

## ACC1 (Ab-80) Antibody [#B0405]

<b>Catalog Number:</b>	B0405
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q13085
<b>Other Names:</b>	ACAC, ACACA, ACC-alpha, ACCA, Acetyl-CoA carboxylase 1, EC 6.3.4.14, EC 6.4.1.2
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-phosphopeptide derived from human ACC1 around the phosphorylation site of serine 79 (S-M-S <sup>P</sup> -G-L).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ACC1 (Ab-80) antibody detects endogenous levels of total ACC1 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
<b>References:</b>	Naomoto Harada, Mol. Cell. Biol., Mar 2007; 27: 1881 - 1888. M. Joachimiak, PNAS, Sep 1997; 94: 9990. D. A. Somers, Plant Physiology, Mar 1993; 101: 1097.