

# Nop30 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Nop30

## Instruction Manual

Catalog Number	PK-AB718-2225
Synonyms	Nop30 Antibody: Nucleolar protein 3, Apoptosis repressor with CARD, ARC
Description	Apoptosis, also known as programmed cell death, plays major roles in development and normal tissue turnover in addition to tumor formation. Apoptosis is regulated by death domain (DD) and/or caspase recruitment domain (CARD) containing molecules and the caspase family of proteases. CARD domain containing cell death regulators include RAIDD, Apaf-1, caspase-9, and caspase-2. A novel CARD domain containing protein was recently identified and designated ARC for apoptosis repressor with CARD. An alternate splicing isoform of ARC was identified as Nop30. While ARC interacts with caspase-2 and -8 and suppresses apoptosis induced by cell death adapters FADD and TRADD and by cell death receptors Fas, TNFR-1 and DR3, Nop30 multimerizes and binds to the splicing factor SRp30c and may act to influence alternative splice site selection in vivo. The Nop30 antibody will not detect ARC protein.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Nop30 antibody was raised against a 14 amino acid peptide from near the carboxy terminus of human Nop30 (GenBank accession no. AAH12798).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
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 INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Application have to be determined individually. Nop30 antibody can be used for detection of Nop30 by Western blot at 0.5 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 0.5 µg/mL.
Images	Available upon request.
References	Jin Z and El Deiry WS. Overview of cell death signaling pathways. Cancer Biol. Ther. 2004; 4:139-63. Koseki T, Inohara N, Chen S, et al. ARC, an inhibitor of apoptosis expressed in skeletal muscle and heart that interacts selectively with caspases. Proc. Natl. Acad. Sci. USA 1998; 95:5156-60. Stoss O, Schwaiger FW, Cooper TA, et al. Alternative splicing determines the intracellular localization of the novel nuclear protein Nop30 and its interaction with the splicing factor SRp30c. J. Biol. Chem. 1999; 274:10951-62
Images	Available upon request.
Related Products	Nop30 Peptide, Cat. No. PK-AB718-2225P Mouse Skeletal Muscle Tissue Lysate, Cat. No. PK-AB718-1407

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