

ARRD1 Antibody [#C12059]

Catalog Number:	C12059
Concentration:	1mg/ml
Swiss-Prot No.:	Q8N5I2
Other Names:	ARRD1; arrestin domain-containing 1
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 peptide derived from internal of human ARRD1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ARRD1 antibody detects endogenous levels of total ARRD1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Humphray S.J., Nature 429:369-374(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004).