

RON

macrophage stimulating 1 receptor

Recombinant Human Active Protein Kinase

HGNC Symbol: MST1R

Synonyms: PTK8, CDw136

Product No.: 0889-0000-1

Lot: 008

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 008, 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: 0.187 µg/µl

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

Biochemical Parameters:

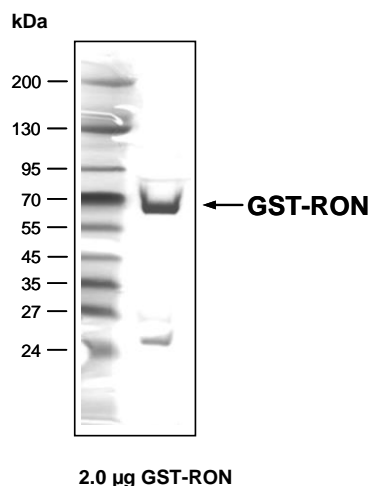
Specific kinase activity (P_i transfer): 47 pmol/µg × min

ATP-K_M: 0.8 µM

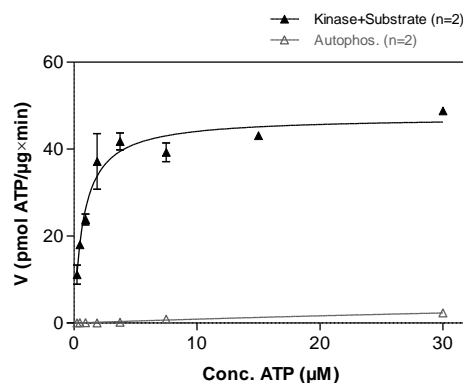
Additional assay technology:

RON Lot 008 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

RON Lot 008:
Coomassie stain



RON Lot 008:
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 80 µg/ml
Kinase: 1 µg/ml
- Filter binding assay
MSPH membrane (Millipore)

RON

Product No.: 0889-0000-1

GST-RON Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG HHHHHG	RDS LEVL FQ C	240
241	PLAMLRKESI	QLRDLSALL	AEVKDVLIPH	ERVVTHSDRV	IGKGHFGVVY	HGEYIDQAQN	300
301	RIQCAIKSL	RITEMQQVEA	FLREGLLMRG	LNHPNVLALI	GIMLPPEGLP	HVLLPYMCHG	360
361	DLLQFIRSPQ	RNPTVKDLIS	FGLQVARGME	YLAEQKRVHR	DLAARNCLMD	ESFTVKVADF	420
421	GLARDILDRE	YYSVQQRHRA	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	DVKYVVPSEFS	AGGLVQAMVT	60
61	YEGDRNESAV	FVAIRNRLHV	LGPDLKSVQS	LATGPAGDPG	QCTCAACGPG	PHGPPGDTDT	120
121	KVLVLDPALP	ALVSCGSSLQ	GRCFLHDLEP	QGTAVHLAAP	ACLFSAHHNR	PDDCPDCVAS	180
181	PLGTRVTVE	QQQASYFYVA	SSLDAVAAS	FSPRSVSIIR	LKADASGFAP	GFVALSVLPK	240
241	HLVSYSTIEYV	HSFHTGAFVY	FLTVQPASVT	DDPSALHTRL	ARLSATEPEL	GDYRELVLDC	300
301	RFAPKRRRRG	APEGGQPYPV	LRVAHSAPVG	AQLATELSIA	EGQEVLFVGF	VTGKDGGPGV	360
361	GPNSVVCAPF	IDLLDTLIDE	GVERCCESPV	HPGLRRGLDF	FQSPSFCPNP	PGLAALSPNT	420
421	SCRHFPLLVS	SSFSRVDLFN	GLLGPVQVTA	LYVTRLDNVT	VAHMGTM DGR	ILQVELVRS	480
481	NYLLVSNFS	LGDSGQPQVR	DVSRLGDHLL	FASGDQVFQV	PIQGPGRHF	LTCGRCLRAW	540
541	HFMGCGWCGN	MCGQQKECPG	SWQQDHCPPK	LTEFHPHSGP	LRGSTRLTLC	GSNFYLHPSG	600
601	LVPEGTHQVT	VGQSPCRPLP	KDSSKLRPVP	RKDFVEEFEC	ELEPLGTQAV	GPTNVSLTVT	660
661	NMPPGKHFRV	DGTSVLRGFS	FMEPVLIIVQ	PLFGPRAGGT	CLTLEGQSL	VGTSRAVLVN	720
721	GTECLLARVS	EGQLLCATPP	GATVASVPLS	LQVGGAVQVG	SWTFQYREDP	VVLSISPNCG	780
781	YINSHITICG	QHLTSAWHLV	LSFHDGLRAV	ESRCERQLPE	QQLCRLPEYV	VRDPQGWVAG	840
841	NLSARGDGAA	GFTLPGFRFL	PPPHPPSANL	VPLKPEEHAI	KFEYIGLGAV	ADCVGINVTV	900
901	GGESCQHEFR	GDMVVCPLPP	SLQLGQDQAP	LQVCVDGECH	ILGRVVRPGP	DGVPQSTLLG	960
961	ILLPLLLLVA	ALATALVFSY	WWRKQLVLP	PNLNDLASLD	QTAGATPLPI	LYSGSDYRSG	1020
1021	LALPAIDGLD	STTCVHGASF	SDSEDESCVP	LLRKESIQLR	DLDSALLAEV	KDVLIPHERV	1080
1081	VTHSDRVIGK	GHFGVVYHGE	YIDQAQNRIO	CAIKSLSRIT	EMQQVEAFLR	EGLLMRGLNH	1140
1141	PNVLALIGIM	LPPEGLPHVL	LPYMCHGDL	QFIRSPQRNP	TVKDLISFGL	QVARGMEYLA	1200
1201	EQKFVHRDLA	ARNCLDES	TVKVADFLA	RDILDREYYS	VQQRHARLP	VKWMMALESIQ	1260
1261	TYRFTTKSDV	WSFGVLLWEL	LTRGAPPYRH	IDPFDLTHFL	AQGRRLPQPE	YCPDSLYQVM	1320
1321	QQCWEADPAV	RPTFRVLVGE	VEQIVSALLG	DHYVQLPATY	MNLGPSTSHE	MNVRPEQPQF	1380
1381	SPMPGNVRRP	RPLSEPPRPT					1440

blue: RON sequence expressed in recombinant protein

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

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.