

Product Information

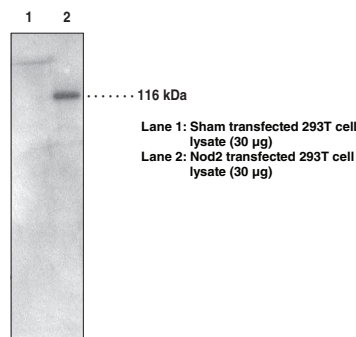


Nod2 Polyclonal Antibody

Item No. 160777

- Contents:** This vial contains 100 µg of lyophilized IgG against human Nod2.
Host: Rabbit
Antigen: The Nod2 polyclonal antibody was raised against residues 28-301 of recombinant human Nod2.^{1,2}
Cross-reactivity: (+) Nod2 transfected 293T cell lysate; (-) SHAM transfected 293T cell lysate
Storage: ≥2 year at -20°C
Application: Reconstitute the lyophilized antibody with 100 µl water containing 0.05% sodium azide. Store the reconstituted antibody at 4°C when not in use. This polyclonal antibody is useful for western blot analysis (at dilutions of 1:500 to 1:1,000) and immunoprecipitation of Nod2 in cells transfected with Nod2-containing plasmids. Nod2 is expressed at low levels in monocytes, and may not be detected using this antibody.

Nod2 is a member of a protein family of apoptosis regulators which includes Apaf-1 and Nod1/CARD4.¹ Nod2 contains two N-terminal caspase recruitment domains (CARDs), a nucleotide binding domain (NBD) and C-terminal leucine rich repeats (LRRs). Nod1/CARD4 is broadly expressed in tissues; however, Nod2 expression is limited to monocytes.^{1,3} Both Nod1 and Nod2 activate NF-κB and act as intracellular receptors for bacterial lipopolysaccharides (LPS).³⁻⁶ The LRRs of both Nod1 and Nod2 are essential for LPS-induced activation of NF-κB.⁵ A frameshift mutation in the Nod2 gene confers susceptibility to Crohn's disease, possibly by causing truncation of the 10th LRR resulting in a protein that is unresponsive to LPS.^{2,7}



References

- Ogura, Y., Inohara, N., Benito, A., *et al.* Nod2, a Nod1/Apaf-1 family member that is restricted to monocytes and activates NF-κB. *J. Biol. Chem.* **276**, 4812-4818 (2001).
- Ogura, Y., Bonen, D.K., Inohara, N., *et al.* A frameshift mutation in NOD2 associated with susceptibility to Crohn's disease. *Nature* **411**, 603-606 (2001).
- Bertin, J., Nir, W.-J., Fischer, C.M., *et al.* Human CARD4 protein is a novel CED-4/Apaf-1 cell death family member that activates NF-κB. *J. Biol. Chem.* **274**, 12955-12958 (1999).
- Inohara, N., Koseki, T., del Peso, L., *et al.* Nod1, an Apaf-1-like activator of caspase-9 and nuclear factor-κB. *J. Biol. Chem.* **274**, 14560-14567 (1999).
- Inohara, N., Ogura, Y., Chen, F.F., *et al.* Human NOD1 confers responsiveness to bacterial lipopolysaccharides. *J. Biol. Chem.* **276**, 2551-2554 (2001).
- Inohara, N., Koseki, T., Lin, J., *et al.* An induced proximity model for NF-κB activation in the Nod1/RICK and RIP signaling pathways. *J. Biol. Chem.* **275**, 27823-27831 (2000).
- Hugot, J.-P., Chamaillard, M., Zouali, H., *et al.* Association of NOD2 leucine-rich repeat variants with susceptibility to Crohn's disease. *Nature* **411**, 599-603 (2001).

Related Products

Caspase-3 (human) Polyclonal Antibody - Item No. 160745 • Apaf-1 Polyclonal Antibody - Item No. 160780 • Caspase-9 Polyclonal Antibody - Item No. 160790 • Nod2 Monoclonal Antibody - Item No. 10004942

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 08/03/2011