

ARFIP1 Antibody [#C14558]

Catalog Number:	C14558
Concentration:	1mg/ml
Swiss-Prot No.:	P53367
Other Names:	ADP-ribosylation factor interacting 1; arfaptin-1; ARFP1
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 ARFIP1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ARFIP1 antibody detects endogenous levels of total ARFIP1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
References:	Kanoh H., J. Biol. Chem. 272:5421-5429(1997). Premont R.T., Submitted (JAN-1999) to the EMBL/GenBank/DDBJ databases. The MGC Project Team; Genome Res. 14:2121-2127(2004).