

## CCKBR Antibody [#G069]

<b>Catalog Number:</b>	G069
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P32239
<b>Other Names:</b>	CCK-B receptor; CCK-BR; CCKRB; Cholecystokinin-2 receptor; GASR; Gastrin/cholecystokinin type B receptor
<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 CCKBR.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CCKBR antibody detects endogenous levels of total CCKBR protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Pisegna J.R., Biochem. Biophys. Res. Commun. 189:296-303(1992). Lee Y.-M., J. Biol. Chem. 268:8164-8169(1993). Ito M., J. Biol. Chem. 268:18300-18305(1993).