

AML1 (Ab-276) Antibody [#B0408]

Catalog Number:	B0408
Concentration:	1mg/ml
Swiss-Prot No.:	Q01196
Other Names:	Acute myeloid leukemia 1 protein, CBF-alpha 2, CBFA2, Core-binding factor, alpha 2 subunit, Oncogene AML-1, PEA2-alpha B, PEBP2-alpha B, Polyomavirus enhancer binding protein 2 alpha B subunit, RUN1, RUNX1, Runt-related transcription factor 1, SL3-3 enhancer factor 1 alpha B subunit, SL3/AKV core-binding factor alpha B subunit
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 non-phosphopeptide derived from human AML1 around the phosphorylation site of 276 (Q-P-S ^P -P-P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AML1 (Ab-276) antibody detects endogenous levels of total AML1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Chaitali Banerjee, PNAS, May 1996; 93: 4968. Jennifer J. Westendorf, Mol. Cell. Biol., Jan 1998; 18: 322 - 333. Ari Melnick, Blood, Dec 2000; 96: 3939 - 3947.