

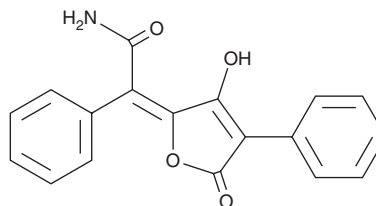
PRODUCT INFORMATION



Pulvinamide

Item No. 44779

CAS Registry No.: 31673-63-9
Formal Name: α -(3-hydroxy-5-oxo-4-phenyl-2(5H)-furanlydene)-benzeneacetamide
MF: $C_{18}H_{13}NO_4$
FW: 307.3
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Pseudocyphellaria* sp.



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

Laboratory Procedures

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

Description

Pulvinamide is a lichen alkaloid pigment that has been found in *P. crocata* and has antibacterial and antifungal activities.¹ It is active against *B. subtilis* and *C. albicans* (MICs = 50 and 200 $\mu\text{g}/\text{ml}$, respectively).

Reference

1. James, P.J.C., Vuong, D., Moggach, S.A., *et al.* Synthesis, characterization, and bioactivity of the lichen pigments pulvinamide, rhizocarpic acid, and epanorin and congeners. *J. Nat. Prod.* **86**(3), 550-556 (2023).

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, 11/24/2025

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