

## ABCA13 Antibody [#C14611]

<b>Catalog Number:</b>	C14611
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q86UQ4
<b>Other Names:</b>	ATP-binding cassette sub-family A member 13
<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 ABCA13.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ABCA13 antibody detects endogenous levels of total ABCA13 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IF: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Prades C., Cytogenet. Genome Res. 98:160-168(2002). Ota T., Nat. Genet. 36:40-45(2004). Schaap F.G., Submitted (APR-2002) to the EMBL/GenBank/DDBJ databases.