

Merlin (Ab-10) Antibody [#B8046]

Catalog Number:	B8046
Concentration:	1mg/ml
Swiss-Prot No.:	P35240
Other Names:	MERL; moesin-ezrin-radixin-like protein; neurofibromatosis type 2 gene; neurofibromatosis-2; Neurofibromin 2; NF-2; NF2; SCH; Schwannomerlin; schwannomerlin; Schwannomin
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 Merlin around the phosphorylation site of serine 10 (R-M-S ^P -F-S).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Merlin (Ab-10) antibody detects endogenous levels of total Merlin protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Trofatter J.A., Cell 72:791-800(1993). Rouleau G.A., Nature 363:515-521(1993). Zucman-Rossi J., Hum. Mol. Genet. 7:2095-2101(1998).