



The Protein Experts

CYTOSKELETON NEWS

NEWS FROM CYTOSKELETON INC.

Helping advance science,
one protein at a time.

December
2019

Phenotypic Profiling: Actin-focused Cancer Therapeutics

Related Publications

Research Tools

Phenotypic Profiling: Actin-focused Cancer Therapeutics

As an integral component of the mammalian cell cytoskeleton, actin is involved in multiple physiological functions, including cell growth, motility, trafficking, and division. These basic functions require remodeling of the actin cytoskeleton, as well as extension and withdrawal of actin-based cellular neurites (e.g., lamellipodia, filopodia), all of which rely upon rapid dynamic cycling between filamentous actin (F-actin) and monomer actin (G-actin)¹. Correspondingly, dysfunctional actin cytoskeletal dynamics are a pathophysiological feature of many human diseases, with perhaps cancer being the prototypical example. For these reasons, actin is a theoretically attractive anti-cancer therapeutic target. However, in practice, actin has proven to be a poor target because of toxic side effects due in large part to the inability of therapeutics to distinguish between actin isoforms^{2,3}. Thus, the actin in cancerous and healthy cells are affected by cancer therapies acting directly upon actin, leading to toxic side effects in different organ systems (e.g., heart and diaphragm)^{2,3}. Recently, drug discovery has shifted from actin to actin-associated structural proteins such as the Arp2/3 complex and tropomyosins (Tpms) as these proteins offer multiple isoforms for selective targeting and the opportunity to avoid toxic side effects²⁻⁴ (Fig. A).

Developing drugs that regulate actin dynamics by targeting structural proteins is not without challenges. Necessary structural knowledge of target proteins is obtained through functional domain mapping and identification of possible drug binding pockets prior to the start of rational drug design³. In the case of actin-associated regulatory proteins such as Arp2/3 and Tpms, these data are lacking (Fig. A). In the absence of complete structural data, researchers screen existing compound libraries against target proteins, searching for drugs that inhibit function³. An attribute of a putative therapeutic target is the existence of multiple isoforms which offer the possibility of selective targeting, especially if only certain isoforms are associated with disease. Moreover, structural isoforms offer more druggable binding pockets which means that different drugs can regulate the same protein, potentially with different mechanisms of action. To take advantage of existing small molecule compound libraries and deal with the lack of structural knowledge of candidate actin-binding proteins, one technique that is employed is phenotypic profiling³. This screening protocol utilizes cellular imaging to measure how compounds affect multiple, complex cellular architectures, including cell size, morphology, and in the case of actin, specific details of the filaments' attributes (number, length, shape, subcellular distribution, signal intensity) by computational and automated image analyses to identify compounds producing relevant biological effects^{3,5-7}. Furthermore, when a relevant compound's mechanism of action is known, this information is used to both further evaluate this compound and predict the mechanisms of action

of other compounds that produce similar phenotypes^{3,7}.

The Arp2/3 complex and Tpms are promising therapeutic targets in cancer drug discovery programs^{3,8-12}. Arp2/3 activity can be inhibited by different drugs through different mechanisms of actions¹¹. Tpm isoforms differentially regulate how susceptible F-actin is to actin-modulating drugs^{3,12}. Tpms consist of over 40 isoforms and display individual spatial and temporal localization patterns during cell/tissue development¹². The actin cytoskeleton consists of both co-polymers of actin and Tpms and polymers of Tpm-free actin¹³; however, a recent study reports that Tpms coat the majority of F-actin in human cells¹⁴. Tpms determine the functional capability of actin filaments in an isoform-dependent manner¹⁵⁻²¹ (Fig. A). One important characteristic of a protein that qualifies it for further study as a therapeutic target is its specific modulation in activity/expression in cancer cell lines. In the case of Tpms, isoform 3.1 is significantly and selectively up-regulated in multiple cancer cell lines; thus, it is a disease-associated isoform that can be specifically targeted^{2,14,22,23} (Fig. A). Importantly, targeting Tpm3.1 also avoids toxic side effects as muscle Tpms do not express this isoform^{22,24,25}. Currently under investigation are three

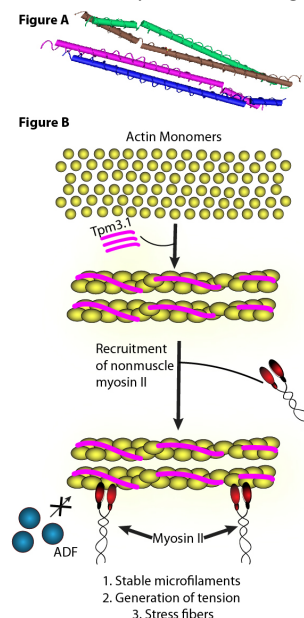


Figure A. Tropomyosin isoform 3.1 (Tpm3.1) is a cancer-associated actin-binding protein. (Top) Crystal structure of an N-terminal fragment of Tpm3.1. PDB ID: 6OTN.

Figure B. (Bottom) Tpm3.1 regulates how F-actin interacts with non-muscle myosin II and ADF/cofilin. Adapted from reference (12).

Meetings

ASCB/EMBO Annual Meeting
December 7-11th
Walter E Washington
Convention Center,
Washington D.C
Booth 701

Cold Spring Harbor
Conference - Development
and 3D Modeling of the
Human Brain
December 9 - 12th
Cold Spring Harbor, NY
Cytoskeleton Supported

Cold Spring Harbor
Conference - Systems
Biology: Global Regulation of
Gene Expression
March 11-14th 2020
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



The Protein
Experts

Helping advance science,
one protein at a time.

Actin PRODUCTS

Continued from Page 1

structurally unique, small molecule compounds that selectively target Tpm3.1 via distinct molecular interactions within Tpm3.1's binding pocket^{3,22,24,25}.

Summary

Investigating actin-associated structural proteins as novel anti-cancer therapeutic targets offers great potential for the development of novel and specific cancer therapies, but there are several caveats and methodological considerations when using phenotypic screening. For instance, 1. candidate cell lines must possess a highly conserved actin cytoskeleton; 2. any actin cytoskeleton stains must be compatible with robotic manipulation; 3. different actin phenotypes must be detectable with the specific screening set-up used and multiple parameters must be measurable (e.g., filament number and length, cell morphology and size, subcellular distribution of F-actin and its distinct structural forms); and 4. actin dynamics within cardiac and skeletal muscle cells must not be affected³. To assist researchers in screening compounds for their interactions with actin and associated proteins, Cytoskeleton offers a variety of purified actin proteins and actin-associated proteins, actin activity and binding assay kits, and live cell imaging probes for F-actin.

Actin Biochem Kits

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

Actin Live Cell Imaging Probes

Product	Ex/Em	Amount	Cat. #
SiR-Actin Kit includes SiR-Actin and Verapamil	630 / 680 nm	50 nmol	CY-SC001
SiR700-Actin Kit Includes SiR700-Actin and Verapamil	690 / 720 nm	35 nmol	CY-SC013
Cytoskeleton Kit Includes SiR-Actin, SiR-Tubulin and Verapamil	630 / 680 nm	50 nmol each	CY-SC006

Labeled Actin Proteins

Labeled Actin	Amount	Cat. #
Biotinylated Actin Protein Rabbit skeletal muscle	5 x 20 µg 20 x 20 µg	AB07-A AB07-C
Pyrene Actin Protein Rabbit skeletal muscle	1 x 1 mg 5 x 1 mg	AP05-A AP05-B
Rhodamine Actin Protein Human platelet, non-muscle	4 x 10 µg 20 x 10 µg	APHR-A APHR-C
Rhodamine Actin Protein Rabbit skeletal muscle	10 x 20 µg 20 x 20 µg	AR05-B AR05-C

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.
- 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.
- Creed S.J. et al. 2008. Tropomyosin isoforms define distinct microfilament populations with different drug susceptibility. *Eur. J. Cell Biol.* **87**, 709-720.
- Feng Y. et al. 2009. Multi-parameter phenotypic profiling: using cellular effects to characterize small-molecule compounds. *Nat. Rev. Drug Discov.* **8**, 567-578.
- Eliceiri K.W. et al. 2012. Biological imaging software tools. *Nat. Methods.* **9**, 697-710.
- Caicedo J.C. et al. 2017. Data-analysis strategies for image-based cell profiling. *Nat. Methods.* **14**, 849-863.
- Jay P. et al. 2000. ARP3beta, the gene encoding a new human actin-related protein, is alternatively spliced and predominantly expressed in brain neuronal cells. *Eur. J. Biochem.* **267**, 2921-2928.
- Millard T.H. et al. 2003. Identification and characterisation of a novel human isoform of Arp2/3 complex subunit p16-ARC/ARPC5. *Cell Motil. Cytoskeleton.* **54**, 81-90.
- Abella J.V. et al. 2016. Isoform diversity in the Arp2/3 complex determines actin filament dynamics. *Nat. Cell Biol.* **18**, 76-86.
- Nolen B.J. et al. 2009. Characterization of two classes of small molecule inhibitors of Arp2/3 complex. *Nature.* **460**, 1031-1034.
- Gunning P.W. et al. 2015. Tropomyosin - master regulator of actin filament function in the cytoskeleton. *J. Cell Sci.* **128**, 2965-2974.
- Lazarides E. 1976. Two general classes of cytoplasmic actin filaments in tissue culture cells: the role of tropomyosin. *J. Supramol. Struct.* **5**, 531-563.
- Meiring J.C.M. et al. 2018. Co-polymers of actin and tropomyosin account for a major fraction of the human actin cytoskeleton. *Curr. Biol.* **28**, 2331-2337.
- Bryce N.S. et al. 2003. Specification of actin filament function and molecular composition by tropomyosin isoforms. *Mol. Biol. Cell* **14**, 1002-1016.
- Tojkander S. et al. 2011. A molecular pathway for myosin II recruitment to stress fibers. *Curr. Biol.* **21**, 539-550.
- Creed S.J. et al. 2011. Tropomyosin isoform 3 promotes the formation of filopodia by regulating the recruitment of actin-binding proteins to actin filaments. *Exp. Cell Res.* **317**, 249-261.
- Janco M. et al. 2016. The impact of tropomyosins on actin filament assembly is isoform specific. *Bioarchitecture.* **6**, 61-75.
- Brayford S. et al. 2016. Tropomyosin promotes lamellipodial persistence by collaborating with Arp2/3 at the leading edge. *Curr. Biol.* **26**, 1312-1318.
- Gateva G. et al. 2017. Tropomyosin isoforms specify functionally distinct actin filament populations *in vitro*. *Curr. Biol.* **27**, 705-713.
- Pathan-Chhatbar S. et al. 2018. Three mammalian tropomyosin isoforms have different regulatory effects on nonmuscle myosin-2B and filamentous b-actin *in vitro*. *J. Biol. Chem.* **293**, 863-875.
- Stehn J.R. et al. 2013. A novel class of anticancer compounds targets the actin cytoskeleton in tumor cells. *Cancer Res.* **73**, 5169-5182.
- Coombes J.D. et al. 2015. Ras transformation overrides a proliferation defect induced by Tpm3.1 knockout. *Cell. Mol. Biol. Lett.* **20**, 626-646.
- Currier M.A. et al. 2017. Identification of cancer-targeted tropomyosin inhibitors and their synergy with microtubule drugs. *Mol. Cancer Ther.* **16**, 1555-1565.
- Kee A.J. et al. 2018. On-target action of anti-tropomyosin drugs regulates glucose metabolism. *Sci. Rep.* **8**, 4604.

Unlabeled Actin Proteins

Unlabeled Actins	Amount	Cat. #
Actin Protein Rabbit skeletal muscle	4 x 250 µg 2 x 1 mg 5 x 1 mg 10 x 1 mg 20 x 1 mg	AKL99-A AKL99-B AKL99-C AKL99-D AKL99-E
Actin Protein Rabbit skeletal muscle	1 x 1 mg 5 x 1 mg	AKL95-B AKL95-C
Actin Protein Bovine cardiac muscle	1 x 1 mg 5 x 1 mg	AD99-A AD99-B
Actin Protein Smooth muscle, chicken gizzard	1 x 1 mg 5 x 1 mg	AS99-A AS99-B
Actin Protein Human platelet, non-muscle	2 x 250 µg 1 x 1 mg 5 x 1 mg	APHL99-A APHL99-C APHL99-E
Pre-formed Actin Filaments Rabbit skeletal muscle	1 x 1 mg 5 x 1 mg	AKF99-A AKF99-B