

Instruction Manual

| | |
|--------------------------------|---|
| Catalog Number | PK-AB718-4547P |
| Quantity | 50 µg |
| Source | 14 amino acids near the amino terminus of human EVER1 |
| Formulation | Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH ₂ PO ₄ , 10 mM Na ₂ HPO ₄ , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide. |
| Reconstitution | During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap. |
| Storage & Stability | Store EVER1 peptide at -20°C, stable for one year. |
| Application | EVER1 peptide is used for blocking the activity of EVER1 antibody. |
| References | Ramoz N, Rueda L-A, Bouadjar B, et al. Mutations in two adjacent novel genes are associated with epidermodysplasia verruciformis. Nat. Genetics 2002; 32:579-81. |

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