

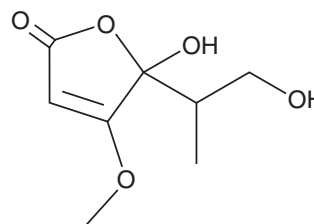
PRODUCT INFORMATION



Dihydro-6-hydroxypenicillic Acid

Item No. 44777

CAS Registry No.: 1661837-02-0
Formal Name: 5-hydroxy-5-(2-hydroxy-1-methylethyl)-4-methoxy-2(5H)-furanone
MF: $C_8H_{12}O_5$
FW: 188.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus//*Penicillium* sp.



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

Laboratory Procedures

Dihydro-6-hydroxypenicillic acid is supplied as a solid. A stock solution may be made by dissolving the dihydro-6-hydroxypenicillic acid in the solvent of choice, which should be purged with an inert gas. Dihydro-6-hydroxypenicillic acid is soluble in DMSO and methanol.

Description

Dihydro-6-hydroxypenicillic acid is a fungal metabolite that has been found in *Talaromyces* and has antiviral activity.¹ It reduces replication of influenza A in a reporter assay using 293T cells ($IC_{50} = 23.8 \mu\text{M}$).

Reference

1. Zhang, R., He, W., Wang, Y., et al. New butyrolactone derivatives from the endophytic Fungus *Talaromyces* sp. CPCC 400783 of *Reynoutria japonica* Houtt. *J. Antibiot. (Tokyo)* **74**(4), 225-232 (2021).

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, 11/25/2025

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