

Rabbit Anti-Human TNF-alpha Converting Enzyme

Instruction Manual

Catalog Number	PK-AB718-2051
Synonyms	ADAM10 Antibody: KUZ
Description	Proinflammatory cytokine tumor necrosis factor- α (TNF- α) contributes to a variety of inflammatory responses and programmed cell death. Notch receptor and its ligand participate in cell fate decisions during vertebrate development and are associated with several human disorders, including a T-cell lymphoma. TNF- α , notch and its ligand delta are all membrane-bound molecules, which are cleaved by proteases to release mature proteins or functional receptor. ADAM10, a metalloprotease-disintegrin in the family of mammalian ADAM (for a disintegrin and metalloprotease), was recently identified to cleave TNF- α , notch and its ligand delta. The genes encoding human, mouse, and bovine ADAM10 were recently cloned and designated ADAM 10, kuzbanian (KUZ), and MADM, respectively. ADAM10 mRNA is expressed in a variety of human and bovine tissues.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	ADAM10 antibody was raised against a peptide corresponding to amino acids 732 to 748 of human ADAM10. This sequence is identical to those of bovine and rat origins and differs from that of mouse ADAM10 by one amino acid (2,4)
Purification Method	Ion exchange chromatography purified.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	
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 INote: 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. ADAM10 can be used for detection of ADAM10 by Western blot at 1:1000 to 1:2000 dilution. An approximately 85 kDa band can be detected, which may represent precursor. This polyclonal antibody can also detect ADAM10 by immunohistochemistry at 2 μ g/mL. For immunofluorescence start at 10 μ g/mL.
Images	Available upon request.
References	Rosendahl MS, Ko SC, Long DL, et al. Identification and characterization of a pro-tumor necrosis factor- α -processing enzyme from the ADAM family of zinc metalloproteases. <i>J Biol Chem</i> 1997;272:24588-93 Pan D, Rubin GM. Kuzbanian controls proteolytic processing of Notch and mediates lateral inhibition during <i>Drosophila</i> and vertebrate neurogenesis. <i>Cell</i> 1997;90:271-80 Qi H, Rand MD, Wu X, Sestan N, Wang W, Rakic P, Xu T, Artavanis-Tsakonas S. Processing of the notch ligand delta by the metalloprotease Kuzbanian. <i>Science</i> 1999;283:91-4
Images	Available upon request.
Related Products	ADAM10 Peptide, Cat. No. PK-AB718-2051P TACE Antibody, Cat. No. PK-AB718-1131 Jurkat Lysate, Cat. No. PK-AB718-1205

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