

PRODUCT INFORMATION



(Z)-9-Hexadecenol

Item No. 30462

CAS Registry No.: 10378-01-5

Synonyms: FOH 16:1, (9Z)-Hexadecen-1-ol,
cis-9-Hexadecenol, Palmitoleyl
Alcohol, (Z)-9-Hexadecen-1-ol,
Z9-16:OH

MF: C₁₆H₃₂O

FW: 240.4

Purity: ≥98%

Supplied as: A liquid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

(Z)-9-Hexadecenol is supplied as a liquid. A stock solution may be made by dissolving the (Z)-9-hexadecenol in the solvent of choice, which should be purged with an inert gas. (Z)-9-Hexadecenol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of (Z)-9-hexadecenol in these solvents is approximately 30 mg/ml.

(Z)-9-Hexadecenol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (Z)-9-hexadecenol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. (Z)-9-Hexadecenol has a solubility of approximately 0.1 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(Z)-9-Hexadecenol is an unsaturated long chain fatty alcohol that has been found in pheromone gland extracts from *H. assulta* and has diverse biological activities.¹⁻³ It is virucidal against herpes simplex virus 2 (HSV-2) when used at a concentration of 10 mM.² (Z)-9-Hexadecenol is active against the bacterium *S. mutans* (MIC = 1.56 µg/ml).³ It decreases flight, upwind, close, landing, and copulatory behavior induced by an *H. assulta* pheromone blend in male *H. assulta* moths in a wind tunnel assay.¹

References

- Xu, M., Guo, H., Hou, C., *et al.* Olfactory perception and behavioral effects of sex pheromone gland components in *Helicoverpa armigera* and *Helicoverpa assulta*. *Sci. Rep.* **6**, 22998 (2016).
- Hilmarsson, H., Kristmundsdóttir, T., and Thormar, H. Virucidal activities of medium- and long-chain fatty alcohols, fatty acids and monoglycerides against herpes simplex virus types 1 and 2: Comparison at different pH levels. *APMIS* **113(1)**, 58-65 (2005).
- Hattori, M., Miyachi, K., Hada, S., *et al.* Effects of long-chain fatty acids and fatty alcohols on the growth of *Streptococcus mutans*. *Chem. Pharm. Bull. (Tokyo)* **35(8)**, 3507-3510 (1987).

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, 02/06/2026

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