

CERTIFICATE of ANALYSIS



iNOS (mouse, recombinant)

Item No. 60864 • Batch No. 0821651 • Expires 13FEB2027

Overview

Formulation:	50 mM HEPES, pH 7.4, with 10% glycerol, 8 μ M BH4, 0.08 mM DTT, 2% protease inhibitor cocktail (EDTA free), and 0.05% nuclease
Concentration:	36.6 mg/mL
Storage/stability:	-80°C/9mo
Specific activity:	19.50 U/mg
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 μ M oxyhemoglobin, 18 μ M tetrahydrobiopterin, and 180 μ M DTT.{1705,1874}
Activity:	713.00 U/ml
Purity:	100,000 x g supernatant

Reviewed and approved by: Beth Hurst

Approved date: 14MAY2026

Elizabeth Hurst

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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