

## TPIP1 Antibody [#C13126]

|                           |                                                                                                                                                               |
|---------------------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Catalog Number:</b>    | C13126                                                                                                                                                        |
| <b>Concentration:</b>     | 1mg/ml                                                                                                                                                        |
| <b>Swiss-Prot No.:</b>    | Q9HCN2                                                                                                                                                        |
| <b>Other Names:</b>       | p53-regulated apoptosis-inducing protein 1, P53aip1, TP53AIP1                                                                                                 |
| <b>Storage/Stability:</b> | Store at -20°C/1 year                                                                                                                                         |
| <b>Form of Antibody:</b>  | Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), 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 TPPI1.                                                                  |
| <b>Purification:</b>      | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.                                         |
| <b>Specificity:</b>       | TPIP1 antibody detects endogenous levels of total TPPI1 protein.                                                                                              |
| <b>Reactivity:</b>        | Human (Identities = 100%, Positives = 100%)                                                                                                                   |
| <b>Applications:</b>      | IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000                                                                                                                |
| <b>References:</b>        | Oda K., Cell 102:849-862(2000).<br>Zhou G., Submitted (MAY-2003) to the EMBL/GenBank/DDBJ databases.<br>The MGC Project Team, Genome Res. 14:2121-2127(2004). |