

UBFD1 Antibody [#C19408]

Catalog Number:	C19408
Concentration:	1mg/ml
Swiss-Prot No.:	O14562
Other Names:	FLJ38870; FLJ42145; ubiquitin family domain containing 1; UBPH
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 UBFD1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	UBFD1 antibody detects endogenous levels of total UBFD1 protein.
Reactivity:	Human, Mouse, Ray
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Loftus B.J., Genomics 60:295-308(1999). Ota T., Nat. Genet. 36:40-45(2004).