

Instruction Manual

Product	Size	Catalog Number
Human Insulin (from yeast)	50 mg	C-52310

Product Description

Insulin is a hormone consisting of a two-chain polypeptide and is produced by the beta-cells of pancreatic islets. It regulates cellular uptake, utilization, and storage of glucose, amino acids, and fatty acids and inhibits the breakdown of glycogen, protein and fat. In serum-free medium insulin is an essential supplement for long-term growth of commonly used mammalian cell lines. Our Human Insulin is a white to off-white powder recombinantly expressed in yeast.

Physical / Chemical Data

Origin: Yeast

Purity: > 98%

Biological Activity: ≥ 25 IU/mg

Zinc Content: $\leq 1\%$

Solubility

10 mg/ml in 0.01 N HCl.

Storage and Stability

Store at -20°C in the dark immediately after arrival. If stored properly, the product is stable until the expiry date stated on the label.

Quality Control

All lots of PromoCell Insulin are subjected to comprehensive quality control tests. Each lot is routinely tested for the use with human primary cells and parameters such as biological activity and endotoxin content. Approved in-house lots are used as a reference.

In addition, all lots have been tested for the absence of microbial contaminants (fungi, bacteria).

Intended Use

The products are for in vitro research use only and not for diagnostic or therapeutic procedures. For safety precautions please see appropriate MSDS.

PromoCell GmbH
Sickingenstr. 63/65
69126 Heidelberg
Germany

info@promocell.com
www.promocell.com

USA/Canada

Phone: 1 - 866 - 251 - 2860 (toll free)
Fax: 1 - 866 - 827 - 9219 (toll free)

Deutschland

Telefon: 0800 - 776 66 23 (gebührenfrei)
Fax: 0800 - 100 83 06 (gebührenfrei)

France

Téléphone: 0800 - 90 93 32 (ligne verte)
Téléfax: 0800 - 90 27 36 (ligne verte)

United Kingdom

Phone: 0800 96 03 33 (toll free)
Fax: 0800 169 85 54 (toll free)

Other Countries

Phone: +49 6221 - 649 34 0
Fax: +49 6221 - 649 34 40

© PromoCell GmbH