

# GroPep Bioreagents

## Human ProIGF-I EB (104-123) antiserum (Rabbit)

### Description

Human ProIGF-I EB (104-123) antiserum was prepared by immunising rabbits with a 21 amino acid peptide derived from human ProIGF-I EB linked to diphtheria toxoid. The peptide is located within the E domain of the 195 amino acid ProIGF-I EB, derived from alternative splicing of the human *igf-1* gene. The antiserum recognises ProIGF-I EB but not ProIGF-I EA, ProIGF-I EC, mature IGF-I or IGF-II. Peptide sequence analysis suggest it may recognise ProIGF-I EB related proteins in a range of non-human primates but not many other species due to alternative splicing of the *igf-1* gene. There is significant homology to similar regions of ProIGF-I variants in bovine, porcine and canine species to suggest it may be effective.

### References:

Rotwein P *et al* (1986) Journal of Biological Chemistry **261**, 4828-4832  
Duguay SJ *et al* (1999) Hormone and Metabolic Research **31**, 43-49

### Immunogen:

Derived from UniProtKB/Swiss-Prot sequence ID P05019 isoform 1 also known as IGF-IB which is recognised as the canonical sequence. The peptide sequence is amino-acids 151-170, derived from ProIGF-I EB, conjugated to diphtheria toxoid.

### Source:

Rabbit antiserum

### Specificity:

Recognises Human ProIGF-I EB (isoform 1, P05019).  
No cross-reactivity against ProIGF-I EA (isoform 2, P05019-2);  
ProIGF-I EC (Q13429) mature IGF-I or IGF-II.

### Appearance:

Lyophilized powder

### Storage/Stability:

At least 2 years at 2 - 4°C (lyophilized).  
After reconstitution store at -20°C or -80°C  
Avoid freeze-thaw cycles

### Reconstitution:

Dissolve the vial contents in 100 µl sterile water and then dilute in a buffer suitable for the application.

### Application and Titre:

Each application and titre should be determined in house. As a guideline the antiserum can be used at a dilution of 1:1000 for Western immunoblotting.

### Product Codes and Pricing

**Human ProIGF-I EB (104-123) antiserum**      **100 µl**      **PAAT1**

### Related Products:

Human ProIGF-I E (71-86) antiserum (Rabbit)  
Human ProIGF-I EA (92-105) antiserum (Rabbit)  
Human ProIGF-I EC (104-110) antiserum (Rabbit)

**\*\*NOT FOR USE IN HUMANS\*\***

**GroPep Bioreagents Pty Ltd**  
51 West Thebarton Road  
Thebarton SA 5031  
Australia

**Postal Address:**  
PO Box 10065  
Adelaide Business Centre  
SA 5000  
Australia

**Email:**  
[info@gropep.com](mailto:info@gropep.com)

**Internet:**  
[www.gropep.com](http://www.gropep.com)

**ABN** 93 147 032 166

**Telephone:** +61 8 7222 1051