

BLCAP Antibody [#C10782]

Catalog Number:	C10782
Concentration:	1mg/ml
Swiss-Prot No.:	P62952
Other Names:	Bladder cancer-associated protein, Bladder cancer 10 kDa protein, Bc10, BLCAP, BC10
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 BLCAP.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BLCAP antibody detects endogenous levels of total BLCAP protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:40000
References:	Olsen S.R., Mol. Endocrinol. 5:1246-1256(1991). Knoell R., Submitted (FEB-2000) to the EMBL/GenBank/DDBJ databases. The MGC Project Team, Genome Res. 14:2121-2127(2004).