quenchers for ROX and Texas Red[®] fluorophores. QXL[®] 610 vinyl sulfone readily reacts with thiol-containing compounds. Product purity is over 95%.

Spectra: Maximum Ex/Em wavelength is 628 nm /none nm.



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.