

Instruction Manual

Catalog Number	PK-AB577-3609P
Quantity	50 µg
Formulation	50 µg (0.2 mg/ml) in phosphate buffered saline (PBS), pH 7.2, containing 0.1% BSA and 0.02% sodium azide.
Storage & Stability	Store at -20°C. Avoid repeated freeze/defrost cycles.
Applications	The peptide is used for blocking the antibody activity of HDAC-9 (PK-AB577-3609). It usually blocks the antibody activity completely in Western blot analysis by incubating the peptide with equal volume of antibody for 30 minutes at 37°C.

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