

Instruction Manual

Catalog Number	PK-AB718-3975P
Quantity	50 µg
Source	19 amino acids near the carboxy terminus of human PDCD4
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 PDCD4 peptide at -20°C, stable for one year.
Application	PDCD4 peptide is used for blocking activity of PDCD4 antibody.
References	Shibihara K, Asano M, Ishida Y, et al. Isolation of a novel mouse gene MA-3 that is induced upon programmed cell death. Gene 1995; 166:297-301.

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