

ATP6V1H Antibody [#C19484]

Catalog Number:	C19484
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UI12
Other Names:	Vacuolar proton pump subunit H; EC 3.6.3.14; V-ATPase subunit H; V-ATPase 50/57 kDa subunits; Vacuolar proton pump subunit SFD; VMA13; Nef-binding protein 1; NBP1; ATP6V1H; CGI-11
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 internal of human ATP6V1H.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ATP6V1H antibody detects endogenous levels of total ATP6V1H protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Liu B., Submitted (DEC-1998) to the EMBL/GenBank/DDBJ databases. Lu X., Immunity 8:647-656(1998). Hu R.-M., Proc. Natl. Acad. Sci. U.S.A. 97:9543-9548(2000). Lai C.-H., Genome Res. 10:703-713(2000).