

CD69 Antibody [#C15597]

Catalog Number:	C15597
Concentration:	1mg/ml
Swiss-Prot No.:	Q07108
Other Names:	activation inducer molecule; AIM; BL-AC/P26; EA1; early activation antigen CD69; early T-cell activation antigen p60; GP32/28; Leu-23; MLR-3
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 CD69.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CD69 antibody detects endogenous levels of total CD69 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 92%); Rat (Identities = 92%, Positives = 92%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Hamann J., J. Immunol. 150:4920-4927(1993). Lopez-Cabrera M., J. Exp. Med. 178:537-547(1993). Ziegler S.F., Eur. J. Immunol. 23:1643-1648(1993).