

## Certificate of Analysis

### IGF-1R, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-465, 14-465-K, 14-465M

Parent Lot # WAB0189

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 IGF-1R residues 959–end, expressed by baculovirus in Sf21 insect cells. Purified using Ni<sup>2+</sup>/NTA agarose. Purity 78% by SDS-PAGE and Coomassie blue staining. MW = 48kDa.

**Specific Activity (Parent lot# WAB0189):** 97U/mg, where one unit of IGF-1R activity is defined as 1nmol phosphate incorporated into 250µM (KKKSPGEYVNIEFG) per minute at 30°C with a final ATP concentration of 100µM.

**Formulation:** 2.03mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 0.2mMMPMSF, 1mM benzamidine, 0.1% 2-mercaptoethanol. Frozen solution.

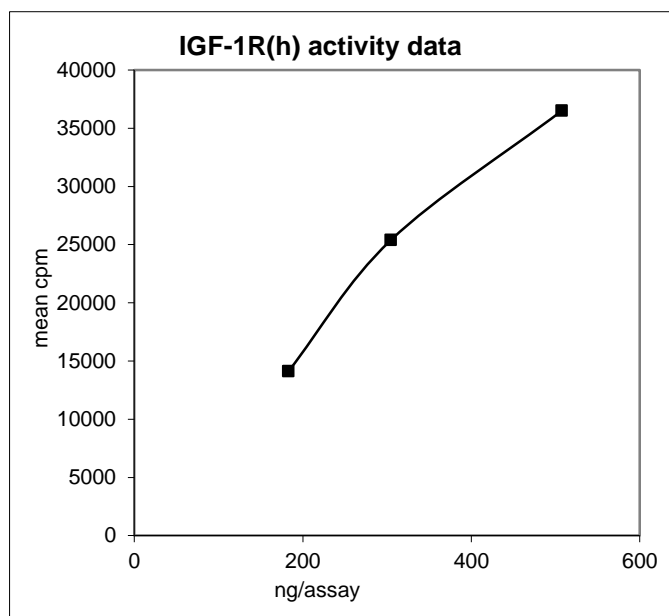
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 6 months from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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

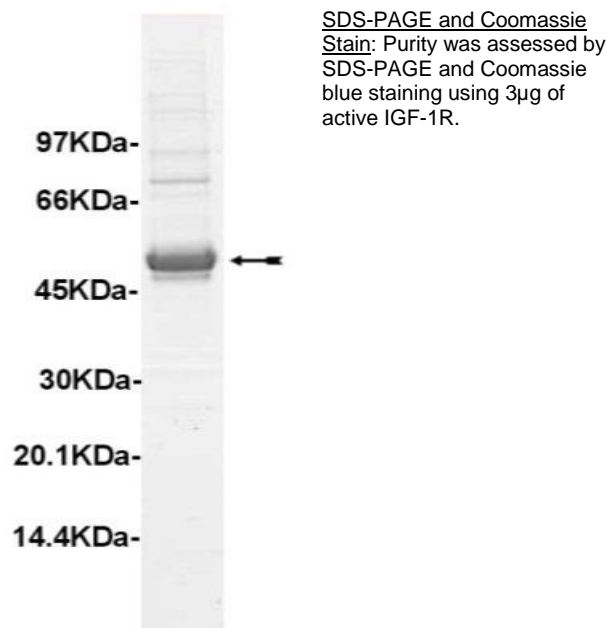
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Kinase Assay:** 183–508ng of this lot of enzyme phosphorylated 250µM (KKKSPGEYVNIEFG) 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 IGF-1R with of the translated sequence listed on page three.



## Certificate of Analysis

## Kinase Assay Protocol

Stock Solutions:

1. **10 x Reaction Buffer:** 500mM Tris/HCl pH7.5, 1mM EGTA, 1mM Na<sub>3</sub>VO<sub>4</sub>, 1% 2-mercaptoethanol.
2. **(KKKSPGEYVNIEFG):** Use at a final assay concentration of 250µM. Prepare a 2.5mM stock and add 2.5µl of stock to reaction mixture.
3. **Manganese Chloride (MnCl<sub>2</sub>):** Use at a final assay concentration of 10mM. Prepare a 200mM stock and add 1.25µl of stock per assay point.
4. **IGF-1R, active:** Dilute with 50mM Tris/HCl pH7.5, 0.1mM EGTA, 0.1mM Na<sub>3</sub>VO<sub>4</sub>, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 183–508ng per assay point.
5. **[γ-<sup>33</sup>P]ATP:** 2.5 x magnesium acetate/[γ-<sup>33</sup>P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [γ-<sup>33</sup>P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

Assay Procedure (96 well plate format):

1. Add 2.5µl of 10 x reaction buffer to wells.
2. Add 2.5µl of **(KKKSPGEYVNIEFG)**.
3. Add 1.25µl of MnCl<sub>2</sub>.
4. Add 6.25µl of dH<sub>2</sub>O.
5. Add **2.5µl (183–508ng) IGF-1R, active**.
6. Add 10µl of diluted [γ-<sup>33</sup>P]ATP mixture.
7. Incubate for 10 minutes at 30°C.
8. Stop the reaction by adding 5µl 3% phosphoric acid to wells.
9. Transfer a 10µl aliquot onto the appropriate area of a **P30 Filtermat**.
10. Wash the filtermat three times for 5 minutes with 75mM phosphoric acid.
11. Wash the filtermat once for 2 minutes with methanol.
12. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
13. 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

## IGF-1R Sequence Information

<b><u>Protein</u></b>	human IGF-1R
<b><u>Tags</u></b>	N-terminal 6His
<b><u>Native sequence</u></b>	H16 of the recombinant protein is equivalent to H959 of human IGF-1R
<b><u>Accession number</u></b>	GenBank X04434

### ***Recombinant IGF-1R amino acid sequence:***

```

1 MAHHHHHHEN LYFQGHRKRN NSRLNGVLY ASVNPEYFSA ADVYVPDEWE VAREKITMSR
61 ELGQGSFGMV YEGVAKGVVK DEPTRVAIK TVNEAASMRE RIEFLNEASV MKEFNCHHV
121 RLLGVVSQGG PTLVIMELMT RGDLSYLRS LRPEMNNPV LAPPSSKMI QMAGEIADGM
181 AYLNANKFVH RDLAARNCMV AEDFTVKIGD FGMTRDIYET DYYRKGGKGL LPVRWMSPE
241 LKDGVTFTYS DVWSFGVVLW EIATLAEQPY QGLSNEQVLR FVMEGGLDK PDNCPDMLFE
301 LMRMCWQYNP KMRPSFLEII SSIKEEMEPG FREVSFYSE ENKLPEPEEL DLEPENMESV
361 PLDPSASSSS LPLDRHSGH KAENGPGPV LVLRSFDER QPYAHMNGGR KNERALPLPQ
421 SSTC
  
```

### ***Recombinant IGF-1R nucleotide sequence:***

```

1 atggcgcatt accatcacca tcatgaaaac ctgtattttc agggccatag aaagagaaat
61 aacagcaggc tggggaaatgg agtgctgtat gcctctgtga acccggagta cttcagcgct
121 gctgatgtgt acgttcctga tgagtgggag gtggctcggg agaagatcac catgagccgg
181 gaacttgggc aggggtcgtt tgggatggct tatgaaggag ttgccaaggg tgtggtgaaa
241 gatgaacctg aaaccagagt ggccattaaa acagtgaacg aggccgcaag catgctgtgag
301 aggattgagt ttcacaacga agcttctgtg atgaaggagt tcaattgtca ccatgtggtg
361 cgattgctgg gtgtggtgtc ccaaggccag ccaacactgg tcatcatgga actgatgaca
421 cggggcgatc tcaaaaagta tctccggtct ctgaggccag aaatggagaa taatccagtc
481 ctgacacctc caagcctgag caagatgatt cagatggccg gagagattgc agacggcatg
541 gcataacctc aagcaataa gttcgtccac agagaccttg ctgcccggaa ttgcatggta
601 gccgaagatt tcacagtcaa aatcggagat tttggtatga cgcgagatat ctatgagaca
661 gactattacc ggaaaggagg gaaaggctg ctgcccgtgc gctggatgtc tcctgagtcc
721 ctcaaggatg gagtctcac cacttactcg gacgtctggt ccttcggggt cgtcctctgg
781 gagatcgcca cactggccga gcagccctac cagggtctgt ccaacgagca agtccttcgc
841 ttgctcatgg agggcggcct tctggacaag ccagacaact gtcctgacat gctgtttgaa
901 ctgatgcgca tgtgctggca gtataacccc aagatgaggc cttccttcct ggagatcatc
961 agcagcatca aagaggagat ggagcctggc ttccgggagg tctccttcta ctacagcgag
1021 gagaacaagc tgcccagacc ggaggagctg gacctggagc cagagaacat ggagagcgtc
1081 cccctggacc cctcggcctc ctctgtctcc ctgccactgc ccgacagaca ctcaggacac
1141 aaggccgaga acggccccgg ccctgggggtg ctggctctcc gcgccagctt cgacgagaga
1201 cagccttacg cccacatgaa cgggggcccgc aagaacgagc gggccttgcc gctgccccag
1261 tcttcgacct gctga
  
```

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