

PRODUCT INFORMATION



Kibdelone A

Item No. 21524

CAS Registry No.: 934464-77-4
Formal Name: (10S,12S,13R)-4-chloro-10,11,12,13-tetrahydro-10,12,13,15-tetrahydroxy-8-methoxy-2-methyl-3-propyl-2H-[1]benzopyrano[2',3':6,7]naphth[2,1-g]isoquinoline-1,5,14,16-tetrone

MF: $C_{29}H_{24}ClNO_{10}$

FW: 582.0

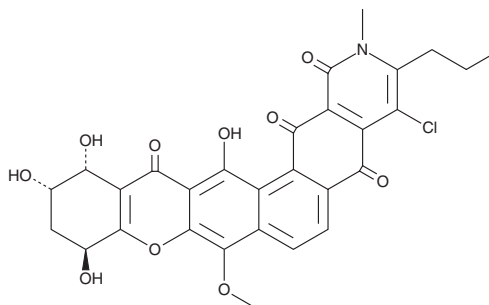
Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Bacterium/*Kibdelosporangium* sp.



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

Laboratory Procedures

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

Description

Kibdelone A is a member of a family of natural heterocyclic polyketides first isolated from a soil actinomycete, *Kibdelosporangium*.¹ Kibdelones have been described as having potent and selective cytotoxicity against a panel of human tumor cell lines.¹ They also display significant antibacterial and nematocidal activity.¹

Reference

1. Ratnayake, R., Lacey, E., Tennant, S., *et al.* Kibdelones: Novel anticancer polyketides from a rare Australian actinomycete. *Chemistry* **13**(5), 1610-1619 (2007).

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, 11/14/2022

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