

Instruction Manual

| | |
|--------------------------------|---|
| Catalog Number | PK-AB718-2289P |
| Quantity | 50 µg |
| Source | 16 amino acids of human Bnip3L |
| 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 Bnip3L peptide at -20°C, stable for one year. |
| Application | Bnip3L peptide is used for blocking the activity of Bnip3L antibody. |
| References | Chen G, Cizeau J, Vande Velde C, et al. Nix and Nip3 form a subfamily of pro-apoptotic mitochondrial proteins. <i>J Biol Chem</i> 1999;274(1):7-10. Daugas E, Susin SA, Zamzami N, et al. Mitochondrio-nuclear translocation of AIF in apoptosis and necrosis. <i>FASEB J</i> 2000;14:729-39 (WD0500) |

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