

PJA2 Antibody [#C15563]

Catalog Number:	C15563
Concentration:	1mg/ml
Swiss-Prot No.:	O43164
Other Names:	E3 ubiquitin-protein ligase Praja2, EC 6.3.2.-; Praja-2; RING finger protein 131
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 PJA2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PJA2 antibody detects endogenous levels of total PJA2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Ishikawa K., DNA Res. 4:307-313(1997). Nakajima D., Submitted (JAN-2004) to the EMBL/GenBank/DDBJ databases. Ota T., Nat. Genet. 36:40-45(2004).