

ProQinase™ p38-delta (non activated)

Mitogen-activated protein kinase 13

Recombinant Human non activated Protein Kinase

HGNC Symbol: MAPK13

Synonyms: PRKM13, SAPK4, p38delta

Product No.: 0941-0000-3

Lot: 004

Description: Human p38-delta, full length, amino acids M₁-L₃₆₅ (as in NCBI/Protein entry NP_002745.1), non activated, untagged, expressed in E.coli

Theoretical MW_{Protein}: 42,501 Da

Expression: E. coli

Activation: No special measures were taken to activate p38-delta Lot 004

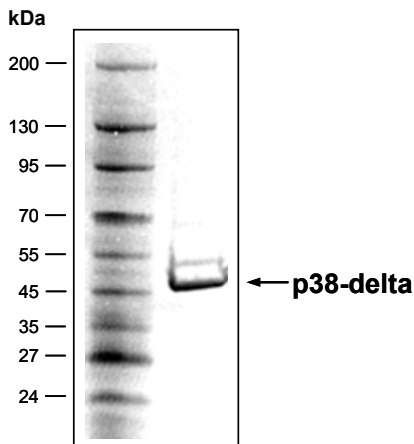
Purification: GST-Affinity Chromatography followed by cleavage of GST-tag

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!

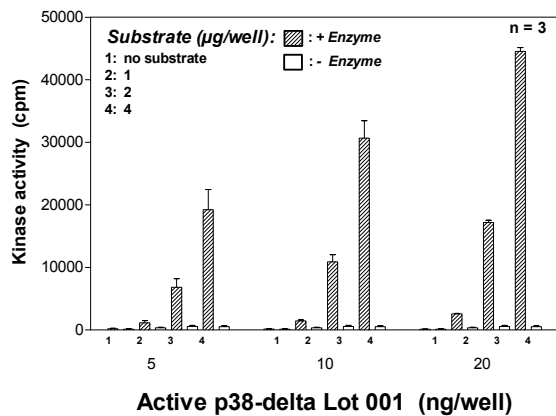
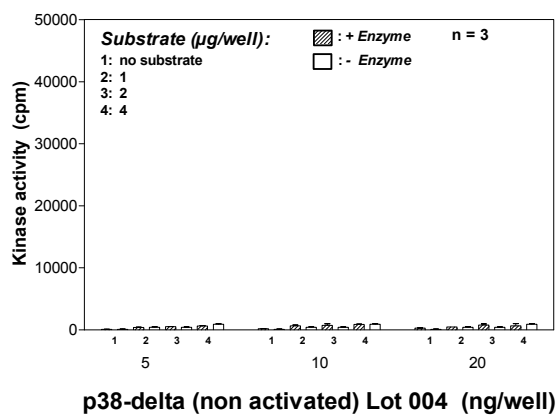
Protein concentration: 0.357 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

p38-delta Lot 004:
Coomassie stain



3.0 µg p38-delta

Kinase activity p38-delta (non activated) vs active p38-delta:



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): ATF2
- Recombinant p38-delta (non activated) or active p38-delta (variable)

Assay: ³³PanQinase® Assay

ProQinase™ p38-delta (non activated)

Product No.: 0941-0000-3

p38-delta Recombinant Fusion Protein Amino Acid Sequence								
1	GPLGS	MSLIR	KKGFYKQDVN	KTAWELPKTY	VSPTHVGSGA	YGSVCSAIDK	RSGEKVAIKK	60
61	LSRPFQSEIF	AKRAYRELLL	LKHMQHENV	GLLDVFTPAS	SLRNFYDFYL	VMPFMQTDLQ		120
121	KINGMEFSEE	KIQYLVYQML	KGLKYIHSAG	VVHRDLKPGN	LAVNEDCELK	ILDFGLARHA		180
181	DAEMTGYVVT	RWYRAPEVIL	SWMHYNQTV	IWSVGCIMAE	MLTGKTLFKG	KDYLDQLTQI		240
241	LKVTGVPGTE	FVQKLNKAA	KSYIQSLPQT	PRKDFTQLFP	RASPQAADLL	EKMLELDVVK		300
301	RLTAAQALH	PFFPEFRDPE	EETEAQQPFD	DSLEHEKLT	DEWKQHIYKE	IVNFSPIARK		360
361	DSRRRSGMKL							420

1-5: remnant of 3C cleavage **blue: p38-delta**

p38-delta wt ¹ amino acid sequence							
1	MSLIRKKGFY	KQDVNKTAW	LPKTYVSPH	VGSGAYGSVC	SAIDKRSGEK	VAIKKLSRPF	60
61	QSEIFAKRAY	RELLLLKHM	HENVIGLLDV	FTPASSLRNF	YDFYLVMPFM	QTDLQKIMGM	120
121	EFSEEKIQYL	VYQMLKGLKY	IHSAGVVHR	LKPGNLAVNE	DCELKILDFG	LARHADAEMT	180
181	GYVTRWYRA	PEVILSWMHY	NQTVDIWSVG	CIMAEMLTGK	TLFKGKDYL	QLTQILKVTG	240
241	VPGTEFVQKL	NDKAAKSYIQ	SLPQTPRKDF	TQLFPRASPQ	AADLLEKMLE	LDVDKRLTAA	300
301	QALHPFFEP	FRDPEEETEA	QQPFDDSL	EKLTVDKWKQ	HIYKEIVNFS	PIARKDSRRR	360
361	SGMKL						420

blue: p38-delta sequence expressed in fusionprotein

¹NCBI/Protein accession number NP_002745.1

Recombinant Proteins