Nitric Oxide (NO2-/NO3-) detection kit

Fast, reliable, easy to use detection kit for separate detection of nitrate and nitrite

The Nitric Oxide (NO_2^-/NO_3^-) , detection kit is a complete kit for the quantitative determination of nitrate and nitrite concentrations. Absorbance is read 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: 27

View Online »

Ordering Information

Order Online »

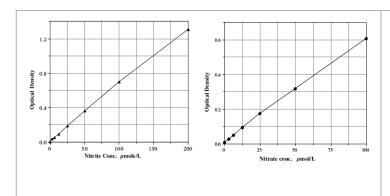
ADI-917-010

2x96 wells

Manuals, SDS & CofA

View Online »

- Homogeneous mix and read assay
- Measures nitrate and nitrite separately
- Fast results in 10 to 40 minutes from up to 39 samples in duplicate



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 nitrite and nitrate

levels in culture supernatants, plasma, serum, saliva, and

urine from any species.

Contents Microtiter plate, Nitrate Reductase, Nitrate Reductase

Storage Buffer, NADH, Nitrate Standard, Nitrite Standard, Reaction buffer concentrate, Griess Reagent I, Griess

Reagent II

Sensitivity (Nitrite) 0.222 µmole/L; (Nitrate) 0.625 µmole/L (range

(Nitrite) 3.125 – 200 µmole/L; (Nitrate) 3.125 – 100

µmole/L)

Species Reactivity Species independent

Last modified: May 29, 2024