

ORCTL-2 Antibody [#C11478]

Catalog Number:	C11478
Concentration:	1mg/ml
Swiss-Prot No.:	Q96BI1
Other Names:	Solute carrier family 22 member 18, Organic cation transporter-like protein 2, ORCTL-2, Imprinted multi-membrane spanning polyspecific transporter-related protein 1, Efflux transporter-like protein, Tumor-suppressing subchromosomal transferable fragment candidate gene 5 protein, Tumor-suppressing STF cDNA 5 protein, Beckwith-Wiedemann syndrome chromosomal region 1 candidate gene A protein, p45-Beckwith-Wiedemann region 1 A, p45-BWR1A, Solute carrier family 22 member 1-like, SLC22A18, BWR1A, BWSCR1A, HET, IMPT1, ITM, ORCTL2, SLC22A1L, TSSC5v
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 C-terminal of human ORCTL-2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ORCTL-2 antibody detects endogenous levels of total ORCTL-2 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Lee M.P., Cancer Res. 58:4155-4159(1998). Morisaki H., DNA Res. 5:235-240(1998). Cooper P.R., Genomics 49:38-51(1998).