TNF-α (human), (recombinant)

Tumor necrosis factor is a multifunctional proinflammatory cytokine, with effects on lipid metabolism, coagulation, insulin resistance, endothelial function, autoimmune disease and allergic responses. TNF- β shares 35% identity and 50% homology in amino acid sequence with the TNF- α and it has been shown that the 2 TNFs share a common receptor on tumor cells.

Ordering Information

Order Online »

BML-SE303-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tumor necrosis factor-α, TNFSF 2

Formulation Sterile, lyophilized powder.

MW 17.5 kDa

Source Produced in *E. coli* as a single, non-glycosylated

polypeptide.

UniProt ID P01375

Last modified: May 29, 2024

