

Recombinant Human NT3

NT-3 is an important factor of the NGF family of proteins like β -NGF and BDNF. NT-3 is thought to promote the survival or differentiation of existing and new neurons in the central nervous system and synapses, and these functions are thought to be communicated by TrkC, a receptor tyrosine kinase thought to induce NT-3 specific signaling.

Materials Provided

Catalog Number:	92-1025
Quantity Supplied:	1 vial(s), 2 μ g per vial
Reconstitution Buffer:	1 mL

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	27200 g/mol
Source:	E. coli
Purity:	\geq 98% by SDS-PAGE and HPLC
Endotoxin Level:	< 0.1 ng/ μ g (<1EU/ μ g).
Formulation:	Sterile filtered through a 0.2 micron filter. Lyophilized with no additives
Storage:	\leq -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	0.1 mg/mL
Reconstitution Volume:	0.020 mL
Reconstitution Solvent:	Reconstitution Buffer

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com