

CA12 Antibody [#C14935]

Catalog Number:	C14935
Concentration:	1mg/ml
Swiss-Prot No.:	O43570
Other Names:	CA-XII; CAH12; carbonate dehydratase XII; carbonic anhydrase 12; CAXII; EC 4.2.1.1
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 CA12.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CA12 antibody detects endogenous levels of total CA12 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%); Rat (Identities = 92%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Tuereci O., Proc. Natl. Acad. Sci. U.S.A. 95:7608-7613(1998). Ivanov S.V., Proc. Natl. Acad. Sci. U.S.A. 95:12596-12601(1998). Kalline N., Submitted (MAY-2003) to the EMBL/GenBank/DDBJ databases.