

NucSpot® 470 Nuclear Stain

Cell membrane-impermeant, nuclear-specific green counterstain.
Suitable for fixed cells or staining dead cells in live cultures.



Product Description

NucSpot® 470 Nuclear Stain is a green cell membrane-impermeant, nuclear-specific counterstain for the FITC channel.

- Nuclear-specific counterstain for fixed cells
- Selectively stains dead cells in live culture
- No wash required
- Suitable for cell cycle profiling by flow cytometry in fixed/permeabilized cells

Note: NucSpot® 470 has limited use for cell viability tracking due to its instability in culture medium during long incubation periods. [NucSpot® 500/515](#) was designed as an improved alternative for NucSpot® 470, offering better stability in culture medium.

Biotium also offers [NucSpot® Nuclear Stains](#) in a wide range of other colors from orange to near-IR. The stains also offer nuclear-specific counterstaining for fixed cells, as well as selective staining of dead cells in live cultures, while avoiding the issue of [photoconversion induced background by DAPI and Hoechst](#).

Learn more about our full selection of [nuclear stains](#) and our stains for [other organelles](#), [viability analysis](#), [membrane staining](#), and more.

Product attributes

Probe cellular localization	Nucleus
For live or fixed cells	For fixed cells
Assay type/options	DNA content/cell cycle profiling by flow cytometry, Live/dead discrimination, No-wash staining, Real-time imaging, Tissue staining
Detection method/readout	Fluorescence microscopy, Flow cytometry
Cell permeability	Membrane impermeant
Apoptosis/viability marker	Dead cell stain
Fixation options	Fix before staining (formaldehyde), Fix before staining (methanol), Permeabilize before staining
Colors	Green
Storage Conditions	Store at 2 to 8 °C