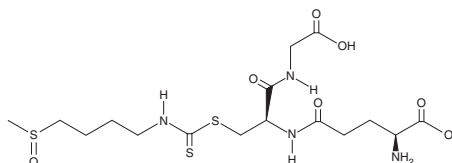


PRODUCT INFORMATION

DL-Sulforaphane Glutathione

Item No. 34731

CAS Registry No.: 289711-21-3
Formal Name: L-γ-glutamyl-S-[[[4-(methylsulfinyl)butyl]amino]thioxomethyl]-L-cysteinyl-glycine
Synonyms: SFN-glutathione, SFN-GSH, Sulforaphane-GSH
MF: C₁₆H₂₈N₄O₇S₃
FW: 484.6
Purity: ≥98%
UV/Vis.: λ_{max}: 253 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

DL-Sulforaphane glutathione is supplied as a solid. A stock solution may be made by dissolving the DL-sulforaphane glutathione in the solvent of choice. DL-Sulforaphane glutathione is soluble in organic solvents such as methanol (warmed), which should be purged with an inert gas. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

DL-Sulforaphane glutathione is a metabolite of the class I and II histone deacetylase (HDAC) inhibitor and anticancer agent sulforaphane (Item No. 10496).¹ It increases mRNA expression of the genes encoding glutathione S-transferase A1 (GSTA1) and the UDP-glucuronosyltransferase (UGT) isoform UGT1A1 in HepG2 cells when used at a concentration of 15 μM.²

References

1. Wang, H., Khor, T.O., Yang, Q., *et al.* Pharmacokinetics and pharmacodynamics of phase II drug metabolizing/antioxidant enzymes gene response by anticancer agent sulforaphane in rat lymphocytes. *Mol. Pharmacol.* **9(10)**, 2819-2827 (2012).
2. Basten, G.P., Bao, Y., and Williamson, G. Sulforaphane and its glutathione conjugate but not sulforaphane nitrile induce UDP-glucuronosyl transferase (UGT1A1) and glutathione transferase (GSTA1) in cultured cells. *Carcinogenesis* **23(8)**, 1399-1404 (2002).

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, 12/08/2022

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