

## NCAM2 Antibody [#C16899]

<b>Catalog Number:</b>	C16899
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O15394
<b>Other Names:</b>	N-CAM 2; NCAM21; Neural cell adhesion molecule 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 peptide derived from internal of human NCAM2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	NCAM2 antibody detects endogenous levels of total NCAM2 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%); Rat (Identities = 92%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Paoloni-Giacobino A., Genomics 43:43-51(1997). Hattori M., Nature 405:311-319(2000). Zhang H., Nat. Biotechnol. 21:660-666(2003).