

## BRD3 Antibody [#C11064]

<b>Catalog Number:</b>	C11064
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q15059
<b>Other Names:</b>	Bromodomain-containing protein 3, KIAA0043, RING3-like protein, RING3L
<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 BRD3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	BRD3 antibody detects endogenous levels of total BRD3 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Nomura N., DNA Res. 1:223-229(1994). Humphray S.J., Nature 429:369-374(2004). Yu L.-R., J. Proteome Res. 6:4150-4162(2007).