

Rabbit Anti-Human/Mouse NADE (Novel Apoptosis Related Protein)

Instruction Manual

Catalog Number	PK-AB718-3359
Synonyms	NADE Antibody:
Description	The p75 neurotrophin receptor (p75NTR) is a member of the tumor necrosis receptor superfamily and can mediate cell death and cell survival in response to nerve growth factor (NGF). The p75NTR-associated cell death executor (NADE) mediates apoptosis by interacting with the cell death domain of p75NTR following the binding of NGF by p75NTR. Recent studies have shown that NADE also interacts with second mitochondria-derived activator of caspase (Smac). Co-expression of NADE and Smac promotes TRAIL-induced apoptosis and inhibits XIAP-mediated Smac ubiquitization. It has been suggested that it is this interaction between NADE and Smac that allows apoptosis to proceed. Finally, although initially discovered as an mRNA expressed in ovarian granulosa cells, NADE has been suggested to play a role in the neuronal death seen in epileptic brain damage.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal NADE antibody was raised against a peptide corresponding to 14 amino acids near the middle of human NADE (GenBank accession no. NP_996798).
Purification Method	Ion exchange chromatography purified.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
Specificity	
Formulation	Antibody is supplied in PBS containing 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	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC I Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually. NADE antibody can be used for detection of NADE by Western blot at 1 µg/mL. Despite its predicted molecular weight, NADE migrates at ~23 kDa in SDS-PAGE. Antibody can also be used for immunohistochemistry starting at 2 µg/mL.
Images	Available upon request.
References	Gentry JJ, Barker PA, and Cater BD. The p75 neurotrophin receptor: multiple interactors and numerous functions. <i>Pro. Brain Res.</i> 2004; 146:25-39. Mukai J, Hachiya T, Shoji-Hoshino S, et al. NADE, a p75NTR-associated cell death executor, is involved in signal transduction mediated by the common neurotrophin receptor p75NTR. <i>J. Biol. Chem.</i> 2000; 275:17566-70. Rapp G, Freudenstein J, Klaudiny J, et al. Characterization of three abundant mRNAs from human ovarian granulosa cells. <i>DNA Cell Biol.</i> 1990; 9:479-85. YI J-S, Lee, S-K, Sato T-A, et al. Co-induction of p75NTR and the associated death executor NADE in degenerating hippocampal neurons after kainate-induced seizures in the rat. <i>Neurosci. Lett.</i> 2003; 347:126-30.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3359P; Human brain Cell Lysate, Cat. No. PK-AB718-1303 TRAIL Antibody, Cat. No. PK-AB718-1113; Smac Antibody (CT), Cat. No. PK-AB718-2409 XIAP Antibody (CT), Cat. No. PK-AB718-3331

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