

# PRODUCT INFORMATION



## iNOS (mouse, recombinant)

Item No. 60864

### Overview and Properties

**Synonyms:** Inducible Nitric Oxide Synthase, NOS II  
**Source:** Active recombinant murine enzyme expressed in *E. coli*  
**Uniprot No.:** P29477  
**Molecular Weight:** 130 kDa/subunit • Homodimer  
**Storage:** -80°C (as supplied)  
**Stability:** ≥6 months  
**Purity:** 100,000 x g supernatant  
**Supplied in:** 50 mM HEPES, pH 7.4, with 10% glycerol, 8 μM BH<sub>4</sub>, 2% protease inhibitor cocktail (EDTA free), and 0.05% nuclease

#### Protein

**Concentration:** *batch specific*  
**Activity:** *batch specific*  
**Specific Activity:** *batch specific*  
**Unit Definition:** One unit of enzyme produces 1 nmol of nitric oxide per minute at 37°C in 50 mM HEPES, pH 7.4, containing 1 mM arginine, 1 mM magnesium acetate, 0.15 mM NADPH, 4.5 μM oxyhemoglobin, 18 μM tetrahydrobiopterin, and 180 μM DTT.<sup>1,2</sup>

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### References

1. Hevel, J.M., and Marletta, M.A. Nitric-oxide synthase assays. *Methods in Enzymology* **233**, 250-258 (1994).
2. Lyons, C.R., Orloff, G.J., and Cunningham, J.M. Molecular cloning and functional expression of an inducible nitric oxide synthase from a murine macrophage cell line. *J. Biol. Chem.* **267**, 6370-6374 (1992).

#### 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, 05/27/2020

#### 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