

# 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 »](#)

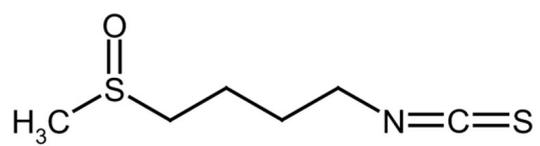
Ordering Information

[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-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)