

(±)4-Hydroxynon-2-enal

Lipid peroxidation product

Cytotoxic product of microsomal lipid peroxidation induced by xenobiotics. Reported to be a specific chemical indicator of lipid peroxidation in rat liver microsomes and in myocardial reperfusion injury. Covalently reacts with LDL and HDL during oxidative modification. Modifies proteins by covalent reaction with histidine and lysine residues. Activates TRPA1.

Citations: 1

[View Online »](#)

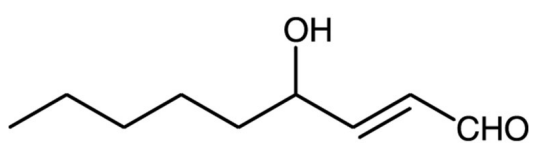
Ordering Information

[Order Online »](#)

BML-HA004-0001	1mg
BML-HA004-0010	10mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Colorless oil.

CAS 75899-68-2

Couple Target TRP channel

Couple Type Activator

Formula $C_9H_{16}O_2$

MW 156.2

Purity ≥97%

Solubility Soluble in 100% ethanol (25mg/ml).



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)