



Helping advance science,  
one protein at a time.

## Microtubules and Motor-Driven Cargo Delivery Related Publications Research Tools

May  
2021

### 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

### Recent Citation Spotlights

- Pro-Inflammatory Stressors Alter The Vesiculome Of MIN6 Pancreatic Beta Cells
- SARS-CoV-2 infection alters the transcriptomic landscape of human airway cells
- A Live Look At Membrane Tension At The Plasma Membrane And Beyond
- RHOA-mediated mechanical force generation through Dectin-1

### Microtubules and Motor-Driven Cargo Delivery

Microtubules, composed of alpha- and beta-tubulin heterodimers, are one of the most recognizable filamentous structural proteins in the dynamic cytoskeleton network. Its function in mitosis and cell division is integral to growth and viability, but its essential role in intracellular organization during interphase is just as important. One core function is its role in cargo transport where microtubule networks have been compared to highways and byways utilized by motor proteins to transport cargo such as vesicles, organelles, protein complexes and many other cell components<sup>(1)</sup>. So what's new in the world of microtubule/motor-driven cargo transport and what's regulating it? Exciting new research is adding to our understanding of the dynamic microtubule/motor cargo system and its effect on disease.

### Microtubule/Motor-Driven Cargo Localization of mRNA and Ribosomes Regulates Cardiac Hypertrophy

When cardiomyocytes are exposed to mechanical or neurohumoral stimuli they activate signaling cascades that upregulate transcription and translation of new sarcomeres to promote hypertrophy. Recent studies suggest that mRNA and ribosomes localize to distinct regions in response to stimuli to promote local sarcomere synthesis<sup>(2)</sup>. Scarborough et al. found that microtubules regulate mRNA and ribosome cargo localization to control local translation and dictate cardiac growth<sup>(3)</sup>. When mice or neonatal rat ventricular myocytes (NRVMs) were exposed to stimulus like phenylephrine (PE) they underwent hypertrophy as expected; this growth could be inhibited with microtubule disrupting drugs like colchicine or nocodazole supporting previous findings that the microtubule system plays an important role in cardiac hypertrophy. Unexpectedly, when the group evaluated changes in the cardiac hypertrophic genes in PE treated NRVMs they found that in the absence or presence of colchicine they saw a similar elevation of these genes. The group determined that the microtubule/motor cargo delivery system was essential to localize mRNA and ribosomes to the correct location for effective translation of sarcomere forming proteins. When proper microtubule function was blocked with microtubule disrupting drugs, mRNAs and ribosomes collapsed around the nucleus leading to rapid protein degradation and suppressed hypertrophy. These data suggest that proper localization of the translation machinery and not just translation levels are critically important for cardiac hypertrophy, and the microtubules/motor-driven cargo delivery system are important players in this process.

### Identifying New Components of The Complex That Transports Mitochondria Cargo

Due to the size of neurons active transport of mitochondria

is necessary; this process is primarily carried out by the microtubule/motor cargo delivery system<sup>(4)</sup>. Impairment of mitochondrial trafficking in neurons results in human neurodegenerative diseases. Previous work has identified Miro and Track as two key components that form a complex used to link mitochondria to kinesin-1 and dynein and transport them in both directions along microtubules<sup>(5)</sup>. However, Miro deletion did not completely block mitochondria movement in a drosophila model, offering the possibility that alternative or additional complexes exist to transport mitochondria. A new study by Zhao et al. identified metaxins, MTX-1 and MTX-2, as new components of the complex for mitochondria transport<sup>(6)</sup>. Their early studies determined that loss of MTX-2 abolished mitochondria distribution. Furthermore, their studies identified MTX-2 and Miro as core adaptor components while interaction with MTX-1 or TRAK-1 specify each complex for interaction with kinesin-1 or dynein respectively. This study identified novel adaptor proteins in the transport complex that is critical for mitochondria distribution in neurons and showed that disruption of this complex and ultimately mitochondrial transport resulted in dendrite degradation. Importantly the group identified unique complexes that mediated cargo interaction with specific motors which adds to our understanding of the regulatory mechanisms of the microtubule/motor cargo delivery system.

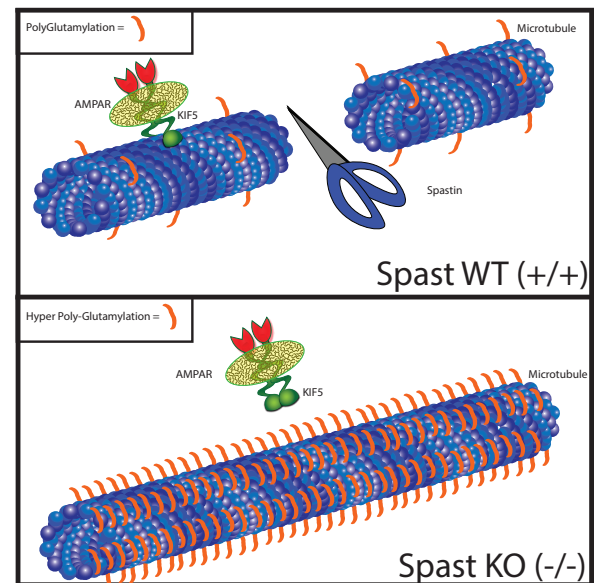


Figure 1: Spastin depletion impairs synaptic cargo transport and cell surface delivery



# Tubulin and Motor PRODUCTS

Helping advance science,  
one protein at a time.

## Continued from Page 1

### Spastin Depletion, Tubulin Polyglutamylation, and Impaired Kinesin Mediated Transport

Mutations in spastin, a microtubule severing protein, is the primary cause of hereditary spastic paraplegia (HSP)<sup>(7)</sup>. Mutations in kinesin family protein (KIF) 5 has also been linked to HSP. Complicated forms of HSP are associated with neurologic impairments and late-onset dementia<sup>(8)</sup>. New findings by Lopes et al. used a spastin knockout mouse to investigate the causes of neurologic deficits identified in complicated HSP; as expected, altered microtubules were observed and played a critical role in a reduction of synapses (in the range of 40%), a measure of neurological dysfunction<sup>(9)</sup>. Neurons from these knockout mice had movement rate defects in microtubule/motor-driven transport of AMPA receptors and had less KIF 5 associated with these microtubules. Additionally, they identified enhanced polyglutamylation in neurons where spastin was knocked out (see figure 1). A link between Spastin function and polyglutamylation on the C-terminal tails of tubulin has been linked previously<sup>(10)</sup>. To define the importance of enhanced polyglutamylation, Lopes et al. utilized shRNA tools to deplete polyglutamylase subunit 1 (PGs1), which catalyzes 80% of the polyglutamylation of neuronal tubulin. Depletion of PGs1 was sufficient to reverse the enhanced polyglutamylation observed in spastin knockout neurons; more importantly, it also restored normal KIF5 interaction, neuronal transport, and synapse numbers. These findings suggests that post-translational modifications (PTMs) that occur on microtubules can have a profound effect on microtubule/motor cargo delivery system, and it will be interesting to further define the interplay between tubulin PTMs and motors.

### Summary

These findings highlight how important proper regulation of the microtubule/motor cargo delivery system is for downstream process that regulate pathologic outcomes like cardiac hypertrophy and neurodegeneration. As our understanding of these biological consequences are uncovered, it will be even more important to identify key players in the transport complexes that may be targets for intervention as a mechanism to properly control this process. It will be important to understand how proteins that regulate the dynamicity and PTM state of tubulin also impact cargo movement and delivery along these microtubule tracks. What an exciting time to be investigating tubulin regulated cargo delivery and Cytoskeleton Inc. is proud to assist researchers with essential tubulin tools to aide in their investigation.

### Purified Tubulins and MAPs

Product	Amount	Cat #
<b>Tubulin protein (97% pure)</b> Source: porcine brain	1 x 4 mg 1 x 40 mg	HTS03-A HTS03-B
<b>Microtubule associated protein rich fraction</b> Source: porcine brain	1 x 100 µg 5 x 100 µg	MAPF-A MAPF-C
<b>Tau protein</b> Source: bovine brain	1 x 50 µg 3 x 50 µg	TA01-A TA01-B
<b>Tubulin protein (&gt;99% pure)</b> Source: porcine brain	1 x 1 mg 5 x 1 mg	T240-A T240-B

### Signal Seeker™ PTM Detection Kits

Product	Assays	Cat. #
<b>Signal-Seeker™ Phosphotyrosine Detection Kit</b>	30	BK160
<b>Signal-Seeker™ Ubiquitination Detection Kit</b>	30	BK161
<b>Signal-Seeker™ SUMOylation 2/3 Detection Kit</b>	30	BK162
<b>Signal-Seeker™ Acetyl-Lysine Detection Kit</b>	30	BK163
<b>Signal-Seeker™ SUMOylation 1 Detection Kit</b>	30	BK165

### Custom Motor Proteins

Product	Amount	Cat. #
<b>Dynein motor protein</b>	1 x 50 µg	CS-DN01-A
<b>Kinesin heavy chain motor domain protein</b> GST tagged: Homo sapiens recombinant	2 x 25 µg 1 x 1 mg	KR01-A KR01-XL
<b>KIF3C kinesin motor domain protein</b> GST tagged: Homo sapiens recombinant	2 x 25µg	KF01-A
<b>KIF3C kinesin motor domain protein</b> GST tagged: Homo sapiens recombinant	2 x 25 µg	KC01-A
<b>Custom purification of kinesin motors</b>	Custom	Inquire for Price

### Tubulin Live Cell Imaging Products

Product	Ex / Em	Amount	Cat #
<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>SiR700-Tubulin Kit</b> 35 nmol SiR700-Tubulin and 1 µmol verapamil	680 / 720 nm	50 nmol	CY-SC014
<b>SPY555-Tubulin Kit</b> Includes SPY555-Tubulin and Verapamil	555 / 580 nm	100 stains	CY-SC203
<b>SPY650-Tubulin</b> Includes SPY650-Tubulin and Verapamil	652 / 674 nm	100 stains	CY-SC503

### Tubulin Kits

Product	Assays	Cat #
<b>Tubulin polymerization HTS assay using &gt;97% pure tubulin</b> OD based - Porcine	24	BK004P
<b>Tubulin polymerization assay using &gt;99% pure tubulin</b> OD based - Porcine	24-30	BK006P
<b>Tubulin polymerization assay using &gt;99% pure tubulin</b> Fluorescence based	96	BK011P
<b>Microtubule Binding Protein Spin-Down Assay Biochem Kit</b>	50-100	BK029
<b>Microtubule/Tubulin In Vivo Assay Biochem Kit</b>	30-100	BK038

### References

- Fourriere L, Jimenez AJ, Perez F, Boncompain G. The role of microtubules in secretory protein transport. *J Cell Sci.* 2020;**133**(2).
- Lewis YE, Moskovitz A, Mutlak M, Heineke J, Caspi LH, Kehat I. Localization of transcripts, translation, and degradation for spatiotemporal sarcomere maintenance. *J Mol Cell Cardiol.* 2018;**116**:16-28.
- Scarborough EA, Uchida K, Vogel M, Erlitzki N, Iyer M, Phyo SA, et al. Microtubules orchestrate local translation to enable cardiac growth. *Nat Commun.* 2021;**12**(1):1547.
- Sheng ZH. Mitochondrial trafficking and anchoring in neurons: New insight and implications. *J Cell Biol.* 2014;**204**(7):1087-98.
- Schwarz TL. Mitochondrial trafficking in neurons. *Cold Spring Harb Perspect Biol.* 2013;**5**(6).
- Zhao Y, Song E, Wang W, Hsieh CH, Wang X, Feng W, et al. Metaxins are core components of mitochondrial transport adaptor complexes. *Nat Commun.* 2021;**12**(1):83.
- Trotta N, Orso G, Rossetto MG, Daga A, Broadie K. The hereditary spastic paraplegia gene, spastin, regulates microtubule stability to modulate synaptic structure and function. *Curr Biol.* 2004;**14**(13):1135-47.
- Murphy S, Gorman G, Beetz C, Byrne P, Dytko M, McMonagle P, et al. Dementia in SPG4 hereditary spastic paraplegia: clinical, genetic, and neuropathologic evidence. *Neurology.* 2009;**73**(5):378-84.
- Lopes AT, Hausrat TJ, Heisler FF, Gromova KV, Lombino FL, Fischer T, et al. Spastin depletion increases tubulin polyglutamylation and impairs kinesin-mediated neuronal transport, leading to working and associative memory deficits. *PLoS Biol.* 2020;**18**(8):e3000820.
- Valenstein ML, Roll-Mecak A. Graded Control of Microtubule Severing by Tubulin Glutamylation. *Cell.* 2016;**164**(5):911-21.