

## PIAS4 Antibody [#C0364]

|                           |  |
|---------------------------|--|
| <b>Catalog Number:</b>    | C0364  |
| <b>Concentration:</b>     | 1mg/ml   |
| <b>Swiss-Prot No.:</b>    | Q8N2W9   |
| <b>Other Names:</b>       | PIAS-gamma, PIASG, PIASy, protein inhibitor of activated STAT protein 4, protein inhibitor of activated STAT protein PIASy, protein inhibitor of activated STAT protein gamma, protein inhibitor of activated STAT, 4                              |
| <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 PIAS4.   |
| <b>Purification:</b>      | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  |
| <b>Specificity:</b>       | PIAS4 antibody detects endogenous levels of total PIAS4 protein.   |
| <b>Reactivity:</b>        | Human, Mouse   |
| <b>Applications:</b>      | WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000  |
| <b>References:</b>        | Zendra E. Zehner, AACR Meeting Abstracts, Apr 2005; 2005: 456.<br>Amador Albor, J. Biol. Chem., Sep 2006; 281: 25850 - 25866.<br>Christina-Maria Hecker, J. Biol. Chem., Jun 2006; 281: 16117 - 16127.<br>Yun Zhao, Stem Cells, Apr 2007; 10.1634. |