

BHLHB3 Antibody [#C11661]

Catalog Number:	C11661
Concentration:	1mg/ml
Swiss-Prot No.:	Q9C0J9
Other Names:	bHLHB3, Differentially expressed in chondrocytes protein 2, hDEC2, Enhancer-of-split and hairy-related protein 1, SHARP-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 BHLHB3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BHLHB3 antibody detects endogenous levels of total BHLHB3 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Fujimoto K., Biochem. Biophys. Res. Commun. 280:164-171(2001). Garriga-Canut M., J. Biol. Chem. 276:14821-14828(2001).