RedDot[™]1 Far-Red Nuclear Stain, 200X in Water

A far-red live cell nuclear stain similar to Draq5™.

Product Description

RedDotTM1 is a far-red cell membrane-permeant nuclear dye similar to Draq5TM. The dye is ideal for specifically staining the nuclei of live cells.

- Ideal for specifically staining the nuclei of live cells for short term imaging
- Highly thermostable and photostable, for convenient handling and demanding imaging applications
- Can be used for DNA content analysis by flow cytometry like Vybrant™ DyeCycle™ Ruby
- Can be used for normalization of In Cell Western® in fixed cells, similar to Drag5TM/Sapphire®700
- $\lambda_{Ex}/\lambda_{Em}$ = 662/694 nm (with DNA), for detection in the Cy®5 channel

RedDotTM1 can be excited at a wide range of wavelengths between 488 nm and 647 nm, including the 488 nm laser for flow cytometry, and can be used for DNA content analysis by flow cytometry. While DAPI and Hoechst dyes show a strong preference for A-T rich regions, RedDotTM1 is relatively insensitive to sequence base composition.

Choose the Right Dye for Your Application

Like Draq5®, RedDot[™]1 is toxic to cells within a few hours, and is recommended for short term staining experiments. For a low-toxicity far-red nuclear stain for live cells, see our NucSpot[™] 650 dye.

RedDotTM1 can be used to stain live bacteria (gram-positive and gram-negative). Staining in live yeast is weak and not nuclear. See our <u>Cellular Stains Table</u> for more information on how our dyes stain various organisms.

RedDotTM1 cannot be used for nuclear-specific staining in fixed cells. Please see RedDotTM2 (cat# 40061), a spectrally similar dye designed for specific nuclear staining of fixed and permeabilized cells or selective nuclear staining of dead cells.

Please note: far-red dyes like RedDotTM1 are not visible to the human eye, but must be imaged with a CCD camera or by confocal microscopy.

Product attributes

Probe cellular localization	Nucleus			
Assay type/options	DNA content/cell cycle profiling by flow cytometry, No-wash staining, Short term staining (<24h)			
Detection method/readout	Fluorescence microscopy, Flow cytometry, Near-IR imager			
Cell permeability	Membrane permeant			
Colors	Far-red			
Excitation/Emission	662 (broad)/694 nm (with DNA)			
Concentration	200X in water			
Storage Conditions	Store at -10 to -35 °C, Protect from light			

Call us: 800-304-5357 Email: btinfo@biotium.com

RedDot™1 and RedDot™2 Far-Red Nuclear Stains

Product	Catalog No.	Cell Permeability	Live/Fixed	Color (Ex/Em)	Applications
RedDot TM 1 Far-Red Nuclear Stain, 200X in Water	40060	Permeant	Live cells only	(Far-red) 662/694 nm	Live cell nuclear stain comparable to Draq5TM For microscopy, cell cycle analysis by flow cytometry Can be used for In-Cell Western normalization
RedDot™2 Far-Red Nuclear Stain, 200X in DMSO	40061	Impermeant	Fixed/permeabilized cells or tissues	(Far-red) 665/695 nm	Fixed cell nuclear stain Greater nuclear specificity than Draq7 TM Selectively stains dead cells in unfixed cultures

Cy Dye is a registered trademark of Cytiva. Draq5 is a trademark of BioStatus Ltd. In Cell Western and Sapphire are registered trademarks of LI-COR Inc. Vybrant and DyeCycle are trademarks of Thermo Fisher Scientific.

References

Download a list of curated RedDot, Not1 references.

This datasheet was generated on November 11, 2025 at 06:53:50 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/reddottm1-far-red-nuclear-stain-200x-in-h2o/