

## Vitamin D3 Receptor (Ab-51) Antibody [#B8205]

<b>Catalog Number:</b>	B8205
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P11473
<b>Other Names:</b>	1,25-dihydroxyvitamin D3 receptor; NR111; vitamin D receptor; vitamin D3 receptor
<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 non-phosphopeptide derived from human Vitamin D3 Receptor around the phosphorylation site of serine 51 (R-R-S <sup>P</sup> -M-K).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Vitamin D3 Receptor (Ab-51) antibody detects endogenous levels of total Vitamin D3 Receptor protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Baker A.R., Proc. Natl. Acad. Sci. U.S.A. 85:3294-3298(1988). Goto H., Biochim. Biophys. Acta 1132:103-108(1992). Miyamoto K., Mol. Endocrinol. 11:1165-1179(1997).