

HDAC-10 peptide

HDAC-10 Antibody Blocking Peptide

Instruction Manual

| | |
|---------------------|---|
| Catalog Number | PK-AB577-3610P |
| Quantity | 50 µg |
| Formulation | 50 µg (0.2 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-10 (PK-AB577-3610). 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.