

MED23 Antibody [#C11881]

| | |
|---------------------------|---|
| Catalog Number: | C11881 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | Q9ULK4 |
| Other Names: | Mediator of RNA polymerase II transcription subunit 23, Mediator complex subunit 23, CRSP complex subunit 3, Cofactor required for Sp1 transcriptional activation subunit 3, Transcriptional coactivator CRSP130, Vitamin D3 receptor-interacting protein complex 130 kDa component, DRIP130, Activator-recruited cofactor 130 kDa component, ARC130, hSur-2, MED23, ARC130, CRSP3, DRIP130, KIAA1216, SUR2 |
| 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 MED23. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | MED23 antibody detects endogenous levels of total MED23 protein. |
| Reactivity: | Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%) |
| Applications: | IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000 |
| References: | Wiemann S., Genome Res. 11:422-435(2001). Mungall A.J., Nature 425:805-811(2003). Akoulitchev S., Nature 407:102-106(2000). |