

Instruction Manual

Catalog Number	PK-AB577-3070
Description	Fas/Apo-1/CD95 is a cell surface type-I membrane glycoprotein that is part of the TNFR family that includes TNFR1, TNFR2, CD27, CD30, and CD40. Binding of Fas-ligand (FasL) to Fas or crosslinking of Fas by anti-Fas antibodies leads to apoptosis in Fas expressing cells. Both Fas and TNF-R contain the 70 amino acids cytoplasmic "death domain" that is responsible for transmitting signal for apoptosis.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide corresponding to residues surrounding amino acid 32 of mouse Fas (Internal ID BV-531)
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	The purified antibody detects a 55 kDa of glycosylated Fas, a 35 kDa native form of Fas and a 29 kDa of cleaved Fas fragment on SDS-PAGE immunoblots.
Formulation	100 µg (0.5 mg/ml) affinity purified polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 0.01% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and freeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	The antibody can be used for Western blotting (0.5-4 µg/ml). However, the optimal conditions should be determined individually. Blocking peptide (PK-AB577-3070P) is available separately.

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