



Product Data Sheet

Product Name: QXL[®] 670 acid, SE

Size: 5 mg

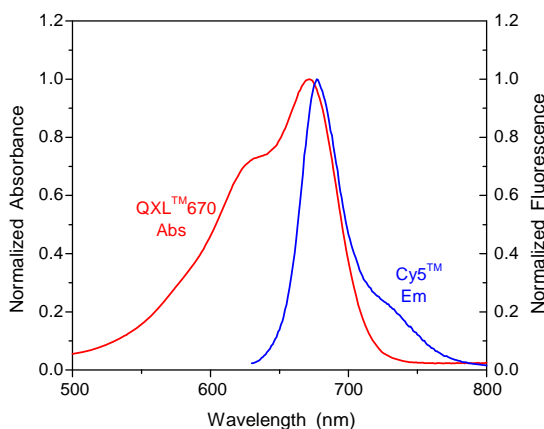
Catalog Number: AS-81841

Molecular Weight: 721.34

Solvent: DMF or DMSO

Product Description: QXL[®] 670 dyes are optimized quenchers for Cy5[®] and Cy5[®]-like fluorophores such as HiLyte Fluor[™] 647. Their absorption spectra perfectly overlap with the fluorescence spectra of Cy5[®]-like dyes. This characteristic makes QXL[®] 670 dyes excellent quenchers for Cy5[®]-like fluorophores. QXL[®] 670 acid, SE readily reacts with amino-containing compounds.

Spectra: Maximum Ex/Em wavelength is 668 nm /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.