ViaFluor® Live Cell Microtubule Stains

Cell-permeant probes for imaging the microtubule cytoskeleton in live cells. Available with blue, green, or far-red fluorescence.



Product Description

ViaFluor® Live Cell Microtubule Stains are cell-permeant probes for imaging the microtubule cytoskeleton in live cells. They are simple, rapid and sensitive stains that can be imaged without a wash step.

- Cell-permeant fluorescent microtubule probes
- Rapid, sensitive, no-wash stains for live cells
- Available with blue, green, or far-red fluorescence
- ViaFluor® 647 dve compatible with SIM or STED

ViaFluor® Live Cell Microtubule Stains bind to polymerized tubulin and may be used to visualize microtubules in live cells for up to 24 hours in immortalized cell types. However, the probes may show cytotoxicity and cell cycle arrest with longer incubation times. The stains are supplied as a 1000X stock solutions in DMSO, and include a vial of 100 mM VRP, an efflux pump inhibitor that may improve probe retention and staining in some cell types.

ViaFluor® Live Cell Microtubule Stains cannot be fixed after staining, and cannot be used with fixed cells or tissues. The probes are not predicted to bind bacterial or yeast microtubules.

Learn more about our other <u>Fluorescent Cellular Stains</u> for nucleus, mitochondria, lysosomes, or see our selection of <u>Microbiology Stains & Kits</u>.

Product attributes

Dye	ViaFluor® 405, ViaFluor® 488, ViaFluor® 647			
Probe cellular localization	cytoskeleton, Microtubules			
Cell permeability	Membrane permeant			
For live or fixed cells	For live/intact cells			
Assay type/options	Long term staining (24-72h)			
Colors	Blue, Green, Far-red			
Concentration	1000X in DMSO			
Storage Conditions	Store at 2 to 8 °C, Protect dye component from light			

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

ViaFluor® Live Cell Microtubule Stains

Dye	Ex/Em	Catalog no.	Size (1000X in DMSO)*	Features
<u>ViaFluor® 405</u> 408/452 nm	408/452 nm	70064-T	10 uL	Blue fluorescence for the DAPI channel
		70064	50 uL	
ViaFluor® 488	500/515 nm	70062-T	10 uL	Green fluorescence for the FITC channel
		70062	50 uL	
ViaFluor® 647	650/675 nm	70063-T	10 uL	Far-red fluorescence for the Cy®5 channel
		70063	50 uL	Compatible with SIM or STED

 $^{^{\}star}$ Supplied with 100 mM VRP in DMSO for optional use.Cy5 is a registered trademark of Cytiva.

This datasheet was generated on November 4, 2025 at 07:10:36 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/viafluor-live-cell-microtubule-stains/