

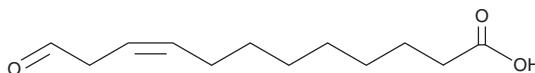
PRODUCT INFORMATION



12-oxo-9(Z)-Dodecenoic Acid

Item No. 37713

CAS Registry No.: 60485-39-4
Formal Name: (9Z)-12-oxo-9-dodecenoic acid
Synonyms: FA 12:2;O, 12-keto-9Z-Dodecenoic Acid,
12-oxo-9-cis-Dodecenoic Acid
MF: $C_{12}H_{20}O_3$
FW: 212.3
Purity: $\geq 95\%$
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥ 2 years



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

Description

12-oxo-9(Z)-Dodecenoic acid is an ω -oxoacid and a metabolite of the ω -6 PUFA linoleic acid (Item Nos. 90150 | 90150.1 | 21909) and the ω -3 PUFA α -linolenic acid (Item Nos. 90210 | 21910).¹ It is formed from these PUFAs via lipoxygenase-produced 13-hydroperoxy intermediates by hydroperoxide lyase in plants. 12-oxo-9(Z)-Dodecenoic acid (100 μM) is active against the plant pathogenic oomycete *P. parasitica*.²

References

1. Vick, B.A. and Zimmerman, D.C. Oxidative systems for modification of fatty acids: The lipoxygenase pathway. *The biochemistry of plants: A comprehensive treatise*. Stumpf, P.K., Elsevier (1987).
2. Prost, I., Dhondt, S., Rothe, G., et al. Evaluation of the antimicrobial activities of plant oxylipins supports their involvement in defense against pathogens. *Plant Physiol.* **139**, 1902-1913 (2005).

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, 01/10/2025

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