

SFT2D3 Antibody [#C19500]

Catalog Number:	C19500
Concentration:	1mg/ml
Swiss-Prot No.:	Q587I9
Other Names:	Vesicle transport protein SFT2C; SFT2 domain-containing protein 3; SFT2D3
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 SFT2D3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SFT2D3 antibody detects endogenous levels of total SFT2D3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Otsuki T., DNA Res. 12:117-126(2005). Hillier L.W., Nature 434:724-731(2005).