

Human IGF-II Polyclonal Antibody (Rabbit)

Description

Human IGF-II polyclonal antibody was prepared by immunising rabbits with a human IGF-II conjugated to ovalbumin. The immunoglobulin fraction was purified by Protein A affinity chromatography. The antibody cross-reacts 100% with human and porcine IGF-II; 96% with chicken IGF-II; 91% with rat IGF-II and 85% with bovine IGF-II. Cross-reactivity to human, chicken and rat IGF-I is < 0.1%. It does not cross react with insulin from a range of species. The antibody is suitable for use in immunoassay, immunoblotting and immunohistochemistry.

References:

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

Immunogen:

Human IGF-II conjugated to ovalbumin

Source:

Rabbit antiserum. The immunoglobulin fraction was purified by Protein A affinity chromatography of serum obtained from a rabbit immunised with a recombinant human IGF-II conjugated to ovalbumin.

Specificity:

Human and porcine IGF-II 100%.
Chicken IGF-II 96%, rat IGF-II 91%, bovine IGF-II 85%.
Human, chicken and rat IGF-I < 0.1%.
No cross-reactivity to insulin

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 400 µl sterile water or buffer solution and stand for 10 minutes at room temperature. Dilute further in a buffer suitable for the application.

Application and Titre:

Suitable for use in immunoassay, immunoblotting and immunochemistry. Each application and titre should be determined in house. As a guideline, following reconstitution a dilution in the range 1:500–1:32,000 should give a suitable absorbance in a conventional immunoassay format.

Product Codes and Pricing

Human IGF-II Polyclonal Antibody

400 µg **PAC1**

Related Products:

IGF-II Polyclonal Antibody (Rabbit)
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

Internet:
www.gropep.com

