

**$\alpha$ -Tumor Necrosis Factor  $\alpha$  (anti-TNF $\alpha$ )**  
**anti-mouse Tumor Necrosis Factor  $\alpha$**   
**rabbit, polyclonal**

Cat. No.	Amount
ABD-025	50 $\mu$ l

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -20°C

**Avoid freeze / thaw cycles**

**Form**

Liquid. Supplied in Serum.

**Description**

Polyclonal antibody anti-mTNF- $\alpha$  is a mouse TNF- $\alpha$  specific rabbit IgG antibody.

**Antigen**

E. coli expressed mouse TNF- $\alpha$

**Applications**

Anti-mTNF- $\alpha$  may be used for Western blotting and ELISA at a dilution of 1:10000-1:50000

**Selected References:**

Hassoun E et al. (2000) Ricin-induced toxicity in the macrophage J744A.1 cells: the role of TNF-alpha and the modulation effects of TNF-alpha polyclonal antibody. *J. Biochem. Mol. Toxicol.* **14**:95.