

Trolox[®]

Cell permeable vitamin E derivative

Cell permeable derivative of vitamin E. Antioxidant. Inhibits peroxynitrite mediated tyrosine nitration and mitochondrial proton permeability. Prevents lipid peroxidation, cytotoxicity and apoptosis induced by cytochrome P450 2E1 in HEP G2 cells.

Citations: 15

[View Online »](#)

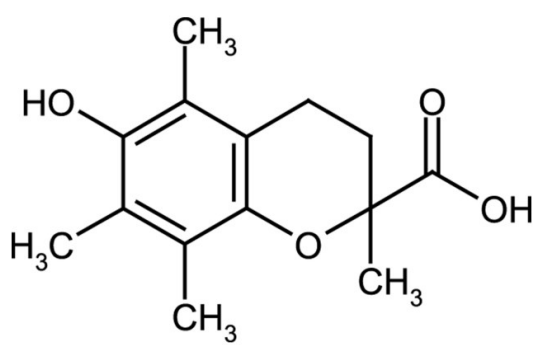
Ordering Information

[Order Online »](#)

ALX-270-267-M100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	6-Hydroxy-2,5,7,8-tetramethylchroman-2-carboxylic acid
Appearance	White to off-white, pale yellow or pale green powder.
CAS	53188-07-1; 56305-04-5
Formula	$C_{14}H_{18}O_4$
Identity	Determined by IR.
MW	250.3
Purity	≥96.5% (Titration (T) NaOH 0.1M)
RTECS	DJ2273000
Solubility	Soluble in 100% ethanol (25mg/ml) or methanol.

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)