

Instruction Manual

Catalog Number	PK-AB718-3631P
Quantity	50 µg
Source	16 amino acids near the carboxy terminus of human PIST
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 PIST peptide at -20°C, stable for one year.
Application	PIST peptide is used for blocking the activity of PIST antibody.
References	<p>Yu Z, Horton A, Bravin M, et al. A novel protein complex linking the delta2 glutamate receptor and autophagy: implications for neurodegeneration in Lurcher mice. <i>Neuron</i> 2002; 35:921-33.</p> <p>Takeuchi T, Misaki A, Chen BK, et al. H-cadherin expression in breast cancer. <i>Histopathology</i> 1999; 35:87-88.</p> <p>Sato M, Mori Y, Sakurada A, et al. The H-cadherin (CDH13) gene is inactivated in human lung cancer. <i>Hum. Gen.</i> 1998; 103:96-101.</p> <p>Shamji AF, Ngheim P, and Schreiber SL. Integration of growth factor and nutrient signaling: implications for cancer biology. <i>Proc. Natl. Acad. Sci. USA</i> 1985; 82:7145-8.</p>

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