



Human ProIGF-II E (89-101) antiserum (Rabbit)

Description

Human ProIGF-II E (89-101) antiserum was prepared by immunising rabbits with a 13 amino acid peptide derived from the E domain of human IGF-II E linked to diphtheria toxoid. The sequence corresponds to IGF-II E₈₉₋₁₀₁ and as a consequence, the antiserum recognises full length IGF-II E and IGF-II E₁₋₁₀₄ but not IGF-II E₁₋₈₈ derived from processing of the human *igf-2* gene and following proteolytic cleavage. It does not recognise mature IGF-II or IGF-I. Sequence analysis suggests it is likely to recognise IGF-II E proteins in non-human primates, dog, guinea-pig, pig and potentially cattle.

References:

Duguay SJ *et al* (1999) *Hormone and Metabolic Research* **31**, 43-49
Khosla S *et al* (2002) *Journal of Clinical Endocrinology and Metabolism* **87**, 3867-3870

Immunogen:

Derived from UniProtKB/Swiss-Prot sequence ID P03144 isoform 1. The peptide sequence is amino-acids 113-125, derived from IGF-II E, conjugated to diphtheria toxoid.

Source:

Rabbit antiserum

Specificity:

Determined by Western blotting. Recognises full length human IGF-II E (IGF-II E₁₋₁₅₆) and proteolytically cleaved variants (IGF-II E₁₋₁₀₄) but not (IGF-II E₁₋₈₇).

No cross-reactivity against mature IGF-II or IGF-I.

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 500 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 1:8000 for Western immunoblotting and 1:4000 for enzyme immunoassay

Product Codes and Pricing

**Human ProIGF-II E (89-101)
antiserum**

100 µl

PAAY1

Related Products:

Human ProIGF-II E (138-156) antiserum (Rabbit)
Human Pro IGF-II E (78-88) antiserum (Rabbit)
Human Pro IGF-II (1-104)
Human Pro IGF-II (1-156)

****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