

Nerve growth factor receptor (human):Fc (human), (recombinant)

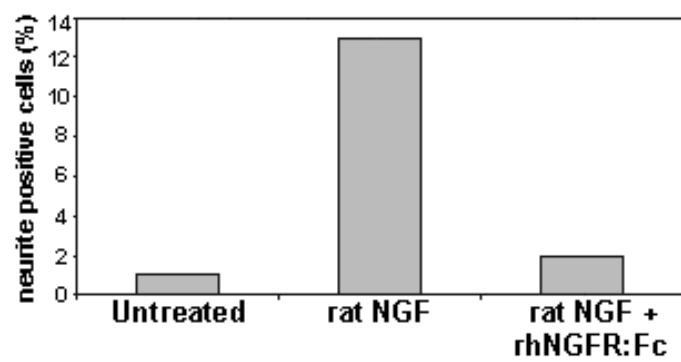
Ordering Information

[Order Online »](#)

ALX-522-044-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C. After reconstitution, it is recommended to make appropriate aliquots and to store them at -20°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	NGFR, TNFRSF 16, Low-affinity nerve growth factor receptor, p75NTR, p75 neurotrophin receptor, CD271
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
MW	~80kDa (under reducing conditions).
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute in 50μl sterile water. Further dilutions can be made with medium containing 5% fetal calf serum.
Source	Produced in HEK 293 cells. The cysteine-rich domain of human NGFR (nerve growth factor receptor; p75 ^{NTR}) (aa 1-250) is fused to the Fc portion of human IgG1.
Specificity	Binds human, mouse and rat NGF (nerve growth factor).
Technical Info / Product Notes	Historical data has shown that NGFR inhibits rat NGF-mediated activation of rat PC12 cells.
UniProt ID	P08138



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)