

Instruction Manual

Catalog Number	PK-AB718-3781P
Quantity	50 µg
Source	16 amino acids near the carboxy terminus of human GAPDH
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 GAPDH peptide at -20°C, stable for one year.
Application	GAPDH peptide is used for blocking the activity of GAPDH antibody.
References	Sirover MA. New nuclear functions of the glycolytic protein, glyceraldehyde-3-phosphate dehydrogenase, in mammalian cells. J. Cell. Biochem. 2005; 95:45-52. Mosialos G, Birkenbach M, Yalamanchili R, et al. The Epstein-Barr virus transforming protein LMP1 engages signaling proteins for the tumor necrosis factor receptor family. Cell 1995; 80:389-99. van Eyndhoven WG, Gamper CJ, Cho E, et al. TRAF-3 mRNA splice-deletion variants encode isoforms that induce NFκB activation. Mol. Immunol. 1999; 36:647-58.

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