



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Light-on LysoView™ 555

Light-On LysoView™ 555 is a red fluorogenic lysosome dye with pH-dependent fluorescence. Light-On LysoView™ dye is unique among commercially available lysosomotropic dyes in that its fluorescence in cells is activated by exposure to UV excitation.



Product Description

Light-On LysoView™ 555 is a red fluorogenic lysosome dye with pH-dependent fluorescence. Light-On LysoView™ dye is unique among commercially available lysosomotropic dyes in that its fluorescence in cells is activated by exposure to UV excitation. In solution, the dye shows pH-dependent fluorescence that does not require UV activation. In cells, the dye initially shows low fluorescence, but brief exposure to UV excitation from a mercury arc lamp through a DAPI filter activates bright red fluorescence localizing to lysosomes.

We hypothesize that the dye assumes a non-fluorescent structure that can be switched to a fluorescent structure by UV excitation. Mercury arc lamp excitation of the dye using other filter sets (FITC, rhodamine) or dye excitation with a 405 nm laser does not activate fluorescence. Lysosomal fluorescence fades several minutes after UV exposure, but can be re-activated in the same cells multiple times by exposure to UV light . Therefore the dye provides a novel tool for UV-activated, reversible fluorescence imaging of lysosomes. See the product protocol for more information.

Also see our red fluorescent [LysoView™ 540](#), and far-red fluorescent [LysoView™ 633](#) lysosome stains, which do not require UV activation.

Product attributes

Probe cellular localization	Lysosomes
For live or fixed cells	For live/intact cells
Assay type/options	Long term staining (24-72h), No-wash staining, Real-time imaging
Cell permeability	Membrane permeant
Colors	Red
Excitation/Emission	554/583 nm

This datasheet was generated on January 21, 2026 at 08:12:25 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/light-on-lysview-555/>