Glowing products for science

NucSpot® Far-Red, 1000X in DMSO

An improved alternative to 7-AAD showing less bleed-through fluorescence in the PE-Texas Red® channel. Ideal for selective detection of dead cells by flow cytometry.



Product Description

NucSpot® Far-Red is an improved alternative to the popular flow cytometry dead cell dye 7-AAD. It has red-shifted fluorescence emission compared to 7-AAD, for less bleed-through fluorescence in the PE-Texas Red® channel.

- Designed as an improved replacement for 7-AAD
- Less bleed into the PE-Texas Red® channel compared to 7-AAD
- Selective detection of dead cells by flow cytometry in the PE-Cy®5 or APC channel
- Perform cell cycle profiling by flow cytometry in fixed/permeabilized cells
- Substitute for 7-AAD in any standard protocol
- $\lambda_{Ex}/\lambda_{Em}$ (with DNA) = 597/667 nm

Like 7-AAD, NucSpot® Far-Red can be excited with the 488 nm laser for detection in the PE-Cy®5 channel, or excited with the 633 nm laser for detection in the APC channel. NucSpot® Far-Red has similar brightness to 7-AAD in the PE-Cy®5 channel, and is much brighter in the APC channel.

NucSpot® Far-Red is excluded by live cells, but stains necrotic and late apoptotic cells with compromised membrane integrity. Like 7-AAD, NucSpot® Far-Red also can be used for DNA content analysis of cell cycle by flow cytometry in fixed and permeabilized cells without the need for RNase-treatment, unlike propidium iodide.

Cy Dye is a registered trademark of Cytiva; Texas Red is a registered trademark of Thermo Fisher Scientific.

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

Product attributes

Probe cellular localization	Nucleus
Assay type/options	DNA content/cell cycle profiling (flow cytometry), Live/dead discrimination, No-wash staining
Detection method/readout	Flow cytometry
Cell permeability	Membrane impermeant
Apoptosis/viability marker	Dead cell stain
Colors	Far-red
Excitation/Emission	597/667 nm (with DNA)
Storage Conditions	Store at -10 to -35 °C

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/nucspot-far-red-1000x-in-dmso/