

F1A α (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat F1A α (Novel Death Receptor Binding Protein)

Instruction Manual

Catalog Number	PK-AB718-2279
Synonyms	F1A alpha Antibody: F1A alpha
Description	Fas and tumor necrosis factor receptor 1 (TNFR1) are two prototype members in the death receptor family. A novel protein that associates with the intracellular domains of Fas and TNFR1 was recently identified and designated F1A α and FEM1 β . F1A α /FEM1 β is the homologue of <i>C. elegans</i> sex determining protein FEM-1. FEM-1/F1A α is cleaved by CED-3 and caspase. FEM-1/F1A α associates with CED-4 and its mammalian homologue Apaf-1. Overexpression of F1A α induces apoptosis. F1A α is therefore a novel member of the death receptor associated protein that mediates apoptosis. F1A α is expressed in a variety of human and mouse tissues.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal F1A α antibody was raised against a peptide (DINYQDQIPRTLEE) corresponding to amino acids 609 to 622 of human F1A α . This sequence is identical to the corresponding sequence of mouse FEM1a.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
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, IHC, 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. F1Ab antibody can be used for detection of F1Ab by Western blot at 1 μ g/mL. An approximately 70 kDa band can be detected. Antibody can also be used for immunohistochemistry starting at 5 μ g/mL. For immunofluorescence start at 20 μ g/mL.
Images	Available upon request.
References	Chan SL, Tan KO, Zhang L, Yee KS , Ronca F, Chan MY, Yu VC. F1A α , a death receptor-binding protein homologous to the <i>Caenorhabditis elegans</i> sex-determining protein, FEM-1, is a caspase substrate that mediates apoptosis. <i>J Biol Chem.</i> 1999;274(45):32461-8. Ventura-Holman T, Maher JF. Sequence, organization, and expression of the human FEM1B gene. <i>Biochem Biophys Res Commun</i> 2000;267(1):317-20 Chan SL, Yee KS , Tan KM, Yu VC. The <i>Caenorhabditis elegans</i> sex determination protein FEM-1 is a CED-3 substrate that associates with CED-4 and mediates apoptosis in mammalian cells. <i>J Biol Chem.</i> 2000;275(24):17925-8. (WD0102)
Images	Available upon request.
Related Products	F1A α Peptide, Cat. No. PK-AB718-2279P Mouse liver Tissue Lysate, Cat. No. PK-AB718-1404

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