

Anti-NOD2 (NT)

(Caspase recruitment domain 15, CARD15)

CATALOG NO.: 54110

BACKGROUND:

Apaf-1 and NOD1 are members of a new family (1), 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 (2). NOD2 interacts with RICK via a homophilic CARD-CARD interaction. NOD2 activates NF- κ B, which is regulated by its C-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 (3, 4) that is a main type of chronic inflammatory bowel disease.

SOURCE & REACTIVITY:

Rabbit anti-NOD2 polyclonal antibody was raised against a synthetic peptide corresponding to 16 amino acids at the N-terminus of human NOD2 (GenBank accession no. Q9HC29). Anti-NOD2 reacts with NOD2 at the molecular weight of approximately 96 kDa on western blot. Species reactivity includes human, while others are not tested. Anti-NOD2 has no cross reactivity towards NOD1.

APPLICATION:

The following concentration ranges are recommended starting points for this product.

Western Blot: 2 to 4 µg/ml.

Positive Control: Jurkat cell lysate



This product is for in vitro research purposes only.

RELATED PRODUCTS:

Anti-NOD2 (CT), Catalog No. **54111** Anti-RICK (NT), Catalog No. **54042**

STORAGE:

The antibody is supplied as immunoaffinity chromatography purified IgG, in 1X PBS containing 0.02% sodium azide. Store at 2-8 °C for up to 1 year. Avoid repeated freeze thaw cycles.

REFERENCES:

- 1. Inohara N. et al. J. Biol. Chem. 274, 14560 (1999).
- 2. Ogura Y. et al. J. Biol. Chem. 276, 4812 (2001).
- 3. Hugot, JP. et al. Nature 411, 599 (2001).
- 4. Ogura, Y. et al. Nature 411, 603 (2001).