

## ARRB1 Antibody [#C12056]

<b>Catalog Number:</b>	C12056
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P49407
<b>Other Names:</b>	ARR1, arrestin 2, arrestin beta 1, beta-arrestin-1
<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 C-terminal of human ARRB1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ARRB1 antibody detects endogenous levels of total ARRB1 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Parruti G., J. Biol. Chem. 268:9753-9761(1993). Yu Q.M., Submitted (AUG-1998) to the EMBL/GenBank/DDBJ databases. MGC Project Team, Genome Res. 14:2121-2127(2004).