

Ebselen

Organoselenium compound. Antioxidant.

Selenium-based inhibitor of protein kinase C, NADPH, 5-lipoxygenase, cyclooxygenase (COX) and NADPH oxidase. Anti-inflammatory antioxidant. Mimics glutathione peroxidase. Inhibits oxidative modifications of low density lipoproteins (LDL). Protects cerebellar granule neurons against 4-hydroxynonenal-induced neuronal death.

Citations: 24

[View Online »](#)

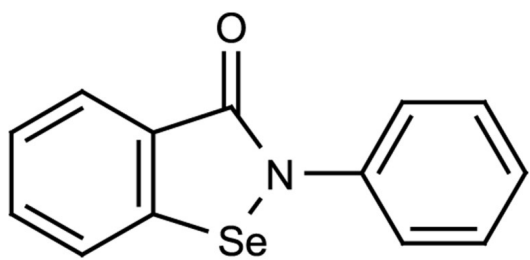
Ordering Information

[Order Online »](#)

ALX-270-097-M025	25mg
------------------	------

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	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White to off-white crystalline solid.
CAS	60940-34-3
Couple Target	COX, Lipoxygenase, NADPH, NADPH oxidase, PKC
Couple Type	Inhibitor
Formula	$C_{13}H_9NOSe$
Identity	Determined by NMR.
MW	274.2
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO or 100% ethanol.

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)