



GroPep Bioreagents



Human Pro IGF-I EC (104-110) Polyclonal Antibody (Rabbit)

Description

Human Pro IGF-I EC (104-110) Polyclonal Antibody was prepared by immunising rabbits with an 8 amino acid peptide derived from the C terminal domain of Pro IGF-I EC linked to diphtheria toxoid. The immunoglobulin fraction was purified by Protein A affinity chromatography. The antibody recognises Pro IGF-I EC but not Pro IGF-I EA, Pro IGF-I EB, mature IGF-I or IGF-II.

References:

Rotwein P *et al* (1986) Journal of Biological Chemistry **261**, 4828-4832
Chew SL *et al* (1995) Endocrinology **136**, 1939-1944
Duguay SJ *et al* (1999) Hormone and Metabolic Research **31**, 43-49

Immunogen:

Derived from UniProtKB/Swiss-Prot sequences ID Q13429 and E9PD02 also known as IGF-IC. The peptide sequence is amino-acids 152-158 (E9PD02) and 133-139 (Q13429) conjugated to diphtheria toxoid.

Source:

Rabbit antiserum

Specificity:

Recognises Human Pro IGF-I EC (sequences E9PD02/Q13429 UniProtKB database)

No cross-reactivity against Pro IGF-I EA (isoform 2, P05019-2); Pro IGF-I EB (isoform P05019); 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 Pro IGF-I EC (104-110)
Polyclonal Antibody**

100 µg

PAAS1

Related Products:

Human Pro IGF-I E (71-86) antiserum (Rabbit)

Human Pro IGF-I EA (92-105) antiserum (Rabbit)

Human Pro IGF-I EB (104-123) 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