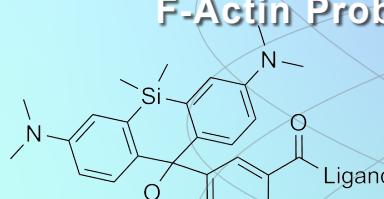
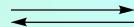
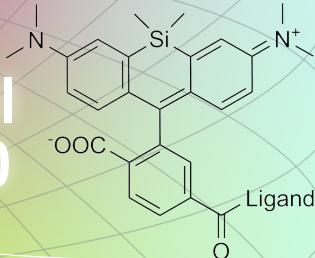


April  
2020



## F-Actin Probes in Living Cells

Related Publications

Research Tools

### Meetings

Cold Spring Harbor  
Conference - Systems  
Biology: Global Regulation of  
Gene Expression  
March 11-14th  
Cold Spring Harbor, NY  
Cytoskeleton Supported

Cold Spring Harbor  
Conference - Neuronal Circuits  
March 18-21st  
Cold Spring Harbor, NY  
Cytoskeleton Supported

Cold Spring Harbor  
Conference - from  
Neuroscience to Artificially  
Intelligent Systems  
March 24-28th  
Cold Spring Harbor, NY  
Cytoskeleton Supported

### Cytoskeleton Products

Actin Proteins  
Activation Assays  
Antibodies  
ECM Proteins  
ELISA Kits  
G-LISA® Kits  
Pull-down Assays  
Motor Proteins  
Small G-Proteins  
Tubulin & FtsZ Proteins

### Contact Us

P: 1 (303) 322.2254  
F: 1 (303) 322.2257  
E: cserve@cytoskeleton.com  
W: cytoskeleton.com

of STED and SiR/SPY-actin probes allows for unparalleled fluorescent visualization of subcellular F-actin structures and their physical characterization in living cells<sup>14-18</sup> (Figs. 1,2). SiR-actin probes have been used to examine F-actin in tissue<sup>19</sup> and a wide variety of cell types, including (but not limited to) human-induced pluripotent stem cell lines, cardiac cells, endothelial cells, epithelial cells, muscle cells, multiple cancer cell lines, and primary neurons<sup>14,16-18,20-22</sup> (Figs. 1,2).

### Fluorescent actin and fluorescent actin-binding domains

The first studies of live cell actin dynamics were performed with fluorescent derivatives of actin protein which were microinjected into cells<sup>50,51</sup>. This was a highly effective procedure but the apparatus took a while to setup. Thus overtime, transfections of GFP-actin conjugates became more popular. Fluorescently labeled actin protein or GFP/eGFP-actins worked well with fluorescence recovery after photobleaching (FRAP) microscopy<sup>11,23-25</sup>, which indicates the dynamic nature of actin cytoskeleton rearrangements. However, GFP/eGFP-actin has several drawbacks<sup>10</sup>. First, the size of GFP (~28 kDa) can impair polymerization<sup>26</sup> and GFP-actin can differentially label F-actin structures<sup>10,24</sup>. Second, some actin-binding proteins (e.g., formin family nucleators) might sterically hinder incorporation of GFP-actin into actin seeds or growing polymers<sup>27,28</sup>. Third, there is a relatively high background signal from non-filamentous fluorescent actin<sup>29</sup>. Finally, expression of eGFP-actin can affect cell behavior<sup>30,31</sup>.

Another method for actin live cell imaging utilizes yeast- or human-derived actin binding domains fused to GFP, eGFP, or m-Cherry fluorophores<sup>10,11,32</sup>. The most common genetically-encoded F-actin probes are Lifeact, utrophin (UtrCH), and F-tractin<sup>10</sup>. Lifeact is a 17 amino acid peptide from yeast Abp140<sup>33,34</sup> used for live cell imaging in mammalian and non-mammalian cells<sup>24,34-36</sup>. Lifeact has several disadvantages, including the possibility of affecting actin dynamics (so-called Lifeact-induced artifacts) and inhibiting the binding of actin-associated proteins such as cofolin<sup>32,37-40</sup>. Although Lifeact-GFP binds strongly to F-actin ( $K_d$ ,  $2.2 \pm 0.3 \mu\text{M}$ ), its binding affinity for G-actin is 10-fold higher<sup>33</sup>, resulting in high background fluorescence. Lifeact does not bind all actin-containing structures<sup>10,38</sup>. Lifeact is introduced into the cell through transfection rather than simply adding it into the medium as is done for the SiR/SiR700/SPY probes. UtrCH is based on the tandem calponin homology domains (CH1 and CH2) of utrophin<sup>41</sup> and consists of the first 261 amino acid residues of human utrophin, an actin binding protein<sup>42</sup>. The CH domains bind to actin with a  $K_d$  of  $\sim 18 \mu\text{M}$ <sup>43</sup>. Utrophin-based probes have been used successfully across a wide range of cell types and species<sup>10,32</sup>. Similar to Lifeact, at high concentrations,

## Continued from Page 1

utrophin-based probes can exert deleterious effects on actin cytoskeletal dynamics<sup>32,37,44</sup>. F-tractin is a 43 amino acid peptide derived from the rat actin-binding inositol 1,4,5-triphosphate 3-kinase A which binds F-actin with a Kd of ~10 μM<sup>45,46</sup>. Due to its larger size (in comparison to other probes), F-tractin might sterically hinder binding of actin-binding proteins that regulate and/or facilitate polymerization<sup>10</sup> and can modify actin-based cellular structures<sup>32</sup>.

### Actin-directed nanobodies and affimer proteins for F-actin

Two new technologies for monitoring actin dynamics in living cells are 1. single-domain antibodies, so-called nanobodies<sup>52</sup>, and 2. actin "affimers" - synthetic, actin-binding proteins isolated from phage library screens<sup>47-49</sup>. If not developed correctly, nanobodies can exhibit a high background signal due to G-actin binding<sup>10</sup>. Recently, three eGFP-fusion actin affimers were described with low micromolar binding affinities for F-actin<sup>47</sup>, but FRAP microscopy suggests that the eGFP-affimers may preferentially bind to a subset of actin filaments and alter actin organization in the cell<sup>47</sup>.

### Summary

Despite multiple options for visualizing F-actin-based structures in living cells, there is no perfect live cell probe (Table 1). Ideally, the best F-actin probe will be sensitive, selective, fluorogenic, produce a very low background signal, non-toxic, and easily introduced into a wide range of cells across multiple species. It is of paramount importance to confirm that changes in actin cytoskeleton dynamics/structural organization are physiologically relevant and not artifacts of the probe itself. To assist researchers in F-actin live cell imaging studies, Cytoskeleton offers the SiR and SPY actin live cell imaging probes.

## References

- Pollard T.D. 2016. What we know and do not know about actin. In: Jockusch B. (eds.) *The Actin Cytoskeleton. Handbook of Experimental Pharmacology*. vol 235. Springer, Cham.
- Dominguez R. and Holmes K.C. 2011. Actin structure and function. *Annu. Rev. Biophys.* **40**, 169-186.
- Verstraelen P. et al. 2018. Image-based profiling of synaptic connectivity in primary neuronal cell culture. *Front. Neurosci.* **12**, 389.
- Hoover B.R. et al. 2010. Tau mislocalization to dendritic spines mediates synaptic dysfunction independently of neurodegeneration. *Neuron*. **68**, 1067-1081.
- Spires-Jones T.L. and Hyman B.T. 2014. The intersection of amyloid beta and tau at synapses in Alzheimer's disease. *Neuron*. **82**, 756-771.
- Dugina V.B. et al. 2019. Biological role of actin isoforms in mammalian cells. *Biochemistry (Mosc)*. **84**, 583-592.
- Hartmann S. et al. 2015. The function of Rho-associated kinases ROCK1 and ROCK2 in the pathogenesis of cardiovascular disease. *Front. Pharmacol.* **6**, 276.
- Bonello T.T. et al. 2009. New approaches to targeting the actin cytoskeleton for chemotherapy. *Future Med. Chem.* **1**, 1311-1331.
- Bryce N.S. et al. 2019. Chemical biology approaches targeting the actin cytoskeleton through phenotypic screening. *Curr. Opin. Chem. Biol.* **51**, 40-47.
- Melak M. et al. 2017. Actin visualization at a glance. *J. Cell Sci.* **130**, 525-530.
- Belin B.J. et al. 2014. Comparative analysis of tools for live cell imaging of actin network architecture. *Bioarchitecture*. **4**, 189-202.
- Lukinavicius G. et al. 2013. A near-infrared fluorophore for live-cell super-resolution microscopy of cellular proteins. *Nat. Chem.* **15**, 132-139.
- Lukinavicius G. et al. 2014. Fluorogenic probes for live-cell imaging of the cytoskeleton. *Nat. Methods*. **11**, 731-733.
- D'Este E. et al. 2015. STED nanoscopy reveals the ubiquity of subcortical cytoskeleton periodicity in living neurons. *Cell Rep.* **10**, 1246-1251.
- Hell S.W. and Wichmann J. 1994. Breaking the diffraction resolution limit by stimulated emission: stimulated-emission-depletion fluorescence microscopy. *Opt. Lett.* **19**, 780-782.
- Bar J. et al. 2016. Periodic F-actin structures shape the neck of dendritic spines. *Sci. Rep.* **6**, 37136.
- Liete S.C. et al. 2016. The actin-binding protein α-adducin is required for maintaining axon diameter. *Cell Rep.* **15**, 490-498.
- Wang L. et al. 2019. Small-molecule fluorescent probes for live-cell super-resolution microscopy. *J. Am. Chem. Soc.* **141**, 2770-2781.
- Klementieva N.V. et al. 2016. Fluorescence imaging of actin fine structure in tumor tissues using SiR-actin staining. *Anticancer Res.* **36**, 5287-5294.
- Kim T.H. et al. 2016. Cancer cells become less deformable and more invasive with activation of β-adrenergic signaling. *J. Cell Sci.* **129**, 4563-4575.
- Hashimoto N. et al. 2017. Reversible differentiation of immortalized human bladder smooth muscle cells accompanied by actin bundle reorganization. *PLoS ONE*. **12**, e0186584.
- Narva E. et al. 2017. A strong contractile actin fence and large adhesions direct human pluripotent colony morphology and adhesion. *Stem Cell Reports*. **9**, 67-76.
- Choidas A. et al. 1998. The suitability and application of a GFP-actin fusion protein for long-term imaging of the organization and dynamics of the cytoskeleton in mammalian cells. *Eur. J. Cell Biol.* **77**, 81-90.
- Lemieux M.G. et al. 2014. Visualization of the actin cytoskeleton: different F-actin-binding probes tell different stories. *Cytoskeleton*. **71**, 157-169.
- Small J. et al. 1999. Visualising the actin cytoskeleton. *Microsc. Res. Tech.* **47**, 3-17.
- Sliogeryte K. et al. 2016. Differential effects of LifeAct-GFP and actin-GFP on cell mechanics assessed using micropipette aspiration. *J. Biomech.* **49**, 310-317.
- Wu J.-Q. and Pollard T.D. 2005. Counting cytokinesis proteins globally and locally in fission yeast. *Science*. **310**, 310-314.
- Chen Q. et al. 2012. Formins filter modified actin subunits during processive elongation. *J. Struct. Biol.* **177**, 32-39.
- Lee C.W. et al. 2013. Dynamic localization of G-actin during membrane protrusion in neuronal motility. *Curr. Biol.* **23**, 1046-1056.
- Aizawa H. et al. 1997. A green fluorescent protein-actin fusion protein dominantly inhibits cytokinesis, cell spreading, and locomotion in Dictyostelium. *Cell Struct. Funct.* **22**, 335-345.
- Nagasaki A. et al. 2017. The position of the GFP tag on actin affects the filament formation in mammalian cells. *Cell Struct. Funct.* **42**, 131-140.
- Patel S. et al. 2017. Functional characterisation of filamentous actin probe expression in neuronal cells. *PLoS ONE*. **12**, e0187979.
- Riedl J. et al. 2008. Lifeact: a versatile marker to visualize F-actin. *Nat. Methods*. **5**, 605-607.
- Riedl J. et al. 2010. Lifeact mice for studying F-actin dynamics. *Nat. Methods*. **7**, 168-169.
- Schachtner H. et al. 2012. Tissue inducible Lifeact expression allows visualization of actin dynamics in vivo and ex vivo. *Eur. J. Cell Biol.* **91**, 923-929.
- Mizoguchi T. et al. 2016. Zebrafish lines expressing UAS-driven red probes for monitoring cytoskeletal dynamics. *Genesis*. **54**, 483-489.
- Du J. et al. 2015. Lifeact and Utr230 induce distinct actin assemblies in cell nuclei. *Cytoskeleton*. **72**, 570-575.
- Munsie L.N. et al. 2009. Lifeact cannot visualize some forms of stress-induced twisted F-actin. *Nat. Methods*. **6**, 317.
- Courtemanche N. et al. 2016. Avoiding artefacts when counting polymerized actin in live cells with LifeAct fused to fluorescent proteins. *Nat. Cell Biol.* **18**, 676-683.
- Flores L.R. et al. 2019. Lifeact-TagGFP2 alters F-actin organization, cellular morphology and biophysical behaviour. *Sci. Rep.* **9**, 3241.
- Burkel B.M. et al. 2007. Versatile fluorescent probes for actin filaments based on the actin-binding domain of utrophin. *Cell Motil. Cytoskelet.* **64**, 822-832.
- Winder S.J. et al. 1995. Calmodulin regulation of utrophin actin binding. *Biochem. Soc. Trans.* **23**, 397S.
- Rybalkova I.N. and Ervasti J.M. 2005. Identification of spectrin-like repeats required for high affinity utrophin-actin interaction. *J. Biol. Chem.* **280**, 23018-23023.
- Spracklen A.J. et al. 2014. The pros and cons of common actin labeling tools for visualizing actin dynamics during Drosophila oogenesis. *Dev. Biol.* **393**, 209-226.
- Schell M.J. et al. 2001. Inositol 1,4,5-triphosphate 3-kinase A associates with F-actin and dendritic spines via its N-terminus. *J. Biol. Chem.* **276**, 37537-37546.
- Brehm M.A. et al. 2004. Identification of the actin-binding domain of Ins(1, 4, 5)P3 3-kinase isoform B (IP3K-B). *Biochem. J.* **382**, 353-362.
- Lopata A. et al. 2018. Affimer proteins for F-actin: novel affinity reagents that label F-actin in live and fixed cells. *Sci. Rep.* **8**, 6572.
- Tiede C. et al. 2014. Adhiron: a stable and versatile peptide display scaffold for molecular recognition applications. *Protein Eng. Des. Sel.* **27**, 145-155.
- Tiede C. et al. 2017. Affimer proteins are versatile and renewable affinity reagents. *Elife* **6**, 24903.
- Shimada Y. et al. 1997. Dynamics of actin in cardiac myofibrils and fibroblast stress fibers. *Cell Structure and Function*, **22**, 59-64.
- Machesky L.M. and Hall A. 1997. Role of actin polymerization and adhesion to extracellular matrix in Rac- and Rho-induced cytoskeletal reorganization. *J. Cell Biol.* **138**, 913-926.
- Rocchetti A. et al. 2014. Fluorescent labelling of the actin cytoskeleton in plants using a camelid antibody. *Plant Methods* **10**, 12. 10.1186/1746-4811-10-12.

## Figures and Table on Page 3

# Figures and Table Appendix

## Figures and Table Referenced in Text

Table 1. Comparision Table of F-actin probes

Table 1. Actin-binding probes for live cell imaging.

	SiR/SPY-actins	Lifeact GFP tagged	Actin fluorescently labeled	Actin GFP tagged	Utrophin GFP tagged	F-Tractin GFP tagged	Nanobody GFP tagged	Affimers GFP tagged
<b>Source</b>	<i>Jaspis johnstoni</i> SiR or SPY probes with desbromodes-methyl-jasplakinolide	<i>Saccharomyces cerevisiae</i> amino-acids 1-17 of ABP140	Skeletal muscle (rabbit) or non-muscle (beta actin human platelet)	Beta-actin fusion protein with GFP	<i>Homo sapiens</i> Amino acids 1-261 of utrophin	<i>Rattus norvegicus</i> Amino acids 10-52 of ITPKA	<i>Vicugna pacos</i> anti-actin-nanobody	Synthetic actin-binding probes isolated from phage library screens
<b>Applications</b>	Live cell imaging of endogenous F-actin including wide field confocal Ref. 14-18 Super-resolution microscopy (e.g., STED, SIM) Ref. 13-18	Live cell imaging Ref. 33,34	FRAP Ref. 50 Live cell imaging Ref. 51	FRAP Ref. 23 Live cell imaging of exogenous and endogenous actin Ref. 24,25	Live cell imaging of endogenous F-actin Ref. 41.	Live cell imaging of endogenous F-actin Ref. 45,46	Live cell imaging of endogenous actin Ref. 52	FRAP Ref. 47-49 Live cell imaging of endogenous actin Ref. 47-49
<b>Advantages</b>	Direct application to cells and tissues Fluorogenic Cell Permeability (No transfection required) Photostability Very Low Background Super-resolution Compatibility (STED, SIM) No Cytotoxicity Multiple Colors (e.g., far-red, red and orange) Binds only F-actin Small organic molecule (more stable)	Multiple Colors (e.g., far-red, red, orange, yellow, green)	Very similar conformation to endogenous actin Small fluorophore size.	Labeled actin is incorporated into endogenous filaments Multiple Colors (e.g., far-red, red, orange, yellow, green)	Does not bind actin monomers (G-actin)	Does not bind actin monomers (G-actin)	Small probe size Low probability of affecting actin dynamics	High nanomolar affinity for F-actin Multiple Colors (e.g., far-red, red, orange, yellow, green) F-actin specific with correct screening protocol.
<b>Disadvantages</b>	Possible effects on actin dynamics at high concentrations Binds G-actin to produce a high background signaling Possible effects on actin dynamics Requires transfection	Large fluorescent reporter GFP Binds G-actin to produce a high background signaling Possible effects on actin dynamics Requires transfection	Requires injection	Large fluorescent reporter GFP Exogenous actin expression Binds G-actin to produce a high background signal Requires transfection	Large size Large fluorescent reporter GFP Possible effects on actin dynamics Requires transfection	Large fluorescent reporter GFP Possible effects on actin dynamics Requires transfection	Large fluorescent reporter GFP Binds G-actin to produce a high background signal Requires transfection	Large fluorescent reporter GFP Requires transfection
<b>Overall rating</b>	+++++	+++	+++	++	+++	+++	+++	+++

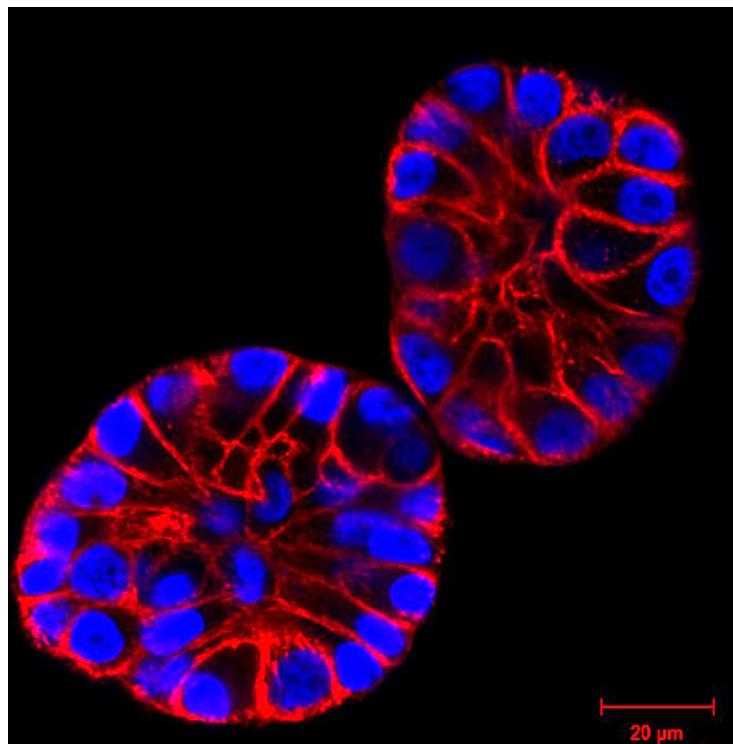


Figure 1. MCF10A cells expressing H2B-GFP (blue) in Matrigel (3D culture) stained with SiR-actin (red). Image taken on an inverted LSM microscope. Courtesy of Christian Conrad and Katharina Jechow, Heidelberg.

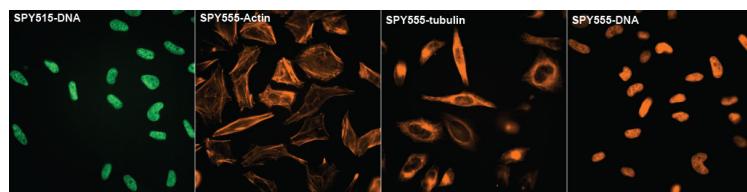


Figure 2. SPY505 and SPY555 staining DNA,Actin,Tubulin, and DNA . Photo comes from front page of Spirochrome's website.

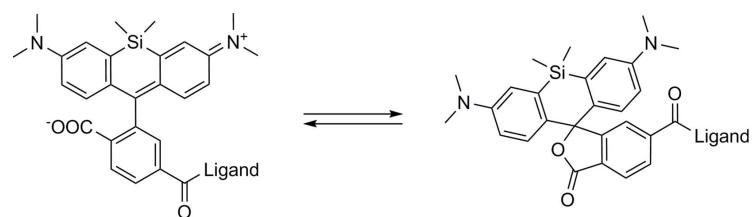


Figure 3. SiR derivatives exist in equilibrium between the fluorescent zwitterionic (open) form (left structure) and the non-fluorescent spiro (closed) form (right structure).

# Live Cell Imaging PRODUCTS

## Cytoskeleton's Live Cell Imaging Tools

### Live Cell Imaging Products

Product	Ex / Em	Amount	Cat #
<b>SPY555-Actin</b> Includes SPY555-Actin	555 / 580 nm	100 stains	CY-SC202
<b>SPY505-DNA</b> Includes SPY505-DNA	512 / 531 nm	100 stains	CY-SC101
<b>SPY555-DNA</b> Includes SPY555-DNA	555 / 580 nm	100 stains	CY-SC201
<b>SPY555-Tubulin</b> Includes SPY555-Tubulin	555 / 580 nm	100 stains	CY-SC203
<b>SPY595-DNA</b> Includes SPY595-DNA	599 / 615 nm	100 stains	CY-SC301
<b>SPY650-DNA</b> Includes SPY650-DNA	652 / 674 nm	100 stains	CY-SC501
<b>SPY650-Tubulin</b> Includes SPY650-Tubulin	652 / 674 nm	100 stains	CY-SC503
<b>SPY700-DNA</b> Includes SPY700-DNA	696 / 718 nm	100 stains	CY-SC601
<b>SiR-Actin™ Kit</b> Includes SiR-Actin and Verapamil	630 / 680 nm	50 nmol	CY-SC001
<b>SiR-Tubulin™ Kit</b> Includes SiR-Tubulin, and Verapamil	630 / 680 nm	50 nmol	CY-SC002
<b>Cytoskeleton Kit</b> Includes SiR-Actin, SiR-tubulin and Verapamil	630 / 680 nm	50 nmol each	CY-SC006
<b>SiR-DNA™ Kit</b> Includes SiR-DNA and Verapamil	630 / 680 nm	50 nmol	CY-SC007
<b>SiR700-Actin Kit</b> Includes SiR700-Actin and Verapamil	690 / 720 nm	35 nmol	CY-SC013
<b>SiR700-Tubulin Kit</b> Includes SiR700-Tubulin and Verapamil	690 / 720 nm	35 nmol	CY-SC014
<b>SiR700-DNA Kit</b> Includes SiR700-DNA and Verapamil	690 / 720 nm	35 nmol	CY-SC015
<b>Flipper-TR™ Kit</b> For fluorescence cell membrane microscopy	480 / 600 nm	50 nmol	CY-SC020

### Featured Papers and Application Notes

- ["Fluorogenic probes for live-cell imaging of the cytoskeleton"](#); G. Lukinavičius, et al. *Nature Methods* **11**, 731–733, 2014.
- ["STED Nanoscopy Reveals the Ubiquity of Subcortical Cytoskeleton Periodicity in Living Neurons"](#); E. D'Este, et al. *Cell Reports*, Volume **10**, Issue **8**, 1246 – 1251, 2015.
- ["A near-infrared fluorophore for live-cell super-resolution microscopy of cellular proteins"](#); G. Lukinavičius, et al.; *Nature Chemistry* **5**, 132–139, 2013.
- ["Dynamic actin filaments control the mechanical behavior of the human red blood cell membrane"](#); D. S. Gokhin, et al.; *Mol. Biol. Cell*; February 25, 2015.
- ["A cleavable cytolsin-neuropeptide Y bioconjugate enables specific drug delivery and demonstrates intracellular mode of action"](#); V. M. Ahrens, et al.; *J. Control. Release*; **209**:170–178, 2015.
- ["Red Si-rhodamine drug conjugates enable imaging in GFP cells"](#); E. Kim, et al.; *Chem. Commun.*, **50**, 4504-4507, 2014.
- ["A marginal band of microtubules transports and organizes mitochondria in retinal bipolar synaptic terminals"](#); M. Graffe, et al.; *J. Gen Physiol.* Vol. **146** No.1: 109-117, 2015.

### Application Notes

- ["A Bright Dye for Live-Cell STED Microscopy"](#); S. Pitsch, I. Köster.

## Explore Cytoskeleton's Novel Kits

### Acti-Stain™ Phalloidins

Product	Amount	Cat #
<b>Acti-stain 488™ phalloidin</b>	300 Slides	PHDG1-A
<b>Acti-stain 555™ phalloidin</b>	300 Slides	PHDH1-A
<b>Acti-stain 670™ phalloidin</b>	300 Slides	PHDN1-A
<b>Rhodamine Phalloidin</b>	1 x 500 µl	PHDR1

### Labeled Actin Proteins

Labeled Actin	Amount	Cat. #
<b>Rhodamine Actin Protein</b> Human platelet, non-muscle	4 x 10 µg 20 x 10 µg	APHR-A APHR-C
<b>Rhodamine Actin Protein</b> Rabbit skeletal muscle	10 x 20 µg 20 x 20 µg	AR05-B AR05-C

### Actin Biochem Kits

Product	Assays	Cat. #
<b>Actin Binding Protein Spin-Down Assay Biochem Kit</b> Rabbit skeletal muscle actin	30-100	BK001
<b>Actin Polymerization Biochem Kit (fluorescence format)</b> Measure actin polymerization <i>in vitro</i> , contains rabbit skeletal muscle actin.	30-100	BK003
<b>Actin Binding Protein Spin-Down Assay Biochem Kit</b> Human platelet actin	30-100	BK013
<b>G-Actin/F-actin In Vivo Assay Biochem Kit</b> Measure the distribution of monomer and polymer actin	30-100	BK037

### G-LISA Activation Assay Kits

Product	Assays	Cat. #
<b>RhoA G-LISA™ Activation Assay</b> (Luminescence format)	96	BK121
<b>RhoA G-LISA™ Activation Assay Kit</b> (Colorimetric format)	96	BK124
<b>Rac1,2,3 G-LISA™ Activation Assay</b> (Colorimetric format)	96	BK125
<b>Rac1 G-LISA™ Activation Assay</b> (Luminescence format)	96	BK126

<b>Rac1 G-LISA™ Activation Assay Kit</b> (Colorimetric Based)	96	BK128
<b>Ras G-LISA™ Activation Assay Kit</b> (Colorimetric Based)	96	BK131

<b>Total RhoA ELISA</b>	96	BK150
-------------------------	----	-------

### Pull Down Activation Assay Kits

Product	Assays	Cat. #
<b>Ras Pull-down Activation Assay Biochem Kit</b> (bead pull-down format)	50	BK008
<b>RhoA / Rac1 / Cdc42 Activation Assay Combo Biochem Kit</b> (bead pull-down format)	3 x 10	BK030
<b>Cdc42 Pull-down Activation Assay Biochem Kit</b> (bead pull-down format)	50	BK034
<b>RhoA Pull-down Activation Assay Biochem Kit</b> (bead pull-down format)	80	BK036