

PVRL4 Antibody [#C17761]

Catalog Number:	C17761
Concentration:	1mg/ml
Swiss-Prot No.:	Q96NY8
Other Names:	Poliovirus receptor-related protein 4; Nectin-4; Ig superfamily receptor LNIR
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 PVRL4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PVRL4 antibody detects endogenous levels of total PVRL4 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Reymond N., J. Biol. Chem. 276:43205-43215(2001). Ota T., Nat. Genet. 36:40-45(2004). Gregory S.G., Nature 441:315-321(2006).