

ERK7 (non activated)

mitogen-activated protein kinase 15

Recombinant Human Active Protein Kinase

HGNC Symbol: MAPK15

Synonyms: ERK8, MAPK15

Product No.: 1106-0000-1

Lot: 001

Description: Human ERK7, N-terminal fragment, amino acids M₁-R₃₅₄ (as in NCBI/Protein entry NP_620590.2), N-terminal 4xFLAG-tag sequence and C-terminal HIS₈-fusion protein, expressed in Sf9 insect cells

Product identity: ERK7 Lot 001, was confirmed as ERK7 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 46,156 Da

Expression: Baculovirus infected Sf9 cells

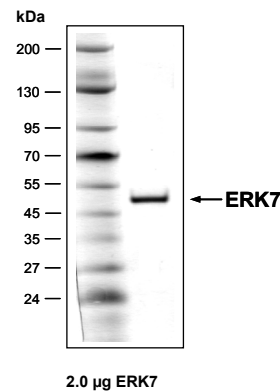
Purification: Immobilized Metal Affinity Chromatography

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

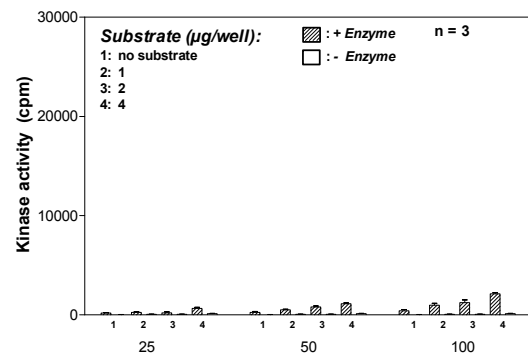
Storage temperature: -80°C
Avoid repeated freeze-thaw cycles!

Protein concentration: 0.563 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

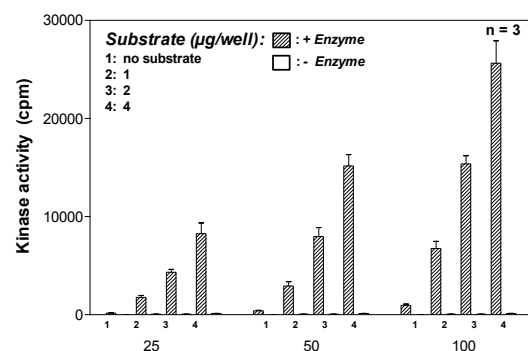
ERK7 Lot 001: Coomassie stain



Kinase activity ERK7 (inactive) vs active ERK7:



ERK7 (non activated) Lot 001 (ng/well)



Active ERK2 (ProQinase Lot 002) (ng/well)

Final assay concentrations:

- 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}
- 1 µM ATP (610,000 cpm ³³P-γ-ATP)
- Substrate (variable): RB-CTF
- Recombinant ERK7 (inactive) or active ERK7 (variable)

Assay: ³³PanQinase® Assay

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.

CoA V1.1 ERK7_Lot001_V1.doc

ERK7 (non activated)

Product No.: 1106-0000-1

ERK7 Recombinant Fusion Protein Amino Acid Sequence							
1	MDYKDDDDKD	YKDDDDKDYK	DDDDKDYKDD	DDKSGGSMC	TVVDPRIVRR	YLLRRQLGQG	60
61	AYGIVWKA	VRTGEVVAIK	KIFDAFRDKT	DAQRTFREIT	LLQEFGDHPN	IISLLDVIRA	120
121	ENDRDIYLVF	EFMDTDLNAV	IRKGGLLDV	HVRSIFYQLL	RATRFLHSGH	VVHRDQKPSN	180
181	VLLDANCTVK	LCDFGLARSL	GDLPEGPEDQ	AVTEYVATRW	YRAPEVLLSS	HRYTLGVDMW	240
241	SLGCILGEML	RGRPLFPGTS	TLHQLELILE	TIPPPSEEDL	LALGSGCRAS	VLHQLGSRPR	300
301	QTLDALPPD	TSPEALDLLR	RLLVFAPDKR	LSATQALQHP	YVQRFHCPSD	EWAREADVPR	360
361	RAHEGVQLSV	PEYRSRVYQM	ILECGSSGT	SRKLL	HHHH	HHHH	420

PINK: 4xFLAG-tag Red: HIS6-tag blue: ERK7 fragment

ERK7 wt ¹ amino acid sequence							
1	MCTVVDPRIV	RRYLLRRQLG	QGAYGIVWKA	VDRRTGEVVA	IKKIFDAFRD	KTDAQRTFRE	60
61	ITLLQEFGDH	PNIISLLDVI	RAENDRDIYL	VFEFMDTDLN	AVIRKGGLLQ	DVHVRSIFYQ	120
121	LLRATRFLHS	GHVVHRDQKP	SNVLLDANCT	VKLCDFGLAR	SLGDLPEGPE	DQAVTEYVAT	180
181	RWYRAPEVLL	SSHRYTLGVD	MWSLGCILGE	MLRGRPLFPG	TSTLHQLELI	LETIPPPSEE	240
241	DLLALGSGCR	ASVLHQLGSR	PRQTLDALLP	PDTSPEALDL	LRLLVFAPD	KRLSATQALQ	300
301	HPYVQRFHCP	SDEWAREADV	RPRAREGVQL	SVPEYRSRVY	QMILECGSS	GTSREKGPPEG	360
361	VSPSQAHLHK	PRADPQLPSR	TPVQGPRPRP	QSSPGHDPAE	HESPRAAKNV	PRQNSAPLLQ	420
421	TALLNGERP	PGAKEAPPLT	LSLVKPSGRG	AAPSLTSQAA	AQVANQALIR	GDWNRGGGVR	480
481	VASVQQVPPR	LPPEARPGRR	MFSTSALQGA	QGGARALLGG	YSQAYGTVCH	SALGHLPLLE	540
541	GHHV						600

blue: ERK7 sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_620590.2