

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-M050	50mg
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

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