

MitoView™ Fix 640

Far-red, mitochondrial stain with excellent tolerance for fixation and permeabilization after staining.



Product Description

MitoView™ Fix 640 is far-red mitochondrial stain. The cell membrane-permeant dye rapidly accumulates in the mitochondria of live cells for no-wash, live cell staining. The dye has a reactive group that can covalently attach to proteins inside the mitochondria, so that staining is well-retained after fixation, permeabilization, and subsequent immunofluorescence staining.

- Stain mitochondria in live cells for subsequent fixation
- Well-retained after PFA or MeOH fixation, permeabilization, and IF staining
- Bright, photostable, and non-toxic for long-term live cell imaging
- Cells retain specific staining in culture for days
- Far-red fluorescence for the Cy@5 channel (Ex/Em 648/670 nm)
- Better specificity after fixation compared to MitoTracker® Deep-Red

MitoView™ Fix is not recommended for staining cells that are already fixed, because staining is non-specific in fixed cells. [MitoView™ Green](#) dye can be used for fixed cell staining, but for best results with fixed cells, we recommend using one of our [mitochondrial marker antibody conjugates](#).

We also offer original [MitoView™ Dyes](#) in a variety of colors, including membrane potential-dependent [MitoView™ 633](#), in addition to a selection of classic mitochondrial dyes. See our [Cellular Stains](#) for more stains for mitochondria, nucleus, cell surface, lysosomes, and other organelles.

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

Product attributes

Probe cellular localization	Mitochondria
Assay type/options	Long term staining (24-72h), No-wash staining
Fixation options	Fix after staining (formaldehyde), Fix after staining (methanol), Permeabilize after staining
Colors	Far-red
Excitation/Emission	648/670 nm
Storage Conditions	Store at 2 to 8 °C or below, Protect from light