

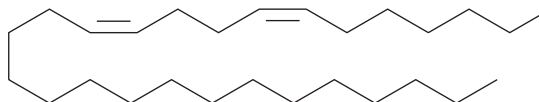
PRODUCT INFORMATION



7(Z),11(Z)-Heptacosadiene

Item No. 10012567

CAS Registry No.: 100462-58-6
Formal Name: heptacos-7Z,11Z-diene
MF: C₂₇H₅₂
FW: 376.7
Purity: ≥98%
Supplied as: A solution in ethanol
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)-Heptacosadiene 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 7(Z),11(Z)-heptacosadiene in these solvents is approximately 20 mg/ml.

Description

The mating and social behaviors of insects are largely orchestrated by a suite of volatile cuticular hydrocarbon pheromones. 7(Z),11(Z)-Heptacosadiene is the predominant female-specific courtship pheromone of the fruit fly *D. melanogaster*. At amounts above 100 ng, 7(Z),11(Z)-heptacosadiene elicits wing vibrations in male *D. melanogaster* flies in a dose-dependent manner.¹

Reference

1. Antony, C., Davis, T.L., Carlson, D.A., *et al.* Compared behavioral responses of male *Drosophila melanogaster* (Canton S) to natural and synthetic aphrodisiacs. *J. Chem. Ecol.* **11(12)**, 1617-1629 (1985).

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, 04/11/2024

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