

Instruction Manual

Catalog Number	PK-AB718-2273
Synonyms	Toso Antibody: Toso, Regulator of Fas-induced apoptosis, Fas apoptotic inhibitory molecule 3, FAIM3
Description	Apoptosis is an important process by which normal tissue homeostasis and function are maintained. One of the major signals that regulate this process is mediated by the activation of the Fas receptor by its ligand. This leads to the formation of a Fas-associated death domain (FADD)-containing death-inducing signaling complex and the activation of caspase-8, which in turn activates downstream effector caspases, such as caspase-3 and -7. Recent experiments have shown that overexpression of Toso, a novel regulator of Fas-induced apoptosis in lymphoid cells, in Jurkat cells as well as transgenic mice render these cells resistant to Fas-induced apoptosis but not to TRAIL-induced apoptosis. Furthermore, Toso was found to associate with FADD, suggesting that Toso functions by disrupting the formation of the death-inducing signaling complex. Despite its predicted molecular weight, Toso often migrates at 60 kDa in SDS-PAGE.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Toso antibody was raised against a 13 amino acid peptide from near the carboxy terminus of human Toso.
Purification Method	Affinity chromatography purified via peptide column.
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 µl 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 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. Toso antibody can be used for detection of Toso by Western blot at 1 µg/mL. Despite its predicted molecular weight, Toso often migrates at 60 kDa in SDS-PAGE.
Images	Available upon request.
References	Curtin JF and Cotter TG. Live and let die: regulatory mechanisms in Fas-mediated apoptosis. Cell Signal. 2003; 15:983-92. Hitoshi Y, Lorens J, Kitada S-I, et al. Toso, a cell surface, specific regulator of Fas-induced apoptosis in T cells. Immunity 1998; 8:461-71. Song Y and Jacob CO. The mouse cell surface protein Toso regulates Fas/Fas ligand-induced apoptosis through its binding to Fas-associated death domain. J
Images	Available upon request.
Related Products	Toso Peptide, Cat. No. PK-AB718-2273P Lung Lysate, Cat. No. PK-AB718-1302

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