

Netrin-1 (human), (recombinant)

The Netrins (whose name means “one who guides” in Sanskrit) are a family of laminin-related secreted protein. In mice, human and rat, netrin-1, netrin-3/NTL2, Netrin-4 and G-Netrins have been cloned. Netrin-1 is predominantly expressed by floor plate cells as commissural axons extend toward the ventral midline. Netrin-1 induces axonal outgrowth, axon orientation and neuronal migration in neuronal development. It is chemoattractive for some neurons and chemorepellent for others. To carry its functions, netrin-1 interacts with specific receptors: DCC/Neogenin and UNC-5. These receptors belong to the family of “dependence receptors” that induce apoptosis in the absence of their respective ligand, and promote survival when bound to their ligand. The Netrin-1 mediated survival signal inhibits p53-induced apoptosis. Netrin-1 is also expressed in organs such as pancreas, lung, bowel, bone and mammary gland. In these tissues, netrin-1 has different roles such as invasion of epithelial cells, leukocytes migration, cell adhesion or angiogenesis. Netrin-1, as a survival factor, also contributes to the control of tumorigenesis. Historical results have shown that netrin-1 induces axon outgrowth and inhibits DCC- and UNC5H2-mediated apoptosis in HEK 293 cells.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|------------------|--------|
| ALX-522-100-C010 | 10µg |
| ALX-522-100-2010 | 2x10µg |

Manuals, SDS & CofA

[View Online »](#)

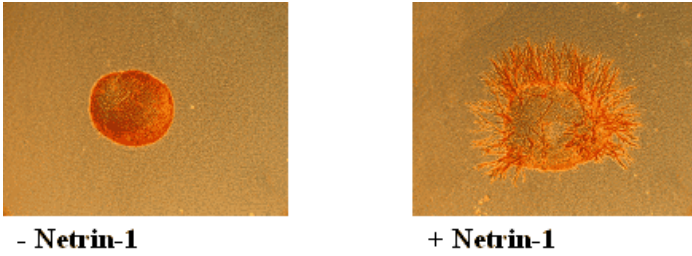
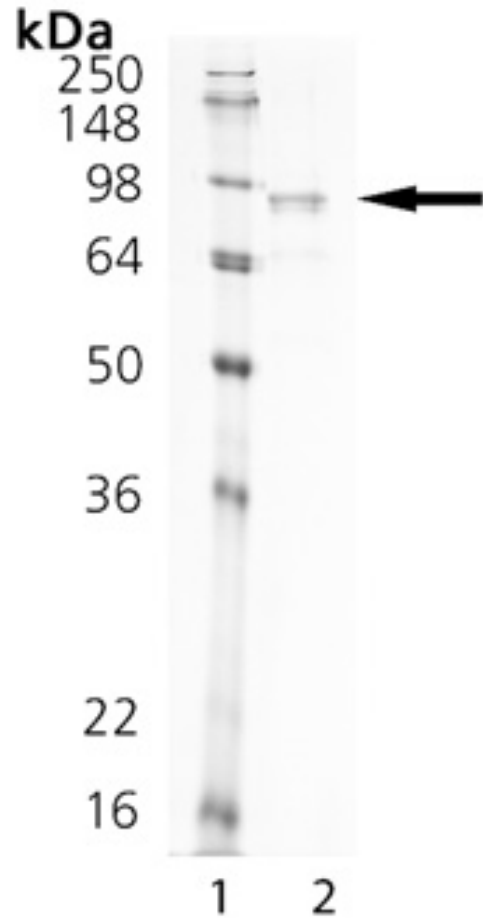
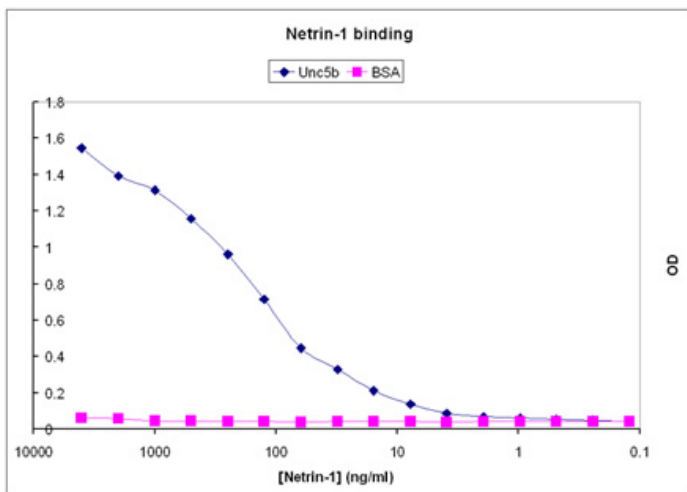


Figure: Netrin-1 (human) (rec.) (Prod. No. [ALX-522-100](#)) mediated commissural axon outgrowth. **Method:** Dorsal spinal cord were dissected out from E13 rat embryos and cultured in collagen matrix in the presence or absence of Netrin-1 (250 ng/ml). Axons were then stained with an anti- β -tubulin antibody. *Picture courtesy of Véronique Corset, Centre Léon Bérard, Lyon.*



SDS-PAGE analysis of Netrin-1 (human), (recombinant): Lane 1: MW Marker, Lane 2: 1 μ g Netrin-1.



ELISA analysis of Netrin-1 (human), (recombinant): Each well is coated with either 500ng Unc5b (Prod. No. ALX-522-095) or 2% BSA. Netrin-1 is titrated in 2-fold series and detected by anti-FLAG antibody.

Handling & Storage

| | |
|--------------------------|--|
| Use/Stability | Stable for at least 6 months after receipt when stored at -20°C. |
| Handling | Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|---------------------------------------|--|
| Application Notes | ELISA: binds to UNC5B at 15-1000 ng/ml. |
| Concentration | 0.1mg/ml after reconstitution. |
| Formulation | Lyophilized. Contains PBS. |
| MW | ~85kDa |
| Purity | ≥90% (SDS-PAGE) |
| Purity Detail | Purified by multi-step chromatography. |
| Reconstitution | Reconstitute with 100µl sterile distilled water. Resuspended protein should be used immediately upon reconstitution to avoid loss in activity due to non-specific binding to the vial. Further dilutions should be made with medium containing 5% fetal calf serum or 0.5% BSA. |
| Source | Produced in HEK 293 cells. The extracellular domain of human netrin-1 (aa 28-604) is fused at the N-terminus to a FLAG [®] -tag. |
| Specificity | Tests positive for binding to human Unc5b by ELISA. Previous lots produced from the same cell line bind to human, mouse and rat receptors UNC5H and DCC. |
| Technical Info / Product Notes | FLAG is a registered trademark of Sigma-Aldrich Co. |
| UniProt ID | O95631 |

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-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)