

hGH

Recombinant human protein expressed in E. coli cells

Catalog # G807-40N

Lot # S346-1

Product Description

Recombinant full-length human GH (Growth Hormone) was expressed in E. coli cells and purified by proprietary chromatographic techniques. The protein accession number is [NP_000506](#).

Gene Aliases

GH; GHN; GH-N; hGH-N; IGHD1B

Endotoxin Level

Less than 0.1 ng/µg (IEU/µg) of hGH.

Formulation

Recombinant protein stored in 1 X PBS, 10% glycerol.

Storage and Stability

Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

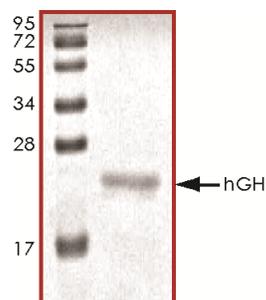
Scientific Background

hGH is a member of the somatotropin/prolactin family of hormones which play an important role in growth control and it is located at the growth hormone locus on chromosome 17 where they are interspersed in the same transcriptional orientation; an arrangement which is thought to have evolved by a series of gene duplications. hGH is synthesized by acidophilic or somatotrophic cells of the anterior pituitary gland (1). The systemic inhibition of GH or IGF1, or both, may have therapeutic potential in preventing some forms of retinopathy (2).

References

1. Niall, H. D. et.al: Sequence of pituitary and placental lactogenic and growth hormones: evolution from a primordial peptide by gene reduplication. Proc. Nat. Acad. Sci. 68: 866-869, 1971.
2. Smith, L. E. H. et.al: Essential role of growth hormone in ischemia-induced retinal neovascularization. Science 276: 1706-1709, 1997.

Purity



The purity was determined to be **>95%** by densitometry analysis in reducing SDS-PAGE Coomassie blue gel. Approx. MW of the major band was **22.3 kDa (22,260 Daltons)** by MW calculation).

hGH

Recombinant human protein expressed in E. coli cells

Catalog Number G807-40N

Specific Lot Number S346-1

Purity	>95%
Concentration	0.2 µg/µl
Stability	1yr At -20°C from date of shipment
Storage & Shipping	Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.