

ProQinase™ p38-beta

Mitogen-activated protein kinase 11

Recombinant Human Active Protein Kinase

Synonyms: MAPK11; PRKM11; SAPK2 ;SAPK2B

Product No.: 0455-0000-3

Lot: 004

Description: Human p38-beta, full length, amino acids M₁-Q₃₆₄ (as in NCBI/Protein entry NP_002742.3), activated, untagged, expressed in E. coli

Product identity: p38-beta, Lot 004, was confirmed as p38-beta by mass spectroscopy LC-ESI-MS/MS (Protagen AG, Germany)

Theoretical MW_{Fusion Protein}: 41,485 Da

Expression: E.coli

Activation: With MKK6 SDTD

Purification: Immobilized Metal Affinity Chromatography

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 3 mM βME, 20% glycerol

Storage temperature: -80°C

For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

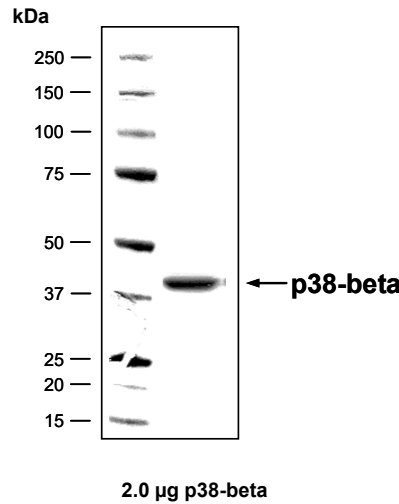
Protein concentration: 0.810 µg/µl (Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Specific activity: 701 pmol/µg×min

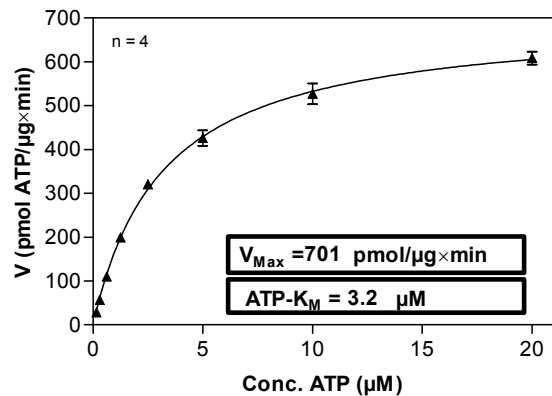
Additional assay technology: p38-beta Lot 004

was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

Coomassie stain:



Determination of K_m value for ATP:



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. ATF2, 200 µg / ml
 - p38-beta: 1.0 µg / ml
- Filter binding assay
 - MSFC membrane (Millipore)

Recombinant Proteins



ProQinase™ p38-beta

Product No.: 0455-0000-3

p38-beta Recombinant Fusion Protein Amino Acid Sequence							
1	GAMSGPRAGF	YRQELNKTVW	EVPQRLQGLR	PVGSGAYGSV	CSAYDARLRQ	KVAVKKLSRP	60
61	FQSLIHARRT	YRELRLKHL	KHENVIGLLD	VFTPATSIED	FSEVYLVTTL	MGADLNNIVK	120
121	CQALSDEHVQ	FLVYQLLRGL	KYIHSAGIIH	RDLKPSNVAV	NEDCELRILD	FGLARQADEE	180
181	MTGYVATRWY	RAPEIMLNWM	HYNQTVDIWS	VGCIMAELLQ	GKALFPGSDY	IDQLKRIMEV	240
241	VGTPSPEVLA	KISSEHARTY	IQSLPPMPQK	DLSSIFRGAN	PLAIDLLGRM	LVLDSQQRVS	300
301	AAEALAHAYF	SQYHDPEDep	EAEPYDESVE	AKERTLEEWK	ELTYQEVLSF	KPPEPKPPG	360
361	SLEIEQ						420

1-2: leagcy from TEV cleavage blue: MAPK11

p38-beta wt ¹ Amino Acid Sequence							
1	MSGPRAGFYR	QELNKTVWEV	PQRLQGLRPV	GSGAYGSVCS	AYDARLRQKV	AVKKLSRPFQ	60
61	SLIHARRTYR	ELRLKHLKH	ENVIGLLDVF	TPATSIEDFS	EVYLVTTIMG	ADLNNIVKCQ	120
121	ALSDEHVQFL	VYQLLRGLKY	IHSAGIIHRD	LKPSNVAVNE	DCELRILD	LARQADEEMT	180
181	GYVATRKYRA	PEIMLNWMHY	NQTVDIWSVG	CIMAELLQK	ALFPGSDYID	QLKRIMEVVG	240
241	TPSPEVLAKI	SSEHARTYIQ	SLPPMPQKDL	SSIFRGANPL	AIDLLGRMLV	LDSDQRVSAA	300
301	EALAHAYFSQ	YHDPEDPEEA	EPYDESVEAK	ERTLEEWKEL	TYQEVLSFKP	PEPPKPPGSL	360
361	EIEQ						420

blue: p38-beta sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_002742.3

Recombinant Proteins