

# β-Nerve growth factor (human), (recombinant)

Human β-NGF is responsible for survival, differentiation and maturation of sympathetic and sensory neurons.

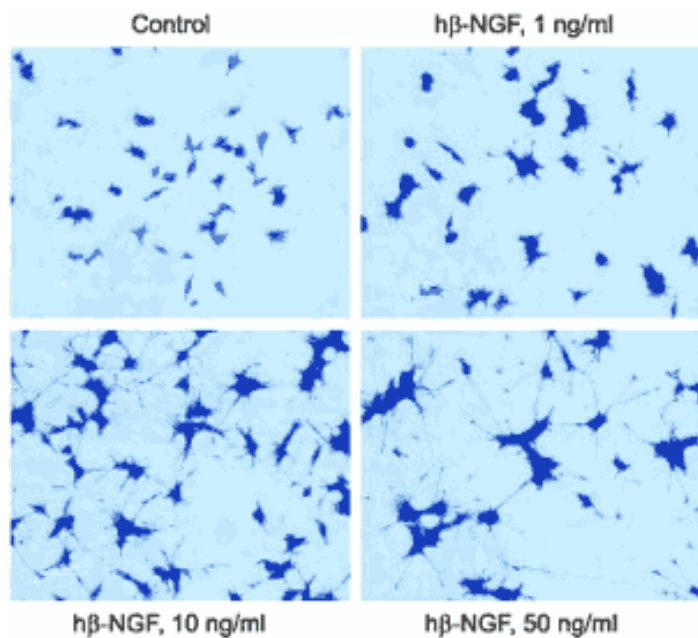
## Ordering Information

[Order Online »](#)

ALX-521-005-C005	5μg
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** Neurite outgrowth in PC12 cells. **Method:** PC12 cells were grown on collagen coated 24-well plates in DMEM medium supplemented with 6% fetal calf serum and 6% horse serum for 24 hours.  $\beta$ -Nerve Growth Factor (human) (rec.) (Prod. No. [ALX-521-005](#)) was added at the indicated concentrations. The culture media was changed every 2-3 days. The development of neurites over a period of 5 days was visualized by methylene blue staining.””” $\beta$ -Nerve growth factor (human)

## Handling & Storage

Use/Stability	Lyophilized protein is stable for 2-3 weeks when stored at room temperature and for 6-9 months when stored at -20°C. Reconstituted protein is stable for up to 1 week when stored at +4°C and for 4-6 weeks when stored at -80°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	β-NGF
Biological Activity	Neurotrophic activity was bioassayed using rat pheochromocytoma PC12 cell cultures over a period of 5-14 days. Activity is in the range of 1-50ng/ml.
CAS	86923-98-0
Formulation	Lyophilized. No additives.
MW	~27kDa (dimer).
Purity	≥98% (HPLC)
Quantity	5µg (determined by amino acid analysis).
Reconstitution	Reconstitute in water or DMEM to a final concentration of at least 10µg/ml and store at -80°C before use.
Source	Produced in <i>E. coli</i> . Dimer of two identical subunits (119 aa) linked together by very strong hydrophobic interactions.
UniProt ID	P01138



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-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@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)

Last modified: May 29, 2024