

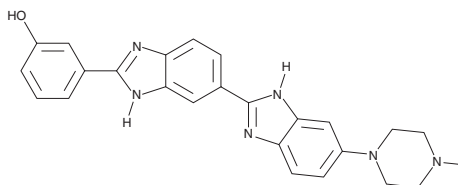
PRODUCT INFORMATION



HOE-S 785026

Item No. 40148

CAS Registry No.: 132869-83-1
Formal Name: 3-[5-(4-methyl-1-piperazinyl)[2,5'-bi-1H-benzimidazol]-2'-yl]-phenol
Synonyms: *m*-Hoechst 33258,
meta-Hoechst 33258,
m-hydroxy Hoechst 33258,
meta-hydroxy Hoechst 33258
MF: C₂₅H₂₄N₆O
FW: 424.5
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-S 785026 is supplied as a solid. A stock solution may be made by dissolving the HOE-S 785026 in the solvent of choice, which should be purged with an inert gas. HOE-S 785026 is soluble in acetonitrile and DMSO.

Description

HOE-S 785026 is a fluorescent DNA probe.¹ It is cytotoxic to L1210 murine skin lymphocytic leukemia cells (IC₅₀ = 670 nM).²

References

1. Breusegem, S.Y., Sadat-Ebrahimi, S.E., Douglas, K.T., *et al.* Increased stability and lifetime of the complex formed between DNA and meta-phenyl-substituted Hoechst dyes as studied by fluorescence titrations and stopped-flow kinetics. *J. Mol. Biol.* **308(40)**, 649-663 (2001).
2. Finlay, 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, 02/15/2024

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