

# Certificate of Analysis

**ICK, active**  
**(Recombinant enzyme expressed in Sf21 insect cells)**  
**Item # 15-023, 15-023-K, 15-023M**  
**Parent Lot # 202962**

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, amino acids 1-312 expressed by baculovirus in Sf21 insect cells. Purified using immobilized metal affinity chromatography. Purity 69% by SDS-PAGE and Coomassie blue staining. MW = 40kDa.

**Specific Activity (Parent lot# 202962):** 520U/mg, where one unit of ICK activity is defined as 1nmol phosphate incorporated into 250µM RRRFRPASPLRGPPK per minute at 30°C with a final ATP concentration of 100µM.

**Formulation:** 0.25mg/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.

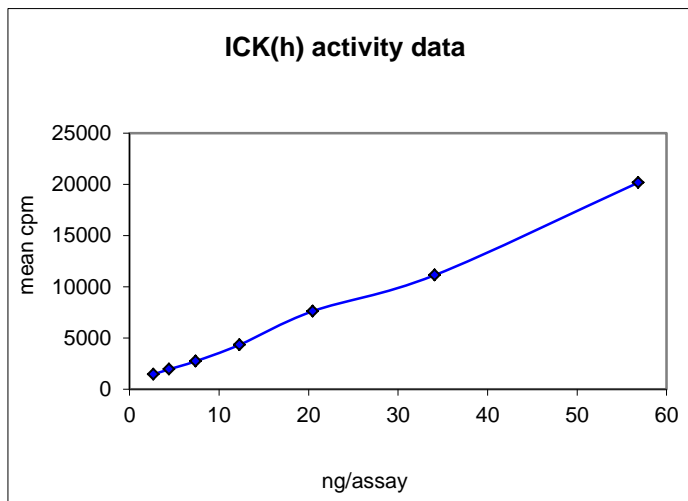
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**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.

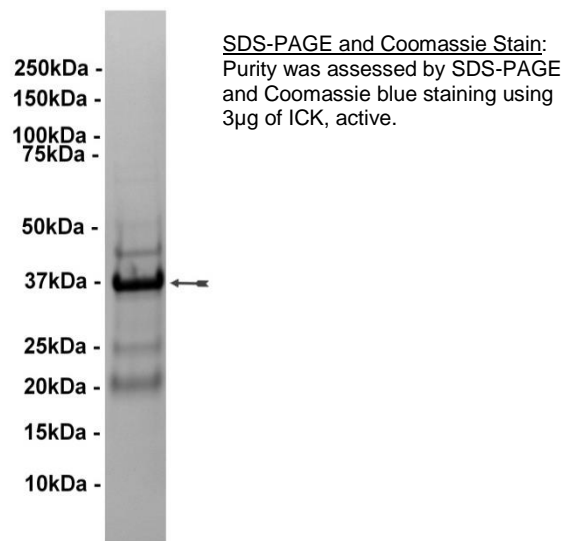
**FOR IN VITRO RESEARCH USE ONLY**  
**NOT FOR USE IN HUMANS OR ANIMALS**

## Quality Control Testing

**Kinase Assay:** 2.7–56.8ng of this lot of enzyme phosphorylated 250µM RRRFRPASPLRGPPK 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 identity as ICK with the translated sequence listed on page three.



# Certificate of Analysis

## Kinase Assay Protocol

### Stock Solutions:

- 1. 5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
- 2. RRRFRPASPLRGPPK:** Use at a final assay concentration of 250 $\mu$ M. Prepare a 2.5mM stock and add 2.5 $\mu$ l of stock per assay point.
- 3. Enzyme:, active** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 2.7–56.8ng per assay point.
- 4. [ $\gamma$ -<sup>33</sup>P]ATP:** 2.5 x MgAc/[ $\gamma$ -<sup>33</sup>P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [ $\gamma$ -<sup>33</sup>P]ATP (specific activity approximately 500 - 800cpm/pmol as required).

### Assay Procedure (96 well plate format):

1. Add 5 $\mu$ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 $\mu$ l of RRRFRPASPLRGPPK.
3. Add **2.5 $\mu$ l (2.7–56.8ng) ICK, active.**
4. Add 5 $\mu$ l of dH<sub>2</sub>O.
5. Add 10 $\mu$ l of diluted [ $\gamma$ -<sup>33</sup>P]ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 $\mu$ l of 3% phosphoric acid.
8. Transfer a 10 $\mu$ 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 dried 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 $\mu$ l of 30% phosphoric acid.

# Certificate of Analysis

## ICK, active Sequence Information

<b>Protein</b>	Human
<b>Tags</b>	N-terminal 6His
<b>Native sequence</b>	M31 of the recombinant protein is equivalent to M1 of human ICK
<b>Accession number</b>	GenBank NM_014920.2

### Recombinant ICK amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMDPEF MNRYYTTIRQL GDGTYGSVLL GRSIESGELI
61 AIKKMKRKFY SWEECMNLRE VKSLKKNLHA NVVKLKEVIR ENDHLYFIFE YMKENLYQLI
121 KERNKLPES AIRNIMYQIL QGLAFIHKHG FFHRDLKPEN LLCMGPELVK IADFGGLAREI
181 RSKPPYTDYV STRWYRAPEV LLRSTNYSSP IDVWAVGCIM AEVYTLRPLF PGASEIDTIF
241 KICQVLGTPK KTDWPEGYQL SSAMNFRWPQ CVPNNLKTLI PNASSEAVQL LRDMLQWDPK
301 KRPTASQALR YPYFQVGHPL GSTTQNLQDS EKPQKGILEK AG

```

### Recombinant ICK nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg gcgccatgga tccggaattc atgaatagat acacaacaat caggcagctc
121 ggggatggaa cctacggttc cgtcctgctg ggaagaagca ttgagtctgg ggagctgac
181 gctattaaaa aaatgaaaag aaaattttat tcctgggagg aatgcatgaa ccttcgggag
241 gttaagtctt taaagaagct caaccatgcc aatgtagtca aattaaaaga agttatcagg
301 gaaaatgata atctttatct tatcttcgag tacatgaagg aaaatcttta ccagctcatt
361 aaagagagaa ataagttggt tcctgagtct gctataagga atatcatgta tcagatatta
421 caaggactcg catttattca caaacacggc ttctttcatc gagacttaaa gcctgagaac
481 tccctctgca tgggaccaga acttgtgaaa attgcagact ttggtttggc ccgagaaata
541 cgatcaaaac ctccatatac agattatgta tctaccagat ggtacagggc tccagaagta
601 ctctgagggt ctaccaacta cagctcccc attgacgtct gggcggtggg ctgcatcatg
661 gcagaagttt acaccctcag gccactcttc cctggagcca gtgaaattga cacaatatc
721 aaaatttgcc aagtgctggg gacacaaaa aagactgact ggctgaagg ctatcaactt
781 tcaagtgcaa tgaacttccg ttggccacag tgtgtacca ataacttaaa gaccttgatt
841 cccaatgcta gcagtgaagc agtccagctc ctgagagaca tgcttcagtg ggatcccaag
901 aaacgaccaa cagctagtca ggcacttcga tctccttact tccaagttgg acaccacta
961 ggcagacca cacaaaacct tcaggattca gaaaaaccac agaaaggcat cctggaaaag
1021 gcaggctaa

```

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.