



Glowing products for science

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

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.



Product attributes

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

Product Description

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

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

Biotium also offers CF® Dye conjugates of [Vitamin D-Binding Protein](#) (Vitamin D-BP, also known as GC Globulin) for visualizing monomeric G-actin in fixed and permeabilized cells. Biotium offers a wide selection of traditional and novel stains for labeling cell structures, organelles, or monitoring viability. Learn more about our [cellular stains](#), or view our [cellular stains selection guides](#). [Conjugates for Annexin, lectins, streptavidin and other biomolecules](#) are also available with bright and photostable [CF® Dyes](#).

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 January 14, 2026 at 01:35:46 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/actinbrite-high-affinity-phalloidin-conjugates/>