

Discovery Services

Certificate of Analysis

WNK2, active

(Recombinant enzyme expressed in Sf21 insect cells) Item # 14-678, 14-678-K, 14-678M Parent Lot # D7BN052U

The data presented in this document apply to the parent lot shown above and to all pack sizes derived from subsequent vialling runs of this parent lot. An alphabetical suffix after the parent lot number is used to denote each vialling run.

Product Description: *N*-terminal 6Histagged, recombinant, human WNK2, amino acids 166–489, expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA agarose. Purity 84% by SDS-PAGE and Coomassie blue staining. MW = 41.2kDa.

Specific Activity (Parent lot# D7BN052U): 37U/mg, where one unit of WNK2, active activity is defined as 1nmol phosphate incorporated into 0.33mg/ml myelin basic protein per minute at 30°C with a final ATP concentration of 100µM. **Formulation: 2.16mg/ml** of enzyme in 50mM Tris/HCl pH7.5, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled micro-centrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C.

FOR IN VITRO RESEARCH USE ONLY NOT FOR USE IN HUMANS OR ANIMALS

Quality Control Testing

<u>Kinase Assay</u>: 43–540ng of this lot of enzyme phosphorylated 0.33mg/ml myelin basic protein in the assay described on page two. Assay background was subtracted from the actual counts to yield the results shown below.







Eurofins Pharma Discovery Services UK Limited Gemini Crescent Dundee Technology Park DUNDEE DD2 1SW United Kingdom T +44 (0)1382 561600 F +44 (0)1382 561601 www.eurofins.com/pharmadiscovery



Discovery Services

Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

- 1. 5 x Reaction Buffer: 40mM MOPS-NaOH pH7.0, 1mM EDTA.
- 2. Myelin Basic Protein (MBP): Use at a final assay concentration of 0.33mg/ml. Make up a 3.3mg/ml stock. Use 2.5µl of stock per assay point.
- 3. WNK2, active: Dilute with 20mM MOPS-NaOH pH 7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 43–540ng per assay point.
- 4. $[\gamma^{-33}P]ATP$: 2.5 x magnesium acetate/ $[\gamma^{-33}P]ATP$ cocktail: 25mM MgAc and 0.25mM ATP to which is added $[\gamma^{-33}P]ATP$ (specific activity approximately 500 800cpm/pmol as required.)

Assay Procedure (96 well plate format):

- 1. Add 5µl of 5 x reaction buffer per assay to wells.
- 2. Add 2.5µl of myelin basic protein (MBP).
- 3. Add 2.5µl (43-540ng) WNK2, active.
- 4. Add 5µl of dH₂O.
- 5. Add 10µl of diluted $[\gamma^{-33}P]$ ATP mixture.
- 6. Incubate for 10 minutes at 30°C.
- 7. Stop the reaction by adding 5µl of 3% phosphoric acid.
- 8. Transfer a 10µl aliquot onto the appropriate area of a **P30 Filtermat**.
- 9. Wash the filtermat three times for 5 minutes with 75mM phosphoric acid.
- 10. Wash the filtermat once for 2 minutes with methanol.
- 11. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
- 12. Read in a scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all assay components plus 1µl of 30% phosphoric acid.



Certificate of Analysis

WNK2 Sequence Information

Protein	human WNK2
<u>Tags</u>	<i>N</i> -term 6His
Native sequence	G31 of recombinant sequence is equivalent to G166 of native human WNK2
Accession number	GenBank NM_006648

Recombinant WNK2 amino acid sequence:

1	MSYYHHHHHH	DYDIPTTENL	YFQGAMDPEF	GRTRRDEPEE	EEDDEDDLKA	VATSLDGRFL
61	KFDIELGRGS	FKTVYKGLDT	ETWVEVAWCE	LQDRKLTKLE	RQRFKEEAEM	LKGLQHPNIV
121	RFYDFWESSA	KGKRCIVLVT	ELMTSGTLKT	YLKRFKVMKP	KVLRSWCRQI	LKGLLFLHTR
181	TPPIIHRDLK	CDNIFITGPT	GSVKIGDLGL	ATLKRASFAK	SVIGTPEFMA	PEMYEEHYDE
241	SVDVYAFGMC	MLEMATSEYP	YSECQNAAQI	YRKVTCGIKP	ASFEKVHDPE	IKEIIGECIC
301	KNKEERYEIK	DLLSHAFFAE	DTGVRVELAE	EDHGRKSTIA	LRLWVEDPKK	LKGK

Recombinant WNK2 nucleotide sequence:

1	atgtcgtact	accatcacca	tcaccatcac	gattacgata	tcccaacgac	cgaaaacctg
61	tattttcagg	gcgccatgga	tccggaattc	gggcgcactc	gccgggacga	gcccgaagag
121	gaggaggacg	acgaggacga	cctcaaggcc	gtggccacct	ctctggacgg	ccgcttcctc
181	aagttcgaca	tcgagctggg	ccgcggttcc	ttcaagacgg	tctacaaggg	gctggacacg
241	gagacctggg	tggaggtggc	ctggtgtgag	ctgcaggacc	ggaagctcac	caagctggag
301	cggcagcggt	tcaaggaaga	ggctgagatg	ctgaaaggcc	tgcagcaccc	caacatcgtg
361	cgcttctacg	acttctggga	gtccagcgcc	aagggcaagc	ggtgcattgt	gctggtgacg
421	gagctgatga	cctcagggac	gctgaagaca	tacctgaagc	ggttcaaggt	gatgaagccc
481	aaggttctcc	gcagctggtg	ccggcagatc	ctgaagggcc	tgctgttcct	gcacacaagg
541	acgccaccca	tcatccaccg	agacctgaaa	tgtgacaata	ttttcatcac	cggaccaact
601	gggtctgtga	agattggcga	cttgggcctg	gccactctga	aaagagcgtc	atttgccaaa
661	agtgtgatag	gtactcccga	gttcatggcg	cccgagatgt	acgaggagca	ctacgatgag
721	tccgtggacg	tctatgcctt	tgggatgtgc	atgctggaga	tggccacctc	ggagtacccc
781	tactcggagt	gccagaatgc	ggcccagatc	taccgcaagg	tcacctgtgg	tatcaagccg
841	gccagctttg	agaaagtgca	cgatcctgaa	atcaaggaga	ttattgggga	gtgtatctgc
901	aaaaacaagg	aggaaaggta	cgagatcaaa	gacctgctga	gccacgcctt	cttcgcagag
961	gacacaggcg	tgagggtgga	gctcgcggag	gaggaccacg	gcaggaagtc	caccatcgcc
1021	ctgaggctct	gggtggaaga	ccccaagaaa	ctgaagggaa	agtaa	

Reviewed and approved by site quality representative.

Unless otherwise stated in our catalogue or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

© 2014 Eurofins Pharma Discovery Services UK Limited is an independent member of Eurofins Discovery Services