

CERTIFICATE of ANALYSIS



eNOS (bovine, recombinant; E. coli expressed)

Item No. 11868 • Batch No. 0818544 • Expires 23FEB2028

Overview

Formulation:	50 mM HEPES, pH 7.8, with 0.5 mM DTT, 0.5 mM EDTA, 300 mM sodium chloride, 2 mM calcium chloride, 10 μ M tetrahydro-L-biopterin (hydrochloride), and 20% glycerol
Concentration:	23.4 mg/mL
Storage/stability:	-80°C/2yr
Specific activity:	3.23 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, with 5 μ M oxyhemoglobin, 1 mM calcium chloride, 20 μ g/ml calmodulin, 0.1 mM NADPH, 50 μ M arginine, 12 μ M tetrahydrobiopterin, and 120 μ M DTT
Activity:	77.80 U/ml
Purity:	Partially pure

Reviewed and approved by: Beth Hurst

Approved date: 23FEB2026

A handwritten signature in dark ink, appearing to read "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