

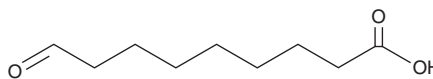
PRODUCT INFORMATION



9-Oxononanoic Acid

Item No. 29882

CAS Registry No.: 2553-17-5
Synonyms: Azelaaldehydic Acid, FA 9:1;O
MF: $C_9H_{16}O_3$
FW: 172.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

9-Oxononanoic acid is supplied as a solid. A stock solution may be made by dissolving the 9-oxononanoic acid in the solvent of choice, which should be purged with an inert gas. 9-Oxononanoic acid is slightly soluble in chloroform and methanol.

Description

9-Oxononanoic acid is an oxidized fatty acid formed via the autoxidation of linoleic acid (Item Nos. 90150 | 90150.1 | 21909).^{1,2} It increases phospholipase A₂ (PLA₂) activity and production of thromboxane B₂ (TXB₂; Item No. 19030) in isolated human plasma.¹ 9-Oxononanoic acid decreases hepatic *de novo* fatty acid synthesis and increases hepatic carnitine palmitoyltransferase activity, a marker of β -oxidation, in rats.²

References

1. Ren, R.F., Hashimoto, T., Mizuno, M., *et al.* A lipid peroxidation product 9-oxononanoic acid induces phospholipase A₂ activity and thromboxane A₂ production in human blood. *J. Clin. Biochem. Nutr.* **52(3)**, 228-233 (2013).
2. Minamoto, S., Kanazawa, K., Ashida, H., *et al.* Effect of orally administered 9-oxononanoic acid on lipogenesis in rat liver. *Biochim. Biophys. Acta* **958(2)**, 199-204 (1988).

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, 02/20/2024

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