

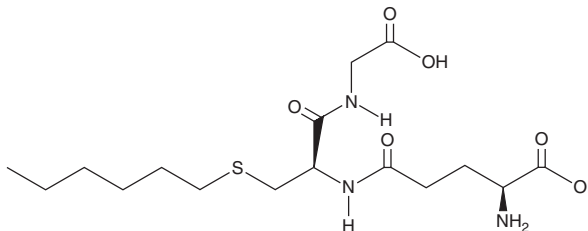
PRODUCT INFORMATION



Hexylglutathione

Item No. 29695

CAS Registry No.: 24425-56-7
Formal Name: L-γ-glutamyl-S-hexyl-L-cysteinyl-glycine
Synonyms: S-Hexylglutathione, NSC 131114
MF: C₁₆H₂₉N₃O₆S
FW: 391.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Hexylglutathione is supplied as a solid. A stock solution may be made by dissolving the RRx-001 in the solvent of choice, which should be purged with an inert gas. RRx-001 is slightly soluble in DMSO.

Hexylglutathione is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, hexylglutathione should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Hexylglutathione has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Hexylglutathione is a mitochondrial membrane-bound microsomal glutathione S-transferase 1 (mtMGST1) inhibitor (IC₅₀ = 0.48 mM).¹ It also inhibits the dopachrome tautomerase activity of macrophage migration inhibitory factor (MIF; IC₅₀ = 3.3 mM).² It has been used as an affinity chromatography ligand to purify glutathione S-transferase (GST) from sheep liver homogenate.³

References

1. Ulziikhishig, E., Lee, K.K., Hossain, Q.S., *et al.* Inhibition of mitochondrial membrane bound-glutathione transferase by mitochondrial permeability transition inhibitors including cyclosporin A. *Life Sci.* **86**(19-20), 726-732 (2010).
2. Swope, M.D., Sun, H.W., Klockow, B., *et al.* Macrophage migration inhibitory factor interactions with glutathione and S-hexylglutathione. *J. Biol. Chem.* **273**(24), 14877-14884 (1998).
3. Reddy, C.C., Burgess, Z.Z., Gong, Z.Z., *et al.* Purification and characterization of the individual glutathione S-transferases from sheep liver. *Arch. Biochem. Biophys.* **224**(1), 87-101 (1983).

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, 12/16/2022

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