

Instruction Manual

Catalog Number	PK-AB718-4627P
Quantity	50 µg
Source	16 amino acids near the amino terminus of human SATB2
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 SATB2 peptide at -20°C, stable for one year.
Application	SATB2 peptide is used for blocking the activity of SATB2 antibody.
References	Szemes M, Gyorgy A, Pawaletz C, et al. Isolation and characterization of SATB2, a novel AT-rich DNA binding protein expressed in development- and cell-specific manner in the rat brain. <i>Neurochem. Res.</i> 2006; 31:237-46.

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