

Instruction Manual

Catalog Number	PK-AB718-4063P
Quantity	50 µg
Source	16 amino acids near the center of human PDL-2
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 PDL-2 peptide at -20°C, stable for one year.
Application	The peptide is used for blocking the activity of PDL-2 antibody.
References	Zhang Y, Chung Y, Bishop C, et al. Regulation of T cell activation and tolerance by PDL2. Proc. Natl. Acad. Sci. USA 2006; 103:11695-700.

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