

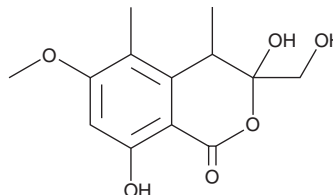
PRODUCT INFORMATION



Banksialactone A

Item No. 30000

CAS Registry No.: 1135775-06-2
Formal Name: 3,4-dihydro-3,8-dihydroxy-3-(hydroxymethyl)-6-methoxy-4,5-dimethyl-1H-2-benzopyran-1-one
MF: C₁₃H₁₆O₆
FW: 268.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years
Item Origin: Fungi/*Aspergillus banksianus*



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

Laboratory Procedures

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

Description

Banksialactone A is a fungal metabolite that has been found in *A. banksianus*.¹

Reference

1. Chaudhary, N.K., Pitt, J.I., Lacey, E., *et al.* Banksialactones and banksiamarins: Isochromanones and isocoumarins from an Australian fungus, *Aspergillus banksianus*. *J. Nat. Prod.* **81(7)**, 1517-1526 (2018).

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, 03/09/2020

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