www.biotium.com

Revised: August 13, 2021

Product Information

LysoView[™] Dyes

Catalog

70066-T

70066

no.

Size

10 ul

50 uL

Ex/Em (nm)**

318.400/464

Detection channel

DAPI, Pacific™

Blue

Product

LysoView[™] 405.

1000X in DMSO

-	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	TRITC, Cy®3, PE		
		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@	Taxaa Dad®		
		70084	50 uL		Texas Reu®		
	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 LysoView^{πi} 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[™] 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). Note: This 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 also can 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 using 1X dye 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[™] 680 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.
- 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.
- Image or detect fluorescence using the appropriate excitation/emission settings or detection channel (see product table). No wash step is required.



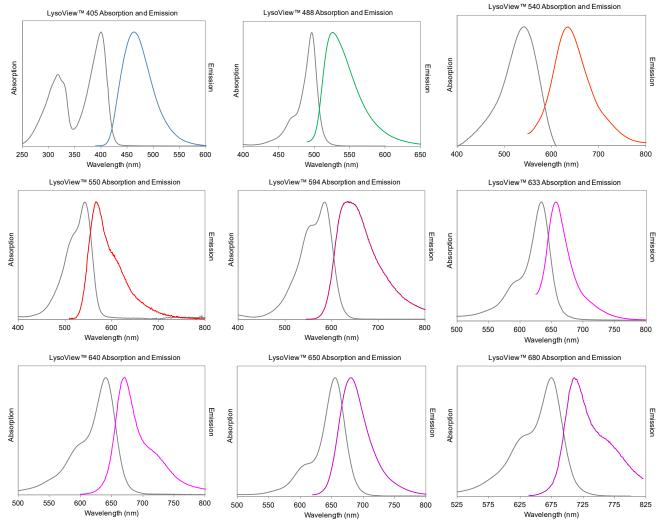


Figure 1. Normalized absorption and emission of LysoView[™] dyes at pH ~5.

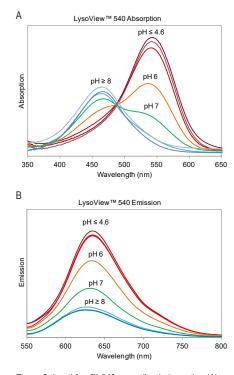
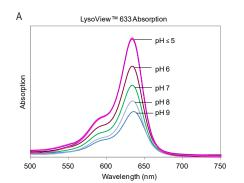


Figure 2. LysoView™ 540 normalized absorption (A) and emission (B) at varying pH.



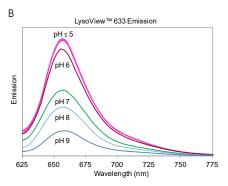


Figure 3. LysoView™ 633 normalized absorption (A) and emission (B) at varying pH.

Related Products

Catalog number	Product	
70060	Light-On LysoView™ 555	
7005270075	MitoView™ Mitochondrial Dyes	
70082	MitoView™ Fix 640	
70065	LipidSpot™ 488 Lipid Droplet Stain	
70069	LipidSpot™ 610 Lipid Droplet Stain	
40081	NucSpot® Live 488 Nuclear Stain	
40082	NucSpot® Live 650 Nuclear Stain	
40060	RedDot™1 far-red nuclear stain for live cells	
40061	RedDot™2 far-red nuclear stain for dead or fixed cells	
70064	ViaFluor® 405 Live Cell Microtubule Stain	
70062	ViaFluor® 488 Live Cell Microtubule Stain	
70063	ViaFluor® 647 Live Cell Microtubule Stain	
30020-30024	CellBrite® Cytoplasmic Membrane Dyes	
3007030079	CellBrite® NIR Cytoplasmic Membrane Dyes	
30088-30090	CellBrite® Fix Membrane Stains	
30092-30104	MemBrite® Fix Cell Surface Staining Kits	
30105-30109 CellBrite® Steady Membrane Staining Kits		
30111-30113	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.

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

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