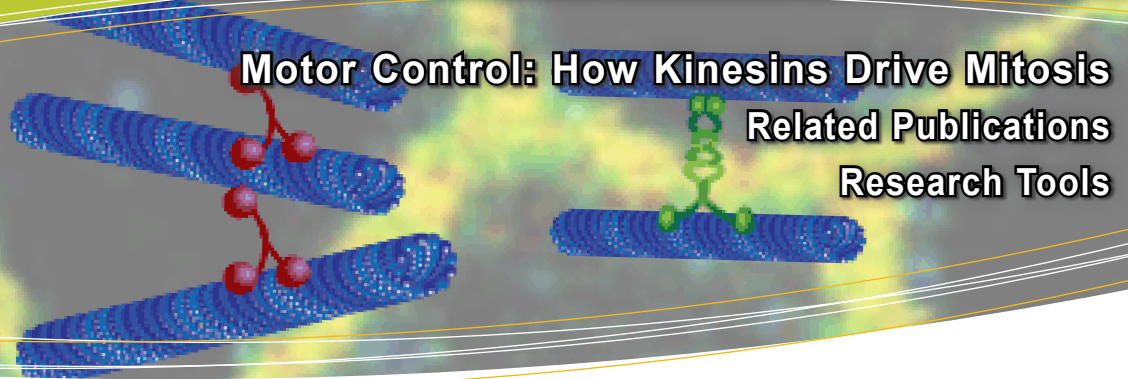




Helping advance science,
one protein at a time.

January 2023



Motor Control: How Kinesins Drive Mitosis Related Publications Research Tools

Sponsored Conferences

Gordon Research Seminar "Fibronectin, Integrins and related Molecules" in Ventura Beach, February 4-5, 2023

2023 Translation Machinery in Health and Disease Gordon Research Conference

Cytoskeleton Products

- Actin Proteins
- Activation Assays
- Antibodies
- ECM Proteins
- ELISA Kits
- G-LISA® Kits
- Live Cell Imaging
- 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

For More News

W: [cytoskeleton.com /blog/](http://cytoskeleton.com/blog/)

Motor Control: How Kinesins Drive Mitosis

One of the main components of the cytoskeleton is microtubules (MTs), which are comprised of heterodimers of α and β -tubulin proteins. Tubulin can polymerize on both ends of MTs at different rates, where polymerization occurs more rapidly on the plus-end and is slower on the minus-end. The molecular motor families, kinesins and dyneins play a role in MT length by regulating the dynamic behavior of growth and shrinkage of MTs, known as dynamic instability¹. These motors work along MTs to provide cellular functions involving organelle transport, vesicle transport, meiosis, and mitosis.

Kinesins were first discovered in 1985 when it was observed that a new class of ATPases could cycle on and off microtubules and induce organelle movement along the axons of squid and vertebrate brain^{2,3}. Since then, 45 kinesin proteins have been identified in humans, with the Kinesins superfamily (KIF) organized into 14 recognized kinesin families⁴. Most kinesins have a plus-end directed motility, but motors in the kinesin-14 subfamily have a C-terminal motor and are minus-end directed. The general architecture of kinesins consists of three domains: 1) a catalytic motor or head domain that is well conserved within each kinesin family that can hydrolyze ATP and binds microtubules, 2) an adjacent neck domain that is involved in coiled-coil interactions and can organize the motor into higher-order oligomers, 3) On the opposite end of the protein is the tail domain, which can be highly divergent within a kinesin family, and is sometimes used to interact with cargo proteins, DNA, or regulatory domains⁵⁻⁸. Below, we will look at recent findings on how kinesins affect the formation of the mitotic spindle and the organization of the chromosomes through the various stages of mitosis⁹.

Kinesins Play a Role in Mitosis

The process of mitosis occurs in somatic cells, where a copy of each chromosome of the parent cell is separated into two daughter cells during cell division. There are 5 phases of mitosis: prophase, prometaphase, metaphase, anaphase, and telophase.

Prophase begins when the duplicated centrosomes move to the opposite sides of the nucleus. Here, the MTs overlap and are orientated antiparallel to each other. The KIF11 motor (aka Eg5), a member of the kinesin-5 subfamily, can crosslink to the MTs and promotes MT sliding outwards to help push the centrosomes apart (Figure 1)^{10,11}. Dynein also assists in moving centrosomes by tethering the MTs to the cell cortex and nuclear envelope.

During prometaphase, the goal is to create bipolar mitotic spindles, form interactions between the chromosome and

MTs, and then promote chromosomal congression of the chromosomes to the spindle equator. The self-organization of the spindle requires KIF11, KIF15, and KIFC1 opposing motor actions to increase the separation of centrosomes and maintain spindle length¹⁰. Once the nuclear envelope breaks down, this allows the MT plus-ends of the spindle access to the macromolecular protein complex known as the kinetochore. Any plus-end MTs that attach to the kinetochore are known as kinetochore-microtubules (kMTs). The collection of all kMTs, typically 10-30 kMTs, within an individual kinetochore makes up a kinetochore-fiber (K-fiber)¹². At the opposite end, the minus-end directed motors of KIFC1 and dynein anchor the minus-ends of MTs to the centrosome to produce focused bipolar spindle poles¹³. Congression of the chromosomes to the spindle equator is aided by dynein and the kinesin-7 subfamily KIF10 motor (aka CENP-E), which promotes lateral sliding of the kinetochore (Figure 1)¹⁴. Kinetochore-mediated pushing and pulling forces need to be generated at the kinetochore by the kinesin motors KIF10, KIF18A, KIF2B, KIF2C, and others for proper chromosomal alignment¹⁵⁻¹⁸. Next, forces must be generated to move the chromosomes from the spindle poles toward the spindle equator. These mechanical forces are known as polar ejection forces, but further studies are needed to understand how much these forces contribute to chromosome congression^{19,20}. It is believed that the chromokinesins KIF22 (aka KID) of the kinesin-10 subfamily, and KIF4 of the kinesins-4 subfamily, interact with chromosomes during mitosis and are the motors associated with generating MT forces that force the chromosomes from the poles (Figure 1)^{20,21}. Depletion of

MOTOR ACTIVITIES IN SPINDLE ASSEMBLY

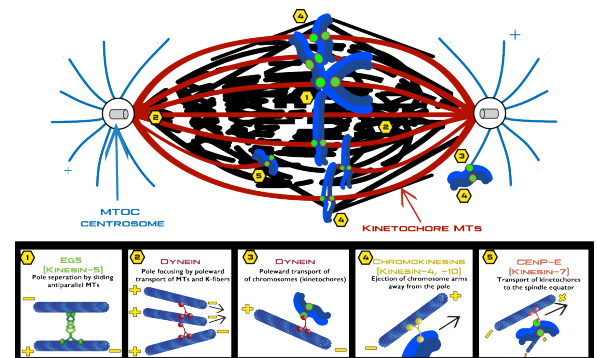


Figure Legend: Schematic representing bipolar spindle formation during mitosis. Microtubules and motor proteins play critical roles in kinetochore orientation and chromosome movement.



The Protein
Experts

Helping advance science,
one protein at a time.

Motor Protein PRODUCTS

Continued from Page 1

KIF22 via RNA interference in HeLa cells led to misaligned chromosomes and defects in spindle length, while KIF4-deficient cells resulted in disorganized chromosomes^{22,23}. These motors play a vital role in chromosomal movement through prometaphase and metaphase.

During metaphase, the sister chromatids must continue to align along the spindle equator and form the metaphase plate. The coordination to synchronously align and separate the sister chromatids is a complex process requiring many factors. The kinesin-13 family members localize to minus-ends of the microtubules at the centrosome, provoking MT depolymerization and generating poleward pulling forces on the chromosomes^{10,24}. KIF18A is also essential for proper metaphase plate alignment and timely anaphase onset²⁵. Studies also point to KIF18A and KIF10's roles in maintaining attachment to kinetochores when MTs are depolymerizing^{10,14,26}.

The final stages of mitosis occur once the chromosomes of the mitotic spindle are correctly aligned to pass the spindle assembly checkpoint. The transition into anaphase leads to the loss of cohesion between sister chromatids, allowing them to be separated and moved to each spindle pole²⁷. Less is known about kinesins and their function in anaphase, but depolymerization of the K-fiber is believed to be largely independent of motor proteins and more reliant on other MT-associated proteins and the intrinsic MT depolymerizing to complete cell division²⁸.

Targeting Kinesins for Drug Therapy

Mitotic drugs have been of great interest as a cancer treatment. For example, Paclitaxel (Taxol) is an anti-mitotic drug that targets MTs to disrupt cell division, and is used as a cancer treatment but is known to have serious side effects as it can also affect normal dividing cells. Screening for potential kinesin inhibitors that target these motor proteins that may be dysregulated in cancer presents an alternative to tubulin-based drugs. Compound screening has identified leads that inhibit kinesins, including KIF11, which have made it to Phase II drug trials²⁹. However, up to this point, most drug trials have been disappointing due to a lack of efficacy and potential side effects³⁰. Nevertheless, other motors like KIF18A appear to be exciting new drug targets, as the viability of cancer cells with chromosomal instability (CIN) depends on KIF18As ability to maintain chromosomal alignment³¹. Targeting motors vital for chromosomal alignment in cancer cells with CIN may make them more vulnerable to anti-mitotic therapies while having minimal effects on normal cells³².

Conclusion

This newsletter offers a glimpse into kinesins' complex mechanistic role in mitosis. Other proteins, microtubule-associated proteins, and types of MTs in the spindle also play significant roles in cell division. Further work is needed to understand how the motors coordinate with each other to generate the type of forces that orient the MTs and align the chromosomes for movement across the cell during the different stages of mitosis. Cytoskeleton Inc. has assisted many drug discovery efforts in this area^{33,34}. With a better understanding, kinesins may present a promising therapeutic target for cancer treatment.

Motor Proteins

Product	Amount	Cat. #
Dynein motor protein	1 x 50 µg	CS-DN01-A
Kinesin heavy chain motor domain protein GST tagged: Homo sapiens recombinant	2 x 25 µg 1 x 1 mg	KR01-A KR01-XL
KIF3C kinesin motor domain protein GST tagged: Homo sapiens recombinant	2 x 25µg	KF01-A
KIF3 kinesin motor domain protein GST tagged: Homo sapiens recombinant	2 x 25 µg	KC01-A
KIF7 Motor Domain Protein (H. Sapien)	1 x 100 µg	CS-KF51

References

- Bringmann, H. et al. A Kinesin-like Motor Inhibits Microtubule Dynamic Instability. *Science* 303, 1519–1522 (2004).
- Vale, R. D., Reese, T. S. & Sheetz, M. P. Identification of a novel force-generating protein, kinesin, involved in microtubule-based motility. *Cell* 42, 39–50 (1985).
- Brady, S. T. A novel brain ATPase with properties expected for the fast axonal transport motor. *Nature* 317, 73–75 (1985).
- Lawrence, C. J. et al. A standardized kinesin nomenclature. *J Cell Biology* 167, 19–22 (2004).
- Wang, W., Cao, L., Wang, C., Gigant, B. & Knossow, M. Kinesin, 30 years later: Recent insights from structural studies: Kinesin, 30 Years Later: Structural Insights. *Protein Sci* 24, 1047–1056 (2015).
- Cochran, J. Kinesin Motor Enzymology: Chemistry, Structure, and Physics of Nanoscale Molecular Machines. *Biophysical Rev* 7, 269–299 (2015).
- Cason, S. E. & Holzbaur, E. L. F. Selective motor activation in organelle transport along axons. *Nat Rev Mol Cell Bio* 23, 699–714 (2022).
- Miki, H., Okada, Y. & Hirokawa, N. Analysis of the kinesin superfamily: insights into structure and function. *Trends Cell Biol* 15, 467–476 (2005).
- Dumont, S. & Mitchison, T. J. Force and Length in the Mitotic Spindle. *Curr Biol* 19, R749–R761 (2009).
- Cross, R. A. & McAnish, A. Prime movers: the mechanochemistry of mitotic kinesins. *Nat Rev Mol Cell Bio* 15, 257–271 (2014).
- Tanenbaum, M. E. & Medema, R. H. Mechanisms of Centrosome Separation and Bipolar Spindle Assembly. *Dev Cell* 19, 797–806 (2010).
- Conway, W. et al. Self-organization of kinetochore-fibers in human mitotic spindles. *Elife* 11, e75458 (2022).
- Merdes, A. & Cleveland, D. W. Pathways of Spindle Pole Formation: Different Mechanisms; Conserved Components. *J Cell Biology* 138, 953–956 (1997).
- Gardner, M. K. CENP-E hangs on at dynamic microtubule ends. *Nat Cell Biol* 15, 1030–1032 (2013).
- Mayr, M. I. et al. The Human Kinesin Kif18A Is a Motile Microtubule Depolymerase Essential for Chromosome Congression. *Curr Biol* 17, 488–498 (2007).
- Stumpff, J., Dassow, G., von, Wagenbach, M., Asbury, C. & Wordeman, L. The Kinesin-8 Motor Kif18A Suppresses Kinetochore Movements to Control Mitotic Chromosome Alignment. *Dev Cell* 14, 252–262 (2008).
- Zhu, C. et al. Functional Analysis of Human Microtubule-based Motor Proteins, the Kinesins and Dyneins, in Mitosis/Cytokinesis Using RNA Interference. *Mol Biol Cell* 16, 3187–3199 (2005).
- Czechanski, A. et al. Kif18a is specifically required for mitotic progression during germ line development. *Dev Biol* 402, 253–262 (2015).
- Brouhard, G. J. & Hunt, A. J. Microtubule movements on the arms of mitotic chromosomes: Polar ejection forces quantified in vitro. *Proc National Acad Sci* 102, 13903–13908 (2005).
- Vukušić, K. & Tolić, I. M. Polar Chromosomes—Challenges of a Risky Path. *Cells* 11, 1531 (2022).
- Almeida, A. C. & Maiato, H. Chromokinesins. *Curr Biol* 28, R1131–R1135 (2018).
- Tokai-Nishizumi, N., Ohsugi, M., Suzuki, E. & Yamamoto, T. The Chromokinesin Kid Is Required for Maintenance of Proper Metaphase Spindle Size. *Mol Biol Cell* 16, 5455–5463 (2005).
- Short, B. KIF4 helps mitotic chromosomes get in shape. *J Cell Biol* 199, 715–715 (2012).
- Diez, S. Multi-talented MCAK: Microtubule depolymerizer with a strong grip. *Nat Cell Biol* 13, 738–740 (2011).
- Orr, B. & Maiato, H. No chromosome left behind: The importance of metaphase alignment for mitotic fidelity. *J Cell Biol* 218, 1086–1088 (2019).
- McIntosh, J. R., Molodtsov, M. I. & Ataullakhanov, F. I. Biophysics of mitosis. *Q Rev Biophys* 45, 147–207 (2012).
- Pandey, H., Popov, M., Goldstein-Levitin, A. & Gheber, L. Mechanisms by Which Kinesin-5 Motors Perform Their Multiple Intracellular Functions. *Int J Mol Sci* 22, 6420 (2021).
- Vukušić, K., Buđa, R. & Tolić, I. M. Force-generating mechanisms of anaphase in human cells. *J Cell Sci* 132, jcs231985 (2019).
- Garcia-Saez, I. & Skoufias, D. A. Eg5 targeting agents: From new anti-mitotic based inhibitor discovery to cancer therapy and resistance. *Biochem Pharmacol* 184, 114364 (2021).
- Yan, V. C. et al. Why Great Mitotic Inhibitors Make Poor Cancer Drugs. *Trends Cancer* 6, 924–941 (2020).
- Tamayo, N. A. et al. Targeting the Mitotic Kinesin KIF18A in Chromosomally Unstable Cancers: Hit Optimization Toward an In Vivo Chemical Probe. *J Med Chem* 65, 4972–4990 (2022).
- Marquis, C. et al. Chromosomally unstable tumor cells specifically require KIF18A for proliferation. *Nat Commun* 12, 1213 (2021).
- Perdomo, D. et al. TbKINX1B: a novel BILBO1 partner and an essential protein in bloodstream form *Trypanosoma brucei*. *Parasite* 29, 14 (2022).
- Fukai, R. et al. Design, synthesis, and evaluation of a novel prodrug, a S-trityl-L-cysteine derivative targeting kinesin spindle protein. *Eur J Med Chem* 215, 113288 (2021).

Preformed Microtubules

Product	Amount	Cat #
Microtubules (Taxol Stabilized and Lyophilized)	4 x 500 µg 1 x 10 mg	MT002-A MT002-XL

Tubulin Proteins

Product	Amount	Cat #
Microtubule associated protein rich fraction Source: porcine brain	1 x 100 µg 5 x 100 µg	MAPF-A MAPF-C
Tubulin protein (>99% pure) Source: porcine brain	1 x 1 mg 5 x 1 mg	T240-A T240-B
Acetylated Tubulin Protein Source: Porcine Brain Source: porcine brain	1 x 500 µg	TAC01