

ProQinase™ ERK2

mitogen-activated protein kinase 1

Recombinant Human Active Protein Kinase

HGNC Symbol: MAPK1

Synonyms: PRKM2, PRKM1, ERK, p41mapk, MAPK2

Product No.: 0634-0000-7

Lot: 011

Description: Human ERK2, full length, amino acids M₁-S₃₆₀ (as in [NCBI/Protein](#) entry NP_002736.3), activated, untagged, expressed in E.coli

Theoretical MW_{Fusion Protein}: 41762 Da

Expression host: E.coli

Purification: GST-Affinity Chromatography

Activation: With MEK1

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

Storage temperature: -80°C

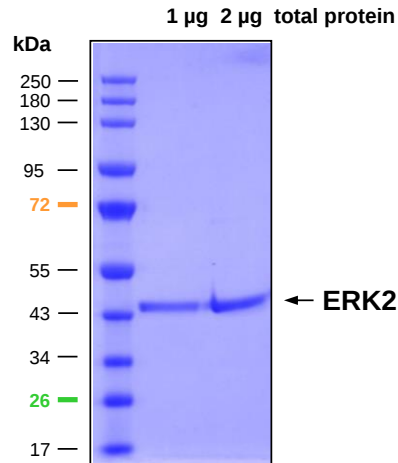
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10 µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.614 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

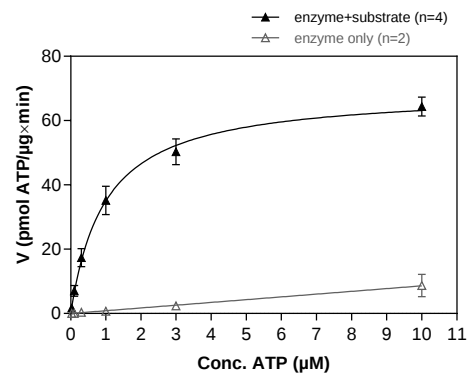
Biochemical Parameters:

Specific kinase activity (P_i transfer): 69 pmol/µg*min
ATP-K_M: 0.98 µM

ERK2 Lot 011:
Coomassie stain



ERK2 Lot 011:
Determination of V_{max} and K_M value for ATP



- 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: RBER-CHKtide 80 µg/ml
Kinase: 0.2 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

Recombinant Proteins

Sequence information

ERK2 Recombinant Fusion Protein Amino Acid Sequence							
1	GPLAMAAAAA	AGAGPEMVRG	QVFDVGPRYT	NLSYIGEGAY	GMVCSAYDNV	NKVRVAIKKI	60
61	SPFEHQTYCQ	RTLREIKILL	RFRHENIIGI	NDIIRAPTIE	QMKDVYIVQD	LMETDLYKLL	120
121	KTQHLSNDHI	CYFLYQILRG	LKYIHSANVL	HRDLKPSNLL	LNTTCDLKIC	DFGLARVADP	180
181	DHDHTGFLTE	YVATRWRYP	EIMLSKGYT	KSIDIWSVGC	ILAEMLSNRP	IFPGKHYLDQ	240
241	LNHILGILGS	PSQEDLNCII	NLKARNYLLS	LPHKNKVPWN	RLFPNADSKA	LDLLDKMLTF	300
301	NPHKRIEVEQ	ALAHPLYEQY	YDPSDEPIAE	APFKFDMELD	DLPKEKLKEL	IFEETARFQP	360
361	GYRS						420

1-4: leagcy from tag cleavage blue: ERK2

ERK2 wt ¹ Amino Acid Sequence							
1	MAAAAAGAG	PEMVRGQVFD	VGPRYTNSLY	IGEGAYGMVC	SAYDNVNKVR	VAIKKISPFE	60
61	HQTYCQRTL	EIKILLRFRH	ENIIGINDII	RAPTIEQMKD	VYIVQDLMET	DLYKLLKTQH	120
121	LSNDHICYFL	YQILRGLKYI	HSANVLHRDL	KPSNLLNTT	CDLKICDFGL	ARVADPDHHD	180
181	TGFLTEYVAT	RWYRAPEIML	NSKGYTKSID	IWSVGCILAE	MLSNRPIFPG	KHYLDQLNHI	240
241	LGILGSPSQE	DLNCIINLKA	RNYLLSLPHK	NKVPWNRLFP	NADSKALDLL	DKMLTFNPHK	300
301	RIEVEQALAH	PYLEQYYDPS	DEPIAEAPFK	FDMELDDLPK	EKLKELIFEE	TARFQPGYRS	360

blue: ERK2 sequence expressed in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/protein/NP_002736.3) accession number NP_002736.3