

Certificate of Analysis

PRAK, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-334, 14-334-K, 14-334M

Parent Lot # 25789U

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 6His-tagged recombinant, human, full-length PRAK expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA agarose. Purity 79% by SDS-PAGE and Coomassie blue staining. MW = 55kDa.

Specific Activity (Parent lot# 25789U): 230U/mg, where one unit of PRAK, active activity is defined as 1nmol phosphate incorporated into 30µM PRAK substrate peptide (KKLRRTLSVA) per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 4.16mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 0.2mM PMSF, 1mM benzamidine, 0.1% 2-mercaptoethanol, 270mM sucrose. Frozen solution.

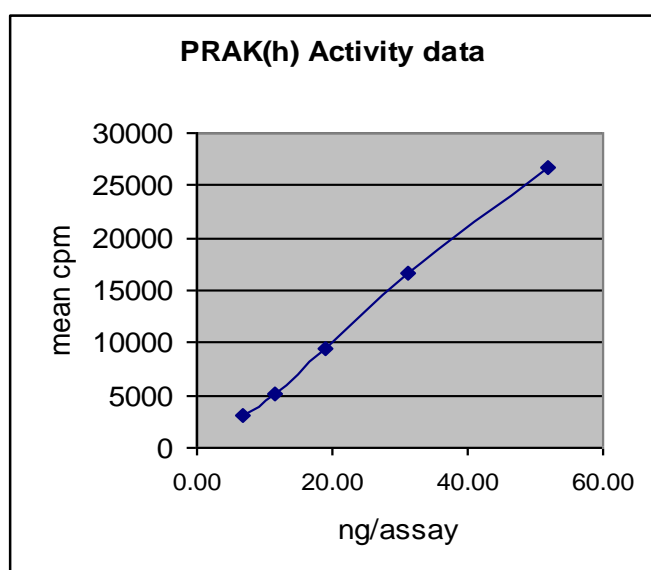
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

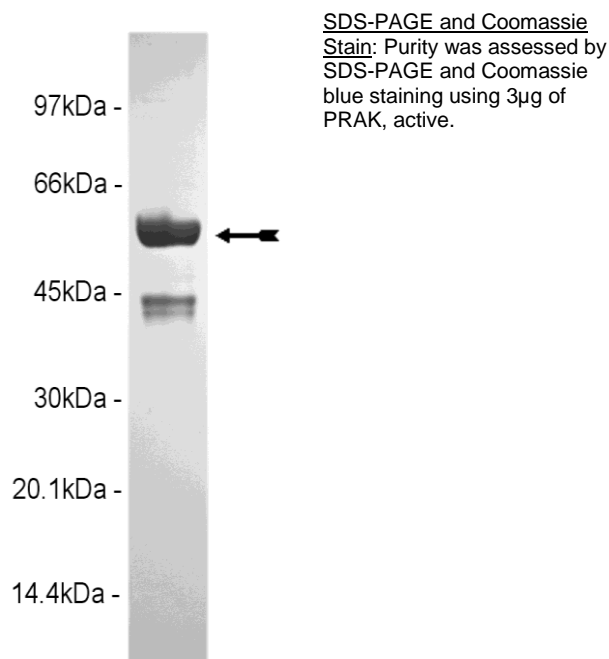
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 7–52ng of this lot of enzyme phosphorylated 30µM (KKLRRTLSVA) in the assay described on page two. Assay background was subtracted from the actual counts to yield the results shown below.



MS Tryptic Fingerprint: Confirmed product identity as PRAK with the translated sequence listed on page three.



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

- 1. 20 x Reaction Buffer:** 1M sodium- β -glycerophosphate pH7.5, 2mM EGTA.
- 2. (KKLRRTLSVA):** Use at a final assay concentration of 30 μ M. Prepare a 300 μ M stock and add 2.5 μ l of stock per assay point.
- 3. PRAK, active:** Dilute with 50mM 1M sodium- β -glycerophosphate pH7.5, 0.1mM EGTA, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 7–52ng per assay point.
- 4. [γ -³³P]ATP:** 2.5 x magnesium acetate/[γ -³³P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [γ -³³P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

Assay Procedure:

1. Add 1.25 μ l of 20 x reaction buffer to a microcentrifuge tube.
2. Add 2.5 μ l of **(KKLRRTLSVA)**, per assay point.
3. Add **2.5 μ l (7–52ng) of PRAK active**.
4. Add 8.75 μ l of dH₂O.
5. Add 10 μ l of diluted [γ -³³P] ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 μ l of 3% phosphoric acid
8. Transfer a 10 μ l 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 1 μ l of 30% phosphoric acid.

Certificate of Analysis

PRAK Sequence Information

<u>Protein</u>	human PRAK
<u>Tags</u>	N-terminal 6His
<u>Native sequence</u>	M8 of the recombinant protein is equivalent to M1 of human PRAK
<u>Accession number</u>	EMBL AF032437

Recombinant PRAK amino acid sequence:

```

1 MHHHHHHMSE ESDMDKAIKE TSILEEYSIN WTQKLGAGIS GPVRVCVKKS TQERFALKIL
61 LDRPKARNEV RLHMMCATHP NIVQIIEVFA NSVQFPHESS PRARLLIVME MMEGGELFHR
121 ISQHRHFTEK QASQVTKQIA LALRHCHLLN IAHRDLKPEN LLFKDNSLDA PVKLCDFGFA
181 KIDQGLMTP QFTPYVYVAPQ VLEAQRHQQK EKSGIIPSTP TPYTYNKSCD LWSLGVIIYV
241 MLCGYPPFYF KHHSTRIPKD MRRKIMTGSF EFPEEWSQI SEMAKDVVRK LLKVKPEERL
301 TIEGVLDPHW LNSTEALDNV LPSAQLMMDK AVVAGIQQAH AEQLANMRIQ DLKVSCLKPLH
361 SVNNPILRKR KLLGTPKPKDS VYIHDHENG AEDSNVALEKL RDVIAQCILP QAGENEDEKL
421 NEVMQEAWKY NRECKLLRDT LQSFWSWNGR FTDKVDRLKL AEIVKQVIEE QTTSHESQ
  
```

Recombinant PRAK nucleotide sequence:

```

1 atgcaccatc accatcacca tatgtcggag gagagcgaca tggacaaagc catcaaggaa
61 acttccattt tagaagaata cagtatcaat tggactcaga agctgggagc tgggaattagt
121 ggtccagtta gactctgtgt aaagaaatct actcaagaac ggtttgcgct gaaaattctt
181 cttgatcgtc caaaagctag aaatgaggta cgtctgcaca tgatgtgtgc cacacacca
241 aacatagttc agattattga agtgtttgct aacagtgtcc agtttcccca tgagtccagc
301 cctagggccc gactcttaat tgtaatggag atgatggaag ggggagagct atttcacaga
361 atcagccagc accggcactt tacagagaag caagccagcc aagtaacaaa gcagatagct
421 ttggctctgc ggcaactgtc cttgttaaac attgctgaca gagacctcaa gcctgaaaaat
481 ctgcttttta aggataactc tttggatgcc ccagtgaagt tgtgtgactt tggatttgcc
541 aagattgacc aaggtgactt gatgacacc cagttcacc cttattatgt agcaccaccag
601 gtactggagg cgcaaagaag gcatcagaag gagaaatctg gcatcatacc tacctcaccg
661 acgcccatac cttacaacaa gagctgtgac ttgtggtccc taggggtgat tatctatgtg
721 atgctgtgcg gataccctcc ttttactcc aacaccaca gccggactat ccaaaggat
781 atgcaagaa agatcatgac aggcagtttt gatttccag aggaagagtg gactcagatc
841 tcagatggg ccaaagatgt tgtgaggaag ctctgaagg tcaaaccgga ggagagactc
901 accatcgagg gagtctgga ccaccctgg ctcaattcca ccgaggccct ggataatgtg
961 ctgcttctg ctcagctgat gatggacaag gcagtgggtg caggaatcca gcaggctcac
1021 gcggaacagt tggccaacat gagaatccag gatctgaaag tcagcctcaa acccctgcac
1081 tcagtgaaca acccattct gcggaagagg aagtacttg gcaccaagcc aaaggacagt
1141 gtctatatcc acgaccatga gaatggagcc gaggattcca atgttgcctt ggaaaaactc
1201 cgagatgtga ttgctcagtg tattctccc caggctggag agaatgaaga tgagaaactg
1261 aatgaagtaa tgcaggaggc ttggaagtat aaccgggaat gcaaactcct aagagatact
1321 ctgcagagct tcagctggaa tggctgtgga ttcacagata aagtagatcg actaaaactg
1381 gcagaaattg tgaagcaggt gatagaagag caaaccacgt cccacgaatc ccaataa
  
```

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