

hGH

Human growth hormone (HGH, GH, somatotropin) is a 191-amino acid peptide pituitary hormone that stimulates growth, cell reproduction, and regeneration. Also regulates body composition, fluids, and sugar and fat metabolism. It is a stress hormone that increases glucose and fatty acid levels and stimulates the production of IGF-1. Can be used medically to treat growth disorders and GH deficiency. Has been abused by competitors for its anabolic properties, and is banned by the IOC and NCAA.

Materials Provided

Catalog Number:	92-1338
Quantity Supplied:	1 vial(s), 100 µg per vial
Reconstitution Buffer:	1 mL

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	22,000 g/mol
Source:	Recombinant
Purity:	≥ 98% by HPLC
Endotoxin Level:	<0.1 EU per 1 µg
Formulation:	N/A
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	100 µg/mL
Reconstitution Volume:	1 mL
Reconstitution Solvent:	Reconstitution Buffer

Handling Instructions:

N/A

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com