

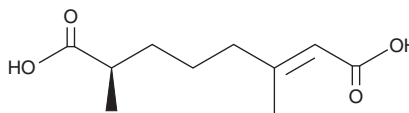
PRODUCT INFORMATION



Callosobruchusic Acid

Item No. 32927

CAS Registry No.: 87172-91-6
Formal Name: 3,7R-dimethyl-2E-octenedioic acid
Synonym: R-(-)-Callosobruchusic Acid
MF: C₁₀H₁₆O₄
FW: 200.2
Purity: ≥95%
Supplied as: A solid
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

Callosobruchusic acid is supplied as a solid. A stock solution may be made by dissolving the callosobruchusic acid in the solvent of choice, which should be purged with an inert gas. Callosobruchusic acid is soluble in methanol and DMSO.

Description

Callosobruchusic acid is an insect pheromone and monoterpene that has been found as a component of the copulation release pheromone erectin in azuki bean weevils (*C. chinensis*).¹ It induces copulation release activity in male azuki bean weevils (EC₅₀ = 6.5 ng/insect).

Reference

1. Mori, K., Ito, T., Tanaka, K., *et al.* Synthesis and biological activity of optically active forms of (E)-3, 7-dimethyl-2-octene-1, 8-dioic acid (callosobruchusic acid): A component of the copulation release pheromone (erectin) of the azuki bean weevil. *Tetrahedron* **39(13)**, 2303-2306 (1983).

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, 11/10/2022

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