

PRODUCT INFORMATION



Bequinostatin A

Item No. 37772

CAS Registry No.: 151013-37-5
Formal Name: (-)-5,6,8,13-tetrahydro-1,6,7,9,11-pentahydroxy-8,13-dioxo-3-pentylbenzo[a]naphthacene-2-carboxylic acid

MF: $C_{28}H_{24}O_9$

FW: 504.5

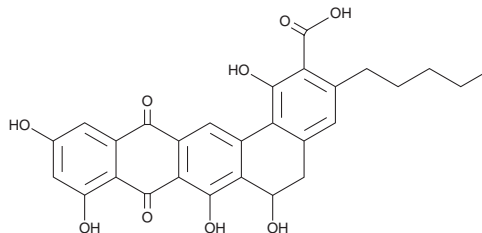
Purity: $\geq 75\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Bacterium/*Streptomyces* sp. MI384-DF12



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

Laboratory Procedures

Bequinostatin A is supplied as a solid. A stock solution may be made by dissolving the bequinostatin A in the solvent of choice, which should be purged with an inert gas. Bequinostatin A is soluble in organic solvents such as methanol, DMSO, and pyridine.

Description

Bequinostatin A is a naphthoquinone originally isolated from *Streptomyces* and a glutathione S-transferase pi 1 (GSTP1) inhibitor ($\text{IC}_{50} = 0.6 \mu\text{g/ml}$ for human GSTP1).¹

Reference

1. Aoyama, T., Kojima, F., Abe, F., *et al.* Bequinostatins A and B new inhibitors of glutathione S-transferase, produced by *Streptomyces* sp. MI384-DF12. Production, isolation, structure determination and biological activities. *J. Antibiot. (Tokyo)* **46(6)**, 914-920 (1993).

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, 02/14/2023

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