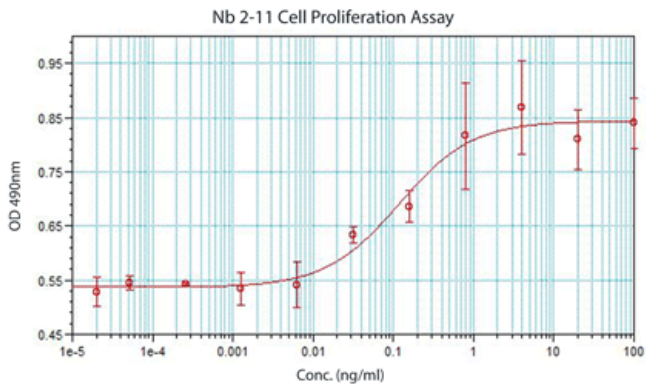


Human growth hormone, (recombinant)

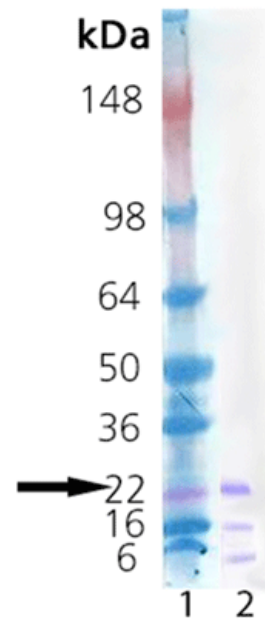
Human growth hormone is a growth factor that has been shown to play an important role in the growth control of cells. It is a member of a group of hormones produced by the pituitary gland that also include prolactin. Human Growth hormone has been shown to stimulate proliferation of Nb2-11 cells in a dose dependent manner. Its ability to promote growth is mediated by the somatomedin family of peptide hormones. It has been shown to enhance glucose transport and facilitates glycogen storage, as well as the ability to stimulate IGF-1 secretion in liver and other tissue.

Manuals, SDS & CofA

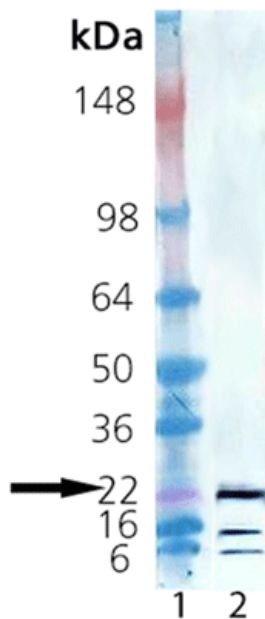
[View Online »](#)



Nb2-11 cell proliferation assay after treatment with Human Growth Hormone: Nb2-11 cells were placed in stationary media at 50,000cells/well in 96 well plates. They were treated with a tittered concentration of Human Growth hormone. The cells were grown at 37°C / 5% CO₂ for 48 hours. Proliferation was measured using an MTT cell viability assay. The ED50 range is around 0.02 – 0.5 ng/ml.



Coomassie stained SDS-PAGE. Lane 1, Molecular weight marker. Lane 2, 1.0µg Human Growth Hormone (ENZ-PRT189).



Western Blot analysis of Human Growth Hormone probed with rabbit anti-FLAG antibody: Lane 1: MWM, Lane 2: 100 ng Human Growth Hormone (ENZ-PRT189) probed with anti-FLAG.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Somatotropin, GH1, HGH
Biological Activity	Induces proliferation of Rat Nb2-11 cells (ED50 range is 0.02ng/ml-0.5ng/ml).
Endotoxin Content	<0.1 EU/μg protein (LAL test; Bio Whittaker)
Formulation	Lyophilized in PBS.
MW	~23 kDa (2 other immunoreactive bands are present at ~16 kDa and ~6 kDa)
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute in 100μl deionized water for a final concentration of 0.1mg/ml.
Source	Produced in HEK cells. The human growth hormone (aa 27-217) is fused at the N-terminus to a FLAG [®] -tag followed by an 8 AA linker.
Technical Info / Product Notes	FLAG is a registered trademark of Sigma-Aldrich Co.
UniProt ID	P01241

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com