## 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 attributes

Call us: 800-304-5357

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

Email: btinfo@biotium.com

## **Product Description**

LysoView<sup>TM</sup> 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<sup>TM</sup> 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<sup>™</sup> 550 is a bright and more photostable choice.

Also see our unique UV-activatable Light-On LysoViewTM 555, which is non-fluorescent until it is reversibly activated by UV 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
LysoView™ 488, 1000X in DMSQ	70066 70067-T	50 uL 10 uL	496/526	GFP, FITC	Green fluorescent dye validated in SIM
LysoView™ 540,	70067 70061-T	50 uL 10 uL	540/634	TRITC, Cy®3, PE	Orange, pH- sensitive fluorescence**
1000X in DMSO	70061	50 uL			
<u>LysoView™ 550,</u> 1000X in DMSO	70083-T 70083	10 uL 50 uL	542/567	TRITC, Cy®3, PE	Bright & photostable orange dye
LysoView™ 594, 1000X in DMSO	70084-T 70084	10 uL 50 uL	585/634	Texas Red®	Bright & photostable red dye
LysoView™ 633 (lyophilized solid)	70058-T 70058	1 vial* 10 vials*	634/657	Cy®5, APC	Far-red, pH-sensitive fluorescence
LysoView™ 640, 1000X in DMSO	70085-T	10 uL	640/671	Cy®5, APC	Bright & photostable far-red dye
LysoView™ 650, 1000X in DMSO	70085 70059-T	50 uL 10 uL	650/675	Cy®5, APC	Photostable far-red dye compatible with SIM, STED
1000X III DIVIOO	70059	50 uL			
LysoView™ 680, 1000X in DMSO	70086-T 70086	10 uL 50 uL	674/711	Cy®5.5	Unique near-IR lysosome stain

This datasheet was generated on November 1, 2025 at 06:00:13 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/lysoview-dyes/

<sup>\*</sup>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.