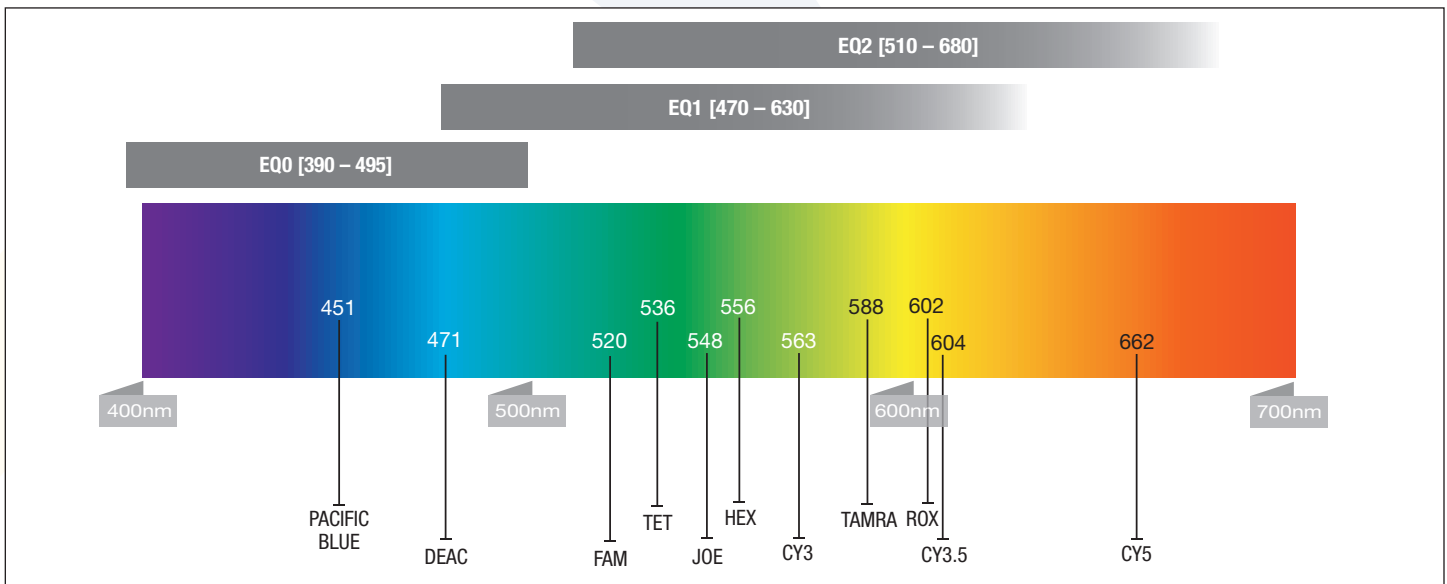




EQ Quencher Dyes for FRET Applications

Amine-Reactive Quencher Dyes

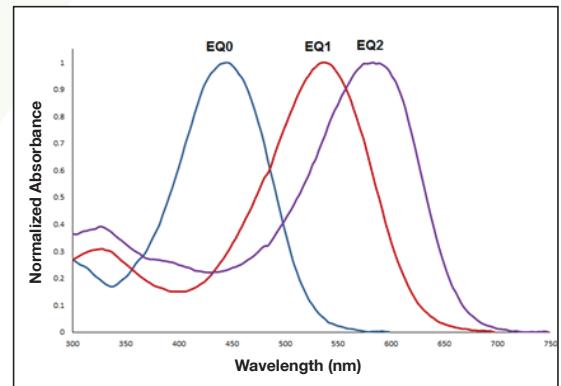
EQ Quencher Dyes are Compatible With a Wide Range of Fluorescent Dyes



EQ quencher dyes are suitable as acceptors in fluorescence resonance energy transfer (FRET) applications due to their broad visible absorption but no detectable fluorescence emission.

- Commonly used in molecular biology FRET applications to quench fluorophores
- NHS ester quencher dye can be conjugated to amine-labeled probes
- Enable multiplexing of probes with the use of different reporter-quencher dye pairs
- Compatible with most common commercially available fluorophores including Alexa Fluor®, ATTO, DyLight®, and BD Horizon dyes

Normalized Absorbance of EQ Dyes



Normalized absorbance of EQ dyes showing their range of quenching through FRET.

RELATED PRODUCTS	
PRODUCT NAME	PRODUCT #
EQ0 NHS ester	ENZ-CHM208
EQ1 NHS ester	ENZ-CHM165
EQ2 NHS ester	ENZ-CHM166

Accelerate Your Research:
www.enzolifesciences.com

GLOBAL HEADQUARTERS
 Enzo Life Sciences Inc.
 10 Executive Boulevard
 Farmingdale, NY 11735
 Phone: 800.942.0430
 Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE
 Enzo Life Sciences (ELS) AG
 Industriestrasse 17
 CH-4415 Lausen, Switzerland
 Phone: +41 61 926 8989
 Fax: +41 61 926 8979
info-eu@enzolifesciences.com