



Glowing products for science

LysoView™ Dyes

LysoView™ fluorescent dyes label lysosomes in live cells without a wash step, and are available in multiple colors with options for super-resolution imaging.



Product Description

LysoView™ dyes are useful for staining lysosomes in live cells.

- No-wash, fluorescent staining of lysosomes in live cells
- Non-toxic for real-time live-cell imaging
- Lower background than LysoTracker® dyes
- Dye options for super-resolution imaging
- Color choices from blue to far-red, plus a unique NIR dye

LysoView™ dyes are fluorescent stains for imaging lysosome localization and morphology in live cells. The dyes accumulate in low pH organelles, resulting in highly specific staining without the need for a wash step. LysoView™ staining is non-toxic, and can imaged over the course of days, even after dye washout for some dye options (see the [Product Information Sheet](#) for more information).

LysoView™ 540 and LysoView™ 633 exhibit pH-sensitive fluorescence, further enhancing the specificity of staining. LysoView™ 488 has been validated in super-resolution imaging by SIM ([Ref. 7](#)), while LysoView™ 650 fluorophore is compatible with super-resolution imaging by SIM and STED.

Note: LysoView™ 540 has limited photostability, and may not be suitable for all microscopy applications. LysoView™ 550 is a bright and more photostable choice.

Also see our unique UV-activatable [Light-On LysoView™ 555](#), which is non-fluorescent until it is reversibly activated by UV light.

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

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	Blue, Green, Orange, Red, Far-red, Near-infrared
Storage Conditions	Store at -10 to -35 °C, Protect from light

LysoView™ Dyes

Product	Catalog number	Size	Abs/Em (nm) (pH ≤ 5)	Detection channel	Features
LysoView™ 405, 1000X in DMSO	70066-T	10 uL	318, 400/ 464	DAPI, Pacific Blue™	Blue fluorescent lysosome stain
70066	50 uL				
LysoView™ 488, 1000X in DMSO	70067-T	10 uL	496/526	GFP, FITC	Green fluorescent dye validated in SIM
70067	50 uL				
LysoView™ 540, 1000X in DMSO	70061-T	10 uL	540/634	TRITC, Cy®3, PE	Orange, pH-sensitive fluorescence**
70061	50 uL				
LysoView™ 550, 1000X in DMSO	70083-T	10 uL	542/567	TRITC, Cy®3, PE	Bright & photostable orange dye
70083	50 uL				
LysoView™ 594, 1000X in DMSO	70084-T	10 uL	585/634	Texas Red®	Bright & photostable red dye
70084	50 uL				
LysoView™ 633 (lyophilized solid)	70058-T	1 vial*	634/657	Cy®5, APC	Far-red, pH-sensitive fluorescence
70058	10 vials*				
LysoView™ 640, 1000X in DMSO	70085-T	10 uL	640/671	Cy®5, APC	Bright & photostable far-red dye
70085	50 uL				
LysoView™ 650, 1000X in DMSO	70059-T	10 uL	650/675	Cy®5, APC	Photostable far-red dye compatible with SIM, STED
70059	50 uL				
LysoView™ 680, 1000X in DMSO	70086-T	10 uL	674/711	Cy®5.5	Unique near-IR lysosome stain
70086	50 uL				

*Each vial of LysoView™ 633 makes 100 uL of 1000X dye after reconstitution.
** LysoView™ 540 has limited photostability and may not be suitable for all microcopy applications.
Cy Dye is a registered trademark of Cytiva; Pacific Blue is a trademark and Texas Red is a registered trademark of Thermo Fisher Scientific.
LysoTracker is a registered trademark of Thermo Fisher Scientific.

