

Product Information

LysoView™ Dyes

Product	Cat. No.	Unit Size	Ex/Em (nm)**	Detection Channel
LysoView™ 405, 1000X in DMSO	70066-T	10 uL	318, 400/464	DAPI, Pacific™
	70066	50 uL		Blue
LysoView™ 488, 1000X in DMSO	70067-T	10 uL	496/526	GFP, FITC
	70067	50 uL		
LysoView™ 540, 1000X in DMSO	70061-T	10 uL	540/634	Texas Red®
	70061	50 uL		
LysoView™ 550, 1000X in DMSO	70083-T	10 uL	542/567	TRITC, Cy®3, PE
	70083	50 uL		
LysoView™ 594, 1000X in DMSO	70084-T	10 uL	585/634	Texas Red®
	70084	50 uL		
LysoView™ 633 (lyophilized solid)	70058-T	1 vial*	634/657	Cy®5, APC
	70058	10 vials*		
LysoView™ 640, 1000X in DMSO	70085-T	10 uL	640/671	Cy®5, APC
	70085	50 uL		
LysoView™ 650, 1000X in DMSO	70059-T	10 uL	650/675	Cy®5, APC
	70059	50 uL		
LysoView™ 680, 1000X in DMSO	70086-T	10 uL	674/711	Cy®5.5
	70086	50 uL		

* One vial of LysoView ™ 633 yields 100 uL of 1000X dye stock solution after reconstitution (see below).

** Spectral properties at $pH \le 5$. See Figure 1.

Storage and Handling

Store at -20°C and protect from light. Product is stable for at least 12 months from date of receipt when stored as recommended. Before use, centrifuge vials briefly to collect all of the liquid at the bottom of the vial.

Reconstitution (LysoView™ 633 only)

To prepare LysoViewTM 633, 1000X stock solution in water: Briefly centrifuge vial to collect any loose material from cap before opening. Add 100 uL dH₂O to one vial of lyophilized dye and vortex to mix. Stock solution is stable for at least 2 weeks when stored at -20°C, protected from light.

Product Description

LysoView[™] Dyes are fluorescent stains for imaging lysosome localization and morphology in live cells. The dyes accumulate in the low pH environment of acidic organelles, resulting in highly specific, no-wash staining. LysoView[™] 540 and LysoView[™] 633 also exhibit pH-sensitive fluorescence (Figures 2-3). We offer a selection of LysoView ${}^{\rm T\!M}$ Dyes for flexibility in multi-color detection:

- LysoView[™] 405: Blue fluorescence for the DAPI or Pacific Blue[™] channel.
- LysoView™ 488: Green fluorescence for the FITC or GFP channel.
- LysoView[™] 540: Orange dye for the Cy®3 channel with pH-dependent fluorescence (Fig. 2). This specific dye has limited photostability and may not be suitable for all microscopy applications.
- LysoView[™] 550: Bright and photostable orange dye for the Cy®3 channel.
- LysoView™ 594: Bright and photostable red dye for the Texas Red® channel.
- LysoView™ 633: Far-red dye with pH-dependent fluorescence (Fig. 3).
- LysoView[™] 640: Bright and photostable far-red dye for the Cy®5 channel.
- LysoView™ 650: Photostable, far-red dye compatible with super-resolution imaging by SIM or STED.
- LysoView™ 680: Bright, photostable, unique near-infrared lysosome dye for the Cy®5.5 channel.

Also see Light-On LysoView $^{\rm TM}$ 555 (Related Products), a unique, UV-activated orange lysosome dye.

Considerations for Staining

- LysoView[™] Dyes can be used to stain adherent cells or cells in suspension.
- We recommend staining in complete medium with serum, but staining can also be done in serum-free medium or other buffer. If buffer is used to stain adherent cells, we recommend using a buffer like HBSS with calcium and magnesium to maintain cell attachment and morphology.
- We recommend 1X dye concentration as a starting point for optimization. Higher or lower concentrations may be optimal for different imaging systems or cell types.
- Do not add 1000X concentrated dye solution directly to cells in culture. This
 will cause transient high dye concentration leading to uneven or non-specific
 staining or toxicity. To add dye to cells without changing the medium, prepare
 an intermediate dilution of 10X dye in medium, then add 1/10 volume to the
 medium on the cells and pipette up and down gently to mix.
- In our tests, cells showed no obvious signs of toxicity after incubation with LysoView[™] Dyes for 3 days or longer, but toxicity may vary by cell type.
- We have observed LysoView[™] 550, LysoView[™] 594, and LysoView[™] 640 to be well retained in stained cells for 3 days after dye washout. Signal is less bright after washout compared to continuous dye incubation. Dye retention after washout may vary between cell types.
- LysoView[™] Dyes are recommended for live cell imaging only. Staining is not well retained after fixation with formaldehyde.

Staining Protocol

- For LysoView[™] 633 only, prepare 1000X stock solution in water as described under Reconstitution. Other LysoView[™] Dyes are supplied as 1000X stock solutions in DMSO.
- 2. Dilute 1000X LysoView[™] stock solution in cell culture medium to a final concentration of 1X.
- 3. Remove the medium from the cells and replace with medium containing dye.
- Incubate live cells with medium containing 1X LysoView[™] for 15-30 minutes, or longer, at 37°C.
- 5. Image or detect fluorescence using the appropriate excitation/emission settings or detection channel (see product table). No wash step is required.



Figure 1. Normalized absorption and emission of LysoView™ Dyes at pH ~5.



Figure 2. LysoView™ 540 normalized absorption (A) and emission (B) at varying pH.





Figure 3. LysoView™ 633 normalized absorption (A) and emission (B) at varying pH.

Related Products

Ca	at. No.	Product			
7	0060	Light-On LysoView™ 555			
7005	270075	MitoView™ Mitochondrial Dyes			
7	0076	Aquaphile™ JC-1			
7	0082	MitoView™ Fix 640			
7	0065	LipidSpot™ 488 Lipid Droplet Stain			
7	0069	LipidSpot [™] 610 Lipid Droplet Stain			
4	0081	NucSpot® Live 488 Nuclear Stain			
4	0082	NucSpot® Live 650 Nuclear Stain			
4008	341038	NucSpot® Nuclear Stains for dead or fixed cells			
4	0060	RedDot™1 Far-Red Nuclear Stain for live cells			
4	0061	RedDot™2 Far-Red Nuclear Stain for dead or fixed cells			
7	0064	ViaFluor® 405 Live Cell Microtubule Stain			
7	0062	ViaFluor® 488 Live Cell Microtubule Stain			
7	0063	ViaFluor® 647 Live Cell Microtubule Stain			
3002	20-30024	CellBrite® Cytoplasmic Membrane Dyes			
3007	030079	CellBrite® NIR Cytoplasmic Membrane Dyes			
3008	38-30090	CellBrite® Fix Membrane Stains			
3009	92-30104	MemBrite® Fix Cell Surface Staining Kits			
3010	05-30109	CellBrite® Steady Membrane Staining Kits			
3011	11-30113	ExoBrite™ EV Membrane Staining Kits			

Please visit our website at www.biotium.com for information on our life science research products, including fluorescent CF® Dye conjugates of transferrin, cholera toxin subunit B, dextrans, lectins, and Annexin V for cellular imaging, plus many more probes and kits for cell biology research.

Cy Dye is a registered trademark of Cytiva; Pacific Blue is a trademark and Texas Red is a registered trademark of Thermo Fisher Scientific.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.