

TRAPPC6A Antibody [#C19149]

| | |
|---------------------------|--|
| Catalog Number: | C19149 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | O75865 |
| Other Names: | Trafficking protein particle complex subunit 6A; TRAPP complex subunit 6A; TRAPPC6A; HSPC289 |
| 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 TRAPPC6A. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | TRAPPC6A antibody detects endogenous levels of total TRAPPC6A protein. |
| Reactivity: | Human (Identities = 100%, Positives = 100%) |
| Applications: | IHC: 1:50~1:100 ELISA: 1:40000 |
| References: | Ye M., Submitted (MAY-1999) to the EMBL/GenBank/DDBJ databases.; Grimwood J., Nature 428:529-535(2004).; The MGC Project Team, Genome Res. 14:2121-2127(2004). |