

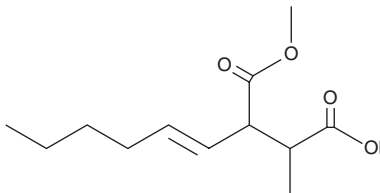
PRODUCT INFORMATION



Methyl Piliformate

Item No. 31757

CAS Registry No.: 898221-37-9
Formal Name: 2-(1E)-1-hexen-1-yl-3-methyl-butanedioic acid, 1-methyl ester
MF: C₁₂H₂₀O₄
FW: 228.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

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

Description

Methyl piliformate is a lichen metabolite that has been found in *Lecanora nipponica*.¹

Reference

1. Takenaka, Y., Hamada, N., and Takao, T. Two aliphatic acid derivatives from the cultured mycobionts of *Lecanora nipponica*. *Zeitschrift fuer Naturforschung, B: Chemical Sciences* **60(12)**, 1324-1326 (2005).

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, 08/11/2023

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