

Instruction Manual

Catalog Number	PK-AB718-3957P
Quantity	50 µg
Source	14 amino acids near the center of human Slug
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 Slug peptide at -20°C, stable for one year.
Application	Slug peptide is used for blocking activity of Slug antibody.
References	Neito MA, Sargent MG, Wilkinson DG, et al. Control of cell behavior during vertebrate development by Slug, a zinc finger gene. Science 1994; 264:835-9

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