# Hoechst 33342 (ultra pure)

Hoechst 33342 is a cell-permeable DNA stain that produces bright blue fluorescence with low cytotoxicity - ideal for dynamic, highcontrast imaging in live-cell applications.

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 »

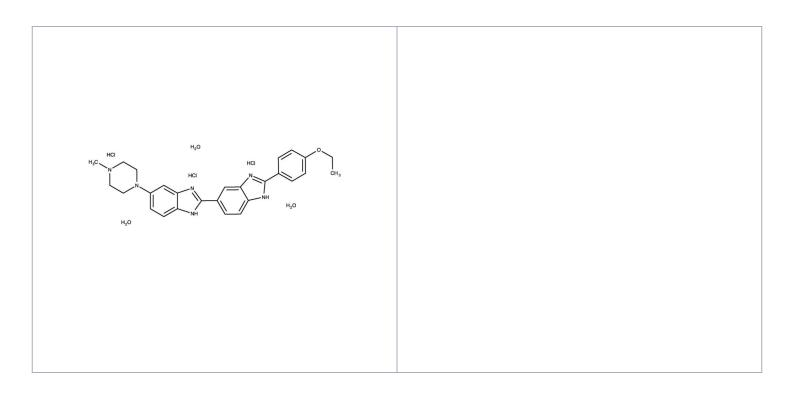
**Ordering Information** 

**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  $C_{27}H_{28}N_6O$  . 3HCl

**MW** 561.9

Purity ≥95% (HPLC)

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



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com