

AATF Antibody [#C11778]

Catalog Number:	C11778
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NY61
Other Names:	CHE-1, CHE1, DED, apoptosis antagonizing transcription factor
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from N-terminal of human AATF.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AATF antibody detects endogenous levels of total AATF protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Lindfors K., Biochem. Biophys. Res. Commun. 276:660-666(2000). Fanciulli M., FASEB J. 14:904-912(2000). Di Padova M., J. Biol. Chem. 278:36496-36504(2003).