

R,S-Sulforaphane

Carcinogenesis inhibitor

R,S-Sulforaphane is a synthetic version of the naturally occurring compound, R-sulforaphane, from Broccoli. It is a racemic mixture of both the R and S isomers.

Potent, selective inducer of phase II detoxification enzymes. Inhibits chemically induced mammary tumor formation in rats. Inhibits LPS-induced HMGB1 (high mobility group box 1) secretion.

Citations: 12

[View Online »](#)

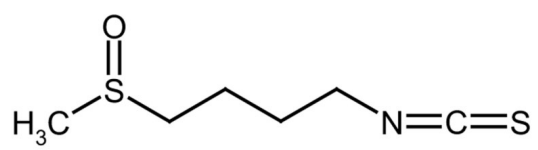
Ordering Information

[Order Online »](#)

ALX-350-232-M025	25mg
------------------	------

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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name D,L-Sulforaphane, 1-Isothiocyanato-4-methylsulfinyl-butane

Appearance Slightly yellowish liquid.

CAS 4478-93-7 (unspecified stereo)

Formula $C_6H_{11}NOS_2$

Identity Identity determined by 1H-NMR and MS.

MI **14:** 8962

MW 177.3

Purity ≥98% (UHPLC)

Solubility Soluble in methanol or DMSO.

Source Synthetic.

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)