



# GroPep Bioreagents

## Barramundi IGF-I (*Lates calcarifer*)

### Description

Barramundi insulin-like growth factor-I (IGF-I) is a single chain 66 amino acid polypeptide that shares structural homology with human IGF-I. Barramundi IGF-I can be used for a variety of *in vivo* and *in vitro* applications including those investigating the actions of growth hormone.

<b>References:</b>	Degger B <i>et al</i> (2000) General and Comparative Endocrinology <b>117</b> , 395-403
<b>Source:</b>	Produced in <i>E.coli</i> .
<b>Molecular Weight:</b>	7309 daltons
<b>Purity:</b>	>95 % (by HPLC analysis)
<b>N-terminal sequence:</b>	Analysis of 5 residues
<b>Biological Activity:</b>	Stimulation of protein synthesis in rat L6 myoblasts (ED <sub>50</sub> < 10 nmol/l)
<b>Endotoxin:</b>	< 0.1 EU/μg
<b>Appearance:</b>	White powder freeze-dried from 0.1M acetic acid and stored under nitrogen at a slight vacuum.
<b>Storage/Stability:</b>	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 and Pricing:

**Barramundi IGF-I**                      **100 μg**                      **YU100**                      **US\$100**

**Related Products:** Barramundi IGF-I antiserum (Rabbit)  
Salmon/Trout IGF-I  
Bream 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](mailto:info@gropep.com)

**Internet:**  
[www.gropep.com](http://www.gropep.com)