

β -Nerve growth factor (mouse), (recombinant)

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

Ordering Information

[Order Online »](#)

ENZ-PRT264-0005	5 μ g
-----------------	-----------

Manuals, SDS & CofA

[View Online »](#)

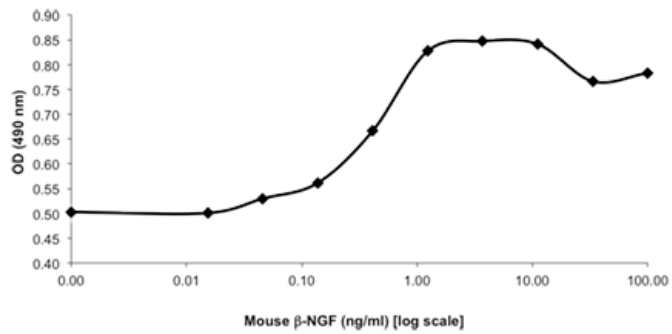


Figure 1: Murine b-NGF induced proliferation of TF-1 cells. Serial dilutions of Mouse β -NGF, starting at 100 ng/mL, were added to TF-1 cells growing in GM-CSF free media. Cell proliferation can be measured and the linear portion of the curve can be used to calculate the ED50.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA solution.
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	β-NGF
Endotoxin Content	≤1 EU/μg protein (LAL test)
Formulation	Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 0.1% Trifluoroacetic Acid (TFA)
MW	~13.6/27.2 kDa (dimer).
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 121/242 amino acids.
UniProt ID	P01139



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