

FOXL1 Antibody [#C10939]

Catalog Number:	C10939
Concentration:	1mg/ml
Swiss-Prot No.:	Q12952
Other Names:	ILF, ILF-1, ILF1, cellular transcription factor ILF-1, forkhead box K2, interleukin enhancer-binding 1
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 FOXL1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	FOXL1 antibody detects endogenous levels of total FOXL1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Fang J., Am. J. Hum. Genet. 67:1382-1388(2000). The MGC Project Team, Genome Res. 14:2121-2127(2004). Pierrou S., EMBO J. 13:5002-5012(1994).