

SLCO1A2 Antibody [#C18884]

Catalog Number:	C18884
Concentration:	1mg/ml
Swiss-Prot No.:	P46721
Other Names:	OATP; OATP-A; OATP1; OATP1A2; organic anion transporting polypeptide A; organic anion-transporting 1; SLC21A3; SO1A2; SO1A4; sodium-independent organic anion transporter; solute carrier family 21 (organic anion transporter), member 3; solute carrier family 21, member 3; solute carrier organic anion transporter family member 1A2
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 SLCO1A2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SLCO1A2 antibody detects endogenous levels of total SLCO1A2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Kullak-Ublick G.-A., Gastroenterology 109:1274-1282(1995). Koenig J., J. Biol. Chem. 275:23161-23168(2000). Lee W., J. Biol. Chem. 280:9610-9617(2005).