

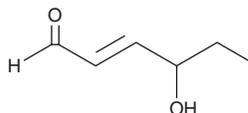
PRODUCT INFORMATION



4-hydroxy Hexenal

Item No. 32060

CAS Registry No.: 17427-21-3
Formal Name: (±)-4-hydroxy-2E-hexenal
Synonym: 4-HHE
MF: C₆H₁₀O₂
FW: 114.1
Purity: ≥98%
UV/Vis.: λ_{max}: 220 nm
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 Hexenal 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 hexenal in these solvents is approximately 50 mg/ml.

4-hydroxy Hexenal is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 4-hydroxy hexenal should be diluted with the aqueous buffer of choice. The solubility of 4-hydroxy hexenal in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

4-hydroxy Hexenal is a lipid peroxidation product derived from oxidized ω-3 fatty acids such as docosahexaenoic acid.^{1,2}

References

1. van Kuijk, F.J.G.M., Holte, L.L., and Dratz, E.A. 4-Hydroxyhexenal: A lipid peroxidation product derived from oxidized docosahexaenoic acid. *Biochim. Biophys. Acta* 1043, 116-118 (1990).
2. van Kuijk, F.J.G.M., Siakotos, A.N., Fong, L.G., *et al.* Quantitative measurement of 4-hydroxyalkenals in oxidized low-density lipoprotein by gas chromatography-mass spectrometry. *Anal. Biochem.* 224, 420-424 (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, 03/07/2024

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