

HECW2 Antibody [#C15541]

Catalog Number:	C15541
Concentration:	1mg/ml
Swiss-Prot No.:	Q9P2P5
Other Names:	E3 ubiquitin-protein ligase HECW2; EC 6.3.2.-; HECT, C2 and WW domain-containing protein 2; NEDD4-like E3 ubiquitin-protein ligase 2
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 HECW2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	HECW2 antibody detects endogenous levels of total HECW2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Nagase T., DNA Res. 7:65-73(2000). Bechtel S., BMC Genomics 8:399-399(2007). Hillier L.W., Nature 434:724-731(2005).