

Certificate of Analysis

PRAK, unactive

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-335, 14-335-K, 14-335M

Parent Lot # 1917434

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 full-length, human PRAK expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA-agarose. Purity 94.4% by SDS-PAGE and Coomassie blue staining. MW = 55kDa.

Formulation: 2.924mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 50% glycerol, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Liquid at -20°C.

Specific Activity (Parent lot# 1917434): As provided, this lot demonstrated <1% of maximum activity. Activated by phosphorylation with SAPK2a, active (cat# 14-251).

Storage and Stability: On receipt of material store at -20°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.

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

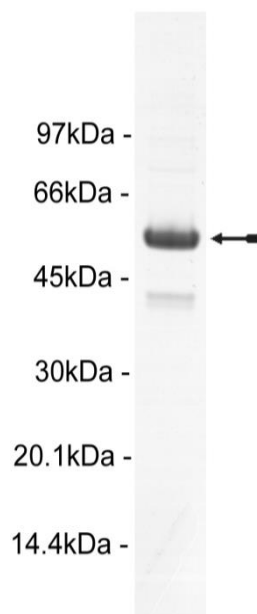
Quality Control Testing

Kinase Assay: 4µM unactive PRAK was activated using SAPK2a (cat# 14-251) at a final concentration of 2U/ml, diluted 10–20 fold and the increased activity against PRAKtide (KKLRRTLSVA) measured. The activation and assay is described on page two. Results of this assay are shown below

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

SDS-PAGE and Coomassie Stain: Purity was assessed by SDS-PAGE and Coomassie blue staining using 3µg of PRAK, unactive.

Active SAPK2a	Unactive PRAK	Mean cpm	Comments
None	5.4µg	5	Background
50mU	5.4µg	6559	Kinase activity



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

1. **5 x Activation Buffer:** 125mM Tris/HCl pH7.5, 0.5mM EGTA, 0.5% 2-mercaptoethanol.
2. **4 x PRAK assay buffer:** 1M Na- β -glycerophosphate pH7.5, 2mM EGTA.
3. **SAPK2a dilution buffer:** 50mM Tris/HCl pH7.5, 0.1mM EGTA, 0.1mM Na₃VO₄, 0.1% 2-mercaptoethanol, 1mg/ml BSA.
4. **PRAK dilution buffer:** 50mM Na- β -glycerophosphate pH7.5, 0.1mM EGTA, 0.1% 2-mercaptoethanol, 1mg/ml BSA.
5. **PRAK, unactive:** Use at a final assay concentration of 4 μ M (0.216mg/ml). Prepare a 1.08mg/ml stock and add 5 μ l of stock per assay point.
6. **SAPK2a, active (Catalogue# 14-251):** Use at a final assay concentration of 2U/ml. Prepare a 500mU stock and add 2.5 μ l of stock per assay point.
7. **5 x Magnesium/ATP Cocktail:** 50mM MgAc, 0.5mM ATP.
8. **[γ -³³P]ATP:** 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.)
9. **PRAKtide (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.

Assay Procedure:

Stage One: *Activation of PRAK by SAPK2a*

1. Add 5 μ l of activation buffer to a microcentrifuge tube.
2. Add 2.5 μ l (50mU) of **SAPK2a, active**.
3. Add **5 μ l of unactive PRAK**.
4. Add 7.5 μ l of dH₂O.
5. Add 5 μ l of Magnesium/ATP cocktail.
6. Incubate for 60 minutes at 30°C.
7. Stop the reaction by diluting 10 and 20-fold in PRAK dilution buffer and store on ice.

Stage Two: *Phosphorylation of PRAKtide by PRAK*

1. Add 6.25 μ l of assay buffer to a microcentrifuge tube.
2. Add 2.5 μ l of **PRAKtide** stock solution.
3. Add **2.5 μ l of Stage One** reaction product.
4. Add 3.75 μ l of dH₂O.
5. Add 10 μ l of the diluted [γ -³³P] ATP.
6. Incubate for 10 minutes at 30°C.
7. Slowly transfer 20 μ l onto the centre of a 2cm x 2cm **P81** paper square.
8. Wash assay squares twice for 5 minutes with 50mM phosphoric acid.
9. Wash assay squares once with acetone for 2 minutes.
10. Transfer assay squares to scintillation vials and add 1ml scintillation cocktail.
11. Read in scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all appropriate 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. Protein is R291E with respect to this accession. This conflict is reported in GenBank AI300977, AL556217 and BI495728.

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 KIDQGDLMTP QFTPYYPVAPQ VLEAQRHQQK EKSGIIPSTP TPYTYNKSCD LWSLGVIIYV
241 MLCGYPPFYS KHHSRTIPKD MRRKIMTGSF EFPEEWSQI SEMAKDVVRK LLKVKPEERL
301 TIEGVLDHPW 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 gagtctgtgt aaagaaatct actcaagaac ggtttgcgct gaaaattctt
181 cttgatcgtc caaaagctag aaatgaggta cgtctgcaca tgatgtgtgc cacacacca
241 aacatagttc agattattga agtgtttgct aacagtgtcc agtttcccca tgagtccagc
301 ctaggggcc gactcttaat tgtaatggag atgatggaag ggggagagct atttcacaga
361 atcagccagc accggcactt tacagagaag caagccagcc aagtaacaaa gcagatagct
421 ttggctctgc ggcactgtca cttgttaaac attgcgcaca gagacctcaa gcctgaaaat
481 ctgcttttta aggataactc tttggatgcc ccagtgaagt tgtgtgactt tggatttgcc
541 aagattgacc aaggtgactt gatgacacc cagttcacc cttattatgt agcaccaccag
601 gtactggagg cgaaagaag gcatcagaag gagaaatctg gcatcatacc tacctcaccg
661 acgccctaca cttacaacaa gagctgtgac ttgtgggtccc taggggtgat tatctatgtg
721 atgctgtgcg gataccctcc tttttactcc aaacaccaca gccggactat cccaaaggat
781 atgcgaagaa agatcatgac aggcagtttt gagttcccag aggaagagtg gagtcagatc
841 tcagagatgg ccaaagatgt tgtgaggaag ctccctgaagg tcaaaccgga ggagagactc
901 accatcgagg gagtgtctgga ccaccctgg ctcaattcca ccgaggccct ggataatgtg
961 ctgccttctg ctgagctgat gatggacaag gcagtgggtg caggaatcca gcaggctcac
1021 gcggaacagt tggccaacat gagaatccag gatctgaaag tcagcctcaa acccctgcac
1081 tcagtgaaca accccattct gcggaagagg aagttacttg gcaccaagcc aaaggacagt
1141 gtctatatcc acgaccatga gaatggagcc gaggattcca atgttgctt ggaaaaactc
1201 cgagatgtga ttgctcagtg tattctcccc caggctggag agaatagaaga tgagaaaactg
1261 aatgaagtaa tgcaggaggc ttggaagtat aaccgggaat gcaaactcct aagagatact
1321 ctgcagagct tcagctggaa tggctcgtgga ttcacagata aagtagatcg actaaaactg
1381 gcagaaattg tgaagcaggt gatagaagag caaacacgt 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.