

Product Information



iNOS Polyclonal Antibody

Catalog No. 160862

Synonyms:	Inducible Nitric Oxide Synthase; NOS II
Contents:	This vial contains 50 µg protein A-purified IgG in 500 µl TBS, pH 7.4, containing 0.5 mg/ml BSA, 50% glycerol, and 0.02% sodium azide
Host:	Rabbit
Antigen:	Purified iNOS from cytokine-induced murine macrophages (RAW 264.7 cells)
Cross-reactivity:	(+) iNOS from most mammalian species and nNOS (~5%); (-) eNOS
Stability:	≥1 year at -20°C
Applications:	Immunoprecipitation, immunocytochemistry Western blot recommended starting dilution: 1:200; 1-3

Laboratory Procedures

This vial contains 50 µg of protein-A purified IgG. For long term storage, we suggest the antibody be stored as supplied at -20°C; it will be stable for at least one year. For immunoblotting experiments, dilute 10 µl of the antibody to 10 ml (1:1,000) with any buffer suitable for your experiment. Higher dilutions of the antibody may be necessary if high background is evident, particularly when using ECL. For other applications the working concentration of the antibody must be determined empirically.

Nitric Oxide Synthase (NOS) catalyzes the biosynthesis of nitric oxide from L-arginine. Constitutively expressed NOS is found in brain (nNOS) and endothelial cells (eNOS). iNOS is a soluble enzyme found in a variety of tissues including macrophages, hepatocytes, vascular smooth muscle cells, and chondrocytes.^{4,5} iNOS expression is increased by a variety of factors including LPS, IFN-γ, IL-1β, and TNF-α, whereas expression is decreased by dexamethasone.⁶⁻¹⁰ The enzyme has been cloned from several species including mouse, rat, and human with homology of at least 80% between these species.⁵ The calculated molecular weight of the protein from the deduced amino acid sequence is 130,000 - 131,000.⁴

References

1. Herencia, F., Ferrándiz, M.L., Ubeda, A., *et al.* Novel anti-inflammatory chalcone derivatives inhibit the induction of nitric oxide synthase and cyclooxygenase-2 in mouse peritoneal macrophages. *FEBS Lett.* **453**, 129-134 (1999).
2. Smith, F.S. and Titheradge, M.A. Detection of NOS isoforms by western-blot analysis. *Methods in Molecular Biology* **100**, 171-180 (1998).
3. Hayes, J.D. and McLellan, L.I. Glutathione and glutathione-dependent enzymes represent a co-ordinately regulated defence against oxidative stress. *Free Rad. Res.* **31**, 273-300 (1999).
4. Knowles, R.G. and Moncada, S. Nitric oxide synthases in mammals. *Biochem. J.* **298**, 249-258 (1994).
5. Michel, T., Xie, Q., Nathan, C. Molecular biological analysis of nitric oxide synthases, Chapter 10, in *Methods in nitric oxide research*. Feelisch, M. and Stamler, J.S. editors. John Wiley & Sons, Chichester, 161-175 (1996).
6. Lorsbach, R.B., Murphy, W.J., Lowenstein, C.J., *et al.* Expression of the nitric oxide synthase gene in mouse macrophages activated for tumor cell killing. Molecular basis for the synergy between interferon-γ and lipopolysaccharide. *J. Biol. Chem.* **268**, 1908-1913 (1993).
7. Balligand, J., Ungureanu-Longrois, D., Simmons, W.W., *et al.* Cytokine-inducible nitric oxide synthase (iNOS) expression in cardiac myocytes. Characterization and regulation of iNOS expression and detection of iNOS activity in single cardiac cyocytes *in vivo*. *J. Biol. Chem.* **269**, 27580-27588 (1994).
8. Schmidt, H.H.H.W., Warner, T.D., Nakane, M., *et al.* Regulation and subcellular location of nitrogen oxide synthases in RAW264.7 macrophages. *Mol. Pharmacol.* **41**, 615-624 (1992).
9. Ogura, T. and Esumi, H. Nitric oxide synthase expression in human neuroblastoma cell line induced by cytokines. *Mol. Neurosci.* **7**, 853-856 (1996).
10. Walker, G., Pfeilschifter, J., and Kunz, D. Mechanism of suppression of inducible nitric-oxide synthase (iNOS) expression in interferon (IFN)-γ-stimulated RAW 264.7 cells by dexamethasone. Evidence for glucocorticoid-induced degradation of iNOS protein by calpain as a key step in post-transcriptional regulation. *J. Biol. Chem.* **271**, 16679-16687 (1996).

Related Products

iNOS (murine macrophage) - Cat. No. 60862 • iNOS Electrophoresis Standard - Cat. No. 360862

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

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 under separate cover to the MSDS supervisor at 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, 06/18/2009

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