

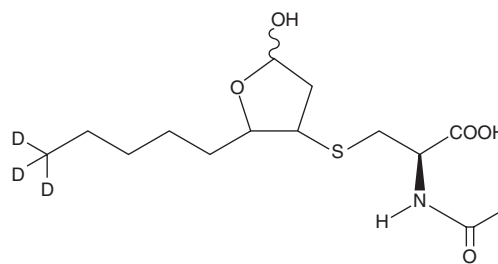
PRODUCT INFORMATION



4-hydroxy Nonenal Mercapturic Acid-d₃

Item No. 9000348

Formal Name: N-acetyl-S-(tetrahydro-5-hydroxy-2-pentyl-3-furanyl)-L-cysteine-11,11,11-d₃
MF: C₁₄H₂₂D₃NO₅S
FW: 322.4
Chemical Purity: ≥98% (4-hydroxy nonenal mercapturic acid)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
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

4-hydroxy Nonenal mercapturic acid-d₃ is intended for use as an internal standard for the quantification of 4-hydroxy nonenal mercapturic acid by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

4-hydroxy Nonenal mercapturic acid-d₃ 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 4-hydroxy nonenal mercapturic acid-d₃ in these solvents is approximately 50 mg/ml.

Description

Peroxidation of polyunsaturated fatty acids in circulating lipid particles and membrane phospholipids leads to transient fatty acid hydroperoxides.¹ Several non-enzymatic routes of decomposition are available to these lipid hydroperoxides, including a β-cleavage reaction which breaks the carbon-carbon bond next to the hydroperoxide and produces alkenals, including 4-hydroxy nonenal (4-HNE). Common ω-6 fatty acids such as linoleic acid, dihomo-γ-linolenic acid, and arachidonic acid can give rise to 4-HNE. 4-HNE is cleared rapidly from the plasma and undergoes enterohepatic circulation as a glutathione conjugate in the rat.² About two thirds of an administered dose of 4-HNE is excreted within 48 hours in the urine, primarily in the form of mercapturic acid conjugates.³ The C-1 aldehyde of 4-HNE is reduced to an alcohol in about half of these metabolites. The remainder are C-1 aldehydes or have been oxidized to C-1 carboxylic acids. These aldehydes and carboxylic acids can also form γ-lactols and γ-lactones, respectively, producing at least four or five end urinary metabolites of 4-HNE *in vivo*.

References

1. Pryor, W.A. and Porter, N.A. Suggested mechanisms for the production of 4-hydroxy-2-nonenal from the autoxidation of polyunsaturated fatty acids. *Free Radic. Biol. Med.* **8**, 541-543 (1990).
2. Laurent, A., Alary, J., Debrauwer, L., *et al.* Analysis in the rat of 4-hydroxynonenal metabolites excreted in bile: Evidence of enterohepatic circulation of these byproducts of lipid peroxidation. *Chem. Res. Toxicol.* **12**, 887-894 (1999).
3. Alary, J., Bravais, F., Cravedi, J.-P., *et al.* Mercapturic acid conjugates as urinary end metabolites of the lipid peroxidation product 4-hydroxy-2-nonenal in the rat. *Chem. Res. Toxicol.* **8**, 34-39 (1995).

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, 02/22/2024

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