

Caspase 2 (Ab-140) Antibody [#B1026]

Catalog Number:	B1026
Concentration:	1mg/ml
Swiss-Prot No.:	P42575
Other Names:	CASP-2; Caspase-2 precursor; EC 3.4.22.-; EC 3.4.22.55; ICH- 1L/1S; ICH-1 protease; ICH1
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 non-phosphopeptide derived from human Caspase 2 around the phosphorylation site of serine 140 (R-L-S ^P -T-D).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Caspase 2 (Ab-140) antibody detects endogenous levels of total Caspase 2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Wang L., Cell 78:739-750(1994). Rieder M.J., Submitted (JAN-2003) to the EMBL/GenBank/DDBJ databases. Hillier L.W., Nature 424:157-164(2003).