

## CD234 Antibody [#G081]

<b>Catalog Number:</b>	G081
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q16570
<b>Other Names:</b>	Duffy antigen/chemokine receptor; Fy glycoprotein; GpFy; Glycoprotein D; Plasmodium vivax receptor; CD234 antigen; DARC; FY, GPD
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human CD234.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CD234 antibody detects endogenous levels of total CD234 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IF: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Chaudhuri A., Proc. Natl. Acad. Sci. U.S.A. 90:10793-10797(1993). Tournamille C., Nat. Genet. 10:224-228(1995). Iwamoto S., Blood 85:622-626(1995).