

Rat IGF-II

Description

Rat insulin-like growth factor-II (IGF-II) is a single chain 67 amino acid polypeptide that shares structural homology with human IGF-II. Only four amino acids are different compared to the sequence of human IGF-II. Rat IGF-II has been shown to stimulate growth in L6 myoblasts and other cell lines. Receptor Grade rat IGF-II is a high quality research reagent particularly for use in studies on cellular and animal growth, for use in specific assays for rat IGF-II and other general research purposes.

References

Dull TJ et al. (1986) Nature **310**, 777-781

Source:

Produced in *E.coli*.

Molecular Weight:

7515 daltons

Purity:

>90 % (by HPLC analysis)

N-terminal sequence:

Analysis of 5 residues

Biological Activity:

IGF type 2 receptor binding assay ($ED_{50} < 5$ ng/ml)
Stimulation of protein synthesis in rat L6 myoblasts
($ED_{50} < 150$ ng/ml)

Endotoxin:

< 0.1 EU/μg

Appearance:

White powder freeze-dried from 0.1M acetic acid and stored under nitrogen at a slight vacuum.

Storage/Stability:

At least 2 years at 2-4°C (as a freeze dried product).

Reconstitution:

Handling of GroPep IGF-I, IGF-II and IGF analogues

Detection:

Procedure for the iodination of IGF peptides
Determination of IGF-I or IGF-II in a range of species by
Radioimmunoassay (RIA)

Product Codes:

Rat IGF-II

100 μg

AAU100

Related Products:

Rat IGF-I

Human IGF-II antiserum (Rabbit)

IGF-II polyclonal antibody

****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