

LRAT Antibody [#C10219]

Catalog Number:	C10219
Concentration:	1mg/ml
Swiss-Prot No.:	O95237
Other Names:	Lecithin retinol acyltransferase, EC 2.3.1.135, Phosphatidylcholine--retinol O-acyltransferase, LRAT
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 LRAT.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LRAT antibody detects endogenous levels of total LRAT protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Ruiz A., J. Biol. Chem. 274:3834-3841(1999). Mondal M.S., Biochemistry 39:5215-5220(2000). Thompson D.A., Nat. Genet. 28:123-124(2001).