

distributed in the US/Canada by:  
**EAGLE BIOSCIENCES, INC.**  
20A NW Blvd, Suite 112 Nashua, NH 03063  
Phone: 617-419-2019 FAX: 617-419-1110  
[www.EagleBio.com](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)



# GroPep Bioreagents

## Human IGF-I monoclonal antibody

### Description

Human IGF-I antibody is a mouse monoclonal antibody (IgG1 kappa) which binds to human IGF-I. The antibody was prepared by immunising mice with IGF-I and synthetic peptides of IGF-I and has 100% cross-reactivity with Des [1-3] IGF-I and LR<sup>3</sup>IGF-I. Cross-reactivity to rat IGF-I is <4%. Cross-reactivity to chicken IGF-I, salmon IGF-I and human IGF-II is <0.4%.

<b>Immunogen:</b>	Human IGF-I and synthetic peptides of IGF-I
<b>Source:</b>	Produced in hybridoma cells derived from the SP2/0 myeloma cell line. Purified from conditioned medium by Protein A affinity chromatography
<b>Specificity:</b>	Cross-reacts 100% with Des [1-3] IGF-I and LR <sup>3</sup> IGF-I. Cross-reactivity to rat IGF-I is <4%. Cross-reactivity to chicken IGF-I, salmon IGF-I and human IGF-II is <0.4%
<b>Appearance:</b>	White powder lyophilized from phosphate buffered saline containing 60 mM trehalose
<b>Storage/Stability:</b>	At least 2 years at 2 - 8°C (lyophilized) After reconstitution store at -20°C or -80°C Avoid freeze-thaw cycles
<b>Reconstitution:</b>	Reconstitute in sterile water as outlined on the Certificate of Analysis
<b>Applications:</b>	GroPep recommends optimal dilutions/concentrations should be determined by the end user. ELISA (1/500 – 1/8000) Western Blot

### Product Codes

<b>Human IGF-I</b>	<b>100 µg</b>	<b>MAMU100</b>
<b>monoclonal antibody</b>	<b>200 µg</b>	<b>MAMU200</b>

<b>Related Products:</b>	Human IGF-I antiserum (Rabbit) IGF-I polyclonal antibody (Rabbit) Human IGF-I (Receptor Grade) Human IGF-I (Media Grade)
--------------------------	---

**\*\*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

WDSP-09.03

**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)