

PureBlu™ Nuclear Staining Dyes



Bio-Rad's PureBlu Nuclear Staining Dyes are high-purity formulations of the traditional 4',6-diamidino-2-phenylindole dihydrochloride (DAPI) and Hoechst 33342 chemistries. PureBlu Dyes are packaged in an easy-to-use format and are the optimal choice to stain the nuclei of cells for cell imaging applications.

Key Benefits of PureBlu Dyes

- No weighing steps
- Single-step dilution upon resuspension
- High-purity formulation

Visit bio-rad.com/web/PureBlu1 for more information.

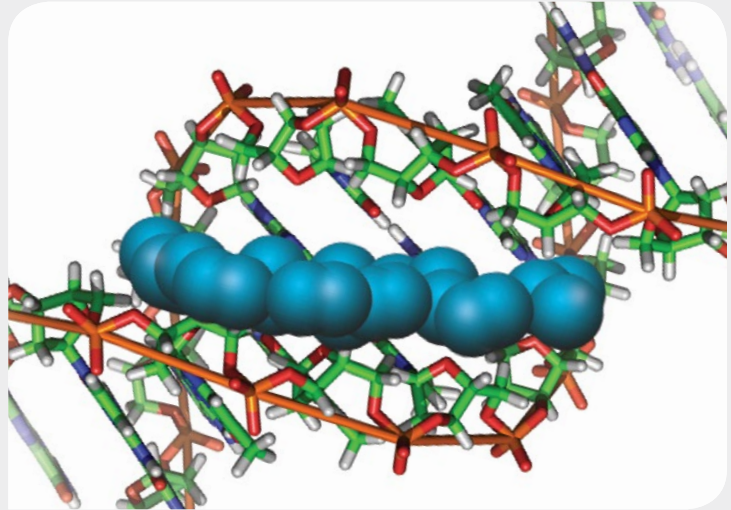
BIO-RAD

Easily Stain the Nuclei of Cells

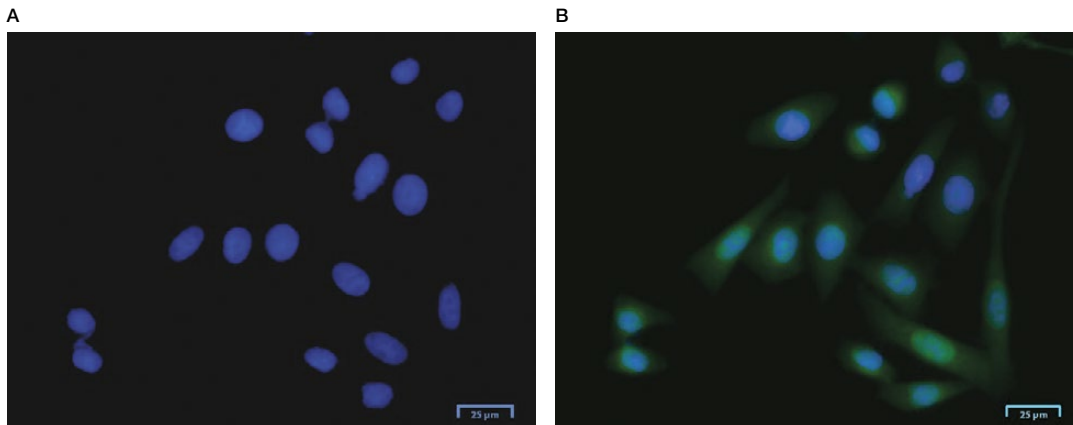
The format of PureBlu Nuclear Staining Dyes ensures that only the amount needed is reconstituted, thus reducing the amount of wasted reagents. Available as 95% pure compounds, PureBlu DAPI is suggested for fixed cells and PureBlu Hoechst 33342 for live cells. Both reagents are compatible with all multicolor cell imaging applications that include the use of an ultraviolet light source.

PureBlu Dye Chemistry

PureBlu Dyes bind with high affinity to the minor groove of AT-rich DNA sequences in eukaryotic and prokaryotic cells. A strong blue fluorescence is then released upon excitation in the blue range of the visible light spectrum, allowing the localization of the nuclei within the eukaryotic cells.



PureBlu Hoechst 33342 Nuclear Staining Dye binding to the minor groove of an AT-rich DNA sequence in a cell.



PureBlu Dyes are the optimal choice for staining the DNA of eukaryotic cells. HeLa cells were transfected with a green fluorescent protein (GFP)-expressing plasmid DNA, stained, and viewed in a six-well dish using a ZOE™ Fluorescent Cell Imager (catalog #145-0031). **A**, cells counterstained with PureBlu DAPI Nuclear Staining Dye (#135-1303); **B**, merged view of cells stained with PureBlu DAPI and GFP.

Ordering Information

Catalog Number	Available Dye	Excitation Maximum, nm	Emission Maximum, nm	Optimal Channel on ZOE™ Fluorescent Cell Imager
135-1303	PureBlu DAPI Nuclear Staining Dye	359	461	Blue
135-1304	PureBlu Hoechst 33342 Nuclear Staining Dye	350	461	Blue



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 3065 7550
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000