

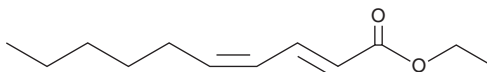
PRODUCT INFORMATION



2(E),4(Z)-Decadienoic Acid ethyl ester

Item No. 30835

CAS Registry No.: 3025-30-7
Formal Name: 2E,4E-decadienoic acid, ethyl ester
Synonyms: Pear ester, SFE 12:2
MF: C₁₂H₂₀O₂
FW: 196.3
Purity: ≥95%
UV/Vis.: λ_{max}: 264 nm
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

2(E),4(Z)-Decadienoic acid ethyl ester is supplied as a neat oil. A stock solution may be made by dissolving the 2(E),4(Z)-decadienoic acid ethyl ester in the solvent of choice, which should be purged with an inert gas. 2(E),4(Z)-Decadienoic acid ethyl ester is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml. 2(E),4(Z)-Decadienoic acid ethyl ester is slightly soluble in dimethyl formamide.

Description

2(E),4(Z)-Decadienoic acid ethyl ester is a volatile compound that has been found in pears and has kairomonal attractant activity.¹ 2(E),4(Z)-Decadienoic acid ethyl ester attracts male and female codling moths (*C. pomonella*), an apple, pear, and walnut pest, to traps placed in orchards when applied at a concentration of 40 mg/trap. It also attracts *C. splendana* males and females, as well as *P. fasciana* males.

Reference

- Schmidt, S., Anfora, G., Ioriatti, C., *et al.* Biological activity of ethyl (e,z)-2,4-decadienoate on different tortricid species: Electrophysiological responses and field tests. *Environ. Entomol.* **36(5)**, 1025-1031 (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, 03/04/2024

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