

AMOTL1 Antibody [#C14473]

Catalog Number:	C14473
Concentration:	1mg/ml
Swiss-Prot No.:	Q8IY63
Other Names:	Angiomotin-like protein 1
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 AMOTL1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AMOTL1 antibody detects endogenous levels of total AMOTL1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Bratt A., Gene 298:69-77(2002). Bratt A., Gene 310:231-231(2003). Yu L.-R., J. Proteome Res. 6:4150-4162(2007).