

HORM1 Antibody [#C11285]

Catalog Number:	C11285
Concentration:	1mg/ml
Swiss-Prot No.:	Q86X24
Other Names:	HORMA domain-containing protein 1, Newborn ovary HORMA protein, Cancer/testis antigen 46, CT46, HORMAD1, NOHMA
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 C-terminal of human HORM1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	HORM1 antibody detects endogenous levels of total HORM1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Pangas S.A., Gene Expr. Patterns 5:257-263(2004). Wiemann S., Genome Res. 11:422-435(2001). Gregory S.G., Nature 441:315-321(2006).