

Apaf-1 antibody (pAb)

Rabbit Anti-
Human/Mouse/Rat/Bovine/Rabbit/Hamster/Monkey/Horse/Guinea Pig

Instruction Manual

Catalog Number	PK-AB577-3018
Description	During apoptosis, Apaf-1 (apoptosis protease activating factor 1) binds to cytochrome c (Apaf-2), a reaction that requires dATP, and this leads to the formation of a large multimeric complex. The complex recruits pro-caspase-9, and activates caspase-9 (Apaf-3). Activated caspase-9 in turn cleaves and activates caspase-3, one of the proteases responsible for the proteolytic cleavage of many key proteins in apoptosis. Apaf-1 is ubiquitously expressed in human tissues.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 1161 of human Apaf-1.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat, bovine, rabbit, hamster, guinea pig, monkey, horse
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.2 mg/ml) immunoaffinity purified, rabbit anti-Apaf-1 polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.05% 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 refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	The antibody can be used for Western blotting (0.5-4 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far.

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