

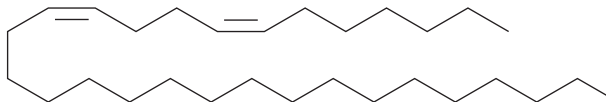
PRODUCT INFORMATION



7(Z),11(Z)-Nonacosadiene

Item No. 9000314

CAS Registry No.: 104410-91-5
Formal Name: (7Z,11Z)-nonacosadiene
MF: C₂₉H₅₆
FW: 404.8
Purity: ≥98%
Supplied as: A solution in hexane
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

7(Z),11(Z)-Nonacosadiene is supplied as a solution in hexane. To change the solvent, simply evaporate the hexane 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 7(Z),11(Z)-nonacosadiene in these solvents is approximately 20 mg/ml.

Description

Unsaturated cuticular hydrocarbons serve as pheromones in insects. In *D. melanogaster*, these hydrocarbons are sexually dimorphic in both their occurrence and their effects. 7(Z),11(Z)-Nonacosadiene is a cuticular pheromone in female fruit flies that stimulates male courtship behavior.¹ The biosynthesis of this C₂₉ diene appears to be mediated by an elongase with female-based expression.²

References

1. Ferveur, J.-F. and Sureau, G. Simultaneous influences on male courtship of stimulatory and inhibitory pheromones produced by live sex-mosaic *Drosophila melanogaster*. *Proc. Biol. Sci.* **263**(1373), 967-973 (1996).
2. Chertemps, T., Duportets, L., Labuer, C., et al. A female-biased expressed elongase involved in long-chain hydrocarbon biosynthesis and courtship behavior in *Drosophila melanogaster*. *Proc. Natl. Acad. Sci. USA* **104**(11), 4273-4278 (2007).

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, 02/15/2024

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