

# R,S-Sulforaphane, High Purity

Phase II enzyme inducer; Nrf2 activator;  
antioxidant with anticarcinogenic properties.

Sulforaphane is a phase II enzyme inducer present in broccoli. It was found to inhibit chemically induced mammary tumor formation in rats. R,S-sulforaphane is a synthetic compound. Inhibits benzo[a]pyrene-DNA and 1,6-dinitropyrene-DNA adducts formation. Inducer of quinone reductase and glutathione S-transferase. Exerts dose-dependent cytotoxicity, inhibits quiescent cells to re-enter the cell cycle, and induces cell cycle arrest and apoptosis in human colon adenocarcinoma HT29 cells. Inhibitor of phase I cytochrome P450 isoenzymes 2E1 and 1A2. Inducer of multidrug resistance associated protein 2 in primary rat and human hepatocytes.

## Ordering Information

[Order Online »](#)

LKT-S8044-M025	25mg
LKT-S8044-M100	100mg
LKT-S8044-M500	500mg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

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

**CAS** 4478-93-7

**Formula**  $C_6H_{11}NOS_2$

**MI** 13: 9051

**MW** 177.29

**Purity** ≥99%

**Solubility** Soluble in ethanol, methanol, or ethyl acetate. Miscible with water and/or DMSO.



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)