

Instruction Manual

Catalog Number	PK-AB577-3576
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 mapping to the N-terminus of mouse Oct-4.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat, bovine, pig
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit anti-Oct-3/4 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), immunoprecipitation (20 µg/ml) and immunohistochemistry (20 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody reacts with 45 kDa Oct-4 (also designated Oct-3A) and in a less extent 33 kDa Oct-3B from human, mouse, and rat samples. Rat kidney tissue lysate can be used as a positive control.

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