

ATP-Citrate Lyase (Ab-454) Antibody [#B0618]

Catalog Number:	B0618
Concentration:	1mg/ml
Swiss-Prot No.:	P53396
Other Names:	ATP-citrate (pro-S-)-lyase, ATP-citrate lyase, Citrate cleavage enzyme, EC 2.3.3.8
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 non-phosphopeptide derived from human ATP-Citrate Lyase around the phosphorylation site of serine 454 (T-A-S ^P -F-S).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ATP-Citrate Lyase (Ab-454) antibody detects endogenous levels of total ATP-Citrate Lyase protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Elshourbagy N.A., Eur. J. Biochem. 204:491-499(1992). Lord K.A., Protein Expr. Purif. 9:133-141(1997). Beausoleil S.A., Proc. Natl. Acad. Sci. U.S.A. 101:12130-12135(2004).