

## 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

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