

Discovery Services

Certificate of Analysis

PKA, catalytic subunit, recombinant

(Recombinant enzyme expressed in *E.coli* cells) Item # 14-440, 14-440-K, 14-440M

Parent Lot # D8MN079U

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: Untagged, recombinant full-length human PKA, catalytic subunit type alpha, expressed in *E.coli* cells. Purified using phosphocellulose P11 followed by gel filtration. Purity 97.8% by SDS-PAGE and Coomassie blue staining. MW = 41kDa.

Specific Activity (Parent lot# D8MN079U): 8580U/mg, where one unit of PKA activity is defined as 1nmol phosphate incorporated into 30µM Kemptide per minute at 30°C with a final ATP concentration of 100µM. **Formulation: 0.432mg/ml** of enzyme in 30mM potassium phosphate pH7.4, 150mM KCI, 1mM EDTA, 1mM DTT, 50% glycerol. 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>: 0.1–1.8ng of this lot of enzyme phosphorylated 30µM Kemptide in the assay described on page two. Assay background was subtracted from the actual counts to yield the results shown below.



<u>MS Tryptic Fingerprint:</u> Confirmed identity as PKA with the translated sequence listed on page three.



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, pH7.0, 1mM EDTA.
- Kemptide: Use at a final concentration of 30µM. Dilute with reaction buffer for a 300µM stock. Use 2.5µl of stock per assay point.
- 3. PKA, active: Dilute with 20mM MOPS, pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2mercaptoethanol, 1mg/ml BSA. Use 0.1–1.8ng 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 **kemptide**.
- 3. Add 2.5µl (0.1–1.8ng) PKA, active.
- 4. Add 5 μ l of dH₂O.
- 5. Add 10µl of diluted [γ -³³P] 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

PKA Sequence Information

<u>Protein</u>	Human PKA, catalytic subunit type alpha				
<u>Tags</u>	untagged				
Native sequence	M1 of the recombinant protein is equivalent to M1 of human PKA				
Accession number	GenBank X07767				

Recombinant PKA amino acid sequence:

1	MGNAAAAKKG	SEQESVKEFL	AKAKEDFLKK	WESPAQNTAH	LDQFERIKTL	GTGSFGRVML
61	VKHKETGNHY	AMKILDKQKV	VKLKQIEHTL	NEKRILQAVN	FPFLVKLEFS	FKDNSNLYMV
121	MEYVPGGEMF	SHLRRIGRFS	EPHARFYAAQ	IVLTFEYLHS	LDLIYRDLKP	ENLLIDQQGY
181	IQVTDFGFAK	RVKGRTWTLC	GTPEYLAPEI	ILSKGYNKAV	DWWALGVLIY	EMAAGYPPFF
241	ADQPIQIYEK	IVSGKVRFPS	HFSSDLKDLL	RNLLQVDLTK	RFGNLKNGVN	DIKNHKWFAT
301	TDWIAIYQRK	VEAPFIPKFK	GPGDTSNFDD	YEEEEIRVSI	NEKCGKEFSE	F

Recombinant PKA nucleotide sequence:

1	atgggcaacg	ccgccgccgc	caagaagggc	agcgagcagg	agagcgtgaa	agaattctta
61	gccaaagcca	aagaagattt	tcttaaaaaa	tgggaaagtc	ccgctcagaa	cacagcccac
121	ttggatcagt	ttgaacgaat	caagaccctc	ggcacgggct	ccttcgggcg	ggtgatgctg
181	gtgaaacaca	aggagaccgg	gaaccactat	gccatgaaga	tcctcgacaa	acagaaggtg
241	gtgaaactga	aacagatcga	acacaccctg	aatgaaaagc	gcatcctgca	agctgtcaac
301	tttccgttcc	tcgtcaaact	cgagttctcc	ttcaaggaca	actcaaactt	atacatggtc
361	atggagtacg	tgcccggcgg	ggagatgttc	tcacacctac	ggcggatcgg	aaggttcagt
421	gagccccatg	cccgtttcta	cgcggcccag	atcgtcctga	cctttgagta	tctgcactcg
481	ctggatctca	tctacaggga	cctgaagccg	gagaatctgc	tcattgacca	gcagggctac
541	attcaggtga	cagacttcgg	tttcgccaag	cgcgtgaagg	gccgcacttg	gaccttgtgc
601	ggcacccctg	agtacctggc	ccctgagatt	atcctgagca	aaggctacaa	caaggccgtg
661	gactggtggg	ccctgggggt	tcttatctat	gaaatggccg	ctggctaccc	gcccttcttc
721	gcagaccagc	ccatccagat	ctatgagaag	atcgtctctg	ggaaggtgcg	cttcccttcc
781	cacttcagct	ctgacttgaa	ggacctgctg	cggaacctcc	tgcaggtaga	tctcaccaag
841	cgctttggga	acctcaagaa	tggggtcaac	gatatcaaga	accacaagtg	gtttgccaca
901	actgactgga	ttgccatcta	ccagaggaag	gtggaagctc	ccttcatacc	aaagtttaaa
961	ggccctgggg	atacgagtaa	ctttgacgac	tatgaggaag	aagaaatccg	ggtctccatc
1021	aatgagaagt	gtggcaagga	gttttctgag	ttttag		

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.