

Decadienal

Induces mitochondrial dysfunction and oxidative stress

Decadienal is a lipid peroxidation product which is an aldehydic bifunctional electrophile. It induces mitochondrial dysfunction and oxidative stress possibly via covalent modification of cytochrome c. It can induce cell proliferation via the p27 pathway in human bronchial epithelial cells.

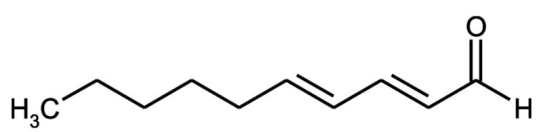
Ordering Information

[Order Online »](#)

BML-HA005-0020	20mg
-----------------------	------

Manuals, SDS & CofA

[View Online »](#)



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.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Pale yellow oil.
CAS	25152-84-5
Formula	$C_{10}H_{16}O$
MW	152.2
Purity	≥95%
Solubility	Soluble at 25mg/ml in DMSO, ethanol, or DMF.



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)