

Instruction Manual

Catalog Number	PK-AB718-4327P
Quantity	50 µg
Source	17 amino acids near the amino terminus of human RAP80
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 RAP80 peptide at -20°C, stable for one year.
Application	RAP80 peptide is used for blocking the activity of RAP80 antibody.
References	Yan Z, Kim YS, and Jetten AM. RAP80, a novel nuclear protein that interacts with the retinoid-related testis-associated receptor. J. Biol. Chem. 2002; 277:32379-88.

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