

VHL Antibody [#B8450]

Catalog Number:	B8450
Concentration:	1mg/ml
Swiss-Prot No.:	P40337
Other Names:	G7 protein; pVHL; Von Hippel-Lindau disease tumor suppressor; von Hippel-Lindau syndrome gene
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 VHL.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	VHL antibody detects endogenous levels of total VHL protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Latif F., Science 260:1317-1320(1993). Richards F.M., Hum. Mol. Genet. 5:639-644(1996). Pause A., Proc. Natl. Acad. Sci. U.S.A. 94:2156-2161(1997).