

## BCAR3 Antibody [#C10193]

<b>Catalog Number:</b>	C10193
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O75815
<b>Other Names:</b>	SH2-containing protein Nsp2, breast cancer antiestrogen resistance 3
<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 BCAR3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	BCAR3 antibody detects endogenous levels of total BCAR3 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	van Agthoven T., EMBO J. 17:2799-2808(1998). Lu Y., J. Biol. Chem. 274:10047-10052(1999). Clark H.F., Genome Res. 13:2265-2270(2003).