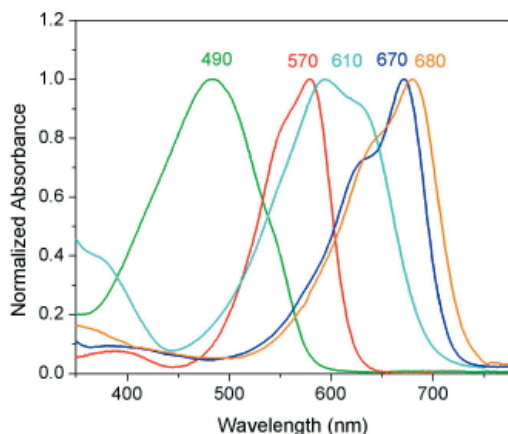


QXL™ Quenchers

* SIMPLY THE BEST *



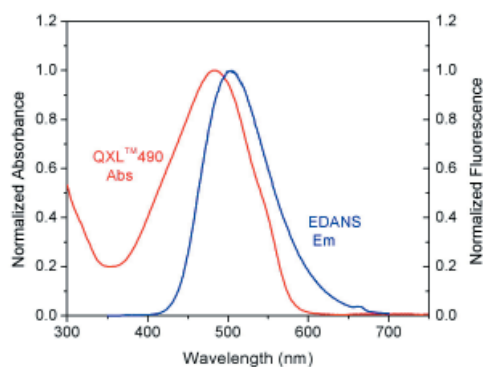
- **Most powerful:** explore new FRET potentials not achievable with other quenchers
- **Versatile reactive forms:** convenient for labeling different biomolecule functionalities
- **Complete set of quenchers:** matching the most commonly used fluorescent donors and covering the full spectrum
- **Enhanced value:** competitive price with the best performance

Product	Size	Catalog #
QXL™ 490 acid	10 mg	81825-10
QXL™ 490 acid, SE	5 mg	81826-5
QXL™ 490 C2 amine	5 mg	81828-5
QXL™ 490 hydrazide	5 mg	81846
QXL™ 490 C2 maleimide	5 mg	81827-5
QXL™ 570 acid	10 mg	81835-10
QXL™ 570 acid, SE	5 mg	81836-5
QXL™ 570 C2 amine	5 mg	81837
QXL™ 570 hydrazide	5 mg	81848
QXL™ 570 C2 maleimide	5 mg	81838
QXL™ 610 acid	10 mg	81815-10
QXL™ 610 acid, SE	5 mg	81816-5
QXL™ 610 C2 amine	5 mg	81818-5
QXL™ 610 hydrazide	10 mg	81849
QXL™ 610 vinyl sulfone *Thiol-Reactive*	25 mg	81817
QXL™ 670 acid	10 mg	81840
QXL™ 670 acid, SE	5 mg	81841
QXL™ 670 C2 amine	5 mg	81842
QXL™ 670 hydrazide	5 mg	81850
QXL™ 670 C2 maleimide	5 mg	81843
QXL™ 680 acid	10 mg	81855
QXL™ 680 acid, SE	5 mg	81851
QXL™ 680 C2 amine	1 mg	81852
QXL™ 680 hydrazide	1 mg	81853
QXL™ 680 C2 maleimide	1 mg	81854

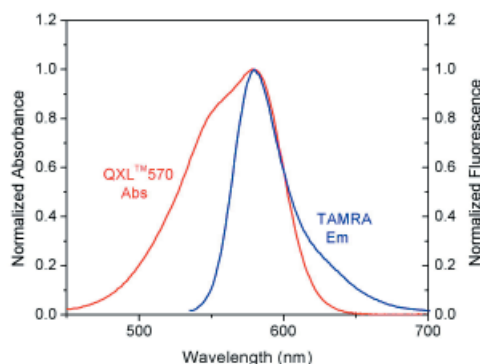
Chemical reactivities and spectral properties of QXL™ Quenchers.

Quencher (Acceptor)	λ_{\max} (in nm)	Amine-Reactive	Thiol-Reactive	Carboxy-Reactive (NH ₂ -Containing)	Recommended FRET Donor
QXL™ 490	488	QXL™ 490 acid; QXL™ 490 acid, SE	QXL™ 490 C2 maleimide	QXL™ 490 C2 amine, QXL™ 490 hydrazide	EDANS, AMCA, HiLyte Fluor™ 430
QXL™ 570	538, 577	QXL™ 570 acid; QXL™ 570 acid, SE	QXL™ 570 C2 maleimide	QXL™ 570 C2 amine, QXL™ 570 hydrazide	Cy3™, TAMRA, ROX, HiLyte Fluor™ TMR
QXL™ 610	594, 628	QXL™ 610 acid; QXL™ 610 acid, SE	QXL™ 610 C2 maleimide	QXL™ 610 C2 amine, QXL™ 610 hydrazide	ROX, Sulforhodamine 101, HiLyte Fluor™ TR
QXL™ 670	668	QXL™ 670 acid; QXL™ 670 acid, SE	QXL™ 670 C2 maleimide	QXL™ 670 C2 amine, QXL™ 670 hydrazide	Cy5™, HiLyte Fluor™ 647
QXL™ 680	679	QXL™ 680 acid; QXL™ 680 acid, SE	QXL™ 680 C2 maleimide	QXL™ 680 C2 amine, QXL™ 680 hydrazide	Cy5™, HiLyte Fluor™ 647

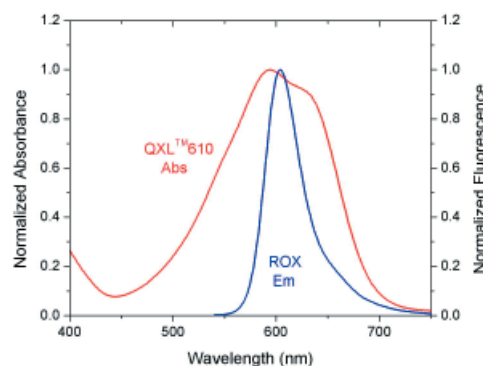
Note: Cy3™ and Cy5™ are trademarks of Amersham Biosciences.



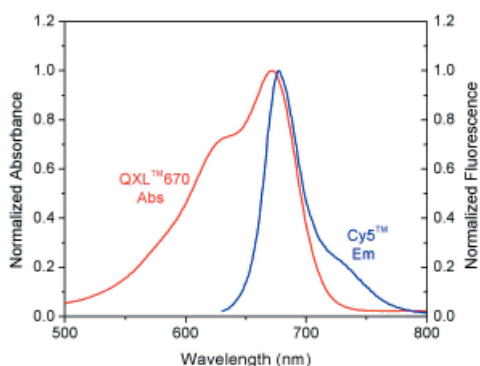
FRET Spectral overlap of EDANS and QXL™ 490.



FRET Spectral overlap of TAMRA and QXL™ 570.



FRET Spectral overlap of ROX and QXL™ 610.



FRET Spectral overlap of Cy5™ and QXL™ 670.