

# NO<sub>2</sub>/NO<sub>3</sub> Assay Kit-FX(Fluorometric) -2,3-Diaminonaphthalene Kit-

## Technical Manual

Technical Manual (Japanese version) is available at <http://www.dojindo.co.jp/manual/nk08.pdf>

### General Information

Nitric oxide (NO), a gaseous paramagnetic radical, is a very important and versatile messenger in biological systems. NO is synthesized from L-arginine by NO synthase (NOS). It has been identified as an endothelial derived relaxation factor (EDRF) and antiplatelet substance. It serves as a neurotransmitter derived from a neutrophil and a cytotoxic substance from an activated macrophage. Although NO's molecular action in the biological system is very versatile, the most important role of NO is the activation of guanylate cyclase.

The Griess assay is one of the most popular and simplest methods to detect the NO concentration. However, the sensitivity of Griess assay is sometimes not sufficient to determine μmol/l level of NO<sub>2</sub><sup>-</sup>. 2,3-Diaminonaphthalene (DAN) reacts with NO<sub>2</sub><sup>-</sup> and generates a fluorescent product as shown in Figure 1.

Therefore, the fluorescence detection method using DAN can be used to determine the NO<sub>2</sub><sup>-</sup> concentration more sensitively.

The reaction of 2,3-Diaminonaphthalene with NO<sub>2</sub><sup>-</sup> proceeds under pH 2 at room temperature within 5 minutes, and the fluorescence of produced naphthalenetriazole

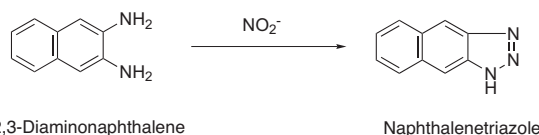


Figure 1. NO<sub>2</sub><sup>-</sup> reaction with 2,3-Diaminonaphthalene

can be determined at pH10 or higher. The NO<sub>2</sub>/NO<sub>3</sub> Assay Kit-FX contains 2,3-diaminonaphthalene, nitrate reductase, enzyme cofactor, buffer solution and NO<sub>2</sub><sup>-</sup>, NO<sub>3</sub><sup>-</sup> solutions as standard. Therefore, NO metabolites, NO<sub>2</sub><sup>-</sup> and NO<sub>3</sub><sup>-</sup>, are easily detectable using this kit. The suitable detection range is from 1 to 10 μmol/l.

The excitation and emission maximum of naphthalenetriazole is 365 nm and 410-425 nm, respectively. However, in order to reduce the fluorescence blank of 2,3-diaminonaphthalene and increase the sensitivity of the NO detection, the use of a 450-465 nm emission filter is recommended.

### Contents of this kit

- Buffer Solution	x 1	- NO <sub>3</sub> Reductase	x 2
- NaNO <sub>2</sub> Standard Solution (200 μmol/l)	x 2	- Enzyme Cofactors	x 2
- NaNO <sub>3</sub> Standard Solution (200 μmol/l)	x 2	- Stop Solution	x 1
- Fluorescence Reagent (DAN) Solution	x 1		

### Storage

Please store this kit at 0-5 °C. The reconstituted solutions of NO<sub>3</sub> Reductase and Enzyme Cofactors should be stored at 0-5 °C, and be used up within 2 weeks. Do not freeze this kit and the reconstituted solutions.

### Precaution

- This kit contains glass vials with an aluminium cap. Please handle carefully.
- Since inside of each vial of NO<sub>3</sub> Reductase and the Enzyme Cofactors is under reduced pressure, add 1.2 ml Buffer Solution into the vial with a syringe in order to avoid the dispersal of the powder. Please do not open the rubber septum at the beginning.

### Required Equipments and Materials

- Fluorescence microplate reader (excitation filter: 360-365 nm, emission filter: 450-465 nm)
- 96 well microplate (black plate or white plate)
- 10 μl, 100-200 μl pipettes, multi-channel pipette

### Preparation of Reagent solutions

- 1) 20 μmol/l NaNO<sub>2</sub> Standard Solution :  
Add 9 ml of Buffer Solution ( or medium ) to a bottle of NaNO<sub>2</sub> Standard Solution.
- 2) 20 μmol/l NaNO<sub>3</sub> Standard Solution :  
Add 9 ml of Buffer Solution ( or medium ) to a bottle of NaNO<sub>3</sub> Standard Solution.
- 3) NO<sub>3</sub> Reductase Solution :  
Add 1.2 ml of Buffer Solution to a bottle of NO<sub>3</sub> Reductase.
- 4) Enzyme Cofactors Solution :  
Add 1.2 ml of Buffer Solution to a bottle of Enzyme Cofactors.

### Preparation of NO<sub>2</sub><sup>-</sup> Standard Curve

- 1) Dilute 80 μl of 20 μmol/l NaNO<sub>2</sub> Standard Solution by serial dilution with 80 μl of Buffer Solution (or medium) in 96-well plate as indicated in Figure 2.
- 2) Add 20 μl of Buffer Solution to each well (total volume is 100 μl/well).
- 3) Add 10 μl of Fluorescence Reagent (DAN) Solution to each well, and mix. Then, leave the plate for 15 minutes at room temperature.
- 4) Add 40 μl of Stop Solution to each well.
- 5) Measure the fluorescence intensity of each well at 450-465 nm (excitation 360-365nm) with a fluorescence microplate reader.

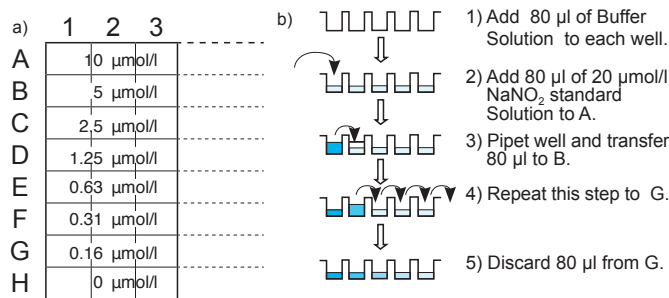


Figure 2. A plate arrangement for preparation of a standard curve.

### General protocol for 96-well micropalate assay

### Preparation of $\text{NO}_3^- + \text{NO}_2^-$ Standard Curve

- 1) Dilute 80  $\mu\text{l}$  of 20  $\mu\text{mol/l}$   $\text{NaNO}_3$  Standard Solution by serial dilution with 80  $\mu\text{l}$  of Buffer Solution (or medium) in 96-well plate as indicated in Figure 2.
- 2) Add 10  $\mu\text{l}$  of  $\text{NO}_3$  Reductase Solution and 10  $\mu\text{l}$  of Enzyme Cofactors Solution to each well.
- 3) Incubate the plate at 37  $^\circ\text{C}$  for 30 minutes.
- 4) Add 10  $\mu\text{l}$  of Fluorescence Reagent (DAN) Solution to each well, and mix. Then, leave the plate for 15 minutes at room temperature for complete reaction.
- 5) Add 40  $\mu\text{l}$  of Stop Solution to each well.
- 6) Measure the fluorescence intensity of each well at 450-465 nm (excitation 360-365 nm) with a fluorescence microplate reader.

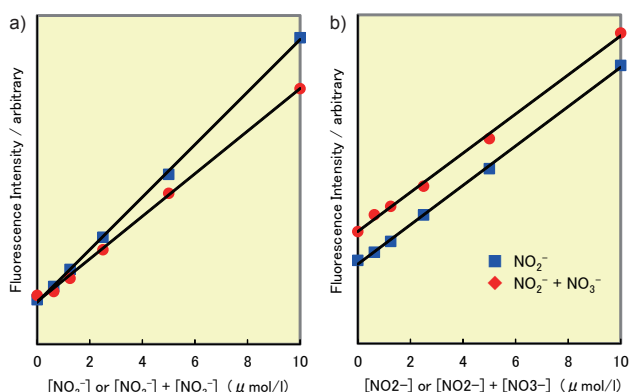


Figure 3. Standard curves of  $\text{NO}_2^-$  and  $\text{NO}_2^- + \text{NO}_3^-$  using  $\text{NO}_2/\text{NO}_3$  Assay Kit-FX. a) in Buffer Solution, b) in a medium (DMEM).

### Preparation of Sample Solution

Cell or tissue culture medium:

Use phenol red free medium. Centrifuge the medium at 1,000  $\times g$  for 15 minutes, and use the supernatant as a sample solution. A cell culture medium that contains  $\text{NO}_3^-$  as a component (such as RPMI1640) cannot be used to detect  $\text{NO}_3^-$ . However,  $\text{NO}_2^-$  in the medium can be determined using this kit.

Serum:

Serum or plasma sample should be treated with Amicon Ultra-4 Centrifugal Filter Unit with Ultracel-10 membrane (7,000  $\times g$ , 4  $^\circ\text{C}$ , 20minutes) to remove hemoglobin and proteins.

### Determination of $\text{NO}_2^-$ Concentration in Sample Solution

- 1) Add 80  $\mu\text{l}$  of a sample solution to a well.
- 2) Add 20  $\mu\text{l}$  of Buffer Solution to each well.
- 3) Add 10  $\mu\text{l}$  of the Fluorescence Reagent (DAN) Solution to each well, and mix. Then, leave the plate for 15 minutes at room temperature for complete reaction.
- 4) Add 40  $\mu\text{l}$  of Stop Solution to each well.
- 5) Measure the fluorescence intensity of each well at 450-465 nm (excitation 360-365 nm) with a fluorescence microplate reader.
- 6) Determine the concentration of  $\text{NO}_2^-$  in the sample solution from the standard curve.

### Determination of $\text{NO}_3^- + \text{NO}_2^-$ Concentration in Sample Solution

- 1) Add 80  $\mu\text{l}$  of a sample solution to a well.
- 2) Add 10  $\mu\text{l}$  of the  $\text{NO}_3$  Reductase Solution and 10  $\mu\text{l}$  of the Enzyme Cofactors Solution to each well.
- 3) Incubate the plate at 37  $^\circ\text{C}$  for 30 minutes.
- 4) Add 10  $\mu\text{l}$  of the Fluorescence Reagent (DAN) Solution to each well, and mix. Then leave the plate for 15 minutes at room temperature for complete reaction.
- 5) Add 40  $\mu\text{l}$  of Stop Solution to each well.
- 6) Measure the fluorescence intensity of each well at 450-465 nm (excitation 360-365 nm) with a fluorescence microplate reader.
- 7) Determine the concentration of  $\text{NO}_3^- + \text{NO}_2^-$  in the sample solution using the standard curve.

### Determination of $\text{NO}_3^-$ Concentration in Sample Solution

$\text{NO}_3^-$  concentration can be obtained by the following equation.

$$[\text{NO}_3^-] = [\text{NO}_2^- + \text{NO}_3^-] - [\text{NO}_2^-]$$

### Notes

1. The  $\text{NO}_3$  Reductase Solution and the Enzyme Cofactors Solution should be kept on an ice bath during the use. Please store these solutions at 0-5  $^\circ\text{C}$ . The solutions should be used up within 2 weeks.
2. This kit can be used to determine  $\text{NO}_3^-$  and  $\text{NO}_2^-$  in the concentration range from 1 to 10  $\mu\text{mol/l}$ . Dilute the sample solution if it contains high concentration of  $\text{NO}_3^-$  and  $\text{NO}_2^-$ . If the sample volume is less than 80  $\mu\text{l}$ , add Buffer Solution to be the total volume of 80  $\mu\text{l}$ .
3. Do not mix the  $\text{NO}_3$  Reductase Solution and the Enzyme Cofactors Solution prior to use.
4.  $\text{NO}_3$  Reductase and the Enzyme Cofactors are packaged under reduced pressure. Please open the caps with caution, or add solvent before opening.
5. The Fluorescent Reagent (DAN) Solution is light sensitive. Please keep it in dark at all times.
6. For fluorophotometric assay, dilute 100  $\mu\text{l}$  of the solution, treated with Stop Solution at the final step with 3.9 ml of ddH<sub>2</sub>O, and measure the fluorescence intensity.

If you need more information, please contact Dojindo technical service.

Dojindo Laboratories

2025-5 Tabaru, Mashiki-machi, Kamimashiki-gun, Kumamoto  
861-2202, Japan Phone: +81-96-286-1515 Fax: +81-96-286-1525  
E-mail: info@dojindo.co.jp Web: www.dojindo.co.jp

Dojindo Molecular Technologies, Inc.

Tel: +1-301-987-2667 Web: http://www.dojindo.com/

Dojindo EU GmbH

Tel: +49-89-3540-4805 Web: http://www.dojindo.eu.com/

Dojindo China Co., Ltd

Tel: +86-21-6427-2302 Web: http://www.dojindo.cn/

NK08 :  $\text{NO}_2/\text{NO}_3$  Assay Kit-FX