

Instruction Manual

Catalog Number	PK-AB577-3444P
Quantity	50 µg
Formulation	50 µg (0.5 mg/ml) in phosphate buffered saline (PBS), pH 7.2, containing 50% glycerol, 1% BSA and 0.02% thimerosal.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	The peptide is used for blocking the antibody activity of Pak4 (Cat. No. PK-AB577-3444). It usually blocks the antibody activity completely in Western blot analysis by incubating the peptide with equal volume of antibody for 30-60 minutes at 37°C.

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