



# GroPep Bioreagents



## IGF-II Polyclonal Antibody (Rabbit)

### Description

This IGF-II antibody was prepared by immunising rabbits with a human IGF-II. The antiserum was then affinity purified. It detects IGF-II in human, mouse, rat and chicken. It cross-reacts up to 20% with human IGF-I. The antibody can be used for a number of applications including immunohistochemistry and Western ligand blotting.

### References:

Owens PC *et al* (1990) Journal of Endocrinology **124**, 269-275

### Immunogen:

Human IGF-II

### Source:

Affinity purified rabbit antiserum

### Specificity:

Characterised for IGF-II primarily by Western ligand blotting and in sections from human, rat and mouse and tissues.  
Cross-reactivity to IGF-I is 20% so appropriate blocking is required

### 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 500µl deionised water and then dilute into an appropriate buffer for the application.

### Application and Titre:

Each application and titre should be determined in house but as a guideline, use at 1:100 – 1:1000 for immunohistochemistry or Western ligand blotting. Examples are shown below.

### Product Codes and Pricing

#### IGF-II

#### Polyclonal Antibody

50 µg

PAAL1

### Related Products:

Human IGF-II antiserum  
Human IGF-II (Receptor grade)  
Chicken IGF-II  
Rat IGF-II

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

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

ABN 93 147 032 166

Telephone: +61 8 7222 1051

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)

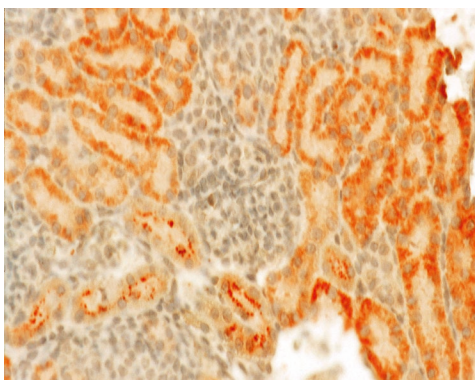


Figure 1 Rat embryo kidney

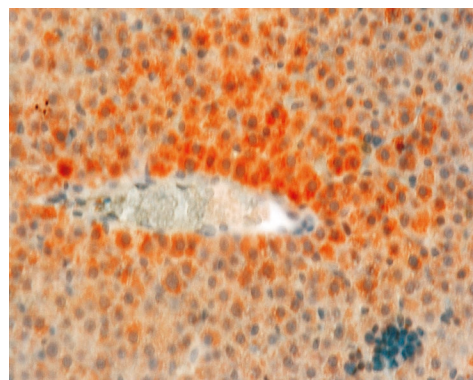


Figure 4 Rat embryo liver

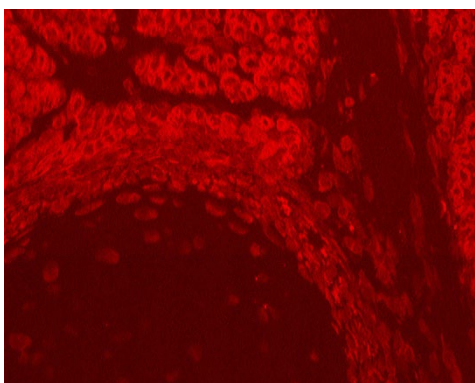


Figure 2 Rat embryo rib

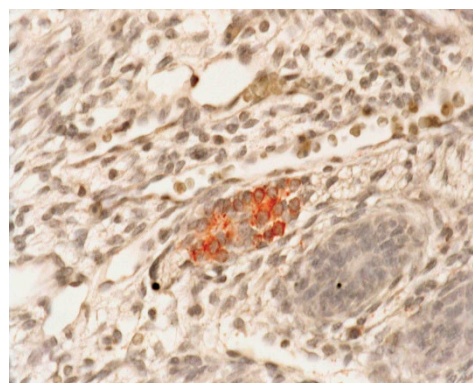


Figure 5 Mouse embryo hair follicle

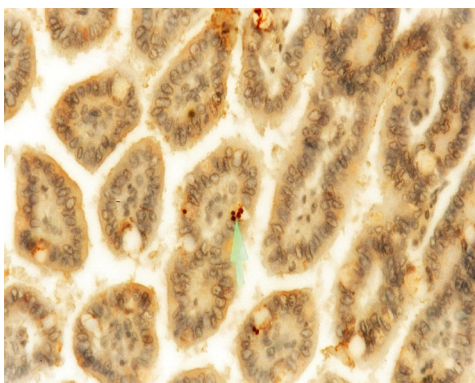


Figure 3 Mouse embryo intestines

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

**ABN** 93 147 032 166

**Telephone:** +61 8 7222 1051

**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)