

Instruction Manual

Catalog Number	PK-AB718-4259P
Quantity	50 µg
Source	17 amino acids near the center of human FABP7
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 FABP7 peptide at -20°C, stable for one year.
Application	FABP7 peptide is used for blocking the activity of FABP7 antibody.
References	Shimizu F, Watanabe TK, Shinomiya H, et al. Isolation and expression of a cDNA for human brain fatty acid-binding protein (B-FABP). <i>Biochim. Biophys. Acta</i> 1997; 1354:24-8.

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