

Instruction Manual

Catalog Number	PK-AB577-3267
Description	Lamins are nuclear membrane structural components that are important in maintaining normal cell functions such as cell cycle control, DNA replication and chromatin organization. Lamin A is specifically cleaved by caspase-6 and therefore serves as a marker for caspase-6 activation. During apoptosis the 70 kDa lamin A is cleaved to a large (40-45 kDa) and small (23 kDa) fragment. The cleavage of lamins results in nuclear disregulation and cell death.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 220 of human Lamin A.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.2 mg/ml) affinity purified rabbit anti-Lamin A 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 blot analysis (0.5-4 µg/ml) and Immunohistochemistry (15-20 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. Cross-reactivity to IκB-alpha of other species has not been determined. The antibody mainly detects full length (~70 kDa) and in a much lesser extent the cleaved fragments (45 kDa and 23 kDa) of Lamin A/C.

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