

Sedanolid

GST activator

Inducer of glutathione S-transferases and inhibitor of chemically induced carcinogenesis. Mosquitocidal, nematocidal and antifungal. Was shown to inhibit cyclooxygenase-1 and -2 (COX-1 and COX-2) as well as topoisomerase I and topoisomerase II.

Citations: 4

[View Online »](#)

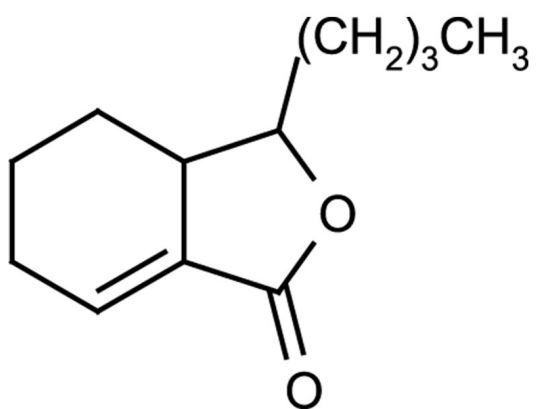
Ordering Information

[Order Online »](#)

ALX-350-229-M100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 3-Butyl-3a,4,5,6-tetrahydro-1(3H)-isobenzofuranone

Appearance White to off-white solid.

CAS 6415-59-4

Couple Target GST

Couple Type Activator

Formula $C_{12}H_{18}O_2$

MW 194.1

Purity ≥98% (UHPLC)

Solubility Soluble in 100% ethanol or ethyl acetate.

Source Isolated from *Apium graveolens* L.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com