

ProQinase™ NDR1

serine/threonine kinase 38

Recombinant Human Active Protein Kinase

HGNC Symbol: STK38

Synonyms: NDR

Product No.: 1775-0000-1

Lot: 001

Description: Human NDR1, full length, amino acids M₁-K₄₆₅ (as in [NCBI/Protein](#) entry NP_009202.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: NDR1 Lot001, was confirmed as NDK1 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 82,280 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: By co-expression of MOB1B

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: 0.266 µg/µl

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

Biochemical Parameters:

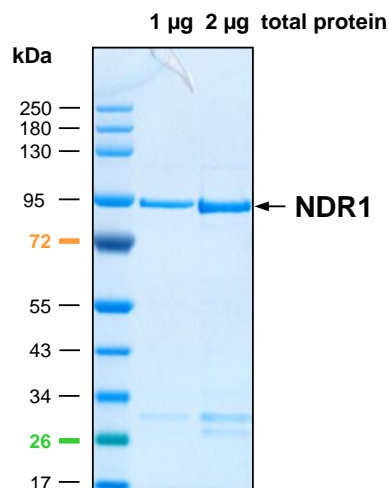
Specific kinase activity (P_i transfer): 2.2 pmol/µg × min

ATP-K_M: 4.5 µM

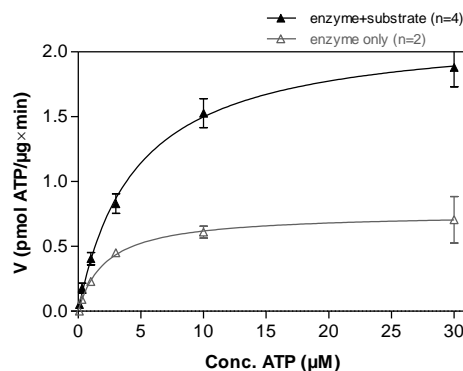
Additional assay technology:

NDR1 Lot 001 was also successfully tested by Reaction Biology 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

NDR1 Lot001: Coomassie stain



NDR1 Lot001: 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: RBER-GSK3 80 µg/ml
Kinase: 4 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

ProQinase™ NDR1

Product No.: 1775-0000-1

GST-NDR1 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFGG	240
241	PMAMTGSTPC	SSMSNHTKER	VTMTKVTLEN	FYSNLIAQHE	EREMRQKLE	KVMEEEGLK	300
301	EKRLRRSAH	ARKETEFLRL	KRTRLGLEDF	ESLKVIGRGA	FGEVRLVQKK	DTGHVYAMKI	360
361	LRKADMLEKE	QVGHIRAERD	ILVEADSLWV	VKMFYSFQDK	LNLYLIMEFL	PGDMMTLLM	420
421	KKDTLTEET	QFYIAETVLA	IDSIHQLGFI	HRDIKPDNLL	LDSKGHVKLS	DFGLCTGLKK	480
481	AHRTEFYRNL	NHSLPSDFTF	QMNNSKRKAE	TWKRNRRLA	FSTVGTDPDI	APEVFMQTGY	540
541	NKLCDWWSLG	VIMYEMLIGY	PPFCSETPQE	TYKVMNWKE	TLTFPPEVPI	SEKAKDLILR	600
601	FCCEWEHRIG	APGVVEIKSN	SFFEGVDWEH	IRERPAAISI	EIKSIDTNS	FDEFPPESDIL	660
661	KPTVATSNHP	ETDYKNKDWV	FINYTYKRFE	GLTARGAIPS	YMKAAG		720

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: NDR1

NDR1 wt ¹ Amino Acid Sequence							
1	MAMTGSTPCS	SMSNHTKERV	TMTKVTLENF	YSNLIAQHEE	REMRQKLEK	VMEEEGLKDE	60
61	EKRLRRSAHA	RKETEFLRLK	RTRLGLEDFE	SLKVIGRGAF	GEVRLVQKKD	TGHVYAMKIL	120
121	RKADMLEKEQ	VGHIRAERDI	LVEADSLWVV	KMFYSFQDKL	NLYLIMEFLP	GGDMMTLLMK	180
181	KDTLTEEETQ	FYIAETVLAI	DSIHQLGFIH	RDIKPDNLLL	DSKGHVKLS	FGLCTGLKKA	240
241	HRTEFYRNLN	HSLPSDFTFQ	NMNSKRKAET	WKRNRRLA	STVGTDPDIA	PEVFMQTGYN	300
301	KLCDWWSLGV	IMYEMLIGYP	PFCSETPQET	YKVMNWKET	LTFPPEVPIS	EKAKDLILRF	360
361	CCEWEHRIGA	PGVVEIKSN	FFEGVDWEHI	RERPAAISIE	IKSIDTNSF	DEFPPESDILK	420
421	PTVATSNHPE	TDYKNKDWVF	INYTYKRFE	LTARGAIPSY	MKAAG		480

blue: NDR1 sequence expressed in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/protein/1009202.1) accession number NP_009202.1