

## ProQinase™ WNK1

WNK lysine deficient protein kinase 1

Recombinant Human Active Protein Kinase

HGNC Symbol: WNK1

**Synonyms:** HSAN2, HSN2, hWNK1, KDP, p65, PRKWNK1, PSK

**Product No.:** 1111-0000-1

**Lot:** 002

**Description:** Human WNK1, internal fragment, amino acids L<sub>171</sub>-E<sub>529</sub> (as in [NCBI/Protein](#) entry NP\_061852.3), N-terminal 4xFLAG, C-terminal HIS<sub>8</sub> fusion protein, expressed in Sf9 insect cells

**Product identity:** WNK1 Lot 002, was confirmed as WNK1 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Fusion Protein</sub>:** 47,013 Da

**Expression host:** Sf9 insect cells

**Purification:** Immobilized Metal Affinity Chromatography

**Activation:** This kinase was not activated by special procedures

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 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!

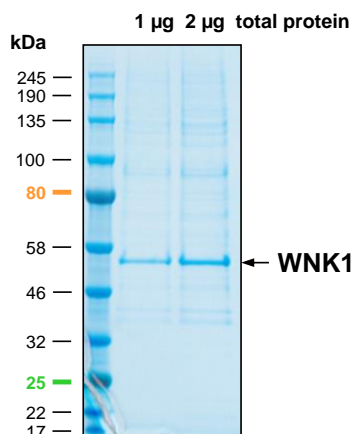
**Protein concentration:** 0.553 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**  
Specific kinase activity (P<sub>i</sub> transfer): 18 pmol/µg × min  
ATP-K<sub>M</sub>: 14 µM

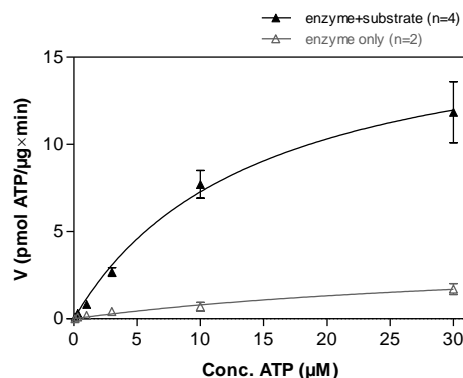
**Additional assay technology:**

WNK1 Lot 002 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

**WNK1 Lot 002:  
Coomassie stain**



**WNK1 Lot 002:  
Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP**



- Assay conditions:  
60 mM HEPES-NaOH, pH 7.5  
3 mM MgCl<sub>2</sub>  
3 mM MnCl<sub>2</sub>  
3 µM Na-orthovanadate  
1.2 mM DTT  
50 µg/ml PEG<sub>20,000</sub>  
ATP (variable)  
Substrate: Bio-RS-Peptide 80 µg/ml  
Kinase: 4 µg/ml
- Filter binding assay  
MSIP membrane (Millipore)

# ProQinase™ WNK1

Product No.: 1111-0000-1

HIS-WNK1 Recombinant Fusion Protein Amino Acid Sequence							
1	<b>MDYKDDDDDKD</b>	<b>YKDDDDDKDYK</b>	<b>DDDDDKDYKDD</b>	<b>DDK</b> SGGGS <b>LV</b>	<b>GSKEEPPPAR</b>	<b>SGSGGSSAKE</b>	60
61	<b>PQEERSQQQD</b>	<b>DIEELETKAV</b>	<b>GMSNDGRFLK</b>	<b>FDIEIGRGSF</b>	<b>KTVYKGLDTE</b>	<b>TTVEVAWCEL</b>	120
121	<b>QDRKLTKSER</b>	<b>QRFKEEAEML</b>	<b>KGLQHPNIVR</b>	<b>FYDSWESTVK</b>	<b>GKKCIVLVTE</b>	<b>LMTSGTLKTY</b>	180
181	<b>LKRKFVKMIK</b>	<b>VLRSWCRQIL</b>	<b>KGLQFLHTRT</b>	<b>PPIIHRDLKC</b>	<b>DNIFITGPTG</b>	<b>SVKIGDLGLA</b>	240
241	<b>TLKRASFAKS</b>	<b>VIGTPEFMAP</b>	<b>EMYEEKYDES</b>	<b>VDVYAFGMCM</b>	<b>LEMATSEYPY</b>	<b>SECQNAAQIY</b>	300
301	<b>RRVTSGVKPA</b>	<b>SFDKVAIPEV</b>	<b>KEIIEGCIRQ</b>	<b>NKDERYSIKD</b>	<b>LLNHAFQEE</b>	<b>TGVRVELAE</b>	360
361	<b>DDGEKIAIKL</b>	<b>WLRIEDIKKL</b>	<b>KGKYKDNEAI</b>	<b>EFSFDLE</b> KLK <b>L</b>	<b>EHHHHHHH</b>		420

**Pink:** 4xFLAG-tag **blue:** WNK1 fragment **Red:** HIS8-tag

WNK1 wt <sup>1</sup> Amino Acid Sequence							
1	MSGGAAEKQS	STPGSLFLSP	PAPAPKNGSS	SDSSVGEKLG	AAAADAVTGR	TEEYRRRRHT	60
61	MDKDSRGAAA	TTTTTEHRFF	RRSVICDSNA	TALELPLGLPL	SLPQPSIPAA	VPQSAPPEPH	120
121	REETVTATAT	SQVAQQPPAA	AAPGEQAVAG	PAPSTVPSST	SKDRPVSQPS	<b>LVGSKEEPPP</b>	180
181	<b>ARSGSGGSA</b>	<b>KEPQEERSQQ</b>	<b>QDDIEELETK</b>	<b>AVGMSNDGRF</b>	<b>LKFDIEIGRG</b>	<b>SFKTVYKGLD</b>	240
241	<b>TETTVEVAWC</b>	<b>ELQDRKLTBS</b>	<b>ERQRFKEEAE</b>	<b>MLKGLQHPNI</b>	<b>VRFYDSWEST</b>	<b>VKGKKCIVLV</b>	300
301	<b>TELMTSGTLK</b>	<b>TYLKRKFVKM</b>	<b>IKVLRSWCRQ</b>	<b>ILKGLQFLHT</b>	<b>RTPPIIHRDL</b>	<b>KCDNIFITGP</b>	360
361	<b>TGSVKIGDLG</b>	<b>LATLKRASFA</b>	<b>KSVIGTPEFM</b>	<b>APEMYEEKYD</b>	<b>ESVDVYAFGM</b>	<b>CMLEMATSEY</b>	420
421	<b>PYSECQNAAQ</b>	<b>IYRRVTSGVK</b>	<b>PASFDKVAIP</b>	<b>EVKEIIEGCI</b>	<b>RQNKDERYSI</b>	<b>KDLLNHAFQ</b>	480
481	<b>EETGVRVELA</b>	<b>EEDDGEKIAI</b>	<b>KLWLRIEDIK</b>	<b>KLKGKYKDNE</b>	<b>AIEFSFDLER</b>	<b>DVPEDVAQEM</b>	540
541	VESGYVCEGD	HKTMAKAIKD	RVSLIKRKRK	QRQLVREEQE	KKKQEESLK	QQVEQSSASQ	600
601	TGIKQLPSAS	TGIPTASTTS	ASVSTQVEPE	EPEADQHQQ	QYQQPSISVL	SDGTVDSSGQ	660
661	SSVFTESRVS	SQQTVSYGSQ	HEQAHSTGTV	PGHIPSTVQA	QSQPHGVYPP	SSVAQGSQSG	720
721	QPSSSLLTGV	SSSQPIQHPQ	QQQGIQQTAP	PQQTQVQYSL	QTSTSSEATT	AQPVSQPQAP	780
781	QVLPQVSAGK	QLPVSQPVPT	IQGEPPQIPVA	TQPSVVPVHS	GAHFLPVGQP	LPTPLLPQYP	840
841	VSQIPISTPH	VSTAQTGFSS	LPITMAAGIT	QPLTLASSA	TAAIIPGVST	VVPSQLPTLL	900
901	QPVTQLPSQV	HPQLLQPAVQ	SMGIPANLQ	AAEVPLSSGD	VLYQGFPPRL	PPQYPGDSNI	960
961	APSSNVASVC	IHSTVLSPPM	PTEVLATPGY	FPTVVQPYVE	SNLLVPMGGV	GGQVQVSQPG	1020
1021	GSLAQAPTTS	SQQAVLESTQ	GVSQVAPAEF	VAVAQTQATQ	PTTLASSVDS	AHSDVASGMS	1080
1081	DGNENVPSSS	GRHEGRTTKR	HYRKSVRSR	RHEKTSRPKL	RILNVSNGD	RVVECQLETH	1140
1141	NRKMVTFKFD	LDGDNPEDIA	TIMVNNDFIL	AIERESFVDQ	VREIIEKADE	MLSEDVSVPE	1200
1201	EGDQGLLESQ	GKDDYGFSGS	QKLEGEFKQP	IPASSMPQQI	GIPTSSLTQV	VHSAGRRFIV	1260
1261	SPVPESRLRE	SKVFPSEITD	TVAASTAQSP	GMNLSHSASS	LSLQAFSEL	RRAQMTGPN	1320
1321	TAPPNFSHTG	PTFPVVPFPL	SSIAGVPTTA	AATAPVPATS	SPPNDISTSV	IQSEVTVPTE	1380
1381	EGIAGVATST	GVVTSGGLPI	PPVSESPVLS	SVVSSITIPA	VVSISTTSPS	LQVPTSTSEI	1440
1441	VVSSTALYPS	VTVSATSASA	GGSTATPGPK	PPAVVSQQA	GSTTVGATLT	SVSTTTSFPS	1500
1501	TASQLCIQLS	SSTSTPTLAE	TVVVSASHLD	KTSHSSTTGL	AFSLSAPSSS	SSPGAGVSSY	1560
1561	ISQPGLHPL	VIPSVIASTP	ILPQAAGPTS	TPLLPQVPSI	PPLVQPVANV	PAVQQTLIHS	1620
1621	QPQALLPNP	PHTHCPEVDS	DTQPKAPGID	DIKTLEELR	SLFSEHSSSG	AQHASVSLET	1680
1681	SLVIESTVTP	GIPTTAVAPS	KLLTSTSTSC	LPPTNLPLGT	VALPVPVVT	PGQVSTPST	1740
1741	TTSGVKPGTA	PSKPPLTKAP	VLPVGTLEPA	GTLPSEQLPP	FPGPSLTQSQ	QPLEDLDAQL	1800
1801	RRTLSPEMIT	VTSAVGPVSM	AAPTAITEAG	TQPQKGVSV	KEGPVLATSS	GAGVFKMGRF	1860
1861	QVSVAADGAQ	KEGKNKSEDA	KSVHFESSTS	ESSVLSSSSP	ESTLVKPEPN	GITIPGISSD	1920
1921	VPESAHKTTA	SEAKSDTGQP	TKVGRFQVTT	TANKVGRFSV	SKTEDKITDT	KKEGPVASPP	1980
1981	FMDLEQAVLP	AVIPKKEKPE	LSEPSHLNGP	SSDPEAAFLS	RDVDDGSGSP	HSPHQLSSKS	2040
2041	LPSQNLSQL	SNFSNYSMS	SDNESDIEDE	DLKLELRLR	DKHLKEIQDL	QSRQKHEIES	2100
2101	LYTKLGVVPP	AVIIPPAAPL	SGRRRRPTKS	KGSKSSRSS	LGNKSPQLSG	NLSGQSAASV	2160
2161	LHPQQLHPP	GNIPESGQNP	LLQPLKPSPS	SDNLYSAFTS	DGAISVPSLS	APGQGTSSTN	2220
2221	TVGATVNSQA	AQAQPPAMTS	SRKGTFTDDL	HKLVDNWARD	AMNLSGRRGS	KGHMNYEGPG	2280
2281	MARKFSAPGQ	LCISMTSNLG	GSAPISAASA	TSLGHFTKSM	CPPQQYGFPA	TPFGAQWSGT	2340
2341	GGPAPQPLGQ	FQPVGTASLQ	NFNISNLQKS	ISNPPGSNLR	TT		2400

**blue:** WNK1 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein) accession number NP\_061852.3

This product was manufactured at Reaction Biology in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals.  
© European Union, 2020. Material may not be reproduced or distributed without written permission from Reaction Biology Europe GmbH