

Amplite™ Fluorimetric NAD/NADH Ratio Assay Kit *Red Fluorescence*

Ordering Information

Product Number: 15263 (250 assays)

Storage Conditions

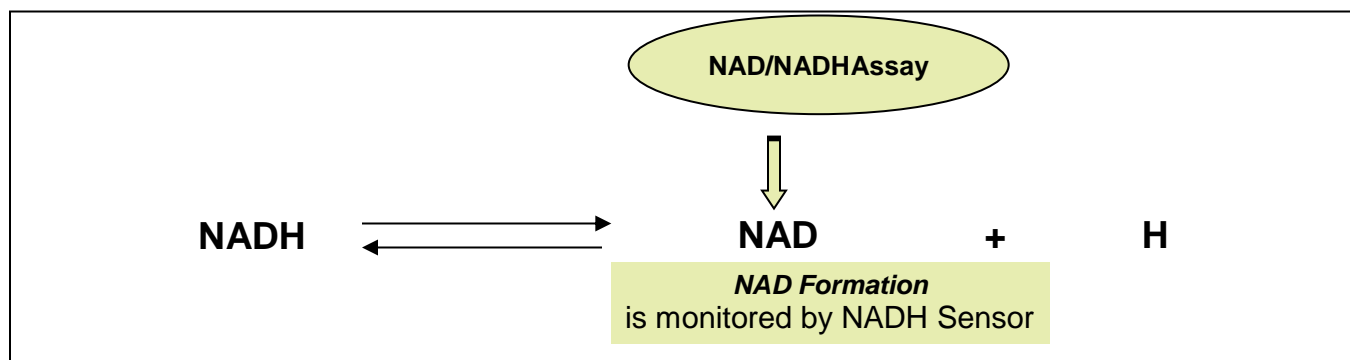
Keep in freezer Avoid exposure to light.

Instrument Platform

Fluorescence microplate readers

Introduction

Nicotinamide adenine dinucleotide (NAD⁺) and nicotinamide adenine dinucleotide phosphate (NADP⁺) are two important cofactors found in cells. NADH is the reduced form of NAD⁺, and NAD⁺ is the oxidized form of NADH. NAD forms NADP with the addition of a phosphate group to the 2' position of the adenylyl nucleotide through an ester linkage. NADP is used in anabolic biological reactions, such as fatty acid and nucleic acid synthesis, which require NADPH as a reducing agent. In chloroplasts, NADP is an oxidizing agent important in the preliminary reactions of photosynthesis. The NADPH produced by photosynthesis is used as reducing power for the biosynthetic reactions in the Calvin cycle of photosynthesis. The traditional NAD/NADH and NADP/NADPH assays are done by monitoring the changes in NADH or NADPH absorption at 340 nm. This method suffers low sensitivity and high interference since the assay is done in the UV range that requires expensive quartz microplates.



This Amplite™ Fluorimetric NAD/NADH Ratio Assay Kit provides a convenient method for sensitive detection of NAD, NADH and their ratio. The enzymes in the system specifically recognize NAD/NADH in an enzyme cycling reaction that significantly increases detection sensitivity. In addition, this assay has very low background since it is run in the red visible range that considerably reduces the interference from biological samples. There is no need to purify NAD/NADH from sample mix. The assay can be performed in a convenient 96-well or 384-well microtiter-plate format. Its signal can be easily read by either a fluorescence microplate reader at Ex/Em = 530 - 570/590 - 600 nm (maximum Ex/Em = 540/590 nm) or an absorbance microplate reader at ~576 nm. This kit provides NAD and NADH extraction buffer, and cell lysis buffer for your convenience. It has been frequently used for determining NAD/NADH from cell lysates.

Kit Components

| Components | Amount |
|--|--------------------------------|
| Component A: NAD/NADH Recycling Enzyme Mixture | 2 bottles (lyophilized powder) |
| Component B: NADH Sensor Buffer | 1 bottle (20 mL) |
| Component C: NADH Standard | 1 vial (142 µg) |
| Component D: NADH Extraction Solution | 1 bottle (10 mL) |
| Component E: NAD Extraction Solution | 1 bottle (10 mL) |
| Component F: NAD/NADH Control Solution | 1 bottle (10 mL) |
| Component G: NAD/NADH Lysis Buffer | 1 bottle (10 mL) |

Assay Protocol for One 96-Well Plate

Brief Summary

Prepare 25 µL of NADH standards and/or test samples → Add 25 µL of NADH or NAD Extraction Solution
→ Incubate at room temperature for 15 minutes → Add 25 µL of NAD or NADH Extraction Solution
→ Add 75 µL of NAD/NADH reaction mixture → Incubate at RT for 15 minutes to 2 hours
→ Monitor fluorescence intensity at Ex/Em = 540/590 nm

Note: Thaw one of each kit component at room temperature before starting the experiment.

1. Prepare NADH stock solution:

Add 200 μL of PBS buffer into the vial of NADH standard (Component C) to have 1 mM (1 nmol/ μL) NADH stock solution.

Note: The unused NADH stock solution should be divided into single use aliquots and stored at -20°C .

2. Prepare NAD/NADH reaction mixture:

Add 10 mL of NADH Sensor Buffer (Component B) to the bottle of NAD/NADH Recycling Enzyme Mixture (Component A), and mix well.

Note: This NAD/NADH reaction mixture is enough for two 96-well plates. The unused NAD/NADH reaction mixture should be divided into single use aliquots and stored at -20°C .

3. Prepare serially diluted NADH standards (0 to 10 μM):

3.1 Add 30 μL of 1 mM NADH stock solution (from Step 1) into 970 μL PBS buffer (pH 7.4) to generate 30 μM (30 pmols/ μL) NADH standard solution.

Note: Diluted NADH standard solution is unstable, and should be used within 4 hours.

3.2 Take 200 μL of 30 μM NADH standard solution (from Step 3.1) to perform 1:3 serial dilutions to get 10, 3, 1, 0.3, 0.1, 0.03 and 0 μM serially diluted NADH standards.

3.3 Add serially diluted NADH standards and/or NAD/NADH containing test samples into a solid black 96-well microplate as described in Tables 1 and 2.

Note: Prepare cells or tissue samples as desired.

Table 1. Layout of NADH standards and test samples in a solid black 96-well microplate

| | | | | | | | | | | | |
|-----|-----|------|------|-----------|-----------|----------|----------|--|--|--|--|
| BL | BL | TS | TS | TS (NADH) | TS (NADH) | TS (NAD) | TS (NAD) | | | | |
| NS1 | NS1 | | | | | | | | | | |
| NS2 | NS2 | | | | | | | | | | |
| NS3 | NS3 | | | | | | | | | | |
| NS4 | NS4 | | | | | | | | | | |
| NS5 | NS5 | | | | | | | | | | |
| NS6 | NS6 | | | | | | | | | | |
| NS7 | NS7 | | | | | | | | | | |

Note: NS= NAD/NADH Standards; BL=Blank Control; TS=Test Samples; TS (NADH) = Test Samples treated with NADH Extraction Solution for 10 to 15 minutes, then neutralized by NAD Extraction Solution; TS (NAD) = Test Samples treated with NAD Extraction Solution for 10 to 15 minutes, then neutralized by NADH Extraction Solution.

Table 2 Reagent composition for each well

| NADH Standard | Blank Control | Test Sample (NAD/NADH) | Test Sample (NADH Extract) | Test Sample (NAD Extract) |
|---|-------------------------------|-------------------------------|-------------------------------|-------------------------------|
| Serial Dilutions*: 25 μL | PBS: 25 μL | Test Sample: 25 μL | Test Sample: 25 μL | Test Sample: 25 μL |
| Component F: 25 μL | Component F: 25 μL | Component F: 25 μL | Component D: 25 μL | Component E: 25 μL |
| Incubate at room temperature for 10 to 15 minutes | | | | |
| Component F: 25 μL | Component F: 25 μL | Component F: 25 μL | Component E: 25 μL | Component D: 25 μL |
| Total: 75 μL | Total: 75 μL | Total: 75 μL | Total: 75 μL | Total: 75 μL |

**Note: Add the serially diluted NADH standards from 0.03 μM to 30 μM into wells from NS1 to NS7 in duplicate. High concentration of NADH (e.g., > 300 μM , final concentration) may cause reduced fluorescence signal due to the over oxidation of NADH sensor (to a non-fluorescent product).*

3.4 **For NADH Extraction (NADH):** Add 25 μL of NADH Extraction Solution (Component D) into the wells of NAD/NADH containing test samples. Incubate at room temperature for 10 to 15 minutes, then add 25 μL of NAD Extraction Solution (Component E) to neutralize the NADH extracts as described in Tables 1 & 2.

For NAD Extraction (NAD): Add 25 μL of NAD Extraction Solution (Component E) into the wells of NAD/NADH containing test samples. Incubate at room temperature for 10 to 15 minutes, then add 25 μL of NADH Extraction Solution (Component D) to neutralize the NAD extracts as described in Tables 1 & 2.

For Total NAD and NADH: Add 25 μL of NAD/NADH Control Solution (Component F) into the wells of NADH standards and NAD/NADH containing test samples. Incubate at room temperature for 10 to 15 minutes, and then add 25 μL of Control Solution (Component F) as described in Tables 1 and 2.

Note1: Prepare cells or tissue samples as desired. NAD/NADH Lysis Buffer (Component G) can be used for lysing the cells (See appendix for details).

Note2: In healthy mammalian cells, there is more NAD compare to NADH, so one can simply use total NAD and NADH minus the NAD to calculate the amount of NADH.

4. Run NAD/NADH assay in supernatants reaction:

4.1 Add 75 μL of NADH reaction mixture (from Step 2) into each well of NADH standard, blank control, and test samples (from Step 3.4) to make the total NADH assay volume of 150 μL /well.

4.2 Incubate the reaction at room temperature for 15 minutes to 2 hours, protected from light.

4.3 Monitor the fluorescence increase with a fluorescence plate reader at Ex/Em = 540/590 nm (cutoff 570 nm).

Note: The contents of the plate can also be transferred to a white clear bottom plate and read by an absorbance microplate reader at the wavelength of 576 ± 5 nm. The absorption detection has lower sensitivity compared to fluorescence reading.

Data Analysis

The fluorescence in blank wells (with the PBS buffer only) is used as a control, and is subtracted from the values for those wells with the NADH reactions. The typical data are shown in Figure 1 (Total NAD and NADH vs. NAD or NADH Extract).

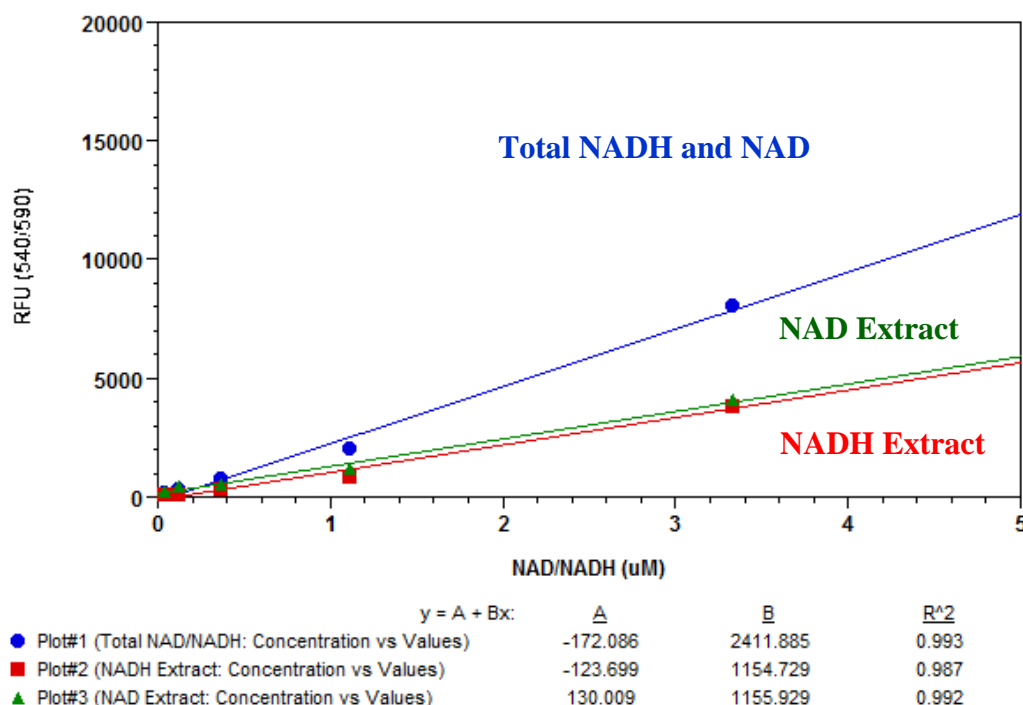


Figure 1. Total NADH and NAD, and their extract dose response were measured with Amplite™ NAD/NADH Ratio Assay Kit in a 96-well black plate using a Gemini microplate reader (Molecular Devices). 25 μL of equal amount of NAD and NADH was treated with or without NADH or NAD extraction solution for 15 minutes, and then neutralized with extraction solutions at room temperature. The signal was acquired at Ex/Em = 540/590 nm (cut off at 570 nm) 30 minutes after adding 75 μL of NADH reaction mixture. The blank signal was subtracted from the values for those wells with the NADH reactions. (*Note: The fluorescence background increases with time, thus it is important to subtract the fluorescence intensity value of the blank wells for each data point.*)

Appendix: Test Sample Preparations Using Component G (NAD/NADH Lysis Buffer)

1. Plant Cell Samples:

Homogenize leaves with the lysis buffer at 200 mg/mL, and centrifuge at 2500 rpm for 5-10 minutes, use the supernatant for tests.

2. Bacterial Cell Samples:

Collect bacterial cells by centrifugation ((10,000 g, 0°C, 15 min). Use about 100 to 10 million cells/mL lysis buffer, keep the treated solution at room temperature for 15 minutes. Centrifuge at 2500 rpm for 5 minutes, and use the supernatant for tests.

3. Mammalian Cell Samples:

Remove medium from plate wells, use about 100 µL lysis buffer per 1-5 million cells (or 50-100 µL/well in a 96-well cell culture plate), and keep the treated solution at room temperature for 15 minutes. Use the cell lysate directly or centrifuge it at 1500 rpm for 5 minutes, use the supernatant for tests.

4. Tissue Samples:

Weigh ~20 mg tissue, wash with cold PBS. Homogenize with 400 µl of lysis buffer in a micro-centrifuge tube. Centrifuge at 2500 rpm for 5-10 minutes, use the supernatant for the assay.

References

- Minglong Shao, Xian Zhang, Zhiming Rao, Meijuan Xu, Taowei Yang, Hui Li, Zhenghong Xu and Shangtian Yang. Efficient testosterone production by engineered *Pichia pastoris* co-expressing human 17β-hydroxysteroid dehydrogenase type 3 and *Saccharomyces cerevisiae* glucose 6-phosphate dehydrogenase with NADPH regeneration. *Green Chem.*, 2016,18, 1774-1784. DOI: 10.1039/C5GC02353J.
- Shi-Qi Ji, Bing Wang, Ming Lu, Fu-Li Li. Direct bioconversion of brown algae into ethanol by thermophilic bacterium *DeFluviitalea phaphyphila*. *Biotechnology for Biofuels*. December 2016, 9:81.
- Tinggang Li, Yu Yan and Jianzhong He. Enhanced direct fermentation of cassava to butanol by *Clostridium* species strain BOH3 in cofactor-mediated medium. *Biotechnology for Biofuels*. 2015,8:166. DOI: 10.1186/s13068-015-0351-7.
- Eugenia Villa-Cuesta, Marissa A. Holmbeck and David M. Rand. Rapamycin increases mitochondrial efficiency by mtDNA-dependent reprogramming of mitochondrial metabolism in *Drosophila*. *Journal of Cell Science* (2014) 127, 2282–2290 doi:10.1242/jcs.142026.
- Tinggang Li, Yu Yan, Jianzhong He. Reducing cofactors contribute to the increase of butanol production by a wild-type *Clostridium* sp. strain BOH3. *Bioresour Technol*. 2014 Mar;155:220-8. doi: 10.1016/j.biortech.2013.12.089. Epub 2014 Jan 2.
- Geng Tian, Jinko Sawashita, Hiroshi Kubo, Shin-ya Nishio, Shigenari Hashimoto, Nobuyoshi Suzuki, Hidekane Yoshimura, Mineko Tsuruoka, Yaoyong Wang, Yingye Liu, Hongming Luo, Zhe Xu, Masayuki Mori, Mitsuki Kitano, Kazunori Hosoe, Toshio Takeda, Shin-ichi Usami, and Keiichi Higuchi. Ubiquinol-10 Supplementation Activates Mitochondria Functions to Decelerate Senescence in Senescence-Accelerated Mice. *Antioxid Redox Signal*. 2014 June 1; 20(16): 2606–2620. doi: 10.1089/ars.2013.5406.
- Bao, Teng; Zhang, Xian; Rao, Zhiming; Zhao, Xiaojing; Zhang, Rongzhen; et al Efficient Whole-Cell Biocatalyst for Acetoin Production with NAD+ Regeneration System through Homologous Co-Expression of 2,3-Butanediol Dehydrogenase and NADH Oxidase in Engineered *Bacillus subtilis*: e102951. *PLoS One* 9.7 (Jul 2014).
- Pengbo Yao, Xiaobo Chen, Yan Yan, Feng Liu, Yuanying Zhang, Xingqi Guo, Baohua Xu. Glutaredoxin 1, glutaredoxin 2, thioredoxin 1, and thioredoxin peroxidase 3 play important roles in antioxidant defense in *Apis cerana cerana*. *Free Radical Biology and Medicine*. Volume 68, March 2014, Pages 335–346.
- Lamonte G, Tang X, Chen JL, Wu J, Ding CK, Keenan MM, Sangokoya C, Kung HN, Ilkayeva O, Boros LG, Newgard CB, Chi JT. Acidosis induces reprogramming of cellular metabolism to mitigate oxidative stress. *Cancer Metab*. 2013 Dec 23;1(1):23. doi: 10.1186/2049-3002-1-23.
- Chao Tong, Alex Morrison, Samantha Mattison, Su Qian, Mark Bryniarski, Bethany Rankin, Jun Wang, D. Paul Thomas, and Ji Li. Impaired SIRT1 nucleocytoplasmic shuttling in the senescent heart during ischemic stress. *FASEB J*, Nov 2013; 27: 4332 - 4342.
- Rubin Tan, Jiansha Li, Xiaochun Peng, Liping Zhu, Lei Cai, Tao Wang, Yuan Su, Kaikobad Irani, and Qinghua Hu. GAPDH is critical for superior efficacy of female bone marrow-derived mesenchymal stem cells on pulmonary hypertension. *Cardiovasc Res*, Oct 2013; 100: 19 - 27.
- Stephen Y. Xue, Valeria Y. Hebert, Danicia M. Hayes, Corie N. Robinson, Mitzi Glover, and Tammy R. Dugas. Nucleoside Reverse Transcriptase Inhibitors Induce a Mitophagy-Associated Endothelial Cytotoxicity That Is Reversed by Coenzyme Q10 Cotreatment. *Toxicol. Sci.*, Aug 2013; 134: 323 - 334.
- Kate J. Roberts, Andrew Cross, Olga Vasieva, Robert J. Moots, and Steven W. Edwards. Inhibition of pre-B cell colony-enhancing factor (PBEF/NAMPT/visfatin) decreases the ability of human neutrophils to generate reactive oxidants but does not impair bacterial killing. *J. Leukoc. Biol.*, Sep 2013; 94: 481 - 492.
- Weijing Cai, Maya Ramdas, Li Zhu, Xue Chen, Gary E. Striker, and Helen Vlassara. Oral advanced glycation endproducts (AGEs) promote insulin resistance and diabetes by depleting the antioxidant defenses AGE receptor-1 and sirtuin 1. *PNAS*, Sep 2012; 109: 15888 - 15893.
- Huang Han-Chang, Xu Ke, Jiang Zhao-Feng, Curcumin-Mediated Neuroprotection Against Amyloid-β-Induced Mitochondrial Dysfunction Involves the Inhibition of GSK-3β. DOI: 10.3233/JAD-2012-120688. *Journal of Alzheimer's Disease*, vol. 32, no. 4, pp. 981-996, 2012.
- Yue Qiu, Claus Tittiger, Claude Wicker-Thomas, Gaëlle Le Goff, Sharon Young, Eric Wajnborg, Thierry Fricaux, Nathalie Taquet, Gary J. Blomquist, and René Feyereisen. An insect-specific P450 oxidative decarboxylase for cuticular hydrocarbon biosynthesis. *PNAS*, Sep 2012; 109: 14858 - 14863.
- Jaime Uribarri, Weijing Cai, Maya Ramdas, Susan Goodman, Renata Pyzik, Xue Chen, Li Zhu, Gary E. Striker, and Helen Vlassara. Restriction of Advanced Glycation End Products Improves Insulin Resistance in Human Type 2 Diabetes: Potential role of AGER1 and SIRT1. *Diabetes Care* 2011; 34: 1610 - 1616.

Warning: This kit shall be only sold to our authorized distributors and end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of the kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at info@aatbio.com if you have any questions.