

MAK (Ab-159) Antibody [#B8128]

Catalog Number:	B8128
Concentration:	1mg/ml
Swiss-Prot No.:	P20794
Other Names:	EC 2.7.11.22; Male germ cell-associated kinase; Serine/threonine-protein kinase MAK
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 MAK.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MAK (Ab-159) antibody detects endogenous levels of total MAK protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Mungall A.J., Nature 425:805-811(2003). Matsushime H., Mol. Cell. Biol. 10:2261-2268(1990). Greenman C., Nature 446:153-158(2007).