



GroPep Bioreagents



Human ProIGF-I E (71-86) antiserum (Rabbit)

Description

Human ProIGF-I E (71-86) antiserum was prepared by immunising rabbits with a 16 amino acid peptide derived from the E domain of human ProIGF-I E linked to diphtheria toxoid. The peptide sequence is located immediately downstream of the cleavage site at the C terminus of mature IGF-I and as a consequence, the antiserum recognises all ProIGF-I E proteins derived from alternative splicing of the human *igf-1* gene. It does not recognise mature IGF-I or IGF-II. It is likely to recognise ProIGF-I E proteins in non-human primates, rabbits and guinea-pigs but not other species 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

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 119-134, derived from ProIGF-I EB, conjugated to diphtheria toxoid.

Source:

Rabbit antiserum

Specificity:

Recognises Human ProIGF-I EB (isoform 1, P05019); ProIGF-I EA (isoform 2, P05019-2); ProIGF-I EC (Q13429).
No cross-reactivity against 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 E (71-86)
antiserum**

100 µl

PAAV1

Related Products:

Human ProIGF-I EA (92-105) 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 7222 1051

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

Email:
info@gropep.com

Internet:
www.gropep.com