

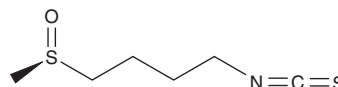
PRODUCT INFORMATION



L-Sulforaphane

Item No. 14797

CAS Registry No.: 142825-10-3
Formal Name: 1-isothiocyanato-4-[(R)-methylsulfinyl]-butane
Synonyms: L-SFN, (R)-SFN, (R)-Sulforaphane
MF: C₆H₁₁NOS₂
FW: 177.3
Purity: ≥98%
UV/Vis.: λ_{max}: 246 nm
Supplied as: A solution in ethanol
Storage: -80°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

L-Sulforaphane is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of L-sulforaphane in these solvents is approximately 16 and 3 mg/ml, respectively.

L-Sulforaphane is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanol solution of L-sulforaphane should be diluted with the aqueous buffer of choice. The solubility of L-Sulforaphane in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

L-Sulforaphane is an isothiocyanate derived from the natural compound glucoraphanin, which is abundant in cruciferous vegetables, including broccoli.¹ It has powerful antioxidant, anti-inflammatory, and anti-carcinogenic effects.^{1,2} L-Sulforaphane, at 40 μM, activates nuclear factor erythroid 2-related factor 2- (Nrf2) mediated gene expression by disrupting its association with Kelch-like ECH-associated protein 1 (Keap1), allowing Nrf2 to enter the nucleus to alter transcription.^{3,4} In this way, sulforaphane induces the expression of phase II detoxification enzymes, which in turn provide diverse beneficial effects.^{1,5}

References

1. James, D., Devaraj, S., Bellur, P., *et al.* Novel concepts of broccoli sulforaphanes and disease: Induction of phase II antioxidant and detoxification enzymes by enhanced-glucoraphanin broccoli. *Nutr. Rev.* **70(11)**, 654-665 (2012).
2. Liang, H. and Yuan, Q. Natural sulforaphane as a functional chemopreventive agent: Including a review of isolation, purification and analysis methods. *Crit. Rev. Biotechnol.* **32(3)**, 218-234 (2012).
3. La Marca, M., Beffy, P., Della Croce, C., *et al.* Structural influence of isothiocyanates on expression of cytochrome P450, phase II enzymes, and activation of Nrf2 in primary rat hepatocytes. *Food Chem. Toxicol.* **50(8)**, 2822-2830 (2012).
4. Dinkova-Kostova, A.T., Holtzclaw, W.D., Cole, R.N., *et al.* Direct evidence that sulfhydryl groups of Keap1 are the sensors regulating induction of phase 2 enzymes that protect against carcinogens and oxidants. *Proc. Natl. Acad. Sci. USA* **99(18)**, 11908-11913 (2012).
5. Munday, R. and Munday, C.M. Induction of phase II detoxification enzymes in rats by plant-derived isothiocyanates: Comparison of allyl isothiocyanate with sulforaphane and related compounds. *J. Agric. Food. Chem.* **52(7)**, 1867-1871 (2004).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/29/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM