

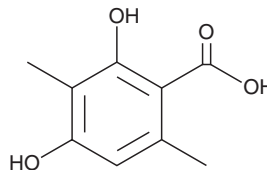
PRODUCT INFORMATION



3-methyl Orsellinic Acid

Item No. 25113

CAS Registry No.: 4707-46-4
Formal Name: 2,4-dihydroxy-3,6-dimethyl-benzoic acid
Synonyms: 2,4-Dihydroxy-3,6-dimethylbenzoic Acid,
 β -Orcinolcarboxylic Acid
MF: $C_9H_{10}O_4$
FW: 182.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*Aspergillus* sp.



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

Laboratory Procedures

3-methyl Orsellinic acid is supplied as a solid. A stock solution may be made by dissolving the 3-methyl orsellinic acid in the solvent of choice, which should be purged with an inert gas. 3-methyl Orsellinic acid is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

3-methyl Orsellinic acid is a fungal metabolite that has been found in *A. terreus*.¹ It has been used as a standard in dereplication of natural products.²

References

1. Takenaka, S., Ojima, N., and Seto, S. The isolation of 2,4-dihydroxy-3,6-dimethylbenzoic acid (3-methylorsellinic acid) from a culture of *Aspergillus terreus*. *J.C.S. Chem. Comm.* **0(7)**, 391-392 (1972).
2. Nielsen, K.F., Månsson, M., Rank, C., *et al.* Dereplication of microbial natural products by LC-DAD-TOFMS. *J. Nat. Prod.* **74(11)**, 2338-2348 (2011).

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, 08/23/2023

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