## Fc (human):TNF-α (soluble) (human), (recombinant)

**Ordering Information** 

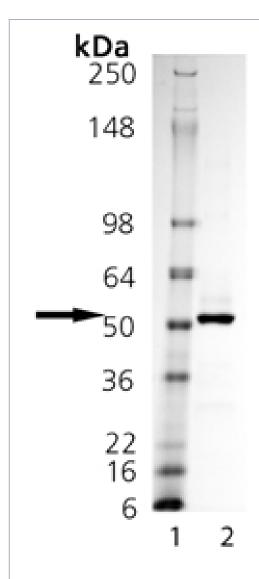
Order Online »

ALX-522-087-C010

10µg

Manuals, SDS & CofA

View Online »



SDS-PAGE analysis of Fc (human):TNF- $\alpha$  (soluble) (human), (recombinant): Lane 1: MW Marker, Lane 2: 1 $\mu$ g TNF- $\alpha$ .

## **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**

Alternative NameTumor necrosis factor α, TNFSF 2

**Concentration** 0.1mg/ml after reconstitution.

Endotoxin Content <0.1EU/µg purified protein (LAL test; Associates of CAPE

COD Inc.).

**Formulation** Lyophilized. Contains PBS.

MW ~48kDa (calculated), ~52kDa (observed)

Purity ≥90% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Reconstitution Reconstitute with 100µl sterile water. Further dilutions

should be made with medium containing 5% fetal calf

serum or a carrier protein.

**Source** Produced in CHO cells. The extracellular domain of human

TNF- $\alpha$  (aa 85-233) is fused at the N-terminus to the Fc portion of human IgG1 and a linker peptide (20 aa).

Technical Info / Product Notes Historical data has shown that TNFα induces killing of

WEHI-164 cells in a concentration range of 0.5-2ng/ml

(does not require an enhancer).

UniProt ID P01375

Last modified: May 29, 2024



info-

eu@enzolifesciences.com