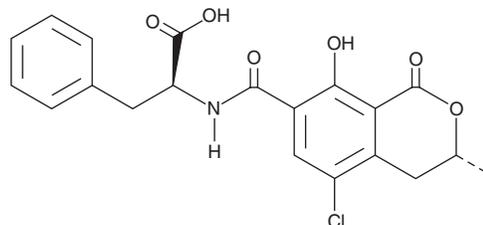


PRODUCT INFORMATION



Obscurolide A₁ Item No. 14439

CAS Registry No.: 144397-99-9
Formal Name: 4-[[tetrahydro-2-(3-hydroxy-1-butenyl)-5-oxo-3-furanyl]amino]-benzoic acid
MF: C₁₅H₁₇NO₅
FW: 291.3
Purity: ≥98%
Supplied as: A powder
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Obscurolide A₁ is supplied as a powder. A stock solution may be made by dissolving the obscurolide A₁ in the solvent of choice. Obscurolide A₁ is soluble in organic solvents such as methanol, DMSO, and acetone.

Description

Obscurolide A₁ is a natural butyrolactone isolated from *S. viridochromogenes*.¹ It inhibits calcium/calmodulin-dependent phosphodiesterase from bovine brain, presumably PDE1 (IC₅₀ = 15 mM).¹

Reference

1. Hoff, H., Drautz, H., Fiedler, H.-P., *et al.* Metabolic products of microorganisms. 261. Obscurolides, a novel class of phosphodiesterase inhibitors from streptomyces. I. Production, isolation, structural elucidation and biological activity of obscurolides A₁ to A₄. *J. Antibiot.* **45(7)**, 1096-1107 (1992).

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, 10/10/2022

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