

GOLT1A Antibody [#C19496]

Catalog Number:	C19496
Concentration:	1mg/ml
Swiss-Prot No.:	C19496
Other Names:	Vesicle transport protein GOT1A; Golgi transport 1 homolog A; hGOT1b; GOLT1A; GOT1B
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 GOLT1A.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GOLT1A antibody detects endogenous levels of total GOLT1A protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:5000
References:	Ota T., Nat. Genet. 36:40-45(2004). Gregory S.G., Nature 441:315-321(2006). Conchon S., EMBO J. 18:3934-3946(1999).