

Recombinant Human TNF α

Tumor necrosis factor alpha (TNF α) is an endogenous pyrogen cytokine that plays a key part in regulating immune cells, immune response to infection, and systemic inflammation. Also regulates cell proliferation, differentiation, apoptosis, lipid metabolism, and coagulation. Is implicated in multiple autoimmune disorders due to its role in inflammation, including inflammatory bowel disease, rheumatoid arthritis, ankylosing spondylitis, and psoriasis. Agonist of TNF receptor 1 and 2.

Materials Provided

Catalog Number:	92-1097
Quantity Supplied:	1 vial(s), 10 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:	17400 g/mol
Source:	E. coli
Purity:	\geq 98% by SDS-PAGE and HPLC
Endotoxin Level:	Endotoxin level is < 0.1 ng/ μ g (<1EU/ μ g).
Formulation:	N/A
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:	25 μ M
Reconstitution Volume:	22.9 μ L
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