

Instruction Manual

| | |
|---------------------|--|
| Catalog Number | PK-AB577-3389 |
| Description | Heat shock proteins (HSPs) are ubiquitously expressed in all organisms. A major function of Hsp90 and other HSPs is to act as molecular chaperones. Hsp90 forms a complex with glucocorticoid receptor (GR), rendering the non ligand-bound receptor transcriptionally inactive. Hsp90 binds the GR as a heterocomplex composed of either Hsp56 or cyclophilin-40, forming an aporeceptor complex. Hsp90 also exists as a dimer with other proteins such as p60/sti1 and p23, forming an aporeceptor complex with estrogen and androgen receptors. |
| Quantity | 100 µg |
| Source / Host | Rabbit |
| Immunogen | Synthetic peptide surrounding amino acid 714 of human Hsp90. |
| 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) 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), in immunoprecipitation (10-20 µg/ml) and immunohistochemistry (frozen sections, 10-20 µg/ml). The optimal concentrations should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody recognizes Hsp90alpha and Hsp90beta of human, mouse, and rat origins. |

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