

PRODUCT INFORMATION

Kibdelone B

Item No. 19444

CAS Registry No.: 934464-78-5
Formal Name: (10R,12R,13S)-rel-(+)-4-chloro-6,7,10,11,12,13-hexahydro-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: C₂₉H₂₆ClNO₁₀

FW: 584.0

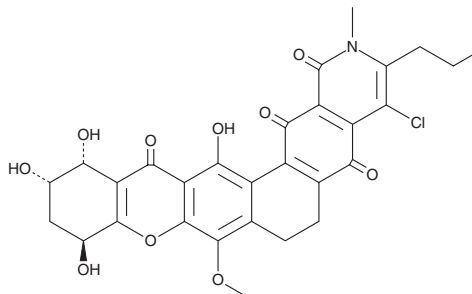
Purity: ≥95%

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 B is supplied as a solid. A stock solution may be made by dissolving the kibdelone B in the solvent of choice, which should be purged with an inert gas. Kibdelone B is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide. Kibdelone B is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Kibdelone B is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Kibdelones are natural aromatic polyketides first isolated from an actinomycete, *Kibdelosporangium*.¹ They exhibit potent and selective cytotoxicity against a panel of human tumor cell lines.¹ Kibdelones also display significant antibacterial and nematocidal activity.¹ Kibdelone B is a member of this group that undergoes facile equilibration to kibdelones A-C under mild conditions.¹ Its mode of action and pharmacology have not been studied.

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, 12/15/2022

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