

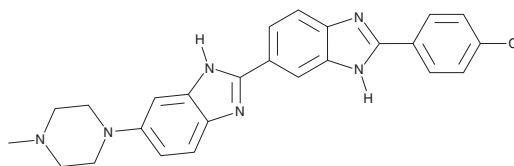
PRODUCT INFORMATION



HOE 32020

Item No. 39910

CAS Registry No.: 23554-99-6
Formal Name: 2'-(4-chlorophenyl)-5-(4-methyl-1-piperazinyl)-2,5'-bi-1H-benzimidazole
Synonym: Hoechst 32020
MF: C₂₅H₂₃ClN₆
FW: 442.9
Purity: ≥95%
Ex./Em. Max: 350/495 nm
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 32020 is supplied as a solid. A stock solution may be made by dissolving the HOE 32020 in the solvent of choice, which should be purged with an inert gas. HOE 32020 is soluble in acetonitrile, methanol, and DMSO.

Description

HOE 32020 is a fluorescent DNA probe.¹ HOE 32020 displays an emission maximum of 495 nm upon excitation at 350 nm when bound to DNA.

Reference

1. Latt, S.A. and Stetten, G. Spectral studies on 33258 Hoechst and related bisbenzimidazole dyes useful for fluorescent detection of deoxyribonucleic acid synthesis. *J. Histochem. Cytochem.* **24(1)**, 24-33 (1976).

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, 03/07/2024

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