

ERK5

Mitogen-activated protein kinase 7

Recombinant Human Active Protein Kinase

HGNC Symbol: MAPK7

Synonyms: BMK1, ERK4, PRKM7

Product No.: 1105-0000-1

Lot: 004

Description: Human ERK5, internal fragment, amino acids L₅-Q₃₉₇ (as in [NCBI/Protein](#) entry NP_002740.2), N-terminal FLAG, C-terminal HIS₈ fusion protein, expressed in Sf9 insect cells

Product identity: ERK5 Lot 004, was confirmed as ERK5 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 50,464 Da

Expression host: Sf9 insect cells

Purification: Immobilized Metal Affinity Chromatography

Activation: This kinase was not activated by special procedures

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 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.298 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

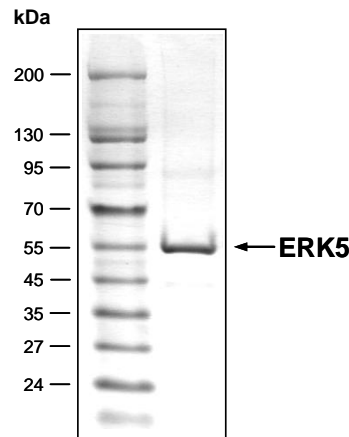
Biochemical Parameters:

Specific kinase activity (P_i transfer): 0.8 pmol/µg × min
ATP-K_M: 3.1 µM

Additional assay technology:

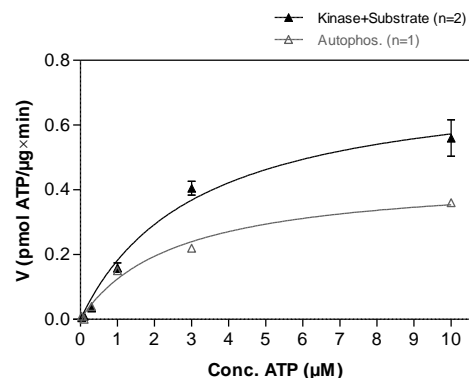
ERK5 Lot 004 was also successfully tested by ProQinase for the use with the ADP-Glo™ Kinase assay from Promega
ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

ERK5 Lot 004: Coomassie stain



2.0 µg ERK5

ERK5 Lot 004: Determination of V_{max} and 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: RBER-CHKtide, 100 µg/ml
 - Kinase: 4 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

ERK5

Product No.: 1105-0000-1

HIS-ERK5 Recombinant Fusion Protein Amino Acid Sequence							
1	MDYKDDDDKD	YKDDDDKDYK	DDDDKDYKDD	DDKSGGSLK	EEDGEDGSAE	PPGPVKAEPA	60
61	HTAASVAAKN	LALLKARSFD	VTFDVGDEYE	IIETIGNGAY	GVVSSARRRL	TGQQVAIKKI	120
121	PNAFDVVTNA	KRTLRELKIL	KHFKHDNIIA	IKDILRPTVP	YGEFKSVYV	LDLMESDLHQ	180
181	IIHSSQPLTL	EHVRYFLYQL	LRGLKYMNSA	QVIHRDLKPS	NLLVNENCEL	KIGDFGMARG	240
241	LCTSPAHEHQY	FMTEYVATRW	YRAPELMLSL	HEYTQAILDW	SVGCIFGEML	ARRQLFPGKN	300
301	YVHQLQLIMM	VLGTPSPAVI	QAVGAERVRA	YIQSLPPRQP	VPWETVYPGA	DRQALSLLGR	360
361	MLRFEPSARI	SAAAALRHPF	LAKYHDPDDE	PDCAPPFDFA	FDREALTRER	IKEAIVAEIE	420
421	DFHARREGIR	QKLLLEHHHHH	HHH				480

1-218: GST **Red**: HIS6-tag **Pink**: FLAG-tag sequence **blue**: ERK5 fragment

ERK5 wt ¹ Amino Acid Sequence							
1	MAEPLKEEDG	EDGSAEPPGP	VKAEPHAHTAA	SVAAKNLALL	KARSFDVTFD	VGDEYEIIET	60
61	IGNGAYGVVS	SARRRLTGQQ	VAIKKIPNAF	DVVTNAKRTL	RELKILKHFK	HDNIIAIKDI	120
121	LRPTVPYGEF	KSVYVVDLM	ESDLHQIHS	SQPLTLEHVR	YFLYQLLRGL	KYMNSAQVIH	180
181	RDLKPSNLLV	NENCELKIGD	FGMARGLCTS	PAEHQYFMTE	YVATRWYRAP	ELMLSLHEYT	240
241	QAIDLWSVGC	IFGEMLARRQ	LFPGKNYVHQ	LQLIMMVLGT	PSPAVIQAVG	AERVRAIQS	300
301	LPPRQVPVWE	TVYPGADRQA	LSLLGRMLRF	EPSARISAAA	ALRHPFLAKY	HDPDDEPDCA	360
361	PPFDFAFDRE	ALTRERIKEA	IVAEIEDFHA	RREGIRQQIR	FQPSLQPVAS	EPGCPDVEMP	420
421	SPWAPSGDCA	MESPPPAPP	CPGPAPDTID	LTLQPPPPVS	EPAPPKKGGA	ISDNTKAALK	480
481	AALLKSLRSR	LRDGPSAPLE	APEPRKPVTA	QERQREREK	RRRRQERAKE	REKRRQERER	540
541	KERGAGASGG	PSTDPLAGLV	LSDNDRSLE	RWTRMARPA	PALTSVPAPA	PAPTPTPTPV	600
601	QPTSPPPGPV	AQPTGPPQS	AGSTSGPVPQ	PACPPPGPAP	HPTGPPGPIP	VPAPPQIATS	660
661	TSLLAQSLV	PPPGLPGSST	PGVLPYFPPG	LPPPDAGGAP	QSSMSESPDV	NLVTQQLSKS	720
721	QVEDPLPPVF	SGTPKGSAG	YGVGFDLEEF	LNQSFDMGVA	DGPQDQGADS	ASLSASLLAD	780
781	WLEGHGMNPA	DIESLQREIQ	MDSPLLADL	PDLQDP			840

blue: ERK5 sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_002740.2

This product was manufactured at ProQinase in Freiburg, Germany, and 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.