

DAPI (ultra pure)

DNA dye

DAPI is a fluorescent dye that binds nucleic acids and demonstrates a high affinity for DNA. It is used extensively in fluorescence microscopy, flow cytometry and microplate assays. DAPI is cell permeable and is readily used for detecting both live, dead and fixed cells. DAPI emits in the blue spectral range and is ideal for multiplexed assays inclusive of green-fluorescent molecules like fluorescein and green fluorescent protein (GFP), or red-fluorescent stains like Texas Red. Aside from labeling cell nuclei, DAPI is also used for the detection of mycoplasma or viral DNA in cell cultures. Wavelength Maxima: Excitation 358nm, Emission 461nm

Citations: 19

[View Online »](#)

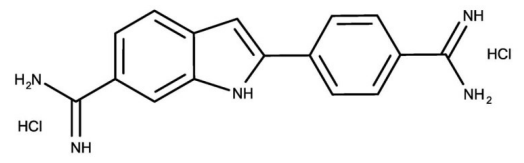
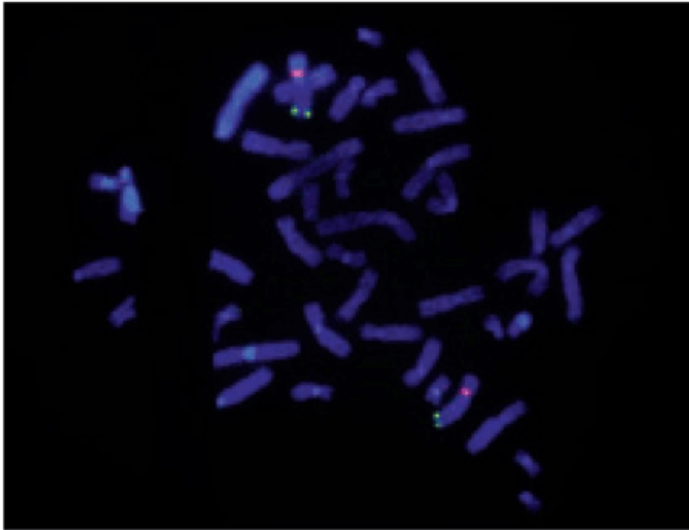
Ordering Information

[Order Online »](#)

ENZ-52404	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',6-Diamidino-2-phenylindole . dihydrochloride
CAS	28718-90-3
Formula	$C_{16}H_{17}Cl_2N_5$
MW	350.3
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.
Technical Info / Product Notes	<p>This product is a member of the CELLESTIAL[®] product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL[®] 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.</p>

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-
be@enzolifesciences.com](mailto:info-be@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)