

Instruction Manual

Catalog Number	PK-AB718-2271P
Quantity	50 µg
Source	16 amino acids near the carboxy terminus of human FLASH
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 FLASH peptide at -20°C, stable for one year.
Application	FLASH peptide is used for blocking the activity of FLASH antibody.
References	Koonin EV, Aravind L, Hofmann K, Tschopp J, Dixit VM Apoptosis. Searching for FLASH domains. Nature 1999;401:662-3.

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