

ARBOR ASSAYS™
Interactive Assay Solutions™



NCal™ International Standard Kit

DetectX®

**Nitric Oxide
Colorimetric Detection Kit**

2 Plate Kit Catalog Number K023-H1

Species Independent

Sample Types Validated:

**Water, Buffers, Serum, Plasma,
Urine, Saliva, and TCM**

Please read this insert completely prior to using the product.
For research use only. Not for use in diagnostic procedures.

www.ArborAssays.com



TABLE OF CONTENTS

Background	3
Assay Principle	4
Related Products	4
Supplied Components	5
Storage Instructions	5
Other Materials Required	6
Precautions	6
Sample Types	7
Sample Preparation	7
Reagent Preparation	7-8
Assay Protocol	9
Calculation of Results	10
Typical Data	11-12
Validation Data Sensitivity, Linearity, etc.	12-14
Warranty & Contact Information	15
Plate Layout Sheet	16

BACKGROUND

Nitric oxide (NO) is a diffusible, transient, reactive molecule that has physiological effects in the picomolar-to-micromolar range. Acting through soluble guanylate cyclase activation, NO is an important physiological regulator of the cardiovascular, nervous, and immunological systems¹. NO is bio-available by two routes. It can be endogenously generated by constitutive or induced enzymes like Nitric Oxide Synthase or it can be orally ingested as nitrates and nitrites for rapid uptake into circulation and subsequent conversion².

The reactive nature of nitric oxide allows it to act as a cytotoxic factor when released during an immune response by cells such as macrophages. It also allows NO to be easily converted to a toxic radical that can produce nitrosative damage to cells, organelles and molecules such as DNA. Nitrosylation is also a regulated post-translational modification in cell signaling³. The balance and dynamics of the regulatory/damage facets of NO are major forces in mitochondrial signaling and dysfunction⁴. NO is linked not only to coronary heart disease, endothelial dysfunctions, erectile dysfunction, and neurological disorders, but also diabetes, chronic periodontitis, autism, cancer, and assorted age-related diseases⁵⁻⁹.

The physical properties of Nitric Oxide make it challenging for direct detection methods. However, colorimetric methods can be applied to measure its stable break-down products, nitrate (NO_3^-) and nitrite (NO_2^-)¹⁰.

1. Moncada, S and EA Higgs. (1991) "Endogenous Nitric Oxide: Physiology, Pathology and Clinical Relevance." Eur. J. Clin. Invest., 21:361-374.
2. Kapil, V. et al. (2010) "Inorganic Nitrate and the Cardiovascular System". Heart, 96:1703-1709.
3. Seth, D and Stamler, JS., (2007) "The SNO-proteome: Causation and Classifications." Curr. Opin. Chem. Biol., 15:1-8.
4. Eursolimsky, JD and Moncada, S. (2007) "Nitric Oxide and Mitochondrial Signaling: From Physiology to Pathophysiology." ATVB 27:2524-2531.
5. Knott, AB and Bossy-Wetzel, E., (2010) "Impact of Nitric Oxide on Metabolism in Health and Age-related Disease." Diab.Obes.Metab. 12(Suppl2):126-133.
6. Van Dyke, K. et al. (2010) "Oxidative/Nitrosative Stresses Trigger Type I Diabetes: Preventable in Streptozotocin Rant and Detectable in Human Disease." Ann. N.Y. Acad. Sci. 1203:138-145.
7. Reher, VSG. et al. (2007) "Nitric Oxide Levels in Saliva Increase with Severity of Chronic Periodontitis." J. Oral Sci. 49(4):271-276.
8. Sogut, S. et al. (2003) "Changes in Nitric Oxide Levels and Antioxidant Enzyme Activities May have a Role in the Pathophysiological Mechanisms Involved in Autism." Clin. Chim. Acta 331:111-117.
9. Balam, E. et al. (2002) "Nitric Oxide Levels and Lipid Peroxidation in Plasma of Patients with Gastric Cancer." Jpn. J. Clin. Oncol. 32(5):162-166.
10. Moshage, H. (1997) "Nitric Oxide Determinations: Much Ado About NO.-Thing?." Clin.Chem. 43(4):553-556.

ASSAY PRINCIPLE

The DetectX® Nitric Oxide Detection Kit is designed to quantitatively measure Nitrate and Nitrite present in a variety of samples. Nitric Oxide content is derived from the sum of Nitrate (NO_3^-) and Nitrite (NO_2^-). Please read the complete kit insert before performing this assay. Both Nitrate and Nitrite standards are provided to generate standard curves for the assay and all samples should be read off the appropriate standard curve. For Nitrite detection, samples are mixed with Color Reagents A and B and incubated at room temperature for 5 minutes. The colored product is read at 540 – 570 nm. The concentration of Nitrite in the sample is calculated, after making a suitable correction for any dilution of the sample, using software available with most plate readers.

Total Nitric Oxide content is measured after the sample is incubated with Nitrate Reductase and NADH. The reductase in combination with NADH reduces Nitrate to Nitrite. After a 20 minute incubation at room temperature, Color Reagents A and B are added and incubated at room temperature for 5 minutes. The colored product is read and calculated as with the Nitrite determination above. The concentration of Nitrate in the sample is calculated by subtracting the measured Nitrite concentration from the Total Nitric Oxide concentration for the sample.

This kit uses Nitrate and Nitrite Standard solutions calibrated to the US National Institute for Science and Technology Standard Reference Materials and ISO/IEC standards.

RELATED PRODUCTS

Kits	Catalog No.
Corticosterone ELISA Kits	K014-H1/H5
Cortisol ELISA Kits	K003-H1/H5
Cortisone ELISA Kits	K017-H1/H5
Cyclic Direct AMP Chemiluminescent ELISA Kits	K019-C1/C5
Cyclic Direct AMP ELISA Kits	K019-H1/H5
Cyclic Direct GMP ELISA Kits (Improved Sensitivity)	K065-H1/H5
Prostaglandin E_2 Multi-Format ELISA Kits	K051-H1/H5
Urinary Creatinine Detection Kits	K002-H1 /H5

SUPPLIED COMPONENTS

Clear Half Area 96 well Plates

Corning CoStar Plate 3694.

2 Plates

Catalog Number X018-2EA

Nitrate Standard

Sodium Nitrate at 2,000 μM in a special stabilizing solution. **Calibrated to NIST Standard Reference Material Lot Number 3185.**

200 μL

Catalog Number C086-200UL

Nitrite Standard

Sodium Nitrite at 2,000 μM in a special stabilizing solution. **Calibrated to ISO/IEC 17025.**

200 μL

Catalog Number C087-200UL

Assay Buffer

A buffer containing detergents and stabilizers.

60 mL

Catalog Number X089-60ML

NADH Concentrate

Reduced β -nicotinamide adenine dinucleotide (NADH) as a stable solution.

1.2 mL

Catalog Number X090-1.2ML

Nitrate Reductase

Nitrate Reductase (NR) as a stable solid stored in a ziploc pouch with dessicant.

1 Vial

Catalog Number C088-1EA

Enzyme Stabilization Buffer

A buffer containing special stabilizers for NR.

1 mL

Catalog Number X091-1ML

Color Reagent A

A solution of Sulfanilamide in acid. **CAUTION: CAUSTIC**

5 mL

Catalog Number X092-5ML

Color Reagent B

A solution of N-(1-Naphthyl)ethylenediamine in acid. **CAUTION: CAUSTIC**

5 mL

Catalog Number X093-5ML

STORAGE INSTRUCTIONS

All components of this kit should be stored at 4°C until the expiration date of the kit.

Once reconstituted, the Nitrate Reductase **must** be stored at -20°C.

OTHER MATERIALS REQUIRED

Distilled or deionized water free of detectable nitrate or nitrite.

10,000 Molecular Weight Cut Off (MWCO) polysulfone filters (Corning Spin-X UF 500, 431478) or similar product.

Repeater pipets using disposable tips for addition of Color Reagents A & B, NADH and Nitrate Reductase.

96 well plate reader capable of reading optical absorption at 540-570 nm.

Software for converting optical density (OD) readings from the plate reader and carrying out four parameter logistic curve (4PLC) fitting. Contact your plate reader manufacturer for details.

PRECAUTIONS

As with all such products, this kit should only be used by qualified personnel who have had laboratory safety instruction. The complete insert should be read and understood before attempting to use the product.

The Color Reagents A and B are both acid solutions and should be handled like any laboratory acid.



ARBOR
ASSAYS

SAMPLE TYPES

NO, Nitrate and Nitrite are identical across species and this kit will measure NO from all sources. We determined NO in human samples and the end user should evaluate recoveries of NO in samples from other species being tested. The kit will measure NO in cell culture medium, however many media contain nitrate salts. Care needs to be taken in the selection of media when NO measurement is to be done.

If samples need to be stored after collection, we recommend storing them at -70°C or lower, preferably after being frozen in liquid nitrogen. This assay has been validated for serum, plasma, urine, and saliva, as well as water and buffer samples. Tris, HEPES, and PBS buffers are compatible at pH 7.2, as is EDTA at ≤ 10 mM. Most cell lysates and tissue homogenates should also be compatible. Detergents such as Triton X-100, Tween 20 and CHAPS are compatible at concentrations of $\leq 0.1\%$. Samples containing these detergents should be diluted at least 1:2 with the Assay Buffer. **Samples containing SDS or azide are not compatible with the assay.** Samples containing visible particulate should be centrifuged prior to filtration and using.

SAMPLE PREPARATION

All samples must be filtered through a 10,000 MWCO spin filter to remove protein.

Serum, plasma, saliva, or urine

Dilute sample with Assay Buffer and filter through a 10,000 MWCO device following the manufacturer's recommendations. Collect the filtrates and either further dilute with Assay Buffer as appropriate or use directly in the assay. For serum and plasma, the recommended final dilution is $\geq 1:4$. For urine and saliva, the recommended final dilution is $\geq 1:8$.

REAGENT PREPARATION

Allow the kit reagents to come to room temperature for 30 minutes. Ensure that all samples have reached room temperature and have been diluted and filtered through a 10,000 MWCO filter prior to running them in the kit.

Nitrate Reductase (NR)

Allow the ziploc pouch to warm to room temperature and remove the vial. Add 550 μ L of Enzyme Stabilization Buffer to the vial. Vortex gently and allow to sit at room temperature for 5 minutes. For extended periods of time (> 2 hours) store reconstituted NR on ice. Store any unused reconstituted NR at -20°C.

Prepare NR for use in the assay by taking one part of reconstituted NR and adding to three parts of Assay Buffer. See Table below.

Nitrate Reductase Dilution Table

	1/2 Plate	One Plate	Two Plates
Reconstituted NR	150 μ L	275 μ L	500 μ L
Assay Buffer	450 μ L	825 μ L	1.5 mL
Total Volume	600 μ L	1.1 mL	2 mL

For extended periods of time (> 2 hours) store reconstituted NR on ice.

REAGENT PREPARATION CONTINUED

NADH Preparation

Prepare NADH by diluting one part of NADH Concentrate with an equal part of Assay Buffer.

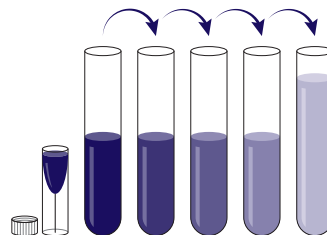
NADH Dilution Table

	1/2 Plate	One Plate	Two Plates
NADH Concentrate	300 µL	550 µL	1 mL
Assay Buffer	300 µL	550 µL	1 mL
Total Reaction Mix Volume	600 µL	1.1 mL	2 mL

Do not store diluted NADH.

Standard Preparation

Nitrate and Nitrite Standards are prepared identically by labeling test tubes as #1 through #7. Briefly vortex to mix and then spin the vial of standard in a microcentrifuge to ensure contents are at bottom of vial. Pipet 360 µL of Assay Buffer into tube #1 and 200 µL into tubes #2 to #7. Carefully add 40 µL of either the NO_2^- or NO_3^- Standard to tube #1 and vortex completely. Take 200 µL of the solution in tube #1 and add it to tube #2 and vortex completely. Repeat this for tubes #3 through #7. The concentration of Nitrate or Nitrite in tubes 1 through 7 will be 200, 100, 50, 25, 12.5, 6.25 and 3.125 µM.



Use all Standards within 2 hours of preparation.

	Std 1	Std 2	Std 3	Std 4	Std 5	Std 6	Std 7
Assay Buffer Vol (µL)	360	200	200	200	200	200	200
Addition	Stock	Std 1	Std 2	Std 3	Std 4	Std 5	Std 6
Vol of Addition (µL)	40	200	200	200	200	200	200
Final Conc (µM)	200	100	50	25	12.5	6.25	3.125

ASSAY PROTOCOL

We recommend that all standards and samples be run in duplicate to allow the end user to accurately determine NO concentrations.

Use the appropriate standards for either Nitrite (NO_2^-) or Nitrate (NO_3^-) determination. All samples should be diluted and filtered through a 10,000 MWCO filter prior to using.

Nitrite Determination Protocol

1. Use the plate layout sheet on the back page to aid in proper sample and standard identification.
2. Pipet 50 μL of samples or Nitrite standards into duplicate wells in the plate.
3. Pipet 50 μL of Assay Buffer into duplicate wells as the Zero standard.
4. Add 25 μL of the Color Reagent A to each well using a repeater pipet.
5. Add 25 μL of the Color Reagent B to each of well using a repeater pipet.
6. Incubate at room temperature for 5 minutes.
7. Read the optical density at 540-570 nm. These readings are for the Nitrite determination.

Total Nitric Oxide Determination Protocol

1. Use the plate layout sheet on the back page to aid in proper sample and standard identification.
2. Pipet 50 μL of samples or Nitrate standards into duplicate wells in the plate.
3. Pipet 50 μL of Assay Buffer into duplicate wells as the Zero standard.
4. Add 10 μL of prepared NADH to each well using a repeater pipet.
5. Add 10 μL of prepared NR to each well using a repeater pipet.
6. Incubate at room temperature for 20 minutes.
7. Add 25 μL of the Color Reagent A to each well using a repeater pipet.
8. Add 25 μL of the Color Reagent B to each of well using a repeater pipet.
9. Incubate at room temperature for 5 minutes.
10. Read the optical density at 540-570 nm. These readings are for the Total Nitric Oxide determination.

CALCULATION OF RESULTS

Average the duplicate optical density readings for each standard and sample. Create a standard curve by reducing the data using the 4PLC fitting routine on the plate reader, after subtracting the mean ODs for the zero standard. The concentrations obtained should be multiplied by the dilution factor to obtain sample values.

Or use the online tool from MyAssays to calculate the data:

www.myassays.com/arbor-assays-nitric-oxide-colorimetric-kit.assay

Nitrite

Nitrite (NO_2) concentrations are calculated from the data obtained from the Nitrite Protocol standard curve data utilizing the curve fitting routine supplied with the plate reader.

Total NO

Total NO concentrations are calculated from the data obtained from the Total Nitric Oxide Protocol (nitrite + nitrate) standard curve data utilizing the curve fitting routine supplied with the plate reader.

Nitrate

Nitrate (NO_3) concentrations are obtained by subtracting the NO_2 concentrations of each sample from the Total NO concentrations. See Below:

$$\text{Nitrate } (\text{NO}_3) = \text{Total NO} - \text{Nitrite } (\text{NO}_2)$$



TYPICAL DATA - NITRITE

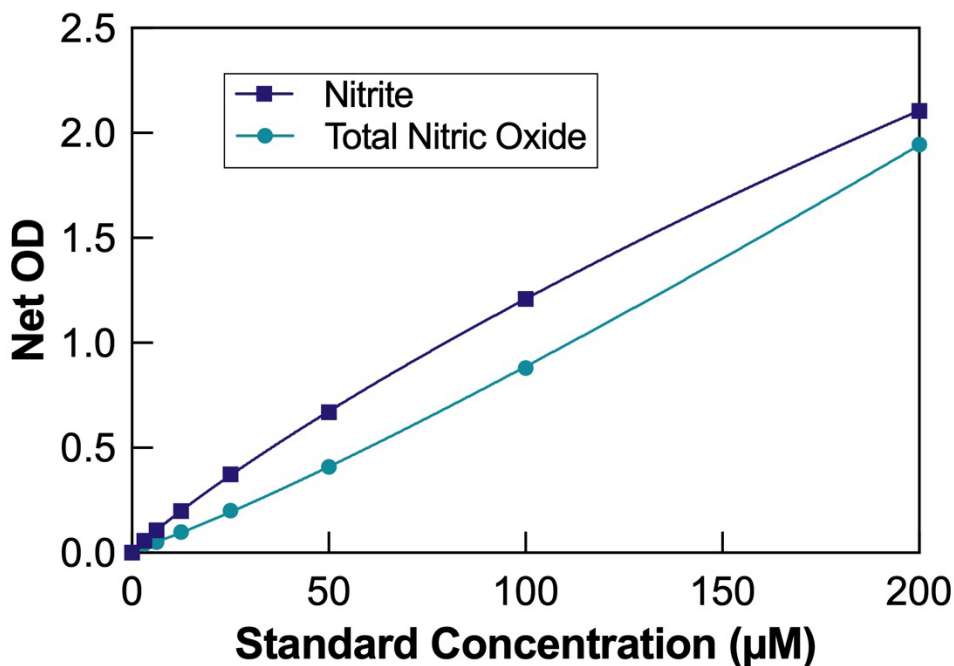
Sample	Mean OD	Net OD	Nitrite Conc. (µM)
Zero	0.038	0	0
Standard 1	2.144	2.106	200
Standard 2	1.248	1.210	100
Standard 3	0.708	0.670	50
Standard 4	0.412	0.374	25
Standard 5	0.236	0.198	12.5
Standard 6	0.145	0.107	6.25
Standard 7	0.095	0.057	3.125
Sample 1	0.639	0.601	43.9
Sample 2	1.484	1.446	124.4

TYPICAL DATA - TOTAL NITRIC OXIDE

Sample	Mean OD	Net OD	Total Nitric Oxide Conc. (µM)
Zero	0.040	0	0
Standard 1	1.984	1.944	200
Standard 2	0.921	0.881	100
Standard 3	0.450	0.410	50
Standard 4	0.240	0.200	25
Standard 5	0.138	0.098	12.5
Standard 6	0.092	0.052	6.25
Standard 7	0.083	0.043	3.125
Sample 1	1.058	1.018	113.1
Sample 2	1.420	1.380	147.9

Always run your own standard curve for calculation of results. Do not use this data.

Typical Standard Curves



Always run your own standard curves for calculation of results. Do not use this data.

VALIDATION DATA

Sensitivity and Limit of Detection

Sensitivity was calculated by comparing the ODs for twenty wells run for each of the zero and standard #7. The detection limit was determined at two (2) standard deviations from the zero along the standard curve. **Sensitivity was determined as 2.63 µM in the Nitrite and 1.02 µM in the Total Nitric Oxide assays.**

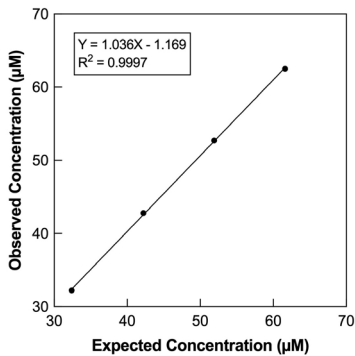
The Limit of Detection was determined in a similar manner by comparing the ODs for twenty wells run for each of the zero and a low concentration human sample. **The Limit of Detection was determined as 0.94 µM in the Nitrite and 3.0 µM in the Total Nitric Oxide assays.**

Linearity

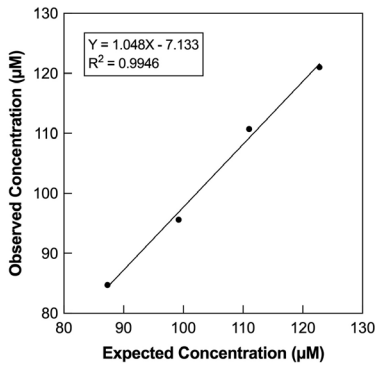
Linearity was determined by taking two diluted human urine samples with known Nitrite and Total NO Concentrations and mixing them in the ratios given below. The measured concentrations were compared to the expected values based on the ratios used.

High Urine Sample	Low Urine Sample	Expected Conc. (µM)		Observed Conc. (µM)		% Recovery	
		Nitrite	Total NO	Nitrite	Total NO	Nitrite	Total NO
80%	20%	61.6	122.8	62.5	121.0	101.5%	98.5%
60%	40%	51.9	111.0	52.7	110.7	101.5%	99.7%
40%	60%	42.2	99.2	42.8	95.6	101.6%	96.4%
20%	80%	32.4	87.3	32.2	84.7	99.5%	97.0%
Mean Recovery						101.0%	97.9%

Nitrite Linearity



Total NO Linearity



Intra Assay Precision

Three samples were diluted in Assay Buffer and run in replicates of 20 in an assay. The mean and precision of the calculated Nitrite or Total NO concentrations were:

Sample	Nitrite Conc. (µM)	%CV	Total NO Conc. (µM)	%CV
1	45.1	4.4	70.5	6.8
2	73.3	9.1	107.4	4.4
3	132.7	1.3	157.8	1.8

Inter Assay Precision

Three samples were diluted in Assay Buffer and run in duplicates in twenty assays run over multiple days by three operators. The mean and precision of the calculated Nitrite or Total NO concentrations were:

Sample	Nitrite Conc. (µM)	%CV	Total NO Conc. (µM)	%CV
1	44.1	3.1	68.8	7.4
2	66.4	4.0	112.1	5.7
3	126.7	6.3	154.4	4.1

LIMITED WARRANTY

Arbor Assays warrants that at the time of shipment this product is free from defects in materials and workmanship. This warranty is in lieu of any other warranty expressed or implied, including but not limited to, any implied warranty of merchantability or fitness for a particular purpose.

We must be notified of any breach of this warranty within 48 hours of receipt of the product. No claim shall be honored if we are not notified within this time period, or if the product has been stored in any way other than outlined in this publication. The sole and exclusive remedy of the customer for any liability based upon this warranty is limited to the replacement of the product, or refund of the invoice price of the goods.

CONTACT INFORMATION

For details concerning this kit or to order any of our products please contact us:

Arbor Assays

1514 Eisenhower Place
Ann Arbor, Michigan 48108 USA

Phone: 734-677-1774

Web: www.ArborAssays.com

Email Addresses:

Info@ArborAssays.com

Orders@ArborAssays.com

Technical@ArborAssays.com



OFFICIAL SUPPLIER TO ISWE

Arbor Assays and the International Society of Wildlife Endocrinology (ISWE) signed an exclusive agreement for Arbor Assays to supply ISWE members with assay kits and reagents for wildlife conservation research.

DetectX[®], ThioStar[®] and the Arbor Assays logo are all registered trademarks.

A	1	2	3	4	5	6	7	8	9	10	11	12
B												
C												
D												
E												
F												
G												
H												



Printed on Forest Stewardship Council certified paper