

API-5 Antibody [#C12052]

Catalog Number:	C12052
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BZZ5
Other Names:	AAC-11, API-5, API5-like 1, API5L1, FIF, XAGL protein, apoptosis inhibitor 5, fibroblast growth factor 2-interacting factor, fibroblast growth factor 2-interacting factor 2, migration-inducing protein MIG8
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 C-terminal of human API-5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	API-5 antibody detects endogenous levels of total API-5 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Tewari M., Cancer Res. 57:4063-4069(1997). Van den Berghe L., Mol. Endocrinol. 14:1709-1724(2000). Olsen J.V., Cell 127:635-648(2006).