

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[®] 570, CAL Fluor[®] Gold 540, CAL Fluor[®] Orange 560, CAL Fluor[®] Red 590, CAL Fluor[®] Red 610, Alexa Fluor[®] 430, Alexa Fluor[®] 488, Alexa Fluor[®] 532, Alexa Fluor[®] 546, Alexa Fluor[®] 555, Alexa Fluor[®] 568, Alexa Fluor[®] 594, DyLight[®] 488, DyLight[®] 550, DyLight[®] 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

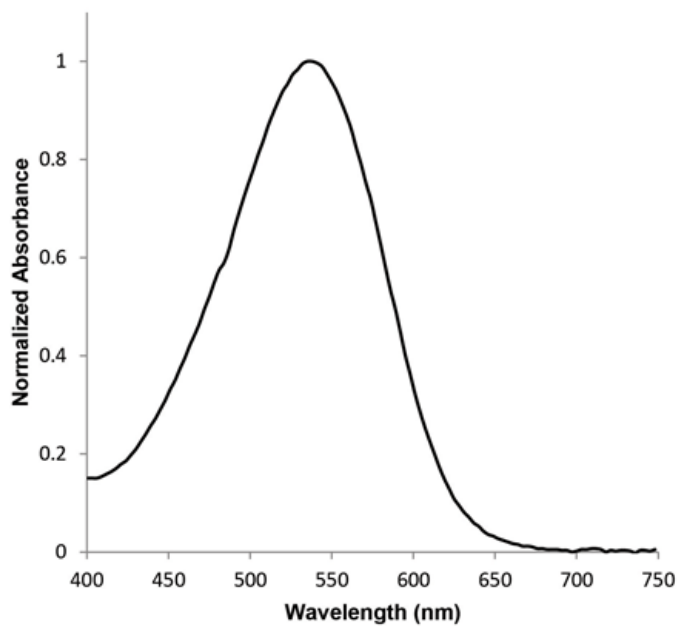
[Order Online »](#)

ENZ-CHM165-0001	1mg
-----------------	-----

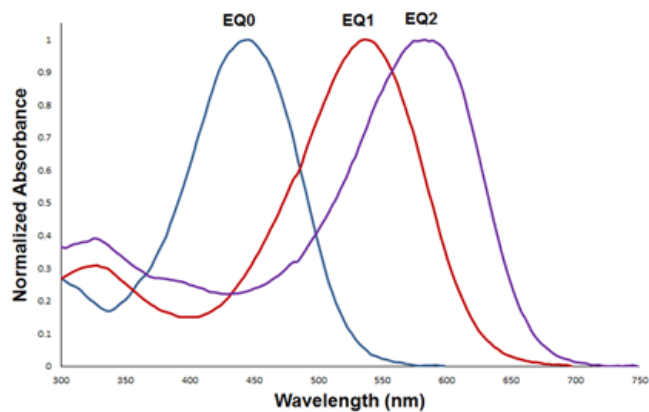
Manuals, SDS & CofA

[View Online »](#)

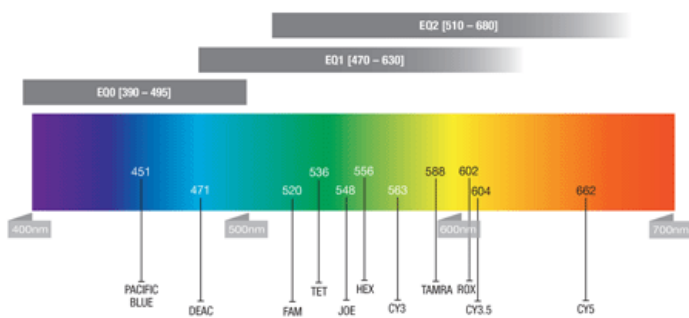
- Commonly used in molecular biology FRET applications to quench fluorophores
- NHS ester quencher dye can be conjugated to amine-labeled probes
- Enable multiplexing of probes with the use of different reporter-quencher dye pairs
- Compatible with most common commercially available fluorophores



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	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](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-
fr@enzolifesciences.com](mailto:info-fr@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)