



Helping advance science,
one protein at a time.

Actin Ring-based Periodic Membrane Skeleton in Neuronal Axons

[Related Publications](#)

[Research Tools](#)

AUGUST 2017

Meetings

**The Triangle
Cytoskeleton Meeting**
September 24th-25th
Saxapahaw, NC

**American Society for Cell
Biology 2017**
Booth # 818
December 2-6,
Philadelphia, PA

**American Society for Cell
Biology 2017**
Booth # 818
December 2-6,
Philadelphia, PA

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

Actin Ring-based Periodic Membrane Skeleton in Neuronal Axons

Actin is an integral part of the neuronal cytoskeleton as it is involved in the regulation of neuronal polarization, cell morphology, the development of neuronal processes (i.e., growth cones with lamellipodial and filopodial extensions and dendritic spines), intracellular trafficking, and synaptic plasticity (dynamic changes in dendritic spine number and/or morphology)¹⁻³. Actin's presence in growth cones and dendritic spines have garnered the attention of scientists for decades; however, actin is also found in neuronal axons, though its presence there has been described as the "black sheep of the neuronal actin family"⁴. This is because the exact details of actin's structure and role in the axon are unknown. Recently, significant advances have been made in unraveling the structure of axonal actin with the discovery of the periodic membrane skeleton (PMS) by nanoscopic microscopy⁵ (Fig. 1). This newsletter discusses the discovery, structure, and possible functions of the PMS in axons.

PMS Discovery and Structure

Discovered in 2013, the PMS is a type of cortical actin and the primary component of the actin cortex, a mixture of F-actin and actin binding proteins which supports eukaryotic cells' plasma membrane and membrane-associated processes such as endo- and exocytosis and cell motility^{4,5}. In neuronal axons, including the initial segment⁶, the PMS consists of short actin filaments bundled into evenly spaced rings that wrap around the circumference of the axon with a periodicity of 180-190 nanometers⁵⁻⁹ (Fig. 1). The short filaments are stabilized by an adducin cap which controls the diameter of actin rings and axons, as well as actin filament growth within the rings^{6,10}. Adjacent actin rings are secured through cross-linkage by spectrin tetramers (β II in the axon proper and β IV in the axon initial segment)^{6,8,11}.

The PMS was described for the first time in fixed mammalian neurons and brain tissue slices using stochastic optical reconstruction microscopy (STORM)⁵. Shortly thereafter, these findings were confirmed by others using STORM, stimulated emission depletion [STED] nanoscopy, and structured illumination microscopy [SIM] in fixed mammalian and non-mammalian neurons^{5,6,8,11,12}, and most importantly, living mammalian neurons using either the F-actin live cell imaging probe SIR-actin^{7,10,13} (silicon rhodamine actin; Fig. 1) or exogenous

expression of fluorescently-labeled β II spectrin⁸.

In cultured mammalian neurons, the PMS is established in the first few days of development in the axon proximal to the cell body before moving distally along the axon as the neuron develops^{5,7,8,11}. In *Drosophila* primary neurons, development of the PMS begins within hours of plating¹¹. Notably, the authors reported differences between the PMS of very young (hours to 2 days old) vs mature (≥ 3 days old) *Drosophila* neurons. For instance, the PMS in young, growing axons, but not older axons, depends upon actin polymerization (i.e., nucleation factors)¹¹.

The actin binding proteins involved in nucleation, assembly, and maintenance of the actin rings are relatively unknown. In *Drosophila* neurons, two nucleation factors, Arp2/3 and formin DAAM [disheveled-associated activator of morphogenesis], participate in PMS formation, likely nucleating the multiple short filaments that comprise the PMS¹¹.

PMS Functions

Neuronal axons are the means by which neurons communicate and transport cargo anterogradely and retrogradely between the cell body and axon terminal. Maintaining healthy axons is necessary for normal brain function as axonal loss is permanent

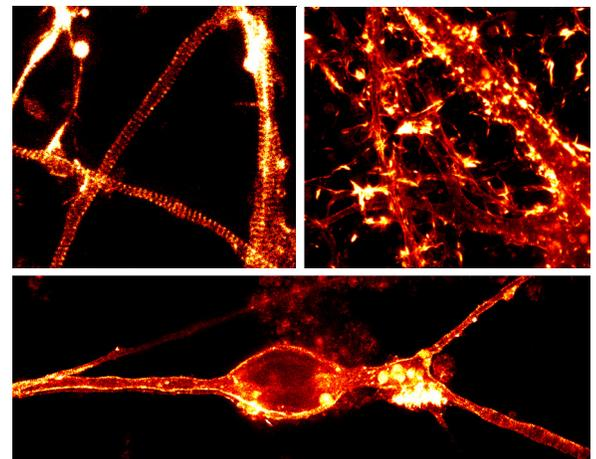


Figure 1. Live cell imaging in neurons with STED microscopy and 100 nM SIR-actin. Upper left: Hippocampal rat neuron, 8 DIV. Upper right: Hippocampal rat neuron, 17 DIV. Bottom image: Rat cerebellar granule neuron, 21 DIV. All images courtesy of Elisa D'Este, MPI Biophysical Chemistry, Göttingen.



Helping advance science,
one protein at a time.

Continued from Page 1

and underlies both normal aging deficits and various neurodegenerative disorders and diseases¹⁴. Functionally, the PMS has been hypothesized to serve as a scaffold for the axonal plasma membrane; spatially organize molecules important for axonal structure and action potential generation, such as ankyrin and sodium channels, respectively, into a periodic distribution in axons^{5,8}; assist in the docking of motor-associated cargo vesicles, especially along the distal part of the axon^{4,9,15}; and organize transmembrane proteins along the axon^{11,15-18}.

Actual functional data are sparse; however, a recent report using *Drosophila* primary neurons fixed and stained with SiR-actin concluded that the PMS is important in maintaining axon integrity by stimulating the polymerization of axonal microtubules (MTs), another primary cytoskeletal component of axons¹¹. Cytochalasin D-induced PMS disassembly resulted in breaks in MT bundles, reduced MT polymerization, and reduced axon numbers¹¹. Other MT-associated functional roles for PMS might involve assembling MTs into bundles¹⁹⁻²¹; serving as anchors for the minus end of MTs^{22,23}; and supporting dynein-mediated sliding of MTs and transport²⁴.

Summary

The recent discovery of actin rings that comprise a sub-membranous lattice in neuronal axons offers an exciting opportunity to better understand the role of axonal actin. Moreover, actin rings have also been described in the nodes of Ranvier in the peripheral nervous system and in at least some dendrites and dendritic spines^{7,25}. Similar to actin rings, actin waves and trails are other actin structures that have received little attention⁴. Without super-resolution microscopy (e.g., STORM, STED, SIM) and the development of live cell imaging probes such as SiR-actin, the existence and composition of the PMS would remain undiscovered. To help scientists further discover and define the composition, assembly, maintenance, and function of actin and other cytoskeletal structures, Cytoskeleton, Inc., provides SiR live cell imaging probes for F-actin, microtubules, DNA, and lysosomes, along with Acti-stain phalloidins for use in fixed cells.

Spirochrome™ Live Cell Imaging Probes

Description	Ex / Em	Cat. #	Amount
SiR700-Actin Kit Includes Verapamil	690 / 720 nm	CY-SC013	35 nmol
SiR-Actin Kit Includes Verapamil	630 / 680 nm	CY-SC001	50 nmol
SiR700-Tubulin Kit Includes Verapamil	690 / 720 nm	CY-SC014	35 nmol
SiR-Tubulin Kit Includes Verapamil	630 / 680 nm	CY-SC002	50 nmol
SiR-Lysosome Kit Includes Verapamil	630 / 680 nm	CY-SC012	50 nmol
SiR700-Lysosome Kit Includes Verapamil	690 / 720 nm	CY-SC016	35 nmol
SiR-DNA Kit includes Verapamil	630 / 680 nm	CY-SC007	50 nmol
SiR700-DNA Kit Includes Verapamil	690 / 720 nm	CY-SC015	35 nmol
Cytoskeleton Kit Includes SiR-Actin, SiR-Tubulin, and Verapamil	630 / 680 nm	CY-SC006	50 nmol each

Actin Proteins

Actin Protein	Cat. #	Amount
Rhodamine Actin Protein Human Platelet, Non-Muscle, >99% Pure	APHR-A APHR-C	4 x 10 µg 20 x 10 µg
Rhodamine Actin Protein Rabbit Skeletal Muscle, >99% Pure	AR05-B AR05-C	10 x 20 µg 20 x 20 µg

References

- Luo L. 2002. Actin cytoskeleton regulation in neuronal morphogenesis and structural plasticity. *Annu. Rev. Cell Dev. Biol.* **18**, 601-635.
- Cingolani L.A. and Goda Y. 2008. Actin in action: The interplay between the actin cytoskeleton and synaptic efficacy. *Nat. Rev. Neurosci.* **9**, 344-356.
- Stieess M. and Bradke F. 2011. Neuronal polarization: The cytoskeleton leads the way. *Dev. Neurobiol.* **71**, 430-444.
- Roy S. 2016. Waves, rings, and trails: The scenic landscape of axonal actin. *J. Cell Biol.* **212**, 131-134.
- Xu K. et al. 2013. Actin, spectrin, and associated proteins form a periodic cytoskeletal structure in axons. *Science.* **339**, 452-456.
- Letierrier C. et al. 2015. Nanoscale architecture of the axon initial segment reveals an organized and robust scaffold. *Cell Rep.* **13**, 2781-2793.
- D'Este E. et al. 2015. STED nanoscopy reveals the ubiquity of subcortical cytoskeleton periodicity in living neurons. *Cell Rep.* **10**, 1246-1251.
- Zhong H. et al. 2014. Developmental mechanism of the periodic membrane skeleton in axons. *eLife.* **3**, e04581.
- Leite S.C. and Sousa M.M. 2016. The neuronal and actin commitment: Why do neurons need rings? *Cytoskeleton.* **73**, 424-434.
- Leite S.C. et al. 2016. The actin-binding protein α -adducin is required for maintaining axon diameter. *Cell Rep.* **15**, 490-498.
- Qu Y. et al. 2017. Periodic actin structures in neuronal axons are required to maintain microtubules. *Mol. Biol. Cell.* **28**, 296-308.
- He J. et al. 2016. Prevalent presence of periodic actin-spectrin-based membrane skeleton in a broad range of neuronal cell types and animal species. *Proc. Natl. Acad. Sci. U.S.A.* **113**, 6029-6034.
- Lukinavicius G. et al. 2014. Fluorogenic probes for live-cell imaging of the cytoskeleton. *Nat. Methods.* **11**, 731-733.
- Adalbert R. and Coleman M.P. 2012. Axon pathology in age-related neurodegenerative disorders. *Neuropathol. Appl. Neurobiol.* **39**, 90-108.
- Gallo G. 2013. More than one ring to bind them all: Recent insights into the structure of the axon. *Dev. Neurobiol.* **73**, 799-805.
- Zhang Y. et al. 2016. Axon membrane skeleton structure is optimized for coordinated sodium propagation. *arXiv.1602.06348*.
- Albrecht D. et al. 2016. Nanoscopic compartmentalization of membrane protein motion at the axon initial segment. *J. Cell Biol.* **215**, 37-46.
- Machnicka B. et al. 2014. Spectrins: a structural platform for stabilization and activation of membrane channels, receptors and transporters. *Biochim. Biophys. Acta.* **1838**, 620-634.
- Prokop A. et al. 2013. Using fly genetics to dissect the cytoskeletal machinery of neurons during axonal growth and maintenance. *J. Cell Sci.* **126**, 2331-2341.
- Sanchez-Soriano N. et al. 2009. Mouse ACF7 and *Drosophila* Short stop modulate filopodia formation and microtubule organization during neuronal growth. *J. Cell Sci.* **122**, 2534-2542.
- Alves-Silva J. et al. 2012. Spectraplakins promote microtubule-mediated axonal growth by functioning as structural microtubule-associated proteins and EB1-dependent +TIPs (tip interaction proteins). *J. Neurosci.* **32**, 9143-9158.
- Nashchekin D. et al. 2016. Patronin/Shot cortical foci assemble the noncentrosomal microtubule array that specifies the *Drosophila* anterior-posterior axis. *Dev. Cell.* **38**, 61-72.
- Ning W. et al. 2016. The CAMSAP3-ACF7 complex couples noncentrosomal microtubules with actin filaments to coordinate their dynamics. *Dev. Cell.* **39**, 61-74.
- Myers K.A. et al. 2006. Microtubule transport in the axon: Re-thinking a potential role for the actin cytoskeleton. *Neuroscientist.* **12**, 107-118.
- Bar J. et al. 2016. Periodic F-actin structures shape the neck of dendritic spines. *Sci. Rep.* **6**, 37136.

Acti-stain Phalloidins

Acti-stain Phalloidin	Cat. #	Amount
Acti-stain™ 488 Phalloidin	PHDG1	500 µl
Acti-stain™ 555 Phalloidin	PHDH1	500 µl
Acti-stain™ 670 Phalloidin	PHDN1	500 µl
Phalloidin (rhodamine)	PHDR1	500 µl

Actin Activation Assay Biochem Kit

Actin Activation Assay	Cat. #	Amount
G-Actin/F-Actin In Vivo Assay Biochem Kit	BK037	30-100 Assays