

PIAS3 Antibody [#C0363]

Catalog Number:	C0363
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y6X2
Other Names:	PROTEIN INHIBITOR OF ACTIVATED STAT3, protein inhibitor of activated STAT 3
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 PIAS3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PIAS3 antibody detects endogenous levels of total PIAS3 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
References:	Jiang-Hong Man, Nucleic Acids Res., Nov 2006; 34: 5552 - 5566. Amir Sonnenblick, Mol. Cell. Biol., Dec 2004; 24: 10584 - 10592. Jianyin Long, PNAS, Jan 2004; 101: 99. Hyun Duk Jang, J. Biol. Chem., Jun 2004; 279: 24873 - 24880.