

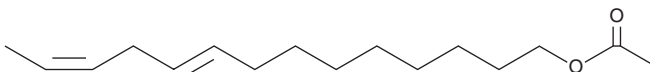
PRODUCT INFORMATION



9(E),12(Z)-Tetradecadienyl Acetate

Item No. 42895

CAS Registry No.: 31654-77-0
Formal Name: 9E,12Z-tetradecadien-1-ol 1-acetate
Synonym: 9E,12Z-TDDA, trans-9,cis-12-Tetradecadienyl Acetate
MF: $C_{16}H_{28}O_2$
FW: 252.4
Purity: $\geq 95\%$
Supplied as: A neat oil
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

9(E),12(Z)-Tetradecadienyl acetate (9E,12Z-TDDA) is supplied as a neat oil. A stock solution may be made by dissolving the 9E,12Z-TDDA in the solvent of choice, which should be purged with an inert gas. 9E,12Z-TDDA is soluble (≥ 10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

9(E),12(Z)-Tetradecadienyl acetate is a derivative of the insect pheromone 9(Z),12(E)-tetradecadienyl acetate.[74089] Formulations containing 9(E),12(Z)-tetradecadienyl acetate have been used to disrupt mating in insects.

References

1. Oyama, M. Mating suppression of *Spodoptera litura* (F.) by (Z, E)-9, 12-tetradecadienyl acetate. *Appl. Ent. Zool.* **12(3)**, 281-286 (1977).

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, 03/11/2025

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