

GroPep Bioreagents

Human [Leu⁶⁰] IGF-I

Description

Human [Leu60] insulin-like Growth Factor-I ([Leu60] IGF-I) is an analogue of human IGF-I which has a Leucine substitution in place of Tyrosine at position 60 of the human IGF-I sequence. [Leu60]IGF-I has normal binding to acid-stripped human serum binding proteins but has reduced binding to the type 1 IGF receptor and to the insulin receptor compared to IGF-I. It is a useful analogue for studying the relative effects of IGF-I through the type 1 receptor and IGF binding proteins.

References

Bayne ML *et al* (1990) Journal of Biological Chemistry **265**, 15648-15652

Source:

Produced in *E.coli*.

Molecular Weight:

7599 daltons

Purity:

>95 % (by HPLC analysis)

N-terminal sequence:

Analysis of 5 residues

Biological Activity:

Stimulation of protein synthesis in rat L6 myoblasts
(ED₅₀ >2000 ng/ml)
IGF type 1 receptor binding assay (ED₅₀ > 100 ng/ml)
IGF binding protein assay: ED₅₀ < 10 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

Product Codes and Pricing

Human [Leu60] IGF-I

100 μg

ABU100

Related Products:

Human [Ala31] IGF-I
Human [Leu24] IGF-I
Human [Ala31] [Leu60] IGF-I
Human [Leu24] [Ala31] IGF-I

****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