

BL-CAM (Ab-807) Antibody [#B8230]

Catalog Number:	B8230
Concentration:	1mg/ml
Swiss-Prot No.:	P20273
Other Names:	B-cell receptor CD22; B-lymphocyte cell adhesion molecule; BL-CAM; Leu-14; Siglec-2
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 BL-CAM around the phosphorylation site of tyrosine 807 (G-D-Y ^P -E-N).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BL-CAM (Ab-807) antibody detects endogenous levels of total BL-CAM protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Wilson G.L., J. Exp. Med. 173:137-146(1991). Wilson G.L., J. Immunol. 150:5013-5024(1993). Stamenkovic I., Nature 345:74-77(1990).