

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.

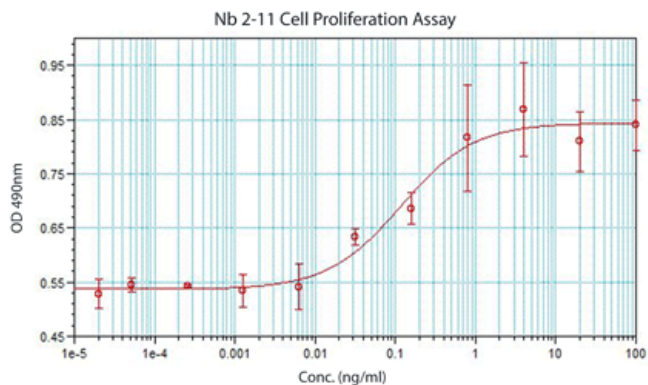
Ordering Information

[Order Online »](#)

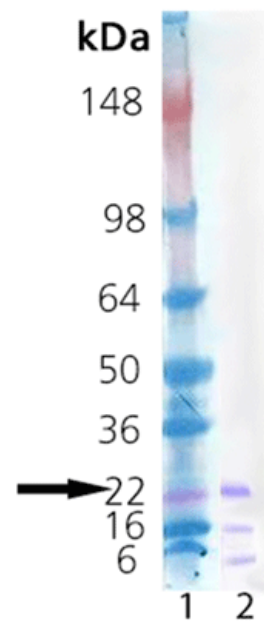
| | |
|-----------------|------|
| ENZ-PRT189-0010 | 10µg |
|-----------------|------|

Manuals, SDS & CofA

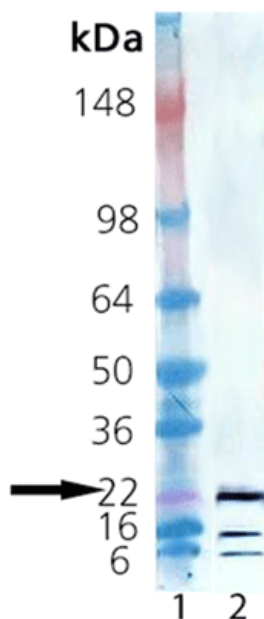
[View Online »](#)



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](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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