

# PRODUCT INFORMATION



## iNOS (mouse recombinant)

Item No. 60864 • Batch No. XXXX

### Overview and Properties

---

**Synonyms:** Inducible Nitric Oxide Synthase, NOS II  
**Source:** Recombinant murine enzyme expressed in *E. coli*  
**Uniprot No.:** P29477

*Batch specific information can be found on the Certificate of Analysis or by contacting Technical Support*

**Molecular Weight:** 130 kDa/subunit • Homodimer  
**Storage:** -80°C (as supplied)  
**Stability:** ≥6 months  
**Purity:** 100,000 x g supernatant  
**Supplied in:** *batch specific*  
**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.

#### 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, 01/18/2016

#### 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