

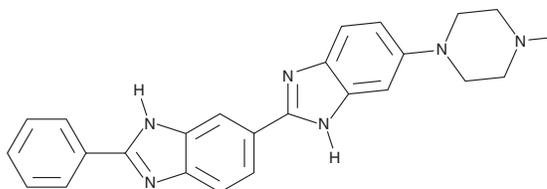
PRODUCT INFORMATION



HOE 33187

Item No. 40147

CAS Registry No.: 23623-08-7
Formal Name: 6-(4-methyl-1-piperazinyl)-2'-phenyl-2,6'-bibenzimidazole
MF: C₂₅H₂₄N₆
FW: 408.5
Purity: ≥98%
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

HOE 33187 is supplied as a solid. A stock solution may be made by dissolving the HOE 33187 in the solvent of choice, which should be purged with an inert gas. HOE 33187 is soluble in organic solvents such as methanol and DMSO. It is slightly soluble in acetonitrile.

HOE 33187 is slightly 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

HOE 33187 is a derivative of the fluorescent DNA probe Hoechst 33258 (Item No. 16756). It is cytotoxic to L1210 murine skin lymphocytic leukemia cells (IC₅₀ = 480 nM).¹

Reference

1. Findlay, G.J. and Baguley, B.C. Potentiation by phenylbisbenzimidazoles of cytotoxicity of anticancer drugs directed against topoisomerase II. *Eur. J. Cancer* **26(5)**, 586-589 (1990).

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, 02/07/2024

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