

FOXB2 Antibody [#C15767]

Catalog Number:	C15767
Concentration:	1mg/ml
Swiss-Prot No.:	Q5VYV0
Other Names:	forkhead box B2
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 FOXB2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	FOXB2 antibody detects endogenous levels of total FOXB2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Humphray S.J., Nature 429:369-374(2004).