

SLC6A16 Antibody [#C17576]

Catalog Number:	C17576
Concentration:	1mg/ml
Swiss-Prot No.:	Q9GZN6
Other Names:	Orphan sodium- and chloride-dependent neurotransmitter transporter NTT5, Solute carrier family 6 member 16
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 SLC6A16.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SLC6A16 antibody detects endogenous levels of total SLC6A16 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Tewari D.S., Biochim. Biophys. Acta 1209:293-295(1994). Hideyama T., Submitted (OCT-1996) to the EMBL/GenBank/DDBJ databases. Yang D., Biochem. Mol. Biol. Int. 38:957-964(1996).