

Property of FGH, Inc.

February 24th, 2022

Project: FGH01.

Contract performed by: Cytoskeleton, Inc.



Cytoskeleton, Inc.

Helping advance science,
one protein at a time.

Contract performed by Cytoskeleton, Inc.
1830 S. Acoma St.
Denver, CO 80223, USA.

Contact information

Telephone	USA-303 - 322 - 2254
Fax	USA-303 - 322 - 2257
Customer Service	cserve@cytoskeleton.com
Technical assistance	USA-303 - 322 - 2254 tservice@cytoskeleton.com
Visit the web page	www.cytoskeleton.com

Contents

	Page
1. Executive Summary (deleted)	3
2. Introduction	4
3. Methods	5-6
4. Results	
a. Summary Table	7
b. Raw activity detail and identification of inhibiting samples	8-20
c. Screenshots of assay plates	21-34
d. Raw Data	35-36

Executive Summary

General Descriptions of the Field of Molecular Motors

Description of the Assays Used in Kinesin and Dynein Targeted Drug Discovery

Eukaryotic kinesin and dynein motor proteins orchestrate a wide range of kinetic events within a cell. They have been shown to move cargoes, such as chromosomes and vesicles, along microtubule tracks (1). They also play a major role in the organization of cytoskeletal architecture as evidenced in the establishment of the microtubule spindle during mitosis (2).

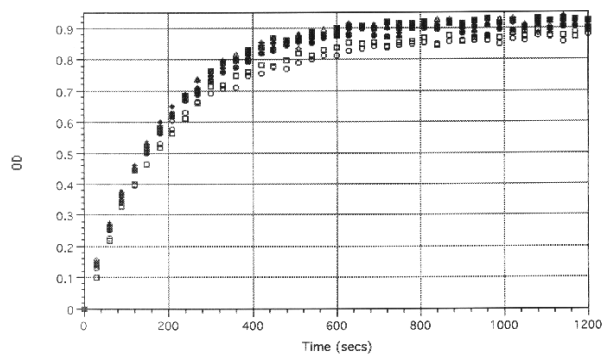
Kinesins operate by utilizing the energy of ATP hydrolysis to move along their microtubule (MT) substrates. Once a kinesin motor binds to its MT track, the ATPase rate of the motor is often enhanced several hundred to several thousand fold (3). MT activated kinesin ATPase is a major parameter in motor function and serves as a powerful method to monitor / study kinesin activity under various experimental conditions.

As part of its Cytoskeleton Motor Werks (CMW) line of research reagents, Cytoskeleton Inc. has developed two assays to measure Microtubule Stimulated Kinesin ATPase. First the kinesin end-point assay is an extremely quick and economical way to measure inorganic phosphate (Pi) generated during the microtubule activated ATPase activity of kinesin motor proteins (Cat. # BK053). Large numbers of assays can be performed simultaneously in a homogenous reaction, making the assay highly suitable for HTS applications. The assay is an adaptation of the method of Kodama, T. et al. J. Biochem. 99: 1465-1472 (1986). Essentially liberated phosphate is detected by a malachite green / molybdate complex which turns a green color on binding to phosphate.

The kinetic kinesin ATPase assay is based on the Kinesin ELIPA (Enzyme Linked Inorganic Phosphate Assay, cat. # BK060) Biochem Kit offered by Cytoskeleton Inc. The assay is an adaptation of a method originally described by Webb for the measurement of glycerol kinase plus D-glyceraldehyde ATPase activity and for actin activated myosin ATPase (4). The assay is based upon an absorbance shift (330 - 360 nm) that occurs when 2-amino-6-mercapto-7-methylpurine ribonucleoside (MESG) is catalytically converted to 2-amino-6-mercapto-7-methyl purine in the presence of inorganic phosphate (Pi). The reaction is catalysed by purine nucleoside phosphorylase (PNP). One molecule of inorganic phosphate will yield one molecule of 2-amino-6-mercapto-7-methyl purine in an essentially irreversible reaction (5). Thus, the absorbance at 360 nm is directly proportional to the amount of Pi generated in the kinesin ATPase reaction. Figure 1 shows a typical set of results from a kinesin heavy chain motor domain protein MT activated ATPase ELIPA.

Figure 1

Kinesin Heavy Chain ELIPA



Method: The reactions were conducted in a 96 well plate format (300 ul reaction volumes). Each reaction contains 94 nM kinesin heavy chain motor domain protein (cat # KR01), 0.66 uM taxol stabilized microtubules (cat # MT001), 0.2 mM MESG, 0.3 U PNP, 15 uM taxol, 15 mM PIPES pH 7, 5 mM MgCl₂, 0.6 mM ATP. Control reactions were carried out in the absence of motor protein and in the absence of MTs, these reactions gave readings of <0.02 (data not shown). Reactions were measured in a SpectraMax 250 (Molecular Devices) set in kinetic mode and 360 nm absorbance wavelength. Readings were taken at room temperature once every 30 seconds for a total reaction time of 20 minutes.

Method: Motor panel screening conditions.

Condition 1, compound preparation.

Prepare compounds by dilution into DMSO at 100x concentration. Pipette 1 μ l of DMSO solution directly into each well.

Condition 2, reaction temperature.

Reactions were run at 22°C for all plates, except plate 3, which was run at 37°C.

Condition 3, reaction's "motor mix".

Mix the following in order at RT to obtain the "motor mix".

2 ml 15 mM Pipes-NaOH pH 7.0, 10 mM MgCl₂, 20 μ M Taxol (Buffer 1).

1 ml 5x MESG (ELIPA 1 reagent, Cat. # BK051).

0.2ml (Plates 1, 5, and 7) or 1 ml (Plates 2, 3, 4, and 6) 2.5 mg/ml MTs (1 x 10mg Cat. # MT002-XL resuspended in 4 ml of Buffer 1).

0.05 ml 100 x PNP (ELIPA 2 reagent, cat. # BK051).

Plate	Motor	Cat. #	Amount per 100 μ l reaction (μ g)
1	Kinesin heavy chain	KR01	0.20
1	CenPE	CP01	0.5
1	Chromokinesin	CR01	4.0
1	Eg5	EG01	2.0
2	MKLP2	CS-MP05	0.5
3	KIF7	CS-KF517	4.0
4	KIFC3	KC01	4.0
4	KIF3C	KF01	4.0
4	MKLP	MP01	1.0
5	KIF22	CS-KF02	2.0
6	KIF3C Repeat	KF01	4.0
6	MCAK	MK01	2.0
7	Kinesin heavy chain Repeat	KR01	0.20

Condition 4, reaction initiation.

Pipette 1.5ml of 3 mM ATP into each motor mix or 3 mM GTP for KIF7.

Pipette 100 μ l into each well.

Condition 5, kinetic protocol.

Start protocol, 5-second rapid circular mixing, 41 readings, 30 seconds apart, 22°C, OD 360nm. For Plate 3, the temperature was increased to 37°C.

Notes:

High DMSO in the dilution of compounds is important to retain solubility.

Plate number	Cat #	Amount per reaction (ug)	Total ug required	Tubes of product required	Lot #
1 (high activity)	KR01	0.2	14.3	1 x 25 ug	038
1 (high activity)	CP01	0.5	12	1 x 25 ug	018
1 (high activity)	CR01	4	96	4 x 25 ug	011
1 (high activity)	EG01	2	48	2 x 25 ug	035
2 (low activity)	CS-MP05	0.5	12	1 x 50 ug	024
3 (low activity)	CS-KIF7 (needs GTP and 37C instead of ATP)	4	96	4 x 25 ug	013
4 (low activity)	KC01	4	96	4 x 25 ug	022
4 (low activity)	KF01	4	96	4 x 25 ug	013
4 (low activity)	MP01	1	24	1 x 25 ug	018
5 (high activity)	CS-KIF02	2	48	1 x 100 ug	013
6 (low activity)	KF01 Repeat	4	96	4 x 25 ug	013
6 (low activity)	MK01	2	48	2 x 25 ug	017
7 (high activity)	KR01 Repeat	0.2	14.3	1 x 25 ug	038

Table 1: Summary of results

Compound ID	Significant inhibition (>20% reduced activity compared to control)	Significant Enhancement (>20% increased activity compared to control)	<20% inhibition or enhancement detected	Notes
A1	Eg5	MKLP2	KHC, CENP-E, Chromokinesin, KIFC3, KIF3C, MCAK, MKLP1, KIF7, KIF22	
A2			KHC, CENP-E, Chromokinesin, Eg5, KIFC3, KIF3C, MCAK, MKLP1, MKLP2, KIF7, KIF22	
A3		Chromokinesin	KHC, CENP-E, Eg5, KIFC3, KIF3C, MCAK, MKLP1, MKLP2, KIF7, KIF22	
Monastrol (100 μ M)	Eg5			Used only for Eg5 control
AMPPNP (1mM)	KHC, CENP-E, Chromokinesin, KIFC3, KIF3C, MCAK, MKLP1, MKLP2, KIF7, KIF22			

Motor: Kinesin HC (Cat# KR01)

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	102.6	106.3		104.4	129.4
	3 μM	114.6	105.9		110.2	136.6
	0.3 μM	100.0	113.1		106.5	132.0
A2	10 μM	96.3	103.9		100.1	124.1
	3 μM	103.4	97.9		100.7	124.8
	0.3 μM	110.0	102.0		106.0	131.4
A3	10 μM	110.8	106.6		108.7	134.7
	3 μM	102.9	99.8		101.4	125.6
	0.3 μM	120.4	115.0		117.7	145.8
AMP-PNP	1 mM	8.7	12.2		10.5	13.0
DMSO	1%	81.1	82.5	78.5	80.7	100.0
No Motor		-1.7			-1.7	0

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Motor: Kinesin HC (Cat# KR01) Repeat

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	121.5	110.6		116.1	116.9
	3 μM	123.4	111.6		117.5	118.4
	0.3 μM	103.0	110.3		106.6	107.4
A2	10 μM	107.9	100.1		104.0	104.8
	3 μM	106.7	106.3		106.5	107.3
	0.3 μM	105.1	107.6		106.3	107.1
A3	10 μM	110.9	105.3		108.1	108.9
	3 μM	111.5	110.2		110.9	111.7
	0.3 μM	109.2	115.8		112.5	113.3
AMP-PNP	1 mM	19.2	17.0		18.1	18.3
DMSO	1%	98.6	91.6	107.5	99.3	100.0
No Motor		6.0			6.0	6.0

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Motor: CENP-E (Cat# CP01)

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	109.9	102.6		106.2	111.3
	3 μM	95.4	94.4		94.9	99.4
	0.3 μM	103.9	97.2		100.6	105.3
A2	10 μM	87.9	97.9		92.9	97.3
	3 μM	108.8	96.5		102.6	107.5
	0.3 μM	114.8	106.5		110.7	115.9
A3	10 μM	89.6	84.7		87.2	91.3
	3 μM	95.4	119.9		107.7	112.8
	0.3 μM	113.4	107.6		110.5	115.7
AMP-PNP	1 mM	15.0	13.9		14.4	15.1
DMSO	1%	97.4	94.4	94.7	95.5	100.0
No Motor		-8.2			-8.2	0

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Motor: Chromokinesin (Cat# CR01)

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	34.0	28.1		31.1	97.0
	3 μM	26.0	28.6		27.3	85.2
	0.3 μM	29.5	28.1		28.8	90.0
A2	10 μM	33.8	32.6		33.2	103.7
	3 μM	26.2	26.8		26.5	82.8
	0.3 μM	35.1	25.2		30.2	94.2
A3	10 μM	55.4	56.4		55.9	174.3
	3 μM	40.5	38.1		39.3	122.7
	0.3 μM	37.3	29.3		33.3	104.0
AMP-PNP	1 mM	15.8	20.7		18.3	57.0
DMSO	1%	27.2	33.7	35.2	32.0	100.0
No Motor		-6.2			-6.2	0

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Motor: Eg5 (Cat# EG01)

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	56.5	61.3		58.9	72.9
	3 μM	70.1	70.8		70.5	87.2
	0.3 μM	96.5	87.7		92.1	114.0
A2	10 μM	82.6	84.5		83.6	103.4
	3 μM	86.5	73.3		79.9	98.9
	0.3 μM	85.8	86.4		86.1	106.6
A3	10 μM	64.9	69.5		67.2	83.2
	3 μM	76.8	83.0		79.9	98.9
	0.3 μM	84.4	80.1		82.3	101.8
Monastrol	100 μM	4.7	5.8		5.3	6.5
DMSO	1%	81.7	83.9	76.7	80.8	100.0
No Motor		-4.7			-4.7	0

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Motor: KIF22 (Cat# KF02)

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	180.6	161.9		171.2	102.7
	3 μM	152.7	157.1		154.9	92.9
	0.3 μM	152.0	157.8		154.9	92.9
A2	10 μM	154.8	166.8		160.8	96.4
	3 μM	164.7	173.2		168.9	101.3
	0.3 μM	167.1	166.0		166.6	99.9
A3	10 μM	170.0	185.6		177.8	106.7
	3 μM	162.7	163.5		163.1	97.8
	0.3 μM	152.8	169.1		161.0	96.6
AMP-PNP	1 mM	73.0	77.8		75.4	45.2
DMSO	1%	160.1	165.8	174.2	166.7	100.0
No Motor		-2.7			-2.7	0

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Motor: KIF3C (Cat# KF01)

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	31.2	28.1		29.6	103.8
	3 μM	29.6	28.8		29.2	102.3
	0.3 μM	27.8	27.3		27.6	96.5
A2	10 μM	37.5	29.8		33.6	117.8
	3 μM	30.0	30.3		30.1	105.5
	0.3 μM	28.2	29.4		28.8	100.9
A3	10 μM	27.6	29.9		28.8	100.8
	3 μM	29.7	28.3		29.0	101.6
	0.3 μM	25.9	27.9		26.9	94.3
AMP-PNP	1 mM	33.4	22.5		28.0	98.0
DMSO	1%	28.6	28.8	28.3	28.5	100.0
No Motor		-5.4			-5.4	0

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Motor: KIF3C (Cat# KF01)

Repeat

Compound	Concentration (μ M)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μ M	24.5	24.0		24.3	100.1
	3 μ M	25.9	24.3		25.1	103.5
	0.3 μ M	24.2	23.7		24.0	98.8
A2	10 μ M	25.7	23.8		24.7	102.0
	3 μ M	24.8	24.1		24.4	100.7
	0.3 μ M	25.1	24.3		24.7	101.8
A3	10 μ M	23.7	23.6		23.6	97.4
	3 μ M	23.8	22.8		23.3	95.9
	0.3 μ M	23.7	24.3		24.0	98.9
AMP-PNP	1 mM	12.6	16.0		14.3	58.9
DMSO	1%	24.0	23.9	24.9	24.3	100.0
No Motor		-1.9			-1.9	0

Significant enhancement (>20% increased activity compared to 1% DMSO control)**Significant inhibition** (>20% reduced activity compared to 1% DMSO control)

Motor: KIFC3 (Cat# KC01)

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	26.4	26.3		26.4	106.6
	3 μM	26.4	29.8		28.1	113.6
	0.3 μM	22.7	23.4		23.0	93.0
A2	10 μM	24.0	28.2		26.1	105.4
	3 μM	26.7	29.9		28.3	114.5
	0.3 μM	23.8	26.6		25.2	101.9
A3	10 μM	25.2	24.7		24.9	100.8
	3 μM	28.3	25.5		26.9	108.7
	0.3 μM	24.8	30.1		27.5	111.0
AMP-PNP	1 mM	4.4	8.7		6.6	26.5
DMSO	1%	26.4	25.0	22.8	24.7	100.0
No Motor		-1.2			-1.2	0

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Motor: MCAK (Cat# MK01)

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	22.8	16.8		19.8	93.8
	3 μM	20.4	22.5		21.4	101.3
	0.3 μM	20.7	24.5		22.6	106.8
A2	10 μM	19.5	25.5		22.5	106.4
	3 μM	19.7	25.4		22.5	106.6
	0.3 μM	27.7	28.3		28.0	132.6
A3	10 μM	24.2	26.0		25.1	118.8
	3 μM	25.7	22.3		24.0	113.3
	0.3 μM	23.9	21.0		22.5	106.4
AMP-PNP	1 mM	12.2	9.1		10.7	50.4
DMSO	1%	21.3	18.7	23.4	21.1	100.0
No Motor		-3.4			-3.4	0

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Motor: MKLP1 (Cat# MP01)

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	130.8	114.5		122.6	95.7
	3 μM	134.9	127.8		131.4	102.6
	0.3 μM	123.0	140.4		131.7	102.8
A2	10 μM	132.9	131.7		132.3	103.2
	3 μM	126.8	131.5		129.1	100.8
	0.3 μM	132.3	126.2		129.3	100.9
A3	10 μM	119.4	116.8		118.1	92.2
	3 μM	132.9	128.9		130.9	102.2
	0.3 μM	115.6	131.8		123.7	96.6
AMP-PNP	1 mM	44.1	54.7		49.4	38.6
DMSO	1%	136.4	133.2	114.6	128.1	100.0
No Motor		3.2			3.2	2.5

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Motor: MKLP2 (Cat# MP05)

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	29.3	31.5		30.4	125.7
	3 μM	30.3	29.0		29.7	122.6
	0.3 μM	35.5	25.3		30.4	125.7
A2	10 μM	30.9	21.6		26.3	108.5
	3 μM	27.3	26.5		26.9	111.0
	0.3 μM	32.1	24.4		28.2	116.6
A3	10 μM	22.5	22.9		22.7	93.7
	3 μM	25.0	26.6		25.8	106.6
	0.3 μM	28.3	28.9		28.6	118.1
AMP-PNP	1 mM		-5.0		-5.0	0
DMSO	1%	24.4	23.7	24.4	24.2	100.0
No Motor		-9.1			-9.1	0

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Motor: KIF7 (Cat# CS-KF51)

Compound	Concentration (μM)	Run 1	Run 2	Run 3	Average (milli units per min)	% Activity
A1	10 μM	36.2	34.9		35.6	98.0
	3 μM	35.0	34.3		34.6	95.5
	0.3 μM	34.0	33.6		33.8	93.2
A2	10 μM	33.6	34.1		33.8	93.2
	3 μM	35.0	35.8		35.4	97.6
	0.3 μM	34.3	34.8		34.6	95.3
A3	10 μM	34.7	33.3		34.0	93.8
	3 μM	36.3	34.3		35.3	97.3
	0.3 μM	33.8	35.3		34.6	95.3
AMP-PNP	1 mM	16.7	35.8		26.2	72.3
DMSO	1%	36.2	35.3	37.4	36.3	100.0
No Motor		-5.2			-5.2	0

Significant enhancement (>20% increased activity compared to 1% DMSO control)

Significant inhibition (>20% reduced activity compared to 1% DMSO control)

Plate 1 Layout

		Compound A1			Compound A2			Compound A3					
		10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M			
		1	2	3	4	5	6	7	8	9	10	11	12
KHC	A	01	1 01	01	01	2 01	01	01	3 01	01	AMPPNP 01	01	KR01 02
	B	01	01	01	01	01	01	01	01	01	01	01	NoMotor 01
CENP-E	C	01	01	01	01	01	01	01	01	01	01	01	CP01 01
	D	01	01	01	01	01	01	01	01	01	01	01	NoMotor 02
Chromokinesin	E	01	01	01	01	01	01	01	01	01	01	01	CR01 01
	F	01	01	01	01	01	01	01	01	01	01	01	NoMotor 03
Eg5	G	01	01	01	01	01	01	01	01	01	Mona 01	01	Eg5 01
	H	01	01	01	01	01	01	01	01	01	01	01	NoMotor 03

Plate1

	1	2	3	4	5	6	7	8	9	10	11	12
A	102.62	114.58	99.966	96.349	103.44	110.03	110.77	102.95	120.36	8.714	81.079	78.471
B	106.27	105.89	113.08	103.86	97.907	101.99	106.61	99.789	115.00	12.228	82.517	-1.737
C	109.90	95.449	103.91	87.898	108.78	114.77	89.628	95.414	113.41	14.965	97.382	94.703
D	102.60	94.353	97.219	97.875	96.498	106.55	84.748	119.92	107.59	13.897	94.365	-8.222
E	34.050	25.965	29.498	33.838	26.224	35.116	55.350	40.500	37.324	15.833	27.197	35.171
F	28.094	28.613	28.145	32.621	26.821	25.226	56.358	38.150	29.316	20.719	33.748	-6.217
G	56.518	70.103	96.462	82.610	86.475	85.844	64.933	76.835	84.413	4.749	81.733	76.724
H	61.296	70.827	87.688	84.531	73.331	86.376	69.474	83.018	80.108	5.827	83.925	-4.745

Settings Information

▲ Kinetic

Time: 20:00

Interval: 00:00:30

Reads: 41

▲ Absorbance

Lm1 360

▲ Plate Type

96 Well Corning Half Area flat cl

Height 14.2 mm

▲ More Settings

Shake before 5s Orb., Med.

Shake between Off

Speed Read On

ReadOrder Row

Read Information

SpectraMax iD5

Plate 2 Layout

		Compound A1			Compound A2			Compound A3					
		10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M			
		1	2	3	4	5	6	7	8	9	10	11	12
MKLP2	A	01	01	01	01	01	MP05 01	01	01	01	AMP-PNP	DMSO	DMSO
	B	01	01	01	01	01	01	01	01	01	AMP-PNP	DMSO	No 01 Motor
	C												
	D												
	E												
	F												
	G												
	H												

Plate 2 MP05

	1	2	3	4	5	6	7	8	9	10	11	12
A	29.324	30.296	35.480	30.933	27.259	32.051	22.454	25.032	28.260	83.932	24.391	24.445
B	31.492	29.033	25.320	21.574	26.454	24.361	22.873	26.555	28.898	-5.049	23.743	-9.050
C	0.238	0.147	0.314	0.205	0.262	-0.229	-0.219	0.252	0.314	-0.281	0.247	0.239
D	0.423	0.364	0.189	-0.255	0.296	0.215	0.175	0.228	0.308	0.350	0.426	-0.250
E	0.200	0.315	0.219	0.431	-0.350	0.378	0.201	0.331	0.201	-0.269	0.332	0.258
F	0.231	0.417	0.244	0.207	0.229	0.260	-0.257	0.253	-0.288	0.210	-0.352	-0.260
G	0.279	-0.210	0.188	0.227	-0.319	0.318	-0.172	0.380	0.281	-0.181	-0.231	0.174
H	0.295	-0.400	0.328	0.346	0.200	-0.243	0.314	0.346	0.207	-0.212	-0.200	0.257

Settings Information

▲ Kinetic

Time: 20:00

Interval: 00:00:30

Reads: 41

▲ Absorbance

Lm1 360

▲ Plate Type

96 Well Corning Half Area flat cl

Height 14.2 mm

▲ More Settings

Shake before 5s Orb., Med.

Shake between Off

Speed Read On

ReadOrder Row

Read Information

SpectraMax iD5

Plate 3 Layout

KIF7

			Compound A1			Compound A2			Compound A3					
			10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M			
			1	2	3	4	5	6	7	8	9	10	11	12
A	03	03	03	03	03	03	03	03	03	03	03	AMP ₂ -PNP	DMSO	DMSO
B	03	03	03	03	03	03	03	03	03	03	03	AMP ₂ -PNP	DMSO	No 03 Motor
C														
D														
E														
F														
G														
H														

Plate 3 Kif7 R

	1	2	3	4	5	6	7	8	9	10	11	12
A	36.166	35.014	34.033	33.594	35.043	34.350	34.679	36.293	33.833	16.709	36.235	37.351
B	34.943	34.276	33.558	34.054	35.804	34.823	33.348	34.274	35.296	35.760	35.255	-5.201
C												
D												
E												
F												
G												
H												

Settings Information

▲ Kinetic

Time: 20:00

Interval: 00:00:30

Reads: 41

▲ Absorbance

Lm1 360

▲ Plate Type

96 Well Corning Half Area flat cl

Height 14.2 mm

▲ More Settings

Shake before 5s Orb., Med.

Shake between Off

Speed Read On

ReadOrder Row

Read Information

SpectraMax iD5

Plate 4 Layout

		Compound A1			Compound A2			Compound A3					
		10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M			
		1	2	3	4	5	6	7	8	9	10	11	12
KIFC3	A	02	02	02	02	02	02 KC01	02	02	02	AMP-PNP	DMSO	DMSO
	B	02	02	02	02	02	02	02	02	02	AMP-PNP	DMSO	No Motor 02
KIF3C	C	02	02	02	02	02	02 KF01	02	02	02	AMP-PNP	DMSO	DMSO
	D	02	02	02	02	02	02	02	02	02	AMP-PNP	DMSO	No Motor 02
	E										AMP-PNP	DMSO	DMSO
	F										AMP-PNP	DMSO	No Motor
MKLP1	G	02	02	02	02	02	02 MP01	02	02	02	AMP-PNP	DMSO	DMSO
	H	02	02	02	02	02	02	02	02	02	AMP-PNP	DMSO	No Motor 02

Plate 4 KC01 KF01 MK01 MP01R

	1	2	3	4	5	6	7	8	9	10	11	12
A	26.429	26.437	22.662	23.979	26.719	23.789	25.236	28.299	24.826	4.387	26.416	22.791
B	26.339	29.753	23.360	28.194	29.942	26.630	24.653	25.465	30.085	8.743	25.016	-1.238
C	31.219	29.608	27.812	37.463	29.962	28.192	27.636	29.665	25.876	33.448	28.563	28.270
D	28.062	28.776	27.297	29.795	30.261	29.379	29.933	28.329	27.947	22.502	28.795	-5.375
E	Masked	Masked	Masked	Masked	Masked	Masked	Masked	Masked	Masked	Masked	Masked	Masked
F	Masked	Masked	Masked	Masked	Masked	Masked	Masked	Masked	Masked	Masked	Masked	Masked
G	130.78	134.93	122.98	132.86	126.76	132.35	119.44	132.88	115.65	44.073	136.42	114.61
H	114.49	127.79	140.42	131.65	131.49	126.24	116.80	128.94	131.78	54.687	133.25	3.173

Settings Information



▲ Kinetic

Time: 20:00

Interval: 00:00:30

Reads: 41

▲ Absorbance

Lm1 360

▲ Plate Type

96 Well Corning Half Area flat c

Height 14.2 mm

▲ More Settings

Shake before 5s Orb., Med.

Shake between Off

Speed Read On

ReadOrder Row

Read Information

SpectraMax iD5

Plate 5 Layout

		Compound A1			Compound A2			Compound A3						
		10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M				
		1	2	3	4	5	6	7	8	9	10	11	12	
A														
B														
C														
D														
E														
F														
G	KIF22	01	01	01	01	01	01	KIF22	01	01	01	AMP-PNP	DMSO	DMSO
H		01	01	01	01	01	01	01	01	01	01	AMP-PNP	DMSO	No Motor 01

Plate 5 KIF22

	1	2	3	4	5	6	7	8	9	10	11	12
A	-0.324	-1.878	-1.147	-5.331	-4.507	-4.635	-5.123	-5.799	-3.614	0.991	-1.069	-0.490
B	-0.767	-3.284	-2.101	-2.343	-1.874	-2.072	1.845	-1.806	-1.888	-2.142	-2.906	5.132
C	0.314	-2.137	-1.255	2.009	2.154	1.437	-2.021	0.910	1.662	4.509	-1.037	-0.499
D	-0.539	-1.728	-0.992	-1.307	1.524	1.593	-1.243	-0.902	-1.007	4.819	-1.456	-0.774
E	-4.363	-3.530	-3.447	-3.783	-3.317	-3.637	-3.848	-3.398	-3.152	-2.557	-3.262	-4.149
F	-3.231	-2.853	-3.782	-2.604	-3.216	-2.776	-3.040	-2.669	-3.763	-2.464	-3.102	-3.778
G	180.61	152.75	152.00	154.76	164.66	167.12	170.00	162.69	152.84	72.987	160.06	174.21
H	161.89	157.14	157.79	166.79	173.15	166.03	185.60	163.50	169.14	77.843	165.82	-2.655

Settings Information

▲ Kinetic

Time: 20:00

Interval: 00:00:30

Reads: 41

▲ Absorbance

Lm1 360

▲ Plate Type

96 Well Corning Half Area flat cl

Height 14.2 mm

▲ More Settings

Shake before 5s Orb., Med.

Shake between Off

Speed Read On

ReadOrder Row

Read Information

SpectraMax iD5

Plate 6 Layout

			Compound A1			Compound A2			Compound A3					
			10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M			
			1	2	3	4	5	6	7	8	9	10	11	12
A														
B														
C														
D														
KIF3C	E	03	03	03	03	03	03	KF01	03	03	03	03	03	03
	F	03	03	03	03	03	03		03	03	03	03	03	03
MCAK	G	04	04	04	04	04	04	MK01	04	04	04	04	04	04
	H	04	04	04	04	04	04		04	04	04	04	04	04

Plate 6 KF01 MK01

	1	2	3	4	5	6	7	8	9	10	11	12
A												
B												
C												
D												
E	24.513	25.920	24.203	25.676	24.752	25.090	23.677	23.769	23.658	12.606	24.000	24.906
F	24.033	24.289	23.721	23.803	24.086	24.319	23.565	22.772	24.309	15.983	23.874	-1.905
G	22.818	20.380	20.702	19.489	19.656	27.723	24.235	25.652	23.948	12.160	21.276	23.419
H	16.837	22.455	24.464	25.515	25.428	28.342	25.995	22.267	21.034	9.148	18.749	-3.447

Settings Information

▲ Kinetic

Time: 20:00

Interval: 00:00:30

Reads: 41

▲ Absorbance

Lm1 360

▲ Plate Type

96 Well Corning Half Area flat

Height 14.2 mm

▲ More Settings

Shake before 5s Orb., Med.

Shake between Off

Speed Read On

ReadOrder Row

Read Information

SpectraMax iD5

Plate 7 Layout

		Compound A1			Compound A2			Compound A3					
		10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M	10 μ M	3 μ M	0.3 μ M			
		1	2	3	4	5	6	7	8	9	10	11	12
A		01	01	01	01	01	01	01	01	01	AMP-PNP	DMSO	DMSO
		01	01	01	01	01	01	01	01	01	AMP-PNP	DMSO	No Motor ₀₁
B													
C													
D													
E													
F													
G													
H													

KHC

KHC

Plate7 KHC repeat

	1	2	3	4	5	6	7	8	9	10	11	12
A	121.51	123.43	102.99	107.86	106.74	105.06	110.89	111.48	109.19	19.234	98.622	107.54
B	110.61	111.59	110.29	100.13	106.32	107.56	105.29	110.23	115.78	17.016	91.601	5.976
C	0.093	0.162	0.179	-0.264	0.150	0.150	0.198	0.167	0.152	0.236	-0.133	0.110
D	0.109	0.112	-0.119	-0.167	0.162	0.200	0.148	0.157	-0.141	0.131	0.155	0.214
E	0.186	-0.136	0.133	-0.209	0.155	0.148	0.143	0.105	0.172	0.131	0.136	-0.138
F	-0.188	0.112	-0.150	-0.167	-0.126	0.136	-0.188	0.157	-0.225	0.138	-0.117	0.095
G	-0.098	0.126	-0.086	0.162	-0.153	-0.127	0.121	-0.171	0.107	-0.117	-0.200	0.136
H	0.153	-0.198	0.241	-0.164	-0.208	-0.169	-0.176	0.143	-0.102	0.119	0.210	0.152

Settings Information

- ▲ Kinetic
 - Time: 20:00
 - Interval: 00:00:30
 - Reads: 41
- ▲ Absorbance
 - Lm1 360
- ▲ Plate Type
 - 96 Well Corning Half Area flat
 - Height 14.2 mm
- ▲ More Settings
 - Shake before 5s Orb., Med.
 - Shake between Off
 - Speed Read On
 - ReadOrder Row

Read Information

SpectraMax iD5

	A	106.8211	106.8567	91.59736	96.91841	107.9953	96.15329	101.9994	114.3823	100.3439	17.73294	106.7717	92.11676	KC01	
	B	106.4587	120.2594	94.41818	113.9561	121.022	107.6352	99.64528	102.9275	121.5989	35.33857	101.1115	-5.00522		
	C	109.3768	103.7313	97.44111	131.2546	104.972	98.77202	96.8239	103.9312	90.65663	117.1855	100.0704	99.04549	KF01	
	D	98.31766	100.8171	95.63467	104.3876	106.0222	102.9316	104.873	99.25255	97.91478	78.83627	100.8842	-18.8327		
	E														
	F														
	G	102.0962	105.3378	96.0116	103.7197	98.96472	103.3237	93.24508	103.7417	90.2848	34.40741	106.4989	89.47473	MP01	
	H	89.38323	99.76496	109.6244	102.7788	102.6523	98.55683	91.18665	100.6593	102.8787	42.69377	104.0264	2.477122		
Plate 5			Cmpd 1		Cmpd 2			Cmpd 3			Inhib	DMSO			
		10 µM	3 µM	0.3 µM	10 µM	3 µM	0.3 µM	10 µM	3 µM	0.3 µM					
	Raw Vmax (slopes)	1	2	3	4	5	6	7	8	9	10	11	12		
	A														
	B														
	C														
	D														
	E														
	F														
	G	180.6062	152.7465	151.9953	154.7564	164.6567	167.1218	170.0034	162.6872	152.8353	72.987	160.0565	174.2105	KIF22 DMSO Avg	
	H	161.8928	157.1397	157.7876	166.7894	173.1542	166.0343	185.5961	163.4955	169.1392	77.84255	165.822	-2.65543		166.6963395
	100*(Vmax/DMSO Avg)	1	2	3	4	5	6	7	8	9	10	11	12		
	A														
	B														
	C														
	D														
	E														
	F														
	G	108.3444	91.63157	91.18096	92.83731	98.77641	100.2553	101.9839	97.59494	91.68486	43.78441	96.0168	104.5077	KIF22	
	H	97.11839	94.26703	94.65569	100.0558	103.8741	99.60285	111.3378	98.07984	101.4655	46.69721	99.47549	-1.59297		
Plate 6			Cmpd 1		Cmpd 2			Cmpd 3			Inhib	DMSO			
		10 µM	3 µM	0.3 µM	10 µM	3 µM	0.3 µM	10 µM	3 µM	0.3 µM					
	Raw Vmax (slopes)	1	2	3	4	5	6	7	8	9	10	11	12		
	A														
	B														
	C														
	D														
	E	24.51281	25.92022	24.20304	25.67625	24.75238	25.08952	23.67718	23.76913	23.65837	12.60581	24.00005	24.90598	KF01 DMSO Avg	
	F	24.03298	24.28856	23.72066	23.80321	24.08576	24.31883	23.56521	22.77152	24.30899	15.98347	23.87416	-1.90462		24.26006544
	G	22.81795	20.37976	20.70179	19.48863	19.65564	27.72327	24.23508	25.65211	23.9482	12.1603	21.27628	23.41925	MK01 DMSO Avg	
	H	16.8372	22.45537	24.46421	25.51515	25.42849	28.34238	25.99499	22.26671	21.034	9.148083	18.74881	-3.44675		21.1481145
	100*(Vmax/DMSO Avg)	1	2	3	4	5	6	7	8	9	10	11	12		
	A														
	B														
	C														
	D														
	E	101.0418	106.8431	99.76492	105.8375	102.0293	103.419	97.59736	97.97637	97.51981	51.96115	98.92823	102.6625	KF01	
	F	99.06396	100.1175	97.77656	98.11684	99.28151	100.2422	97.1358	93.86422	100.2017	65.88385	98.4093	-7.85085		
	G	107.8959	96.36681	97.88951	92.15303	92.94276	131.091	114.5969	121.2974	113.2404	57.50066	100.6061	110.7392	MK01	
	H	79.61561	106.1814	115.6803	120.6497	120.24	134.0185	122.9187	105.2893	99.46038	43.2572	88.65477	-16.2981		
Plate 7			Cmpd 1		Cmpd 2			Cmpd 3			Inhib	DMSO			
		10 µM	3 µM	0.3 µM	10 µM	3 µM	0.3 µM	10 µM	3 µM	0.3 µM					
	Raw Vmax (slopes)	1	2	3	4	5	6	7	8	9	10	11	12		
	A	121.5148	123.4279	102.9913	107.8579	106.7395	105.0574	110.8931	111.4778	109.1938	19.23434	98.62177	107.5374	KHC DMSO Avg	
	B	110.6069	111.5873	110.2949	100.126	106.3202	107.5606	105.2899	110.2252	115.7846	17.01553	91.60052	5.976454		99.25323224
	C														
	D														
	E														
	F														
	G														
	H														
	100*(Vmax/DMSO Avg)	1	2	3	4	5	6	7	8	9	10	11	12		
	A	150.5967	152.9677	127.6401	133.6714	132.2852	130.2007	137.433	138.1576	135.327	23.83767	122.2248	133.2742		
	B	137.0783	138.2933	136.6916	124.089	131.7657	133.3029	130.4888	136.6052	143.4951	21.08783	113.5231	7.406789		
	C														
	D														
	E														
	F														
	G														
	H														