

Rabbit Anti-Human Nucleotide-binding Oligomerization Domain Protein 2 (New CARD Containing Protein)

Instruction Manual

Catalog Number	PK-AB718-2513
Synonyms	NOD2 Antibody: Caspase recruitment domain 15, CARD15
Description	Apaf-1 and NOD1 are members of a new family, which are involved in the regulation of apoptosis and immune response. Each of them contains a caspase recruitment domain (CARD) and a nucleotide-binding oligomerization domain (NOD). A third member in this family was recently identified and designated NOD2. NOD2 interacts with RICK via a homophilic CARD-CARD interaction. NOD2 activates NF- κ B, which is regulated by its carboxy-terminal leucine-rich repeat domain that acts as an intracellular receptor for components of bacteria. The variants of NOD2, either a frameshift or a missense, were associated with Crohn's disease that is a main type of chronic inflammatory bowel disease.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit NOD2 polyclonal antibody was raised against a synthetic peptide corresponding to 14 amino acids at the carboxy terminus of human NOD2 (GenBank accession no. Q9HC29).
Purification Method	Immunoaffinity chromatography purified IgG
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	NOD2 has no cross-reaction with NOD1.
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 μ l 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, ICC, IF INote: 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. NOD2 antibody can be used for detection of Noxa by Western blot at 2 to 4 μ g/mL. Antibody can also be used for immunocytochemistry starting at 10 μ g/mL. For immunofluorescence start at 20 μ g/mL.
Images	Available upon request.
References	Inohara N, Koseki T, del Peso L, et al. Nod1, an Apaf-1-like activator of caspase-9 and nuclear factor- κ B. J. Biol. Chem. 1999; 274:14560-7. Ogura Y, Inohara N, Benito A, et al. Nod2, a Nod1/Apaf-1 family member that is restricted to monocytes and activates NF- κ B. J. Biol. Chem. 2001; 276:4812-8. Hugot JP, Chamaillard M, et al. Association of NOD2 leucine-rich repeat variants with susceptibility to Crohn's disease. Nature 2001; 411:599-603. Ogura Y, Bonen DK, Inohara N, et al. A frameshift mutation in NOD2 associated with susceptibility to Crohn's disease. Nature 2001; 411:603-6.
Images	Available upon request.
Related Products	HeLa Lysate, Cat. No. PK-AB718-1201 NOD2 Antibody, Cat. No. PK-AB718-2511 NOD2 Peptide, Cat. No. PK-AB718-2513P RICK Antibody, Cat. No. PK-AB718-2075

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