

eNOS (human), (recombinant)

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-853-0010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	eNOS is relatively unstable, and loses activity on ice within 30 minutes. Prepare aliquots and snap-freeze in liquid nitrogen until ready for use.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Nitric oxide synthase (endothelial), NOS III
Formulation	Liquid. In 50 mM TRIS, pH 7.4, containing 0.1mM EDTA, 10% glycerol, 100mM NaCl, 0.1mM DTT, 10 µM tetrahydrobiopterin.
MW	~135kDa/subunit; homodimer.
Purity	≥90% (SDS-PAGE, Western Blot)
Source	Produced in Sf9 insect cells. Untagged human eNOS. Produced in a baculovirus expression system.
Specific Activity	≥120 U/mg. 1 U is defined as the amount of enzyme required to produce 1 nmol of NO (nitric oxide) per minute at 37°C. NO synthesis is measured by the conversion of oxyhemoglobin to methemoglobin.
UniProt ID	P29474

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com