

## Instruction Manual

Catalog Number	PK-AB577-3342
Description	Members of Oct family of transcription factors specifically interact with Octamer motif ATGCAAAT, a regulatory element important for tissue- and cell-specific transcription as well as for transcription of a number of housekeeping genes. All of the members of the Oct family contain two highly conserved domains which are separated by 14-26 variable amino acids. These include the POU homeodomain and the POU-specific domain. Both are required for DNA binding and are involved in protein-protein interactions. Evidences indicate that regulation of Oct family transcription factors occurs at the level of phosphorylation.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 727 of human Oct-1.
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 anti-Oct-1 polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 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). Based on researchers feed back, it also works in immunohistochemistry (10-20 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far.

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