

# EQ0 NHS ester

## Amine-reactive quencher dye

EQ0 NHS ester is an amine-reactive quencher dye. It has a broad visible absorption but no detectable fluorescence emission, making it useful as an acceptor in fluorescence resonance energy transfer (FRET) applications.

It can be used to quench reporter dyes: AMCA, AMCA-X, DEAC, DEAC-X, MCA, MCC, CF™ 350, CF™405S, CF™405M, Pacific Blue™, Cascade Blue®, Marina Blue®, Alexa Fluor® 350, Alexa Fluor® 405, DyLight® 350, DyLight® 405, ATTO 390, ATTO 425, ATTO 465, iFluor™350, iFluor™405, BD Horizon V450, BUV395, eFluor®450.

The EQ dyes are a family of non-fluorescent dyes that are designed for fluorescence quenching in FRET applications.

### Ordering Information

[Order Online »](#)

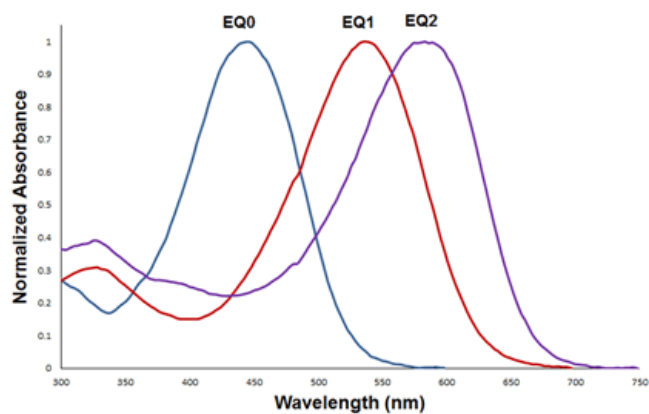
ENZ-CHM208-0001	1mg
-----------------	-----

### Manuals, SDS & CofA

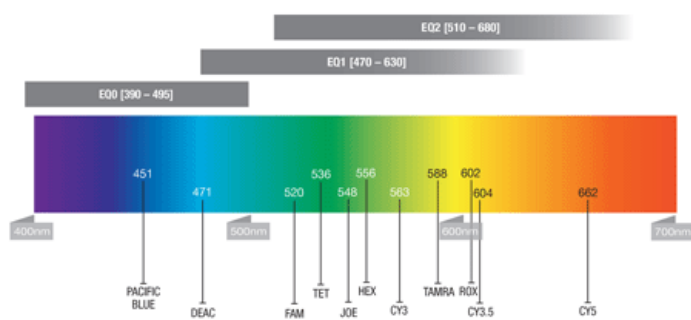
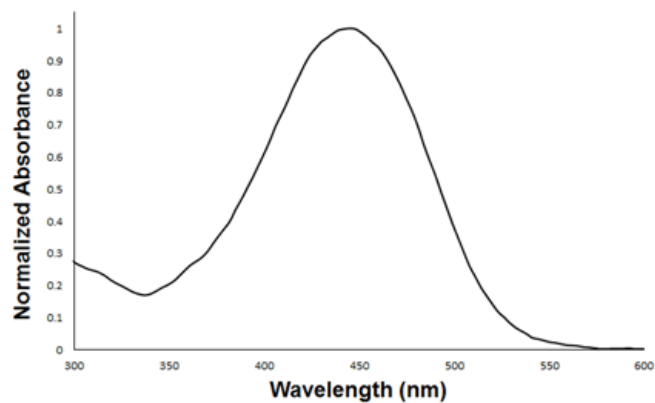
[View Online »](#)

- 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

## Normalized Absorbance of EQ Dyes



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



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Protect from moisture. Protect from light.
Long Term Storage	-20°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	Red brown solid.
Application	Fluorescent detection, FRET
Correction Factor (260nm)	0.48
Correction Factor (280nm)	0.42
Emission Maximum	none
Excitation Maximum	445 nm
Extinction Coefficient	23,600 M-1cm-1
MW	439.5
Purity	≥80% (HPLC)
Solubility	Soluble in DMSO or DMF.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)