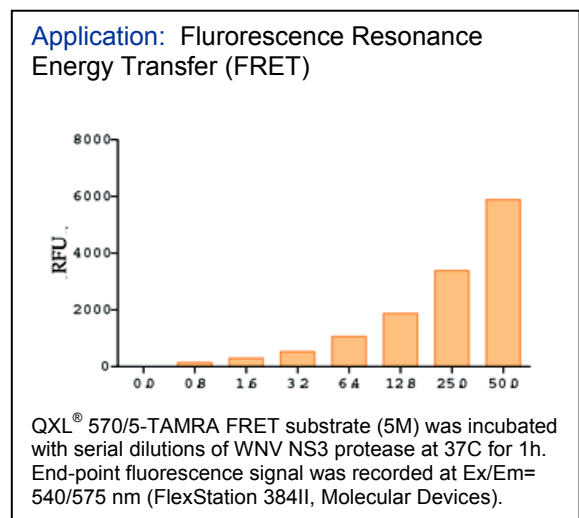
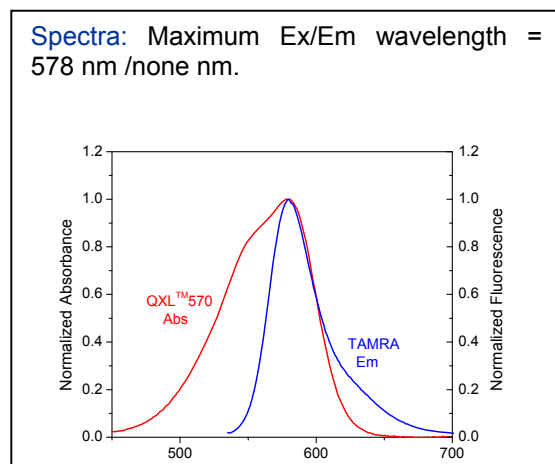




Product Data Sheet

Product Name: QXL[®] 570 acid, SE
Size: 5 mg
Catalog Number: AS-81836-5
Molecular Weight: 775.75
Solvent: DMF or DMSO

Product Description: QXL[®] 570 dyes are optimized quenchers for rhodamines (such as TAMRA, sulforhodamine B, ROX) and Cy3[®] fluorophores. Their absorption spectra perfectly overlap with the fluorescence spectra of TAMRA, sulforhodamine B, ROX and Cy3[®]. This characteristic makes QXL[®] 570 dyes excellent quenchers for rhodamines (such as TAMRA, sulforhodamine B, ROX) and Cy3[®] fluorophores. QXL[®] 570 acid, SE readily reacts with amino-containing compounds. Product purity is over 95%.



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