

## CIB3 Antibody [#C11530]

<b>Catalog Number:</b>	C11530
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q96Q77
<b>Other Names:</b>	Calcium and integrin-binding family member 3, Kinase-interacting protein 3, KIP 3, CIB3, KIP3
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from C-terminal of human CIB3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CIB3 antibody detects endogenous levels of total CIB3 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:20000
<b>References:</b>	Hayashi A., Submitted (NOV-2000) to the EMBL/GenBank/DDBJ databases. The MGC Project Team, Genome Res. 14:2121-2127(2004).