

# Hoechst 33258 (ultra pure)

## DNA dye

Hoechst 33258 is a fluorescent probe useful for detecting DNA by fluorescence microscopy and flow cytometry. Hoechst 33258 may be used on live or fixed cells and is also applicable for cell cycle analysis and monitoring DNA condensation by flow cytometry. Hoechst 33358 is less cell permeable than other derivatives, however this dye also allows for quantitative measurements when plotted in a standard emission-to-content curve. Wavelength Maxima: Excitation 352nm, Emission 461nm

Citations: 11

[View Online »](#)

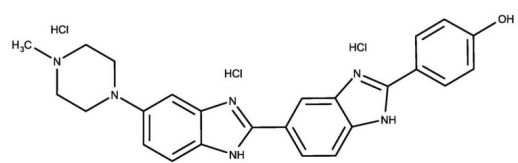
## Ordering Information

[Order Online »](#)

ENZ-52402	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	4-[5-(4-Methyl-1-piperazinyl)][2,5'-bi-1H-benzimidazol]-2'-yl]-phenol . trihydrochloride
Appearance	Yellow solid.
CAS	23491-45-4
Formula	$C_{25}H_{24}N_6O \cdot 3HCl$
MW	533.9 (trihydrochloride)
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.

## 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.

Last modified: May 29, 2024



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)