

CAD Antibody [#B8308]

Catalog Number:	B8308
Concentration:	1mg/ml
Swiss-Prot No.:	P27708
Other Names:	CAD protein; EC 2.1.3.2; EC 3.5.2.3; EC 6.3.5.5; glutamine-dependent carbamoyl-phosphate synthase; PYR1
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 CAD.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CAD antibody detects endogenous levels of total CAD protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 91%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Iwahana H., Biochem. Biophys. Res. Commun. 219:249-255(1996). Davidson J.N., DNA Cell Biol. 9:667-676(1990). Olsen J.V., Cell 127:635-648(2006).