

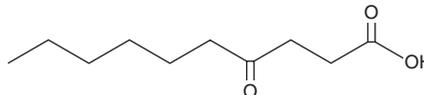
PRODUCT INFORMATION



4-oxo Decanoic Acid

Item No. 44815

CAS Registry No.: 4144-54-1
Synonym: FA 10:0;O, 4-ODA, 4-oxo Capric Acid
MF: $C_{10}H_{18}O_3$
FW: 186.3
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

4-oxo Decanoic acid (4-ODA) is supplied as a solid. A stock solution may be made by dissolving the 4-ODA in the solvent of choice, which should be purged with an inert gas. 4-ODA is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

4-ODA is a sex pheromone produced by virgin female Asian hornets (*V. velutina*).¹ It is produced at a 0.78 ratio of 4-oxo octanoic acid to 4-ODA. 4-ODA, alone or in the natural ratio with 4-oxo octanoic acid, elicits antennal responses from male Asian hornets in bioassays.

Reference

1. Wen, P., Cheng, Y.-N., Dong, S.-H., *et al.* The sex pheromone of a globally invasive honey bee predator, the Asian eusocial hornet, *Vespa velutina*. *Sci. Rep.* **7(1)**, 12956 (2017).

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, 11/14/2025

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