

## Retinoic Acid Receptor $\alpha$ (Ab-77) Antibody [#B8217]

<b>Catalog Number:</b>	B8217
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P10276
<b>Other Names:</b>	NR1B1; RAR-alpha; Retinoic acid receptor alpha; retinoic acid receptor, alpha; RRA
<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 non-phosphopeptide derived from human Retinoic Acid Receptor $\alpha$ around the phosphorylation site of serine 77 (P-P-S <sup>P</sup> -P-P).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Retinoic Acid Receptor $\alpha$ (Ab-77) antibody detects endogenous levels of total Retinoic Acid Receptor $\alpha$ 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: 1:10000
<b>References:</b>	Giguere V., Nature 330:624-629(1987). Chen Z., Proc. Natl. Acad. Sci. U.S.A. 91:1178-1182(1994). Hjalt T.A.H., Mamm. Genome 10:528-529(1999).