

Instruction Manual

Catalog Number	PK-AB718-4413P
Quantity	50 µg
Source	14 amino acids near the center of human GTSP1
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 GSTP1 peptide at -20°C, stable for one year.
Application	GSTP1 peptide is used for blocking the activity of GTSP1 antibody.
References	Pearson WR. Phylogenies of glutathione transferase families. <i>Methods Enzymol.</i> 2005; 401:186-204.

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