

MAPKAPK5

(MAP kinase activated protein kinase 5)

Recombinant Human Active Protein Kinase

Synonyms: PRAK

Product No.: 0712-0000-1

Lot: 005

Description: Human MAPKAPK5, full length, amino acids M₁-Q₄₇₁ (as in GenBank entry NM_003668.2), activated, untagged, expressed in E.coli

Product identity: MAPKAPK5, Lot 005, was confirmed as MAPKAPK5 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW: 54,161 Da

Expression: E.coli

Activation: with MKK6-SDTD / p38-alpha

Purification: Immobilized Metal Affinity Chromatography

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 20% glycerol

Storage temperature: -80°C
Avoid repeated freeze-thaw cycles!

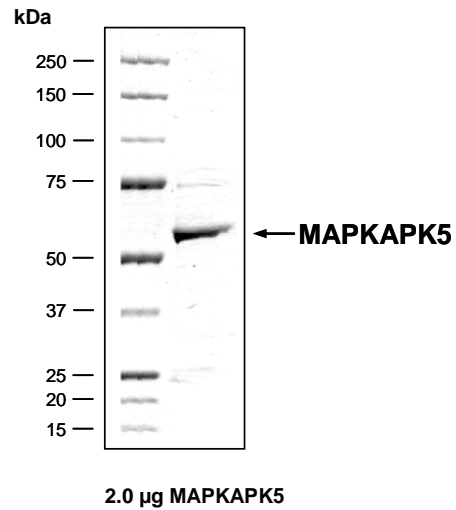
Protein concentration: 0.772 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Determination of K_m value & Specific activity:

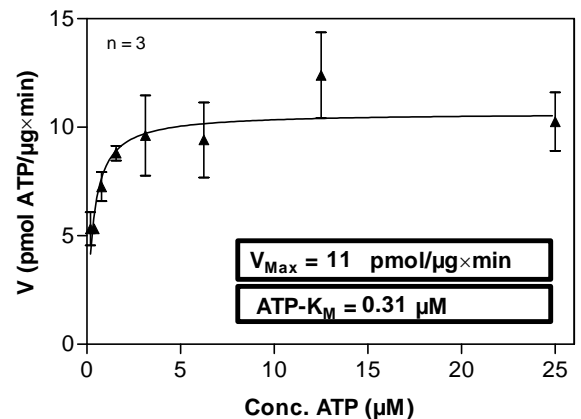
- Assay conditions:
 - 60 mM HEPES-NaOH, pH 7.5
 - 3 mM MgCl₂
 - 3 mM MnCl₂
 - 3 µM Na-orthovanadate
 - 1.2 mM DTT
 - 50 µg / ml PEG_{20,000}
 - ATP (variable)
 - Substrate: rec. RBER-CHKtide 40 µg / ml
 - MAPKAPK5: 400 ng / ml
- Filter binding assay
 - MSFC membrane (Millipore)

Specific activity: 11 pmol/µg×min

Coomassie stain:



Determination of K_m value for ATP:



This product is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.

MAPKAPK5

Product No.: 0712-0000-1

MAPKAPK5 Recombinant Fusion Protein Amino Acid Sequence							
1	GAMSEESDMD	KAIKETSILE	EYSINWTQKL	GAGISGPVRV	CVKKSTQERF	ALKILLDRPK	60
61	ARNEVRLHMM	CATHPNIVQI	IEVFANSVQF	PHSSPRARL	LIVMEMMEGG	ELFHRIHQHR	120
121	HFTEKQASQV	TKQIALALRH	CHLLNIAHRD	LKPENLLFKD	NSLDAPVKLC	DFGFAKIDQG	180
181	DLMTPQFTPY	YVAPQVLEAQ	RRHQKEKSGI	IPTSPTPYTY	NKSCDLWSLG	VIIYVMLCGY	240
241	PPFYSKHHSR	TIPKDMRRI	MTGSFEFPEE	EQISEMAK	DVVRKLLKVK	PEERLTIEGV	300
301	LDHPWLNSTE	ALDNVLPASQ	LMMDKAVVAG	IQQAHAEQLA	NMRIQDLKVS	LKPLHSVNNP	360
361	ILRKRLLGT	KPKDSVYIHD	HENGAEDSNV	ALEKLRDVIA	QCILPQAGEN	EDEKLNEVMQ	420
421	EAWKYNRECK	LLRDTLQSF	WNGRGFTDKV	DRLKLAIEIVK	QVIEEQTTSH	ESQ	480

blue:MAPKAPK5

MAPKAPK5 wt ¹ amino acid sequence							
1	MSEESDMDKA	IKETSILEEY	SINWTQKLGA	GISGPVRVCV	KKSTQERFAL	KILLDRPKAR	60
61	NEVRLHMMCA	THPNIVQIIE	VFANSVQFPH	ESSPRARLLI	VMEMMEGGEL	FHRISQHRHF	120
121	TEKQASQVTK	QIALALRHCH	LLNIAHRDLK	PENLLFKDNS	LDAPVKLCDF	GFAKIDQGDL	180
181	MTPQFTPYV	APQVLEAQR	HQKEKSGIIP	TSPTPYTYNK	SCDLWSLGI	IYVMLCGYPP	240
241	FYSKHHSRTI	PKDMRRKIMT	GSFEFPEEEW	SQISEMAKDV	VRKLLKVKPE	ERLTIEGVLD	300
301	HPWLNSTEAL	DNVLPASQLM	MDKAVVAGIQ	QAHAEQLANM	RIQDLKVSLK	PLHSVNNPIL	360
361	RKRLLGTGTP	KDSVYIHDHE	NGAEDSNVAL	EKLRDVIAQC	ILPQAGENED	EKLNEVMQEA	420
421	WKYNRECKLL	RDTLQSF	WNGRGFTDKVDR	LKLAIEIVKQV	IEEQTTSHES	Q	480

blue: MAPKAPK5 sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_003659.2

This product is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.