

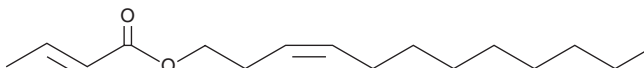
# PRODUCT INFORMATION



## Z3-Dodecenyl E2-butenate

Item No. 19787

CAS Registry No.: 104086-73-9  
Formal Name: 2E-butenic acid,  
3Z-dodecen-1-yl ester  
MF:  $C_{16}H_{28}O_2$   
FW: 252.4  
Purity:  $\geq 98\%$   
Supplied as: A solution in ethanol  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 2$  years



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

### Laboratory Procedures

Z3-Dodecenyl E2-butenate 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 Z3-dodecenyl E2-butenate in these solvents is approximately 30 mg/ml.

### Description

Z3-Dodecenyl E2-butenate is a female sex pheromone produced by the sweet potato weevil, *C. formicarius*. It has been used to bait and trap male sweet potato weevils to manage population counts in agriculture fields with reduced use of insecticides.<sup>1}</sup>

### Reference

1. Reddy, G.V.P., Gadi, N., and Taianao, A.J. Efficient sex pheromone trapping: Catching the sweetpotato weevil, *Cylas formicarius*. *J. Chem. Ecol.* **38**(7), 846-853 (2012).

#### 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, 04/02/2024

#### 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