

AMPD1 Antibody [#C14443]

Catalog Number:	C14443
Concentration:	1mg/ml
Swiss-Prot No.:	P23109
Other Names:	AMP deaminase 1; EC 3.5.4.6; Myoadenylate deaminase; AMP deaminase isoform M
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 AMPD1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AMPD1 antibody detects endogenous levels of total AMPD1 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:	Gregory S.G., Nature 441:315-321(2006). Morisaki H., Hum. Mutat. 16:467-472(2000). Sjoebloom T., Science 314:268-274(2006).