

PARK7 Antibody [#C13096]

Catalog Number:	C13096
Concentration:	1mg/ml
Swiss-Prot No.:	Q99497
Other Names:	BK215D11.1; PARK7
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 PARK7.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PARK7 antibody detects endogenous levels of total PARK7 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Nagakubo D., Biochem. Biophys. Res. Commun. 231:509-513(1997). Gregory S.G., Nature 441:315-321(2006). The MGC Project Team; Genome Res. 14:2121-2127(2004).