## ActinBrite™ High Affinity Phalloidin Conjugates

Novel phalloidin conjugates of fluorescent dyes designed to preserve strong F-actin binding for bright, specific staining that remains stable for over a month.

# Hothum

## **Product Description**

ActinBrite<sup>TM</sup> High Affinity Phalloidin Conjugates are novel phalloidin conjugates of fluorescent dyes that were designed to preserve high phalloidin affinity for F-actin, resulting in bright staining that can be imaged more than one month after staining with minimal loss of signal or specificity.

- Novel fluorescent phalloidin conjugates designed to preserve F-actin binding
- Allows stained samples to be stored for over a month or longer depending on the conjugate
- Available in 7 colors from green to near-IR for seamless multiplexing
- Serve as direct replacements for any phalloidin conjugate

### Reliable F-actin Staining, No Fading

Fluorescent phalloidins are popular tools for staining F-actin in fixed cells and tissues, but traditional dye conjugation can reduce their binding affinity—leading to weaker, less stable staining compared to other probes like labeled antibodies. As a result, phalloidin-stained samples typically lose signal within a week, especially when used with standard fluorescence mounting media.

ActinBrite<sup>TM</sup> High Affinity Phalloidin Conjugates solve this problem by preserving strong F-actin binding, delivering bright, reliable staining that lasts. With ActinBrite<sup>TM</sup>, samples can be imaged after for a month or more (depending on the conjugate and mounting method)—making delayed imaging easier and more dependable.

#### **Product attributes**

Call us: 800-304-5357

Probe cellular localization	Cytoskeleton, F-Actin			
For live or fixed cells	For fixed cells			
Assay type/options	Tissue staining			
Detection method/readout	Fluorescence microscopy			
Cell permeability	Membrane impermeant			
Fixation options	Fix before staining (formaldehyde), Permeabilize before staining			
Toxin	Phalloidin			
Colors	Green, Orange, Red, Far-red, Near-infrared			

Email: btinfo@biotium.com

Conjugation	Ex/Em	Excitation Laser Line	<b>Detection Channel</b>	Size	Catalog No.
ActinBrite™ 488/505 488/	488/505 nm	488 nm	FITC	50 U	<u>00095-T</u>
				300 U	<u>00095</u>
ActinBrite™ 530/555	532/558 nm	512 nm, 532 nm	Alexa Fluor® 532	50 U	<u>00096-T</u>
				300 U	00096
<u>ActinBrite™ 550/565</u> 547/565	547/565 nm	7/565 nm 555 nm, 561 nm	Cy®3, Rhodamine	50 U	<u>00097-T</u>
				300 U	<u>00097</u>
ActinBrite™ 610/630	608/629 nm	594 nm	Texas Red®	50 U	<u>00098-T</u>
				300 U	00098
<u>ActinBrite™ 645/665</u> 644/665 nr	644/665 nm	640 nm	Cy®5	50 U	<u>00101-T</u>
				300 U	<u>00101</u>
<u>ActinBrite™ 665/690</u> 669/685	669/685 nm	685 nm 640 nm	Alexa Fluor® 660	50 U	<u>00099-T</u>
				300 U	00099
<u>ActinBrite™ 750/770</u> 75	750/775 nm	730 nm	Alexa Fluor® 750	50 U	<u>00100-T</u>
				300 U	00100

Biotium offers a wide selection of traditional and novel stains for labeling cell structures, organelles, or monitoring viability. Learn more about our <u>cellular stains</u>, or view our <u>cellular stains</u> selection guides. Conjugates for Annexin, lectins, streptavidin and other biomolecules are also available with bright and photostable <u>CF® Dyes</u>.

CF is a registered trademark of Biotium, Inc. CY DYE is a registered trademark of Cytiva; Alexa Fluor and Texas Red are registered trademarks of Thermo Fisher Scientific.

This datasheet was generated on November 11, 2025 at 07:22:40 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/actinbrite-high-affinity-phalloidin-conjugates/">https://biotium.com/product/actinbrite-high-affinity-phalloidin-conjugates/</a>