

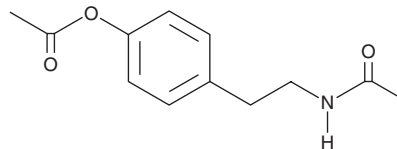
PRODUCT INFORMATION



N,O-Diacetyltyramine

Item No. 28738

CAS Registry No.: 14383-56-3
Formal Name: N-[2-[4-(acetyloxy)phenyl]ethyl]-acetamide
Synonyms: 4-(2-acetamidoethyl) Phenyl Acetate
MF: C₁₂H₁₅NO₃
FW: 221.3
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/Actinoplanes sp.



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

Laboratory Procedures

N,O-Diacetyltyramine is supplied as a solid. A stock solution may be made by dissolving the N,O-diacetyltyramine in the solvent of choice, which should be purged with an inert gas. N,O-Diacetyltyramine is soluble in ethanol, DMSO, methanol, and dichloromethane.

Description

N,O-Diacetyltyramine is a fungal metabolite that has been found in *P. endophytica* and has diverse biological activities.¹ It is active against a panel of 13 bacteria (MICs = 8-128 µg/ml), as well as a panel of nine fungi (MICs = 64-256 µg/ml). N,O-Diacetyltyramine is cytotoxic to MDA-MB-231, HeLa, MCF-7, and OAW42 cancer cells when used at concentrations ranging from 10 to 5,000 nM. It is larvicidal against *A. aegypti*, *A. albimanus*, and *C. quinquefasciatus* mosquitos (LC_{50s} = 724.8, 403.1, and 506.7 ppm, respectively).²

References

1. Mangamuri, U.K., Vijayalakshmi, M., Poda, S., *et al.* Bioactive metabolites produced by pseudonocardia endophytica VUK-10 from mangrove sediments: Isolation, chemical structure determination and bioactivity. *J. Microbiol. Biotechnol.* **25(5)**, 629-636 (2015).
2. Quevedo, R., Nunez-Dallos, N., and Quinones, M.L. Larvicidal activity of single and macrocyclic tyrosine derivatives against three important vector mosquitoes. *Res. Chem. Intermed.* **41(8)**, 5283-5292 (2014).

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, 12/05/2022

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