



Helping advance science,
one protein at a time.

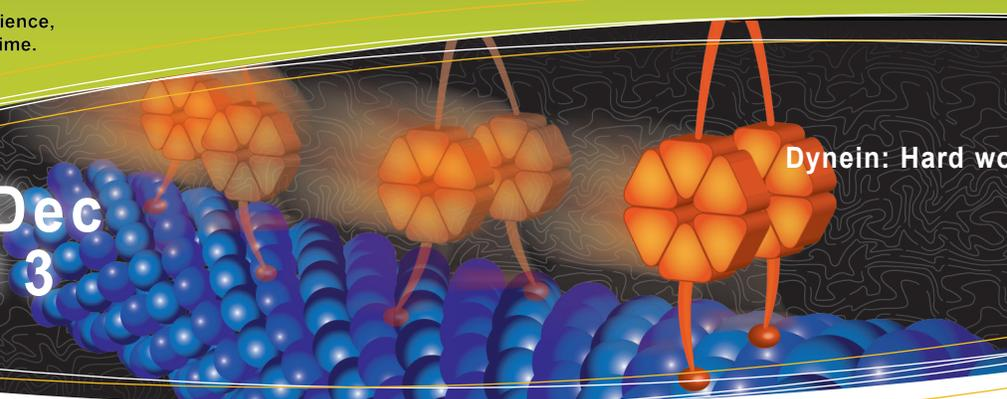
Nov/Dec
2013

this issue

Dynein: Hard working, but team oriented

Related Publications

Research Tools



Meetings

Neuroscience 2013

Nov 9-13, San Diego, CA
Booth # 412

ASCB 2013

Dec 14-18, New Orleans, LA
Booth # 416

Rocky Mountain

Cytoskeleton Meeting

2014

May 23, Fort Collins, CO
Sponsor

Muscle & Molecular Motors GRC 2014

July 6-11, West Dover, VT
Sponsor

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

Distributors

www.cytoskeleton.com/distributors/

Dynein: Hard working, but team oriented

In eukaryote cells, cytoplasmic dynein complex (CDC) and kinesin-14 family members are the only two types of motors known to move cargo to the minus-end of microtubules (MTs)¹. This function is critical for cell metabolism, structure, and movement. The majority of these processes utilize CDC because of its high fidelity and steady processive movement. CDC's wide range of roles is impressive and the roles are further defined by ancillary proteins which select cargo or an intracellular location. Examples of the diversity of these processes are: a) transporting cargo into neuronal dendrites², b) mitotic chromosome movement³, c) a pre-caspase pathway to apoptosis⁴, d) mitochondrial movement and fission⁵, and e) amyloid precursor protein vesicle transport⁶. Likewise, ancillary proteins for these functions are varied and include Presenilin, LIS1, NUDEL, NuMA, p150 Glued in dynactin, Miro, Milton, BimL, and BimEL, to name but a few.

Interestingly, the structure of CDC's motor domain is much like a torsion gear design⁷ (Figure 1), which might represent its actual mechanism. Each dynein molecule is composed of a six domain ring motor head attached to a 20 nm amino terminal coiled-coil tail which binds the cargo and a 9 nm coiled-coil neck with a small MT binding domain at its end. The six domain ring binds up to four ATPs, with one domain hydrolyzing ATP in the power cycle and the other three poised to act as regulators or switches. Domains 5 and 6 do not contain the adenine binding motif (P-loop), which therefore does not allow ATP to bind. The torsion gear mechanism may occur due to the six domain ring being expanded at low loads and being compressed at high loads. At higher loads, the CDC may lie down on the MT surface and help create a catch-bond which stops movement in the plus-end direction.

There are several methods utilized by motors to move along MTs; these are single head one step, double head step over step, stutter step, and linear diffusion⁸⁻¹⁰. The most well-known movement is the step over step method utilized by double head motor proteins like kinesin heavy chain (KHC, kinesin-1 sub-family) and cytoplasmic myosin. These motors move without speed or gear control and more frequently fall off MTs compared to CDC under similar loads. In contrast, CDC is a different case; for example, *in vitro*, it moves in steps

as large as 24-36 nm, which is equivalent to the length of 3 to 4 tubulin heterodimers¹¹. This is surprising because CDC exists as a dimer with a spacer distance (motor to MT) of 9 nm each, but the motor domains are not bound closely to each other as they are in myosin and kinesin, and there is additional spacer length from the cargo binding spacer which is 15 nm long (Figure 1). *In vitro*, there is not much resistance placed on the motor in the form of cargo dragging through a cellular matrix. *In vivo*, the CDC movement is very different, with step length reduced to 8 nm when under load, which corresponds to the length of a tubulin heterodimer¹². Thus, CDC can change its step size and maintain MT contact under different loads.

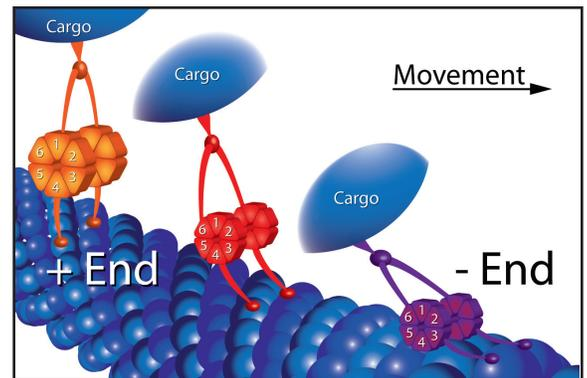


Figure 1 legend: Schematic representation of proposed torsion gear mechanism and microtubule (MT) catch-bonding. The six member ring of CDC's motor domain is shown in a relaxed state in orange and during high load conditions the ring is compressed as shown by the red model. During very high load conditions, CDC's microtubule binding domain clamps down onto the MT as shown by the purple model.

In recent years, some exquisite live cell microscopic work by several groups¹¹⁻¹⁵ has revealed how dynein moves and maintains control of movement *in vivo*. The groups improved on earlier work by a) utilizing laser-captured vesicles to measure force, b) by limiting their data to only straight tracking (<5° angles) particles, c) by more accurately defining what was a stall and a run in one direction, d) measuring the width of a stalled vesicle as an indication of opposing motors



The Protein
Experts

Helping advance science,
one protein at a time.

Motor Protein PRODUCTS

Continued from Page 1

working against each other, and e) utilizing late endosome vesicles, which show more uniform motor protein components, i.e., KHC and CDC only.

When MT minus-end-directed vesicles are observed *in vivo*, they move in plus and minus-end directions and also stall; in effect the vesicles are changing from KHC to CDC driven motion and back again. This scenario is non-intuitive and we might expect only the presence of CDC. Mallik's group studied this movement in more detail, initially utilizing a laser trap to calculate the force generated on a moving vesicle¹². In the majority of cases, the authors measured a force of 5-6 piconewtons (pN) in the plus direction (KHC) and a 5-8 pN force in the minus direction (CDC). From *in vitro* measurements, it is known that these forces reflect the presence of one KHC dimer and 5 to 10 CDC complexes. The length of motion after starting a run was also measured, yielding a ratio of 5:1 (6 nm CDC/1.2 nm KHC), which is highly in favor of minus-end movement. Furthermore, the speed and step size after stall were measured and both were in favor of CDC in the minus-end direction by a factor of 1.5 to 2.0 fold. All of these aspects combined to produce overall motion in the minus-end direction. The authors summarized their findings by concluding that dynein is a tenacious team worker where each molecule contributes 1 pN in force¹², extending others' findings that dynein catch-bonds MTs and prohibits movement in the plus direction¹⁶.

In summary, dynein represents an evolutionarily unique mechanochemical-gear motor which applies strong forces by combining the effects of multiple motors. Dynein is tenacious and catch-bonds tightly to MTs to oppose KHC motor force. At Cytoskeleton, we are developing load and non-load MT stimulated CDC ATPase assays which can be used to identify inhibitors and enhancers such as that reported by Firestone et al.¹⁷. Dynein protein is also available as a custom protein preparation (see New proteins... table), contact tbservice@cytoskeleton.com for more information.

New proteins made to order...

Protein Name & Disease Relevance	Source	Application	Amount	Module #
Dynein (cytoplasmic) Neurodegeneration, vesicle, and organelle transport	Porcine brain	Microtubule stimulated ATPase assay	1 x 50 µg 1 x 1 mg	CS-DN01
MKLP2 motor domain Tumor biology	Human recomb.	Microtubule stimulated ATPase assay	1 x 50 µg	CS-MP05
KIF7 motor domain Development and basal cell carcinoma	Human recomb.	Microtubule stimulated ATPase assay	1 x 100 µg	CS-KF51
Myosin S1 fragment (cardiac) Myocardial weakness	Bovine cardiac	Actin or thin filament stimulated ATPase assay	1 x 250 µg 1 x 1 mg	CS-MYS03
Myosin S1 fragment (skeletal)	Rabbit skeletal	F-actin stimulated ATPase assay	1 x 250 µg	CS-MYS04
Myosin S1 fragment (smooth)	Chicken gizzard	F-actin stimulated ATPase assay	1 x 250 µg	CS-MYS05
Myosin S1 fragment (non-muscle)	Bovine spleen	F-actin stimulated ATPase assay	1 x 250 µg	CS-MYS06
Actin Thin Filaments Pre-assembled complex of Tropomyosin α/β, Troponins C,I,T, and F-actin	Bovine cardiac	Calcium sensitive F-actin stimulated myosin ATPase assay	1 x 1 mg	CS-TFC01
Tropomyosin / Tropomodulin complex Complex of Tropomyosin α/β, Troponins C,I,T	Bovine cardiac	Actin thin filament preparation	1 x 1 mg 5 x 1 mg	CS-TT05

References

- Paschal B.M. & Vallee R.B. 1987. Isolated flagellar outer arm dynein translocates brain microtubules *in vitro*. *Nature*. **330**, 672-674.
- Sasaki S. et al. 2000. A LIS1/NUDEL/cytoplasmic dynein heavy chain complex in the developing and adult nervous system. *Neuron*. **28**, 681-696.
- Merdes A. et al. 1996. A complex of NuMA and cytoplasmic dynein is essential for mitotic spindle assembly. *Cell*. **87**, 447-458.
- Puthalakath H. et al. 1999. The proapoptotic activity of the Bcl-2 family member Bim is regulated by interaction with dynein motor complex. *Mol. Cell*. **3**, 287-296.
- Schwarz T.L. 2013. Mitochondrial trafficking in neurons. *Cold Spring Harb. Perspect. Biol.* doi 10.1101/cshperspect.a011304.
- Gunawardena S. et al. 2013. Presenilin controls kinesin-1 and dynein function during APP-vesicle transport *in vivo*. *Hum. Mol. Genet.* **22**, 3828-2843.
- Vallee R.B. et al. 1988. Microtubule-associated protein 1C from brain is a two-headed cytosolic dynein. *Nature*. **332**, 561-563.
- Hirokawa N. 1998. Kinesin and dynein superfamily proteins and the mechanism of organelle transport. *Science*. **279**, 519-526.
- Kull F.J. & Endow S.A. 2013. Force generation by kinesin and myosin cytoskeletal motor proteins. *J. Cell Sci.* **126**, 9-19.
- Vallee R.B. et al. 2012. Multiple modes of cytoplasmic dynein regulation. *Nat. Cell Biol.* **14**, 224-230.
- DeWitt M.A. et al. 2012. Cytoplasmic dynein moves through uncoordinated stepping of AAA+ ring domains. *Science*. **335**, 221-225.
- Rai A.K. et al. 2013. Molecular adaptations allow dynein to generate large collective forces inside cells. *Cell*. **152**, 172-182.
- Soppina V. et al. 2009. Tug-of-war between dissimilar teams of microtubule motors regulates transport and fission of endosomes. *Proc. Natl. Acad. Sci. U.S.A.* **106**, 19381-19386.
- Qiu W. et al. 2012. Dynein achieves processive motion using both stochastic and coordinated stepping. *Nat. Struct. Mol. Biol.* **19**, 193-200.
- Dufrene Y.F. et al. 2011. Five challenges to bringing single-molecule force microscopy into live cells. *Nat. Methods*. **8**, 123-127.
- Kunwar A. et al. 2011. Mechanical stochastic tug-of-war models cannot explain bidirectional lipid droplet transport. *Proc. Natl. Acad. Sci. U.S.A.* **108**, 18960-18965.
- Firestone A.J. et al. 2012. Small molecule inhibitors of AAA+ ATPase motor cytoplasmic dynein. *Nature*. **484**, 125-129.

Motor Proteins and Kits

Name	Source	Purity	Cat. #	Amount
CENP-E Motor Domain Protein	<i>H. sapiens</i>	>85%	CP01-A	2 x 25 µg
Chromokinesin Motor Domain Protein	<i>H. sapiens</i>	>85%	CR01-A	2 x 25 µg
Eg5 Motor Domain Protein	<i>H. sapiens</i>	>85%	EG01-A	2 x 25 µg
Eg5 Homolog BimC Motor Dom. Protein	<i>A. nidulans</i>	>85%	BM01-A	2 x 25 µg
Eg5 Homolog BimC Motor Dom. Protein	<i>A. fumigatus</i>	>85%	EG02-A	2 x 25 µg
KIF3 Motor Domain Protein	<i>H. sapiens</i>	>85%	KC01-A	2 x 15 µg
KIF3C Motor Domain Protein	<i>H. sapiens</i>	>85%	KF01-A	2 x 25 µg
Kinesin Heavy Chain Motor Dom. Protein	<i>H. sapiens</i>	>85%	KR01-A	2 x 25 µg
MCAK Motor Domain Protein	<i>H. sapiens</i>	>85%	MK01-A	2 x 25 µg
MKLP1 Motor Domain Protein	<i>H. sapiens</i>	>85%	MP01-A	2 x 25 µg
Kinesin ELIPA Biochem Kit	<i>na</i>	<i>na</i>	BK060	96 assays
Kinesin ATPase Endpoint Assay	<i>na</i>	<i>na</i>	BK053	1000 asy
Microtubules (pre-formed)	<i>porcine</i>	>99%	MT002	4 x 500 µg
Actin Filaments (pre-formed)	<i>rabbit</i>	>99%	AKF99	1 x 1 mg