

Instruction Manual

Catalog Number	PK-AB718-3035P
Quantity	50 µg
Source	16 amino amino terminus of human EndoG
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 EndoG peptide at -20°C, stable for one year.
Application	EndoG peptide is used for blocking the activity of EndoG antibody.
References	<p>Tiranti V, Rossi E, Ruiz-Carrillo A, Rossi G, Rocchi M, DiDonato S, Zuffardi O, Zeviani M. Chromosomal localization of mitochondrial transcription factor A (TCF6), single-stranded DNA-binding protein (SSBP), and endonuclease G (ENDOG), three human housekeeping genes involved in mitochondrial biogenesis. <i>Genomics</i>. 1995;25(2):559-64.</p> <p>Gupta A and Tsai LH. Cyclin-dependent kinase 5 and neuronal migration in the neocortex. <i>Neurosignals</i> 2003; 12:173-9.</p> <p>Liang Y, Yu W, Li Y, et al. Nudel functions in membrane traffic mainly through association with Lis1 and cytoplasmic dynein. <i>J. Cell Biol.</i> 2004; 164:557-66.</p>

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