

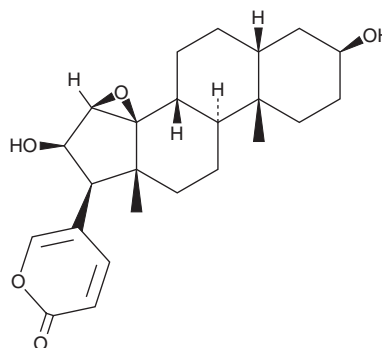
PRODUCT INFORMATION



Desacetylcinobufagin

Item No. 39822

CAS Registry No.: 4026-95-3
Formal Name: (5 β)-14,15 β -epoxy-3 β ,16 β -dihydroxy-
bufa-20,22-dienolide
Synonyms: Deacetylcinobufagin, NSC 234203
MF: C₂₄H₃₂O₅
FW: 400.5
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Bufonis Venenum*



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

Laboratory Procedures

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

Description

Desacetylcinobufagin is an active metabolite of the cardiotoxic steroid cinobufagin (Item No. 19844).¹ It is cytotoxic to HepG2, K562, and HL-60 cells (IC₅₀s = 51.2, 50.1, and 7.5 μ M, respectively).

Reference

1. He, X., Hu, H., Wu, Y., *et al.* Urinary metabolites of cinobufagin in rats and their antiproliferative activities. *Nat. Prod. Res.* **26(6)**, 489-499 (2012).

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, 11/21/2023

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