

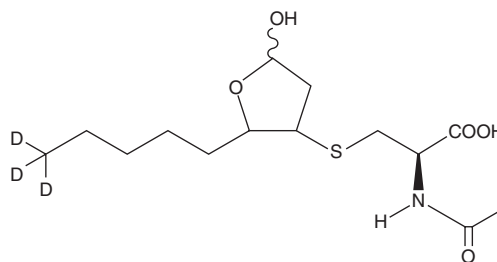
# Product Information



## 4-hydroxy Nonenal Mercapturic Acid-d<sub>3</sub>

Item No. 9000348

<b>Formal Name:</b>	N-acetyl-S-(tetrahydro-5-hydroxy-2-pentyl-3-furanyl)-L-cysteine-11,11,11-d <sub>3</sub>
<b>MF:</b>	C <sub>14</sub> H <sub>22</sub> D <sub>3</sub> NO <sub>5</sub> S
<b>FW:</b>	322.4
<b>Chemical Purity:</b>	≥98% 4-hydroxy nonenal mercapturic acid
<b>Deuterium Incorporation:</b>	≤1% d <sub>0</sub>
<b>Stability:</b>	≥1 year at -80°C
<b>Supplied as:</b>	A solution in ethanol



### Laboratory Procedures

4-hydroxy Nonenal mercapturic acid-d<sub>3</sub> contains three deuterium atoms at the terminal methyl position. It is intended for use as an internal standard for the quantification of 4-hydroxy nonenal mercapturic acid by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that 4-hydroxy nonenal mercapturic acid-d<sub>3</sub> be stored as supplied at -80°C. It should be stable for at least one year.

4-hydroxy Nonenal mercapturic acid-d<sub>3</sub> 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-d<sub>3</sub> in these solvents is approximately 50 mg/ml.

4-hydroxy Nonenal mercapturic acid-d<sub>3</sub> is used as an internal standard for the quantification of 4-hydroxy nonenal mercapturic acid by stable isotope dilution 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).

Peroxidation of polyunsaturated fatty acids in circulating lipid particles and membrane phospholipids leads to transient fatty acid hydroperoxides.<sup>1</sup> 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.<sup>2</sup> 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.<sup>3</sup> 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).

### Related Products

4-hydroxy Nonenal - Item No. 32100 • 4-hydroxy Nonenal Mercapturic Acid - Item No. 32110 • 4-hydroxy Nonenal-d<sub>3</sub> - Item No. 332101 • 4-hydroxy Nonenal - Item No. 9000876 • 4-oxo 2-Nonenal-d<sub>3</sub> - Item No. 10004174 • 4-hydroperoxy 2-Nonenal - Item No. 10004413

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/23/2010

### Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
www.caymanchem.com