

### Instruction Manual

|                     |  |
|---------------------|--|
| Catalog Number      | PK-AB577-3601P   |
| 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 HDAC-1 (PK-AB577-3601). 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.