

## Instruction Manual

Catalog Number	PK-AB577-3624
Description	The nucleosome is made up of four core histone proteins (H2A, H2B, H3 and H4) and is the primary building block of chromatin. The N-terminal tail of core histones undergoes different posttranscriptional modification including acetylation, phosphorylation and methylation. These modifications occur in response to cell signal stimuli and have a direct effect on gene expression.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide corresponding to the N-terminus of human Histone H4.
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-Histone H4 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), in immunoprecipitation (20-40 µg/ml) and Immunohistochemistry (20-40 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody detects ~11 kDa histone H4 protein. It does not cross-react with other histones.

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