# **EQ1 NHS ester**

### Amine-reactive quencher dye

EQ1 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: FAM, Cy2, TET, JOE, VIC, HEX, Cy3, NED, TAMRA, Cy3.5, ROX, Texas Red, Quasar<sup>®</sup> 570, CAL Fluor<sup>®</sup> Gold 540, CAL Fluor<sup>®</sup> Orange 560, CAL Fluor<sup>®</sup> Red 590, CAL Fluor<sup>®</sup> Red 610, Alexa Fluor<sup>®</sup> 430, Alexa Fluor<sup>®</sup> 488, Alexa Fluor<sup>®</sup> 532, Alexa Fluor<sup>®</sup> 546, Alexa Fluor<sup>®</sup> 555, Alexa Fluor<sup>®</sup> 568, Alexa Fluor<sup>®</sup> 594, DyLight<sup>®</sup> 488, DyLight<sup>®</sup> 550, DyLight<sup>®</sup> 594, ATTO 488, ATTO 520, ATTO 532, ATTO Rho6G, ATTO 550, ATTO 565, ATTO Rho11, ATTO 590, ATTO 594.

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

## **Ordering Information**

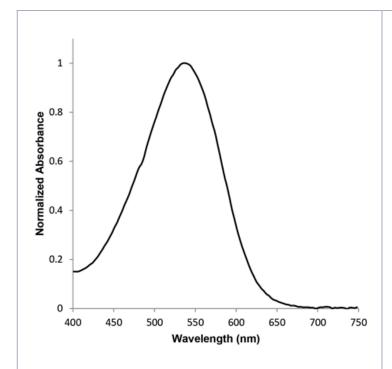
Order Online »

<b>ENZ-CHM165-0001</b> 1mg	
----------------------------	--

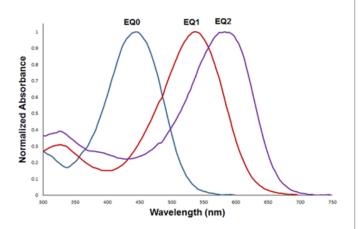
Manuals, SDS & CofA

**View Online »** 

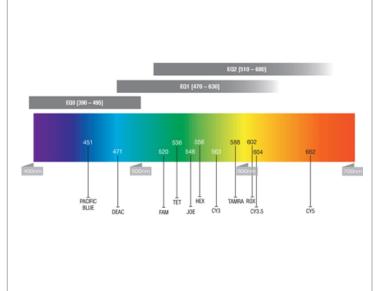
- Commonly used in molecular biology FRET applications to quench fluorophores
- NHS ester quencher dye can be conjugated to amine-labeled probes
- Enable multiplexing of probes with the use of different reporterquencher dye pairs
- Compatible with most common commercially available fluorophores



# 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** Dark red solid.

**Application** Fluorescent detection, FRET

Correction Factor (260nm) 0.34

Correction Factor (280nm) 0.27

Emission Maximum none

Excitation Maximum 535 nm

Extinction Coefficient 39,200 M-1cm-1

**MW** 535.5

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

ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com

European Sales Office

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com