

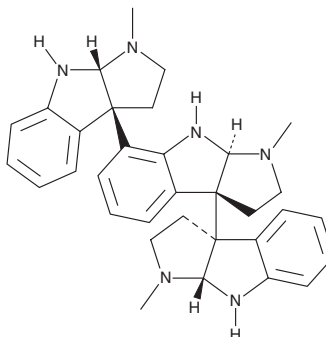
# PRODUCT INFORMATION



## Hodgkinsine B

Item No. 31421

**CAS Registry No.:** 586955-76-2  
**Formal Name:** 2,2',2'',3,3',3'',8,8',8'',8aR,8'aS,8''aR-dodecahydro-1,1',1''-trimethyl-3aR,3'aS(1H,1'H):7',3''aR(1''H)-terpyrrolo[2,3-b]indole  
**Synonym:** (-)-Hodgkinsine B  
**MF:** C<sub>33</sub>H<sub>38</sub>N<sub>6</sub>  
**FW:** 518.7  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

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

### Description

Hodgkinsine B is an alkaloid and isomer of hodgkinsine (Item No. 31413) that has analgesic activity.<sup>1</sup> It increases the latency to tail withdrawal in the tail-flick test in mice when administered at a dose of 10 mg/kg.

### Reference

1. Kodanko, J.J., Hiebert, S., Peterson, E.A., *et al.* Synthesis of all low-energy stereoisomers of the tris(pyrrolidinoindoline) alkaloid hodgkinsine and preliminary assessment of their antinociceptive activity. *J. Org. Chem.* **72(21)**, 7909-7914 (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, 11/09/2022

#### 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](http://WWW.CAYMANCHEM.COM)