

PRODUCT INFORMATION



eNOS Polyclonal Antiserum

Item No. 160880

Overview and Properties

| | |
|----------------------------|---|
| Contents: | This vial contains polyclonal antiserum. |
| Synonyms: | Endothelial NOS, NOS III, ecNOS |
| Immunogen: | Synthetic peptide from the C-terminal region of human eNOS |
| Cross Reactivity: | (-) iNOS and nNOS |
| Species Reactivity: | (+) Human and bovine; other species not tested |
| Uniprot No.: | P29474 |
| Form: | Lyophilized |
| Storage: | -20°C (as supplied) |
| Stability: | ≥1 year |
| Storage Buffer: | Polyclonal antiserum when reconstituted in 100 µl of deionized water |
| Host: | Rabbit |
| Application: | WB; the recommended starting dilution 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

Description

eNOS catalyzes the formation of nitric oxide (NO) from L-arginine in many cell types including vascular endothelium, bronchiolar epithelium, cardiac myocytes, spleen, and kidney.¹⁻³ Although eNOS was initially characterized as a constitutive enzyme, expression of eNOS mRNA is altered by many factors including shear stress, fibroblast growth factor, hypoxia, and LPS.⁴⁻⁶ eNOS is myristoylated and palmitoylated at the N-terminus, modifications which are required for localization to the plasmallemmal caveolae of endothelial cells.⁷⁻¹¹ Additionally, eNOS is phosphorylated on Ser¹¹⁷⁹ by protein kinase Akt resulting in an increase in NOS activity.¹²⁻¹⁴

References

1. Janssens, S.P., Shimouchi, A., Quertermous, T., *et al.* *J. Biol. Chem.* **267**(21), 14519-14522 (1992).
2. Shaul, P.W., North, A.J., Wu, L.C., *et al.* *J. Clin. Invest.* **94**(6), 2231-2236 (1994).
3. Balligand, J.L., Kobzik, L., Han, X., *et al.* *J. Biol. Chem.* **270**(24), 14582-14586 (1995).
4. Xiao, Z., Zhang, Z., and Diamond, S.L. *J. Cell. Physiol.* **171**(2), 205-211 (1997).
5. Arnett, U.A., McMillan, A., Dinerman, J.L., *et al.* *J. Biol. Chem.* **(251)271**, 15069-15073 (1996).
6. Lu, J.L., Schmiede, L.M., III, Kuo, L., *et al.* *Biochem. Biophys. Res. Commun.* **225**(1), 1-5 (1996).
7. Busconi, L. and Michel, T. *J. Biol. Chem.* **268**(12), 8410-8413 (1993).
8. Robinson, L.J. and Michel, T. *Proc. Nat. Acad. Sci. USA* **92**(25), 11776-11780 (1995).
9. Feron, O., Saldana, F., Michel, J.B., *et al.* *J. Biol. Chem.* **273**(6), 3125-3128 (1998).
10. García-Cardena, G., Oh, P., Liu, J., *et al.* *Proc. Nat. Acad. Sci. USA* **93**(13), 6448-6453 (1996).
11. Feron, O., Belhassen, L., Kobzik, L., *et al.* *J. Biol. Chem.* **271**(37), 22810-22814 (1996).
12. Corson, M.A., James, N.L., Latta, S.E., *et al.* *Circ. Res.* **79**(5), 984-991 (1996).
13. Dimmeler, S., Fleming, I., Fisslthaler, B., *et al.* *Nature* **399**(6736), 601-605 (1999).
14. Fulton, D., Gratton, J.P., McCabe, T.J., *et al.* *Nature* **399**(6736), 597-601 (1999).

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, 05/06/2019

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