

PRODUCT INFORMATION



Hodgkinsine trimethyl (iodide)

Item No. 31414

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

MF: C₃₆H₄₇N₆ • 3I

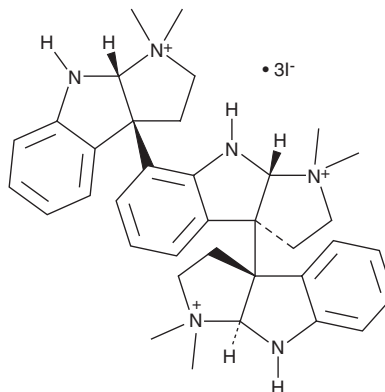
FW: 944.5

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

Description

Hodgkinsine trimethyl is a trimethylated form of hodgekinsine (Item No. 31413).

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