

Instruction Manual

| | |
|--------------------------------|--|
| Catalog Number | PK-AB718-1131 |
| Synonyms | TACE Antibody: adam17 |
| Description | Tumor-necrosis factor- α is a proinflammatory cytokine and contributes to a variety of inflammatory disease responses and programmed cell death. TNF- α is synthesized as a 26K type II membrane-bound precursor that is cleaved by a convertase to generate secreted 17K mature TNF- α . TNF- α converting enzyme (TACE) protein was recently purified and the human and mouse TACE cDNAs were cloned by several groups separately. TACE is a membrane-bound metalloprotease-disintegrin in the family of mammalian ADAM (for a disintegrin and metalloprotease). TACE also processes other cell surface proteins, including TNF receptor, TGF α , the L-selectin adhesion molecule, and alpha-cleavage of amyloid protein precursor (APP). TACE mRNA is expressed in a variety of human and murine tissues. TACE was selected as one of the few targets in cytokine activation by the Eighth International Conference of the Inflammation Research Association. |
| Quantity | 100 μ g |
| Source / Host | Rabbit |
| Immunogen | Rabbit polyclonal TACE antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human TACE. This sequence differs from those of mouse and rat TACE by one amino acid. |
| Purification Method | Affinity chromatography purified via peptide column. |
| Clone / IgG Subtype | Polyclonal antibody |
| Species Reactivity | Human, Mouse, Rat |
| Specificity | 80 to 130 kDa bands can be detected, which may represent mature protein, precursor, and glycosylated TACE. |
| Formulation | Antibody is supplied in PBS containing 0.02% sodium azide. |
| Reconstitution | During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 μ l or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap. |
| Storage & Stability | Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures. |
| Applications | E, WB, ICC, IF, FACS !Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually. |
| Images | Available upon request. |
| References | Black RA, Rauch CT, Kozlosky CJ, et al. A metalloproteinase disintegrin that releases tumour-necrosis factor- α from cells. <i>Nature</i> 1997;385:729-733 Moss ML, Jin SL, Milla ME, et al. Cloning of a disintegrin metalloproteinase that processes precursor tumour-necrosis factor- α . <i>Nature</i> 1997;385:733-736 Mizui Y, Yamazaki K, Sagane K, Tanaka I. cDNA cloning of mouse tumor necrosis factor- α converting enzyme (TACE) and partial analysis of its promoter. <i>Gene</i> 1999;233:67-74 Peschon JJ, Slack JL, Reddy P, et al. An essential role for ectodomain shedding in mammalian development. <i>Science</i> 1998;282:1281-4 |
| Images | Available upon request. |
| Related Products | Blocking Peptide, Cat. No. PK-AB718-1131P HeLa Lysate, Cat. No. PK-AB718-1201 APP Antibody (CT), Cat. No. PK-AB718-2136 ADAM10 (CT) Antibody, Cat. No. PK-AB718-2051 |

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.