

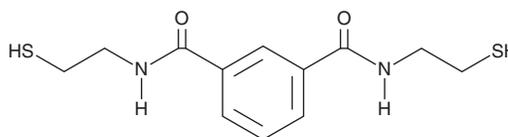
PRODUCT INFORMATION



Emeramide

Item No. 32889

CAS Registry No.: 351994-94-0
Formal Name: N¹,N³-bis(2-mercaptoethyl)-1,3-benzenedicarboxamide
Synonyms: BDTH₂, NBMI
MF: C₁₂H₁₆N₂O₂S₂
FW: 284.4
Purity: ≥98%
Supplied as: A crystalline 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

Emeramide is supplied as a crystalline solid. A stock solution may be made by dissolving the emeraimide in the solvent of choice, which should be purged with an inert gas. Emeramide is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of emeraimide in these solvents is approximately 30 mg/ml.

Emeramide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, emeraimide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Emeramide has a solubility of approximately 0.50 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Emeramide is an antioxidant heavy metal chelator.¹ It prevents methylmercury-induced glutathione (GSH) loss, and cytotoxicity to, isolated mouse aortic endothelial cells when used at a concentration of 50 μM. Emeramide (50 μM) also prevents reactive oxygen species (ROS) production and cytotoxicity induced by bleomycin (Item No. 13877) in primary bovine pulmonary artery endothelial cells.² It increases survival of rats in a model of mercury chloride-induced toxicity when administered at a dose of 2 mg/kg.³

References

1. Secor, J.D., Kotha, S.R., Gurney, T.O., *et al.* Novel lipid-soluble thiol-redox antioxidant and heavy metal chelator, N,N'-bis(2-mercaptoethyl)isophthalamide (NBMI) and phospholipase D-specific inhibitor, 5-fluoro-2-indolyl des-chlorohalopemide (FIPI) attenuate mercury-induced lipid signaling leading to protection against cytotoxicity in aortic endothelial cells. *Int. J. Toxicol.* **30(6)**, 619-638 (2011).
2. Patel, R.B., Kotha, S.R., Sauers, L.A., *et al.* Thiol-redox antioxidants protect against lung vascular endothelial cytoskeletal alterations caused by pulmonary fibrosis inducer, bleomycin: Comparison between classical thiol-protectant, N-acetyl-L-cysteine, and novel thiol antioxidant, N,N'-bis-2-mercaptoethyl isophthalamide. *Toxicol. Mech. Methods* **22(5)**, 383-396 (2012).
3. Clarke, D., Buchanan, R., Gupta, N., *et al.* Amelioration of acute mercury toxicity by a novel, non-toxic lipid soluble chelator N,N'-bis-(2-mercaptoethyl)isophthalamide: Effect on animal survival, health, mercury excretion and organ accumulation. *Toxicol. Environ. Chem.* **94(3)**, 616-640 (2012).

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, 12/13/2022

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