

PRODUCT INFORMATION



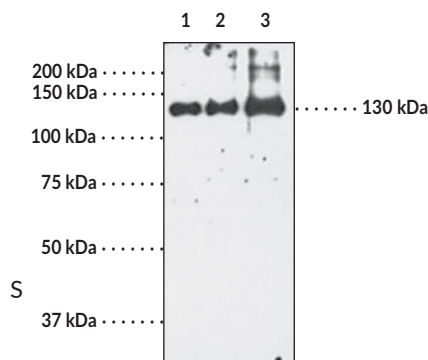
iNOS Polyclonal Antibody

Item No. 160862

Overview and Properties

Contents: This vial contains 500 µl of protein A-purified polyclonal antibody.
Synonyms: Inducible Nitric Oxide Synthase, NOS II
Immunogen: Purified iNOS from cytokine-induced mouse macrophages (RAW 264.7) cells.
Cross Reactivity: (+) nNOS (5%); (-) eNOS
Species Reactivity: (+) Mouse; other species not tested
Uniprot No.: P29477
Form: Liquid
Storage: -20°C (as supplied)
Stability: ≥3 years
Storage Buffer: PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host: Rabbit
Applications: Immunohistochemistry (IHC), Immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution is 1:200.¹⁻³ Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: iNOS electrophoresis standard (50 ng)
Lane 2: iNOS electrophoresis standard (100 ng)
Lane 3: iNOS electrophoresis standard (200 ng)

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/17/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

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-1 β , 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. *FEBS Lett.* **453**, 129-134 (1999).
2. Smith, F.S., and Titheradge, M.A. *Methods Mol. Biol.* **100**, 171-180 (1998).
3. Hayes, J.D., and McLellan, L.I. *Free Rad. Res.* **31**, 273-300 (1999).
4. Knowles, R.G., and Moncada, S. *Biochem. J.* **298**, 249-258 (1994).
5. Michel, T., Xie, Q.W., and Nathan, C. Feilisch, 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. *J. Biol. Chem.* **268**, 1908-1913 (1993).
7. Balligand, J., Ungureanu-Longrois, D., Simmons, W.W., et al. *J. Biol. Chem.* **269**, 27580-27588 (1994).
8. Schmidt, H.H.H.W., Warner, T.D., Nakane, M., et al. *Mol. Pharmacol.* **41**, 615-624 (1992).
9. Ogura, T., and Esumi, H. *Mol. Neurosci.* **7**, 853-856 (1996).
10. Walker, G., Pfeilschifter, J., and Kunz, D. *J. Biol. Chem.* **271**, 16679-16687 (1996).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM