

# HDAC-5 peptide

## HDAC-5 Antibody Blocking Peptide

### Instruction Manual

Catalog Number	PK-AB577-3605P
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. Avoid repeated freeze/defrost cycles.
Applications	The peptide is used for blocking the antibody activity of HDAC-5 (PK-AB577-3605). 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.