

Instruction Manual

Catalog Number	PK-AB577-3447
Description	Five ionomycin-inducible complementary cDNAs, designated PAR1, 2, 3, 4 and 5, have been isolated from the prostate cancer cell line AT-3. Nucleotide sequencing identified PAR1 as the rat homolog of monkeyP-1, PAR2 as the injury-inducible gene HB-EGF, and PAR3 as the serum-induced gene cyr-61. PAR4 and PAR5 sequences were not found to correspond to any previously described proteins. PAR4 (prostate apoptosis response-4) is specifically expressed by cells entering apoptosis and is not inducible by growth factor stimulation, oxidative stress and necrosis, or growth arrest. The PAR4 gene encodes a 38 kDa protein with a putative nuclear localization signal and carboxy terminal leucine zipper.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 324 of human Par4.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat, dog, horse, zebrafish
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) 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), immunoprecipitation, immunohistochemistry. The optimal concentrations 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.