

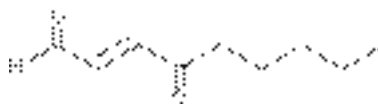
Product Information



4-oxo-2-Nonenal

Item No. 10185

CAS Registry No.: 103560-62-9
Formal Name: 4-oxo-2E-nonenal
Synonym: 4-ONE
MF: C₉H₁₄O₂
FW: 154.2
Purity: ≥98%
Stability: ≥6 months at -80°C
Supplied as: A solution in methyl acetate



Laboratory Procedures

For long term storage, we suggest that 4-oxo-2-nonenal be stored as supplied at -80°C. It should be stable for at least six months.

4-oxo-2-Nonenal is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 4-oxo-2-nonenal in these solvents is approximately 50 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of 4-oxo-2-nonenal is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 4-oxo-2-nonenal in PBS (pH 7.2) is approximately 0.5 mg/ml. Store aqueous solutions of 4-oxo-2-nonenal on ice and use within 12 hours of preparation. Although the aqueous solutions of 4-oxo-2-nonenal may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

4-hydroxy Nonenal is a lipid peroxidation product derived from oxidized ω-6 polyunsaturated fatty acids such as arachidonic acid and linoleic acid.^{1,2} It exhibits various biological activities such as cytotoxicity, growth inhibiting activity, genotoxicity, and chemotactic activity and has been widely used as a marker of lipid peroxidation.¹⁻³ 4-oxo-2-Nonenal is a more recently identified product of lipid peroxidation.⁴⁻⁶ It actively modifies histidine and lysine residues on proteins and causes protein cross-linking.^{7,8} 4-oxo-2-Nonenal also modifies 2'-deoxyguanosine, further implicating lipid peroxidation in mutagenesis and carcinogenesis.⁴

References

1. Pryor, W.A. and Porter, N.A. *Free Radic. Biol. Med.* **8**, 541-543 (1990).
2. Esterbauer, H., Schaur, R.J., and Zollner, H. *Free Radic. Biol. Med.* **11**, 81-128 (1991).
3. Sodom, R.S. and Chung, F.-L. *Cancer Res.* **48**, 320-323 (1988).
4. Rindgen, D., Nakajima, M., Wehrli, S., *et al. Chem. Res. Toxicol.* **12**, 1195-1204 (1999).
5. Lee, S.H. and Blair, I.A. *Chem. Res. Toxicol.* **13**, 698-702 (2000).
6. Spiteller, P., Kern, W., Reiner, J., *et al. Biochim. Biophys. Acta* **1531**, 188-208 (2001).
7. Liu, Z., Minkler, P.E., and Sayre, L.M. *Chem. Res. Toxicol.* **16**, 901-911 (2003).
8. Zhang, W.-H., Liu, J., Xu, G., *et al. Chem. Res. Toxicol.* **16**, 512-523 (2003).

Related Products

4-hydroxy Nonenal - Item No. 32100 • 4-hydroxy Nonenal Mercapturic Acid- Item No. 32110 • 4-hydroxy Nonenal-d₃ - Item No. 332101

WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG 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 under separate cover to the MSDS supervisor at 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, 10/03/2011

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