



The Protein Experts

CYTOSKELETON NEWS

NEWS FROM CYTOSKELETON INC.

Helping advance science, one protein at a time.

NOV/DEC
2011

this issue

Rho GTPase control of neurite extension

Neuroscience Publications

Neuroscience Products

Rho GTPase control of neurite extension

The Rho family of guanine nucleotide triphosphatases (GTPases) governs neurite extension by regulating the interplay of actin and microtubule (MT) cytoskeletal dynamics. During neurite extension, GTPases are activated by growth factors and guidance cue proteins through receptor binding on the membrane of the growth cone, the motile tip of the extending axonal neurite. The most studied GTPases are RhoA, Rac1 and Cdc42 (1-3).

Neurite extension is a process essential for nervous system development, synaptic plasticity and treatment and repair of brain diseases and injuries. Neurite extension or growth can be considered a three step process (4).

Step 1: Protrusion: Rac1 and Cdc42 induce formation of actin filaments via activation of the downstream effector PAK which activates the WAVE (Rac) or N-WASP (Cdc42) protein that complexes with Arp2/3 to produce actin-rich lamellipodia or filopodia, respectively (1-3). These actin-rich protrusions form and extend from the peripheral domain of the growth cone and act as environmental sensors for axon growth and guidance (Fig. 1). There is also increasing evidence that actin filaments and MTs

Step 2: Engorgement: MTs and organelles move into the protruding regions with MTs providing a scaffold for anterograde and retrograde transport and a structural connection between the stabilized axon shaft and the motile central and peripheral domains of the growth cone (5).

Step 3: Consolidation: RhoA and its downstream effector Rho Kinase (ROCK) activate the myosin II motor protein which mediates actomyosin contraction and the formation of stress fibers and focal adhesions. Activated RhoA/myosin II induces F-actin and MT bundling and re-distribution of F-actin and MTs away from the growth cone periphery, causing the membrane to shrink around MT bundles, consolidating a segment of the developing axon shaft (5,6,11-18).

GTPase signaling is much more complex than initially believed and is influenced by many factors that affect the balance of GTPase activation and inactivation (Fig 2). For instance, RhoA signaling via mDia promotes axon extension and stabilization of MTs (2,3), Rac1 activity is sometimes required for growth cone inhibition (19) and over-expression of the Rac GEF Vav2 inhibits neurite extension (20). Both Rac1 and RhoA/ROCK activate LIM kinases which inhibit the actin depolymerizing factor cofilin, leading to neurite stabilization and retraction (3,21). These paradoxical findings suggest that GTPase control of neurite extension results from a balancing of multiple factors, including GEF/GAP activity, the specific downstream pathways activated (20,22) and the degree to which actin filaments and MTs dynamically interact (Fig. 2).

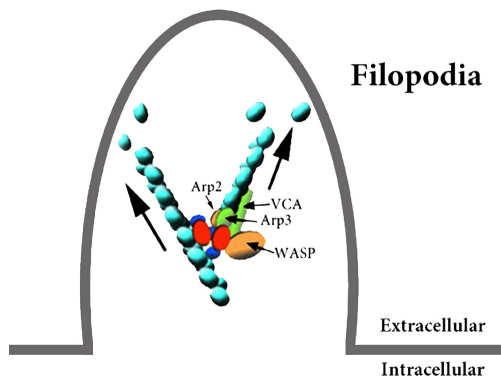


Figure 1. Actin polymerization mediated by Arp2/3/VCA/WASP proteins in filopodia.

interact during neurite extension. MTs defined as dynamic (vs stable), based on specific post-translational modifications, are found in the peripheral domain of the growth cone where MTs can interact with F-actin via proteins that can bind or regulate both cytoskeletal elements (5,6). For instance, Rac1 regulates interplay between actin and MT dynamics during protrusion (6-10).

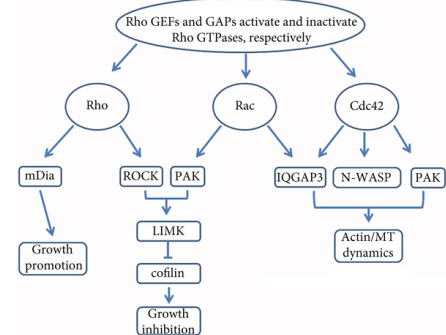


Figure 2. Rho GTPase pathways controlling neurite extension. Adapted from ref. 3

Upcoming Meetings

- Neuroscience Nov 12-16
- AACR-NCI-EORTC Nov 12-16
- ASCB Dec 3-7

Cytoskeleton Products

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

Contact Us

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

Distributors

www.cytoskeleton.com/distributors/

Neuroscience news

Neuroscience publications

Neuroscience tools



Helping advance science,
one protein at a time.

NEUROSCIENCE PRODUCTS

Visit cytoskeleton.com for more information.

Rho GTPase control of neurite extension (Cont'd from Pg 1)

For more information on how GTPases regulate neurite extension, please see the references and reviews below.

References

- Samuel and Hynds, 2010. *Mol. Neurobiol.* v 42, p 133.
- Govek et al., 2005. *Genes Dev.* v 19, p 1.
- Hall and Lalli, 2010. *Cold Spring Harb. Perspect. Biol.* 2:a001818.
- Dent et al., 2010. *Cold Spring Harb. Perspect. Biol.* 3:a001800.
- Fukushima, 2011. *Cytoskeleton of the Nervous System*. R.A. Nixon, A. Yuan (eds.). New York: Springer.
- Conde and Caceres, 2009. *Nat. Rev. Neurosci.* v 10, p 319
- Gallo, 1998. *J. Neurobiol.* v 35, p 121.
- Grabham et al., 2003. *J. Cell Sci.* v 116, p 3739.
- Rochlin et al., 1999. *Mol. Biol. Cell.* v 10, p 2309.
- Rodriguez et al., 2003. *Nat. Cell Biol.* v 5, p 599.
- Fritz and VanBerkum, 2002. *Dev. Biol.* v 252, p 46.
- Loudon et al., 2006. *J. Neurobiol.* v 66, p 847.
- Murray et al., 2010. *Neural Develop.* v 5, p 16.
- Lin et al., 1996. *Neuron.* v 16, p 769.
- Medeiros et al., 2006. *Nat. Cell Biol.* v 8, p 215.
- Fukushima and Morita, 2006. *Brain Res.* v 1094, p 65.
- Burnette et al., 2008. *Dev. Cell.* v 15, p 163.
- Schaefer et al., 2008. *Dev. Cell.* v 15, p 146.
- Jin and Strittmatter, 1997. *J. Neurosci.* v 17, p 6256.
- Moon and Gomez, 2010. *Mol. Cell Neurosci.* v 44, p 118.
- Amano et al., 2010. *Cytoskeleton.* v 67, p 545.
- Woo and Gomez, 2006. *J. Neurosci.* v 26, p 1418.

Pull-down Activation Assays

| Pull-down Activation Assays | Effector Protein | Cat. # | # of Assays |
|--|------------------------|--------|-------------|
| Cdc42 Activation Assay Biochem Kit™ | p21 activated kinase 1 | BK034 | 50 |
| Rac1 Activation Assay Biochem Kit™ | p21 activated kinase 1 | BK035 | 50 |
| Ras Activation Assay Biochem Kit™ | Kinase Raf1 | BK008 | 50 |
| RalA Activation Assay Biochem Kit™ | Ral-BP1 | BK040 | 50 |
| RhoA Activation Assay Biochem Kit™ | Rhotekin | BK036 | 80 |

Fluorescent Probes

| Product | Excitation | Emission | Cat. # | Amount |
|-----------------------------------|------------|----------|---------|------------|
| Acti-stain™ 488 phalloidin | 480 nm | 535 nm | PHDG1-A | 300 Slides |
| Acti-stain™ 555 phalloidin | 535 nm | 585 nm | PHDH1-A | 300 Slides |
| Acti-stain™ 670 phalloidin | 630 nm | 680 nm | PHDN1-A | 300 Slides |
| Rhodamine phalloidin | 535 nm | 585 nm | PHDR1 | 1 x 500 µl |

Neuroscience Tools

G-LISA® Activation Assays

| Product | Cat # | # of Assays |
|--|-------|-------------|
| RhoA G-LISA® Activation Assay, colorimetric | BK124 | 96 |
| RhoA G-LISA® Activation Assay, luminescence | BK121 | 96 |
| Cdc42 G-LISA® Activation Assay, colorimetric | BK127 | 96 |
| RalA G-LISA® Activation Assay, colorimetric | BK129 | 96 |
| Rac1,2,3 G-LISA® Activation Assay, colorimetric | BK125 | 96 |
| Rac1 G-LISA® Activation Assay, colorimetric | BK128 | 96 |

Cytoskeletal Proteins

| Labeled Actins | Source | Purity | Cat. # | Amount |
|---|----------------------------|--------|----------------------|--------------------------|
| Rhodamine Actin Protein | Human platelet, non-muscle | >99% | APHR-A APHR-C | 4 x 10 µg 20 x 10 µg |
| Rhodamine Actin Protein | Rabbit skeletal muscle | >99% | AR05-B AR05-C | 10 x 20 µg 20 x 20 µg |
| Arp2/3 Protein Complex | Bovine Brain | >95% | RP01-A RP01-B | 2 x 50 µg 6 x 50 µg |
| WASP VCA domain GST Fusion Protein | Recombinant human | >95% | VCG03-A VCG03-B | 1 x 500 µg 5 x 500 µg |
| AMCA Labeled Tubulin | Porcine Brain | >99% | TL440M-A TL440M-B | 5 x 20 µg 20 x 20 µg |
| HiLyte Fluor™ 488 Labeled Tubulin | Porcine Brain | >99% | TL488M-A TL488M-B | 5 x 20 µg 20 x 20 µg |
| TRITC Rhodamine Labeled Tubulin | Porcine Brain | >99% | TL590M-A TL590M-B | 5 x 20 µg 20 x 20 µg |
| X-Rhodamine Labeled Tubulin | Bovine Brain | >99% | TL620M-A TL620M-B | 5 x 20 µg 20 x 20 µg |
| HiLyte Fluor™ 647 Labeled Tubulin | Porcine Brain | >99% | TL670M-A TL670M-B | 5 x 20 µg 20 x 20 µg |

G-switch™ Activators and Inhibitors

| G-Protein Modulator | Cell Entry Mechanism | Rho Family Protein Modulation | Cat. # | Amount |
|--|----------------------|-------------------------------|------------------|-------------------------------|
| Rho Activator II Deamidation of Rho Gln-63 | Cell permeable | Direct | CN03-A CN03-B | 1 x 20 µg 5 x 20 µg |
| Rho Inhibitor I ADP ribosylation of Rho Asn-41 | Cell permeable | Direct | CT04-A CT04-B | 1 x 20 µg 5 x 20 µg |
| Rho/Rac/Cdc42 Activator I Deamidation of Rho Gln-63 & Rac/Cdc42 Gln-61 | Cell permeable | Direct | CN04-A CN04-B | 1 x 20 µg 5 x 20 µg |
| Rho Pathway Inhibitor I Rho kinase (ROCK) inhibitor Y-27632 | Cell permeable | Direct | CN06-A CN06-B | 5 x 10 units 20 x 10 units |
| Rho Activator I SHP-2 phosphatase mediated Rho activation | Cell permeable | Indirect | CN01-A CN01-B | 5 x 10 units 20 x 10 units |
| Rac/Cdc42 Activator II EGF receptor mediated Rac/Cdc42 activation | Receptor mediated | Indirect | CN02-A CN02-B | 5 x 10 units 20 x 10 units |

www.cytoskeleton.com