



GroPep Bioreagents



Human ProIGF-I EA (92-105) antiserum (Rabbit)

Description

Human ProIGF-I EA (92-105) antiserum was prepared by immunising rabbits with a 15 amino acid peptide derived from the E domain of human ProIGF-I EA linked to diphtheria toxoid. The peptide is located at the C terminus of the 153 amino acid ProIGF-I EA derived from alternative splicing of the human *igf-1* gene. The antiserum recognises ProIGF-I EA but not ProIGF-I EB, ProIGF-I EC, mature IGF-I or IGF-II. Sequence analysis suggests it may recognise ProIGF-I EA related proteins in a range of other species including non-human primates, rabbit, guinea-pig, dog, horse, sheep and goat but not rodents, due to alternative splicing of the *igf-1* gene

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
Pfeffer LA *et al* (2009) Molecular Biology of the Cell **20**, 3810-3817

Immunogen:

Derived from UniProtKB/Swiss-Prot sequence ID P05019-2 isoform 2 also known as IGF-IA. The peptide sequence is amino-acids 139-153, derived from ProIGF-I EA, conjugated to diphtheria toxoid.

Source:

Rabbit antiserum

Specificity:

Recognises Human ProIGF-I EA (isoform 2, P05019-2).
No cross-reactivity against ProIGF-I EB (isoform 1, P05019); 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 EA (92-105)
antiserum**

100 µl

PAAU1

Related Products:

Human ProIGF-I E (71-86) antiserum (Rabbit)
Human ProIGF-I EB (104-123) 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

ABN 93 147 032 166

Telephone: +61 8 8152 9360

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

Email:
info@gropep.com

Internet:
www.gropep.com