

## UPF1 Antibody [#C18329]

|                           |                                                                                                                                                                    |
|---------------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Catalog Number:</b>    | C18329                                                                                                                                                             |
| <b>Concentration:</b>     | 1mg/ml                                                                                                                                                             |
| <b>Swiss-Prot No.:</b>    | Q92900                                                                                                                                                             |
| <b>Other Names:</b>       | EC 3.6.1.-; ATP-dependent helicase RENT1; Nonsense mRNA reducing factor 1; NORF1; Up-frameshift suppressor 1 homolog; hUpf1                                        |
| <b>Storage/Stability:</b> | Store at -20°C/1 year                                                                                                                                              |
| <b>Form of Antibody:</b>  | Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.                               |
| <b>Immunogen:</b>         | The antiserum was produced against synthesized peptide derived from internal of human UPF1.                                                                        |
| <b>Purification:</b>      | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.                                              |
| <b>Specificity:</b>       | UPF1 antibody detects endogenous levels of total UPF1 protein.                                                                                                     |
| <b>Reactivity:</b>        | Human (Identities = 100%, Positives = 100%);<br>Mouse (Identities = 100%, Positives = 100%)                                                                        |
| <b>Applications:</b>      | WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000                                                                                                                    |
| <b>References:</b>        | Perlick H.A., Proc. Natl. Acad. Sci. U.S.A. 93:10928-10932(1996).<br>Applequist S.E., Nucleic Acids Res. 25:814-821(1997).<br>Nagase T., DNA Res. 3:321-329(1996). |