

Nitric Oxide (total), detection kit

**Sensitive, efficient, convenient method for
detection of total Nitric oxide in just 40min**

The Nitric Oxide (total), detection kit is based on the enzymatic conversion of nitrate to nitrite by the enzyme nitrate reductase, followed by the Griess reaction to form a colored azo dye product. Quantification is performed by measuring absorption at 540-570 nm. Only 0.050 ml diluted sample needed per well saves precious sample.

Nitric oxide (NO) is a major mammalian secretory product that initiates host defense, homeostatic and developmental functions by either direct effect or intercellular signaling. NO is the product of a five-electron oxidation of the amino acid L-arginine mediated by nitric oxide synthase. As a direct effector, NO is thought to activate regulatory proteins, kinases and proteases that are directed by reactive oxygen intermediates. Activation of the immune system is associated with an increase in macrophage NO production. Soluble NO exerts a variety of homeostatic influences as an activator of soluble guanylyl cyclase, a neuronal potentiator, a peripheral nervous system neurotransmitter, and a contraction regulator of both smooth muscle and vascular tissue. In addition, NO has been linked to the formation of olfactory and synaptic memories and remodeling.

Citations: 56

[View Online »](#)

Ordering Information

[Order Online »](#)

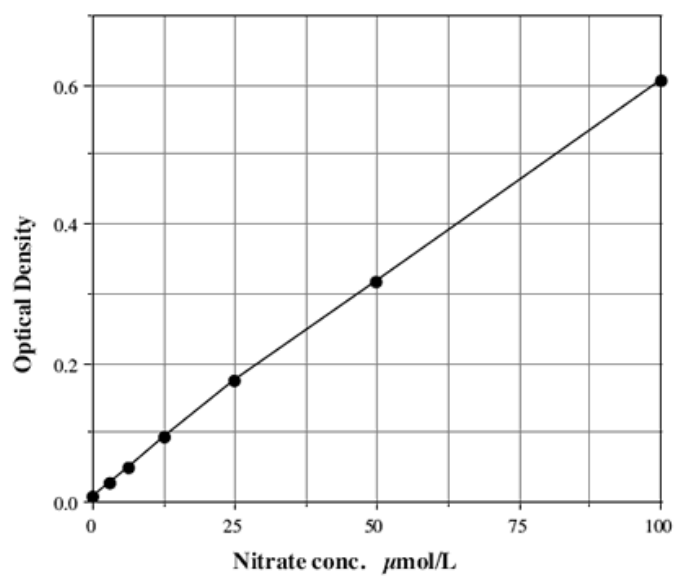
ADI-917-020

2x96 wells

Manuals, SDS & CofA

[View Online »](#)

- Efficient - only 50 µl diluted sample needed, saving your precious sample
- Sensitive - detect as little as 0.625 µmol/L
- Trusted - widely cited in peer-reviewed publications
- Fast - results from up to 80 samples in duplicate in just 40 minutes



Handling & Storage

Use/Stability Store all components at 4°C, except nitrate reductase at -20°C.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

Application Notes For the quantitative determination of total Nitric Oxide in culture supernatants, plasma, serum, saliva, and urine from any species. Cited sample type includes tissue (ref).

Contents Microtiter plate, Nitrate Reductase, Nitrate Reductase Storage Buffer, NADH, Nitrate Standard, Reaction buffer concentrate, Griess Reagent I, Griess Reagent II

Sensitivity 0.625 µmol/L (range 3.125 – 100 µmol/L)

Species Reactivity Species independent

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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