

ARRDC3 Antibody [#C12061]

Catalog Number:	C12061
Concentration:	1mg/ml
Swiss-Prot No.:	Q96B67
Other Names:	Arrestin domain-containing protein 3, ARRDC3, KIAA1376
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 C-terminal of human ARRDC3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ARRDC3 antibody detects endogenous levels of total ARRDC3 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Nagase T., DNA Res. 7:65-73(2000). The MGC Project Team, Genome Res. 14:2121-2127(2004).