



EAGLE  
BIOSCIENCES

# **Biotinylated Rat IGFBP-4 Recombinant Protein**

Catalog Number:

**BP420-G2.5** (1 x 96 wells)

For Research Use Only. Not for use in diagnostic procedures.

v. 1.0

---

**Eagle Biosciences, Inc.**

20A Northwest Blvd., Suite 112, Nashua, NH 03063

Phone: 617-419-2019 Fax: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com)



## **Biotinylated Human Insulin-like Growth Factor 1 (IGF-1) •**

### **Description:**

Biotinylated rat IGFBP-4 has been prepared from receptor grade recombinant rat IGFBP-4 (product code BP4CU020) using the IBT biotinylation technology. The product was purified using chromatographic techniques.

- **Appearance:**  
Lyophilized with no additives.
- **Reconstitution** Reconstitute 2.5 µg aliquot with 0.1 ml double distilled water or buffer of your choice.
- **Storage**  
The product is shipped at room temperature, for long term storage store at – 20 °C. Reconstituted samples can be stored for at least one week at 2-8 °C.

### **Application Notes**

Biotinylated IGF's and analogs have found a wide range of applications in-vitro and in-vivo (see application note IGF004) and are a safe and stable alternative to 125I-labeled IGF's. As 125I-labeled IGFBP's have also been used for methods as e.g. proteolysis assays, binding studies and in-vivo studies, there should be some potential for biotinylated IGFBP's, too. A literature search on the use of biotinylated IGFBP's resulted in a small number of reports, which are limited to IGFBP-1(1-3), IGFBP-2 (4,5,17), IGFBP-3 (6-14, 18) and IGFBP-4 (15). Until today we found no scientific papers on the use of biotinylated IGFBP-5 and IGFBP-6.

Though there is small number of papers, the authors have used the biotinylated IGFBP's in a broad range of techniques to study IGF-dependent and IGF-independent actions of IGFBP's. Biotinylated IGFBP's have been used in immunoassays (1, 6, 11, 15), proteolysis studies (2,3), ligand blot (4,10, 12), cross-linking studies (9, 12) and binding studies (1, 4, 5, 7, 8, 13). Biotinylated IGFBP-3 has been used to demonstrate the nuclear appearance of IGFBP-3 in vivo (14) and as a substrate for transglutaminase (13).

### **Biotinylation Technology:**

We have developed our own biotinylation technology for IGF's that is different from the published methods. Our technology permits biotinylation without loss of binding capacity to IGFBP's or IGF antibodies.

### **Studies with Biotinylated IGFBP's from IBT**

To prove, that biotinylated IGFBP's from IBT are useful tools to demonstrate IGF dependent or independent actions of IGFBP's we used IGFBP-3 as an example. IGFBP-3 was biotinylated using the same technology as for our biotinylated IGF's (as described in notes below).



## **Stability**

Though our biotinylated IGF's are very stable in solution, we have shipped them in the past on dry ice to ensure maximum activity for our customers. This has increased shipping cost so we have decided to lyophilize our IGF's. In this form we have kept them for three months at temperatures up to 35 °C without loss of activity. So they can be shipped now without dry ice at reduced freight cost.

For long term storage, we recommend to store the IGF's refrigerated after arrival. The reconstituted biotinylated IGF's are stable for at least one year at a concentration of 20 µg/ml, if kept frozen at -20 °C, and at least for one month at 4 °C. We recommend to aliquot the reconstituted biotinylated IGF's and to store them at a temperature of - 20 °C.

## **Potential Uses for Biotinylated IGF's**

### **Western-Ligand Blotting**

We have tested our biotinylated IGF's and analogs in western-ligand blotting.

Chemiluminescent substrates as well as colorimetric ones have successfully been used. We also have developed our own buffer system, with buffers, that are stable for at least six months and give an excellent signal-to background ratio. Kits are available for human (bovine, sheep, donkey, pig, guinea pig, goat) IGFBP's and mouse/rat IGFBP's.

### **Western-Ligand blots with biotinylated IGF-I and analogs**

Biotinylated IGF-I binds to all six IGFBP's, but binds to IGFBP-6 very weakly. For all IGF-I analogs at least weak signals have been obtained with IGFBP-1 to -5, but no signals were observed with IGFBP-6.

### **ELISA**

We have tested our biotinylated human IGF-I and IGF-II in a competitive ELISA format using our polyclonal antisera PAA1 and PAC1. Binding to the antibodies was not affected by biotinylation.

### **Immunoprecipitation/Chromatography of IGFBP's**

We have used our biotinylated human IGF-I and IGF-II in immunoprecipitation experiments. Binding of biotinylated IGF's to Streptavidin may also be used to prepare media for affinity chromatography in a column format.

### **Cross-linking Experiments**

We have started to develop methods using our biotinylated IGF's for cross-linking experiments. The method is not fully developed, e.g. the washing procedure needs to be improved and the extraction procedure for proteins needs improvement, too. Preliminary results are available.

### **Immunohistochemistry**

We have started experiments with our biotinylated IGF's for IHC with human cells fixed on slides. The results of histochemistry have not been examined systematically, but staining has been observed on the surface of the cells.

**Flow Cytometry:**

The use of Biotin-Des (1-3) IGF-I in flow cytometry has been reported by Xu et al. (1995), Immunology 85, 394-399.

**References**

Khawaja, X. Z.: Development of a scintillation proximity assay for human insulin-like growth factor-binding protein 4 compatible with inhibitor high-throughput screening. Analytical Biochemistry, Volume 366, Issue 1, 1 July 2007, Pages 80-86



## **Warranty Information**

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at [info@eaglebio.com](mailto:info@eaglebio.com) or at 866-411-8023.