DAPI (ultra pure)

DAPI is a fluorescent dye with strong DNAbinding affinity, commonly used in fluorescence microscopy, flow cytometry, and microplate assays for reliable nuclear visualization.

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

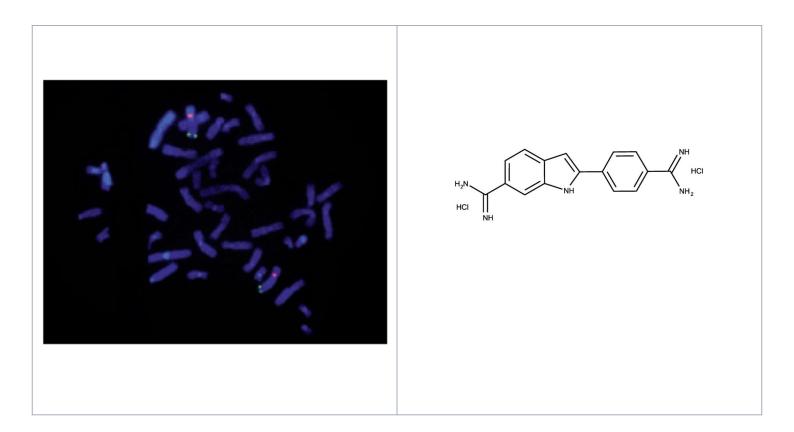
Order Online »

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 NotesThis 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.

Last modified: May 29, 2024

