

## NR1I3 Antibody [#C11053]

<b>Catalog Number:</b>	C11053
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q14994
<b>Other Names:</b>	CAR, Constitutive androstane receptor, Orphan nuclear receptor MB67, Orphan nuclear receptor NR1I3
<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 peptide derived from internal of human NR1I3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	NR1I3 antibody detects endogenous levels of total NR1I3 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Baes M., Mol. Cell. Biol. 14:1544-1552(1994). Choi H.S., J. Steroid Biochem. Mol. Biol. 56:23-30(1996). Xu R.X., Mol. Cell 16:919-928(2004).