

TNF- α (human), (recombinant)

Tumor necrosis factor-alpha (TNF α), also known as cachectin, is a 17.5 kDa, 157 amino acid member of the TNF superfamily of cytokines that is a potent lymphoid factor with effects on a wide range of target cells. Active TNF α is produced as both soluble and membrane-anchored trimers by macrophages, NK cells, and T- and B-lymphocytes. TNF exerts proinflammatory signals via binding and inducing trimerization of TNF-receptor 1 (TNFR1) expressed on most normal and transformed cells, or to TNF-receptor 2 (TNFR2), expressed on endothelial and most immune cells. TNF signaling regulates hematopoiesis differentiation, endothelial cell activation, apoptosis, lipid metabolism, tumor progression, and immune surveillance, and dysregulation of TNF or its receptors is implicated in numerous disease states including cancer, osteoporosis, autoimmune diseases, diabetes, and atherosclerosis.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-908-066-010	10 μ g
-----------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Upon reconstitution TNF- α should be stored at 4°C between 2-7 days and for future use below -20°C. For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA). Please prevent freeze-thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Tumor necrosis factor- α , TNFSF 2
Application Notes	Western blot control.
Formulation	Lyophilized. Sterile filtered liquid, pH 7.2, containing 20mM PB and 100mM sodium chloride
Purity	≥95% (SDS-PAGE; RP-HPLC)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> .
UniProt ID	P01375



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com