

PRS6A Antibody [#C14013]

| | |
|---------------------------|---|
| Catalog Number: | C14013 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | P17980 |
| Other Names: | 26S protease regulatory subunit 6A; proteasome subunit P50; PRS6A; TAT-binding protein 1; TBP-1; TBP1 |
| 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 PRS6A. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | PRS6A antibody detects endogenous levels of total PRS6A protein. |
| Reactivity: | Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%) |
| Applications: | WB: 1:500~1:1000 ELISA: 1:40000 |
| References: | Leung D.W., Front. Biosci. 6:D944-D953(2001). Nagamine K., Submitted (MAR-2000) to the EMBL/GenBank/DDBJ databases. Clark H.F., Genome Res. 13:2265-2270(2003). |