

ProQinase™ RON

macrophage stimulating 1 receptor

Recombinant Human Active Protein Kinase

HGNC Symbol: MST1R

Synonyms: PTK8, CDw136, SEA, CD136

Product No.: 0889-0000-1

Lot: 009

Description: Human RON, internal fragment, amino acids L₁₀₅₂-E₁₃₇₀ (as in [NCBI/Protein](#) entry NP_002438.2), activated, N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: RON Lot 009, was confirmed as RON by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 64,998 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: in vitro auto activation

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 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: 0.374 µg/µl

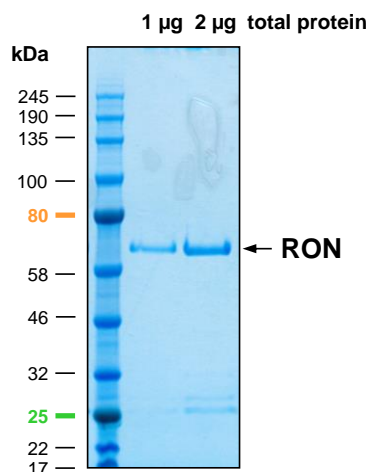
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

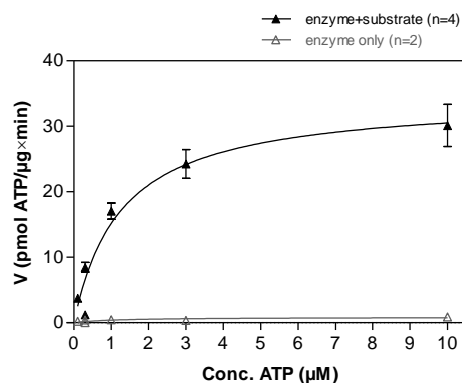
Specific kinase activity (P_i transfer): 34 pmol/µg × min

ATP-K_M: 1.3 µM

RON Lot 009: Coomassie stain



RON Lot 009: 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: TRK-C derived peptide 5 µg/ml
 - Kinase: 1.0 µg/ml
- Filter binding assay
 - MSIP membrane (Millipore)

ProQinase™ RON

Product No.: 0889-0000-1

GST-RON Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPE	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG HHHHHG	RDS LEVLFGQ	240
241	PLAMLRKESI	QLRDLDSALL	AEVKDVLIPH	ERVVTHSDRV	IGKGHFGVVY	HGEYIDQAQN	300
301	RIQCAIKSL	RITEMQQVEA	FLREGLLMRG	LNHPNVLALI	GIMLPPEGLP	HVLLPYMCHG	360
361	DLLQFIRSPQ	RNPTVKDLIS	FGLQVARGME	YLAEQKRVHR	DLAARNCLMD	ESFTVKVADF	420
421	GLARDILDRE	YYSVQQRHHA	RLPVKWMMALE	SLQTYRFTTK	SDVWSFGVLL	WELLTRGAPP	480
481	YRHIDPFDLT	HFLAQGRRLP	QPEYCPDSLY	QVMQQCWEAD	PAVRPTFRVL	VGEVEQIVSA	540
541	LLGDHYVQLP	ATYMNLPST	SHE				600

1-218: GST **Red:** HIS6-tag **Green:** 3C cleavage site **blue:** RON fragment

RON wt ¹ Amino Acid Sequence							
1	MELLPLPQS	FLLLLLLPAK	PAAGEDWQCP	RTPYAASRDF	DVKYVVPSPS	AGGLVQAMVT	60
61	YEGDRNESAV	FVAIRNRLHV	LGPDLKSVQS	LATGPAGDPG	CQCAACGPG	PHGPPGDTDT	120
121	KVLVLDPALP	ALVSCGSSLQ	GRCFLHDLEP	QGTAVHLAAP	ACLFSAHHR	PDDCPDCVAS	180
181	PLGTRVTVE	QGQASYFYVA	SSLDAVAAS	FSPRSVSIRR	LKADASGFAP	GFVALSVLPK	240
241	HLVSYIEYV	HSFHTGAFVY	FLTVPASVT	DDPSALHTRL	ARLSATEPEL	GDYRELVLDC	300
301	RFAPKRRRRG	APEGGQPYPV	LRVAHSAPVG	AQLATELSIA	EGQEVLFGVF	VTGKDGPGV	360
361	GPNSVCAFP	IDLLDTLIDE	GVERCCESPV	HPGLRRGLDF	FQSPSFCPNP	PGLEALSPT	420
421	SCRHFPLLVS	SSFSRVDLFN	GLLGPVQVTA	LYVTRLNVT	VAHMGTMGR	ILQVELVRS	480
481	NYLLYVS NFS	LGDSGQPQVR	DVSRLGDHLL	FASGDQVFQV	PIQGPGRHF	LTCGRCLRAW	540
541	HFMGCGWCGN	MCGQQKECPG	SWQQDHCPPK	LTEFHPSGP	LRGSTRLLC	GSNFYLHPSG	600
601	LVPEGTHQVT	VGQSPCRPLP	KDSSKLRPVP	RKDFVEEFEC	ELEPLGTQAV	GPTNVSLTVT	660
661	NMPPGKHFRV	DGTSVLRGFS	FMEPVLIQV	PLFGPRAGGT	CLTLEGQSL	VGTSRAVLN	720
721	GTECLLARVS	EGQLLCATPP	GATVASVPLS	LQVGAQVPG	SWTFQYREDP	VVLSISPNCG	780
781	YINSHITICG	QHLTSAWHLV	LSFHDGLRAV	ESRCERQLPE	QQLCRLPEYV	VRDPQGWVAG	840
841	NLSARGDGAA	GFTLPGFRFL	PPPHPPSANL	VPLKPEEHAI	KFEYIGLGAV	ADCVGINVT	900
901	GGESCQHEFR	GDMVVCPLPP	SLQLGQDGAP	LQVCVDGECH	ILGRVVRPGP	DGVPQSTLLG	960
961	ILLPLLLLVA	ALATALVFSY	WWRKQLVLP	PNLNDLASLD	QTAGATPLPI	LYSGSDYRSG	1020
1021	LALPAIDGLD	STTCVHGASF	SDSEDESCVP	LLRKESIQLR	DLDSALLAEV	KDVLIPHERV	1080
1081	VTHSDRVIGK	GHFGVVYHGE	YIDQAQRNIQ	CAIKSLSRIT	EMQQVEAFLR	EGLLMRGLNH	1140
1141	PNVLALIGIM	LPPEGLPHVL	LPYMCHGDL	QFIRSPQRNP	TVKDLISFGL	QVARGMEYLA	1200
1201	EQKFVHRDLA	ARNCLDES	TVKVADFGLA	RDILDREYYS	VQQRHARLP	VKWMALESIQ	1260
1261	TYRFTTKSDV	WSFGVLLWEL	LTRGAPPYRH	IDPFDLTHFL	AQGRRLPQPE	YCPDSLYQVM	1320
1321	QQWEADPAV	RPTFRVLVGE	VEQIVSALLG	DHYVQLPATY	MNLGPST	SHE	1380
1381	SPMPGNVRRP	RPLSEPPRPT					1440

blue: RON sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_002438.2