

Instruction Manual

| | |
|---------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Catalog Number | PK-AB577-3636 |
| Description | PAX genes encode nuclear transcription factors that may function as major controllers of developmental processes in both vertebrates and invertebrates. Mutations in murine PAX genes underlie three natural mouse alleles and several corresponding human syndromes (aniridia, foveal hypoplasia and Peters' anomaly). Murine PAX genes have been shown to be proto-oncogenes. Furthermore, human PAX genes have recently been demonstrated to play an influential part in some common human cancers such as brain tumors and lymphomas. All PAX genes encode a DNA-binding domain termed the paired domain and in addition some also encode a second binding domain--the paired type homeobox. PAX6 is involved in the early development of the optical vesicle and has been shown to interact with Six3, another important visual development protein. |
| Quantity | 100 µg |
| Source / Host | Rabbit |
| Immunogen | Synthetic peptide surrounding amino acid 303 of rat PAX6. |
| Clone / IgG Subtype | Rabbit IgG |
| Species Reactivity | human, mouse, rat |
| Specificity | See Applications or Species Reactivity. |
| Formulation | 100 µg (0.5 mg/ml) affinity purified rabbit polyclonal antibody in phosphate-buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, and 0.01% thimerosal. |
| Reconstitution | Can be diluted in other aqueous buffers at the concentrations determined for the respective application. |
| Storage & Stability | Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles. |
| Applications | Western blotting (0.5-4 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody recognizes ~32 kDa and 47 kDa isoforms of PAX6 in samples from human, mouse and rat origins. Reactivity to other species has not been tested. Blocking peptide (Cat.# PK-AB577-3636P) is available separately. |

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.