

PRODUCT INFORMATION



Hodgkinsine

Item No. 31413

CAS Registry No.: 18210-71-4

Formal Name: (3aS,3'aR,3''aR,8aS,8'aR,8''aR)-2,2',2'',3,3',3'',8,8',8'',8a,8'a,8''a-dodecahydro-1,1',1''-trimethyl-,3a,3'a(1H,1'H):7',3''a(1''H)-terpyrrolo[2,3-b]indole

Synonym: (-)-Hodgkinsine

MF: C₃₃H₃₈N₆

FW: 518.7

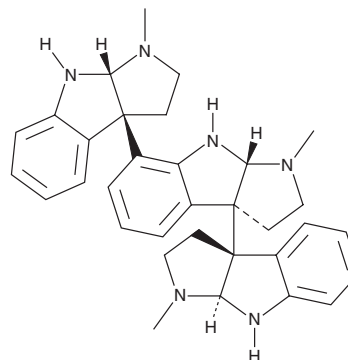
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Hodgkinsonia frutescens*



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

Laboratory Procedures

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

Description

Hodgkinsine is a pyrrolidinoindoline alkaloid that has been found in *P. colorata* and has analgesic activity.¹ It increases latency to paw licking in the hot plate test and latency to tail withdrawal in the tail-flick test in mice when administered at a dose of 5 mg/kg.

Reference

1. Amador, T.A., Verotta, L., Nunes, D.S., *et al.* Antinociceptive profile of hodgekinsine. *Planta Med.* **66**(8), 770-772 (2000).

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, 11/09/2022

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