

Nerve growth factor 2.5S (mouse), (grade II)

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-521-0070	25µg
ALX-521-007-C100	100µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 2-3 weeks when stored at room temperature. After reconstitution, stable for up to 1 week when stored at +4°C and for 4-6 weeks at -80°C.
Handling	Do not reuse medium containing mNGF. Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	NGF 2.5S
Biological Activity	Neurotrophic activity was bioassayed using rat pheochromocytoma cell PC12 cultures over a period of 7-14 days. Activity is in the range of 0.5-50ng/ml. Recommended concentration to be used <i>in vitro</i> for maintenance of sympathetic and sensory nerve cultures is 50ng/ml medium.
CAS	93928-24-6
Formulation	Lyophilized.
MW	~26kDa; homodimer.
Purity	≥95% (HPLC)
Quantity	Concentration is estimated spectroscopically from its extinction coefficient and amino acid analysis.
Reconstitution	Dilute in sterile water or DMEM to a concentration of at least 10µg/ml and store at -80°C before use.
Source	Isolated from mouse submaxillary glands by a modification of the procedure described by T. Darling, et al.; Cell Culture Meth. Mol. Cell Biol. 4 , 79 (1984).

Last modified: March 10, 2025



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)