



The
Protein
Experts



Product Uses

- Measurement of Microtubule-activated ATPase assays
- Identification/characterization of proteins or small molecules that affect motor ATPase activity
- Identification/characterization of proteins or small molecules that affect motor/microtubule interactions.
- Characterization of proteins that have microtubule depolymerization activity

References

- 1 Hackney, D and Jiang, W. (2001) *Methods in Molecular Biology* (Humana Press) 164:65-71
2. Hunter et al. 2003. *Molecular Cell* 11:445-457.
3. Wordeman, L and Mitchison, T.J. 1995. *J. Cell Biol.* 128:95-125.

Related Products and Uses

- Kinesin ELIPA Assay Kit Cat.# BK060
- Kinesin ATPase End Point Assay Kit Cat.# BK053
- Microtubule Spin Down Assay Kit Cat.# BK029
- MCAK Microtubule Depolymerization Assay Kit Cat.# CDS69
- Microtubules Cat.# MT001

Hs MCAK Motor Domain

Cat. # MK01

Lot# 015

Store at 4°C (desiccated) or at -70°C

Material

The conserved motor domain of human MCAK was expressed in a prokaryotic system. The recombinant protein contains a GST-Tag at the amino terminal end and has a combined molecular weight of 79 kD. The protein has been determined to be biologically active in a microtubule-activated ATPase activity test (see below). The protein is supplied as a lyophilized powder.

<u>MK01 size</u>	<u>Minimum amt. per tube</u>	<u>Actual amt. per tube</u>	<u>Vol. of buffer for 5 mg/ml resuspension</u>	<u>Minimum* ATPase (Vmax) (nmol/min/mg)</u>	<u>Minimum* ATPase (Endpoint) (nmol/min/mg)</u>
MK01-A,B	25 ug	25 ug	5 ul	160	120
MK01-XL	1 mg	1 mg	200 ul	160	120

*Guaranteed minimum ATPase activity. For actual activity of this Lot, see p. 3.

Storage and Reconstitution

The lyophilized protein is stable for at least 1 year when stored at 4°C with a desiccant (humidity <10%). Alternatively, the lyophilized protein can be stored at -70°C and is stable at least 1 year. The protein should be reconstituted to 5 mg/ml with distilled water or CMW Buffer 1 (100 mM PIPES pH 7, 200 mM KCl, 2 mM MgCl₂, 1 mM DTT, 20 uM ATP). See above table for resuspension volumes. The protein can be aliquoted into experiment sized tubes and snap frozen in liquid nitrogen. When reconstituted and stored at -70°C, the protein will be stable for at least 4 months. For working concentrations the MCAK should be diluted in CMW Buffer 1. NOTE: Kinesins do not respond well to repeated freeze/thaws and for storage at -70°C the protein concentration should not be less than 5 mg/ml. Kinesin diluted below 5 mg/ml should not be re-frozen as it will lose activity.

Purity

Protein purity is estimated by scanning densitometry of a coomassie-stained SDS-PAGE gradient gel. Figure 1 shows 10 ug of MK01 protein and purity was determined to be >80%. The total protein in each tube will therefore be approximately 20% greater than the amount shown on the tube. The major contaminant at approximately 30 kD is GST protein. The microtubule-activated ATPase activity of the MCAK motor is not inhibited by this contaminant.

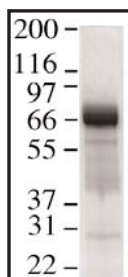


Figure 1. MCAK Motor Domain protein gel. A 10 ug sample of recombinant MCAK Motor Domain protein (GST-tagged) was separated on a 4-20% SDS-PAGE gradient gel, along with Mark12 molecular weight markers (Invitrogen). The fusion protein runs at 70 kD on the polyacrylamide gradient gel. Protein quantitation was determined using Advanced Protein assay (cat.# ADV01).

MICROTUBULE ACTIVATED ATPase ASSAY

MCAK ATPase activity was measured by monitoring real time free phosphate generation using the Kinesin ELIPA Assay Kit (cat.# BK060). The assay is based upon an absorbance shift (330 nm - 360 nm) that occurs when 2-amino-6-mercapto-7-methylpurine ribonucleoside (MESG) is catalytically converted to 2-amino-6-mercapto-7-methylpurine in the presence of inorganic phosphate (Pi). One molecule of Pi will yield one molecule of 2-amino-6-mercapto-7-methylpurine in an essentially irreversible reaction. Hence, the absorbance at 360 nm is directly proportional to the amount of Pi generated in the kinesin ATPase reaction. Under the conditions outlined below, the Vmax for MCAK microtubule-activated ATPase activity for this Lot was 195 nmoles ATP generated per minute per mg of MK01 (Figure 2). The ATPase rate for this Lot using a 10 minute endpoint assay (Kinesin ATPase End Point Assay Kit, cat.# BK053) was 170 nmoles ATP per minute per mg of MK01. Both of these values are above the guaranteed minimum.

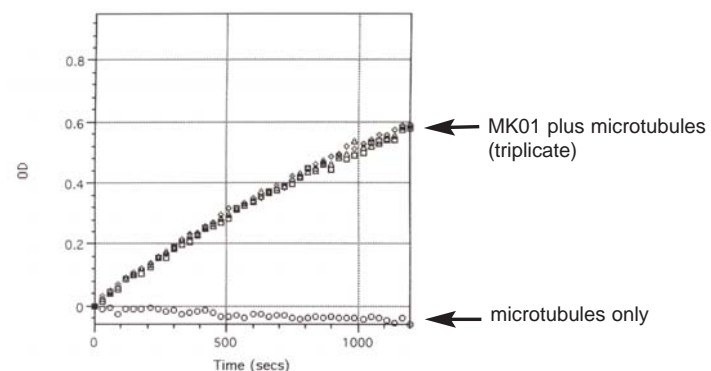


Figure 2. MCAK microtubule-activated ATPase activity using the Kinesin ELIPA Assay Kit (cat.# BK060).

Reagents

- 1) Kinesin ELIPA Assay Kit (cat.# BK060)

Equipment

- 1) Monochromatic spectrophotometer (set to 360 nm) or a filter based spectrophotometer with a 360 nm filter and bandwidth of <10 nm.

Method (ELIPA ATPase assay)

The reactions were conducted in 96 well plates (300 ul reaction volumes). Each reaction contains 10 ug of MCAK protein (MK01), 0.7 uM taxol stabilized Microtubules (cat# MT001), 0.2 mM MESG, 0.3U 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 MK01. These reactions gave readings of <0.1. Reactions were measured in a SpectraMax 250 (Molecular Devices) set in kinetic mode and 360nm absorbance wavelength. Readings were taken at room temperature once every 30 seconds for a total reaction time of 20 minutes. The nmoles of ATP generated in a given time was determined by the use of a phosphate standard curve (not shown).