

# Hoechst 33342 (ultra pure)

Cell permeable DNA dye

Hoechst 33342 is fluorescent probe useful for detecting DNA by fluorescence microscopy and flow cytometry. Hoechst 33342 may be used on live or fixed cells and is also applicable for cell cycle analysis and monitoring DNA condensation by flow cytometry. An additional ethyl group renders Hoechst 33342 more lipophilic, and thus more readily cell permeable than Hoechst 33258 . Exclusion of Hoechst 33342 dye is a characteristic common to stem cells, as well as chemotherapy-resistant cancer cells. Wavelength Maxima: Excitation 350nm, Emission 461nm

Citations: 36

[View Online »](#)

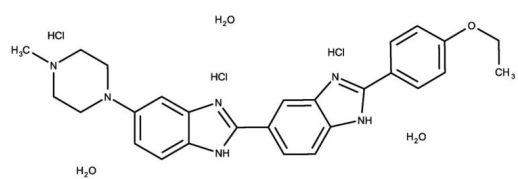
## Ordering Information

[Order Online »](#)

ENZ-52401	100mg
-----------	-------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended.
Handling	Protect from light. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	2'-(4-Ethoxyphenyl)-5-(4-methyl-1-piperazinyl)-2,5'-bi-1H-benzimidazole
Appearance	Yellow solid.
CAS	23491-52-3
Formula	$C_{27}H_{28}N_6O \cdot 3HCl$
MW	561.9
Purity	≥95% (HPLC)
Solubility	Soluble in water (50mg/ml) or DMSO (10mg/ml).

## Technical Info / Product Notes

This product is a member of the CELLESTIAL<sup>®</sup> product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL<sup>®</sup> reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.

### Application Note:

[Monitor Mitochondrial Membrane Potential in Cancer Cell Lines with a Dual-emission Fluorescent Dye](#)



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

Last modified: May 29, 2024