

ABCB10 Antibody [#C14619]

Catalog Number:	C14619
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NRK6
Other Names:	ABC transporter 10; ABCBA; ATP-binding cassette sub-family B member 10, mitochondrial; ATP-binding cassette transporter 10; M-ABC2; mitochondrial ATP-binding cassette 2; MTABC2
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 ABCB10.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ABCB10 antibody detects endogenous levels of total ABCB10 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Zhang F., FEBS Lett. 478:89-94(2000). Ito K., Submitted (APR-1998) to the EMBL/GenBank/DDBJ databases. Gregory S.G., Nature 441:315-321(2006).