

Vitamin D Receptor (Ab-208) Antibody [#B1245]

Catalog Number:	B1245
Concentration:	1mg/ml
Swiss-Prot No.:	P11473
Other Names:	1,25-dihydroxyvitamin D3 receptor, NR1H1, vitamin D receptor, vitamin D3 receptor
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human Vitamin D Receptor around the phosphorylation site of serine 208 (D-L-S ^P -E-E).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Vitamin D Receptor (Ab-208) antibody detects endogenous levels of total Vitamin D Receptor protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Goto H., Biochim. Biophys. Acta 1132:103-108(1992). Miyamoto K., Mol. Endocrinol. 11:1165-1179(1997). Yu X.-P., J. Biol. Chem. 266:7588-7595(1991).