

FOXB1/2 Antibody [#C11579]

Catalog Number:	C11579
Concentration:	1mg/ml
Swiss-Prot No.:	Q99853/Q5VYV0
Other Names:	Forkhead box protein B1/2, Transcription factor FKH-5, FOXB1/2, FKH5
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 FOXB1/2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	FOXB1/2 antibody detects endogenous levels of total FOXB1/2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Bingle C.D., Submitted (JUN-1998) to the EMBL/GenBank/DDBJ databases. Cascio S.M., Submitted (MAR-1998) to the EMBL/GenBank/DDBJ databases. Bingle C.D., Submitted (JAN-1997) to the EMBL/GenBank/DDBJ databases.