

Amplite® Colorimetric NADH and NADPH Assay Kit

Catalog number: 15272
Unit size: 400 Tests

Component	Storage	Amount (Cat No. 15272)
Component A: NADH/NADPH Probe	Freeze (< -15 °C), Minimize light exposure	1 bottle (4 mL)
Component B: Assay Buffer	Freeze (< -15 °C)	1 bottle (16 mL)
Component C: NADPH Standard	Freeze (< -15 °C), Minimize light exposure	1 vial (167 µg, lyophilized)
Component D: Lysis Buffer	Freeze (< -15 °C), Minimize light exposure	1 bottle (10 mL)

OVERVIEW

Nicotinamide adenine dinucleotide (NAD⁺) and nicotinamide adenine dinucleotide phosphate (NADP⁺) are two important cofactors found in cells. NADH is the reduced form of NAD⁺. NAD forms NADP with the addition of a phosphate group to the 2' position of the adenyl nucleotide through an ester linkage. The traditional NAD/NADH and NADP/NADPH assays are based on monitoring the changes in NADH or NADPH absorption at 340 nm. The short UV wavelength of NAD/NADH and NADP/NADPH assays makes the traditional methods suffer low sensitivity and high interference. Due to the weak absorption of NAD and NADH, the UV absorption method requires large sample sizes, making NAD and NADH measurements impractical for limited sample size. AAT Bioquest's Amplite® Colorimetric NADH and NADPH Assay Kit provides a convenient method for the detection of NADPH. The NADPH probe is a chromogenic sensor that has its maximum absorbance at ~460 nm upon NADH reduction. The absorbance increase at ~460 nm is directly proportional to the concentration of NADPH in the solution. The NADPH probe can recognize NADPH in an enzyme-free reaction, and the signal can be easily read by an absorbance microplate reader at ~ 460 nm. The Amplite® Colorimetric NADH and NADPH Assay Kit provides a sensitive assay to detect as little as 3 µM NADPH in a 100 µL assay volume. The assay can be performed in a convenient 96-well or 384-well microtiter-plate format.

AT A GLANCE

Protocol Summary

1. Prepare NADPH working solution (50 µL)
2. Add NADPH standards or test samples (50 µL)
3. Incubate at RT for 15 minutes to 2 hours
4. Monitor Absorbance at 460 nm

Important

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

KEY PARAMETERS

Absorbance microplate reader

Absorbance 460 nm
Recommended plate Clear bottom

CELL PREPARATION

For guidelines on cell sample preparation, please visit:

<https://www.aatbio.com/resources/guides/cell-sample-preparation.html>

PREPARATION OF STOCK SOLUTIONS

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated

freeze-thaw cycles

NADPH standard solution (1 mM)

1. Add 200 µL of PBS buffer to the vial of NADPH standard (Component C) to make a 1 mM (1 nmol/µL) NADPH stock solution.

PREPARATION OF STANDARD SOLUTIONS

For convenience, use the Serial Dilution Planner:
<https://www.aatbio.com/tools/serial-dilution/15272>

NADPH standard

Add 100 µL of NADPH standard solution into 400 µL PBS buffer (pH 7.4) to generate 200 µM (200 pmol/µL) NADPH standard solution (NS7). Take 200 µL of 200 µM NADPH standard solution to perform 1:2 serial dilutions to get remaining serial dilutions of NADPH standard (NS6 - NS1). Note: Diluted NADPH standard solution is unstable, and should be used within 4 hours.

PREPARATION OF WORKING SOLUTION

NADH/NADPH Probe Working Solution

1. Add 1 mL of NADH/NADPH Probe (Component A) to 4 mL of NADPH Assay Buffer (Component B) and mix well.

Note: 5 mL of NADPH working solution is enough for one 96-well plate. The working solution is not stable. It should be used promptly and avoid direct exposure to light.

SAMPLE EXPERIMENTAL PROTOCOL

Table 1. Layout of NADPH standards and test samples in a white/clear bottom 96-well microplate. NS = NADPH standard (NS1 - NS7, 3.13 to 200 µM); BL = blank control; TS = test sample.

BL	BL	TS	TS
NS1	NS1
NS2	NS2
NS3	NS3		
NS4	NS4		
NS5	NS5		
NS6	NS6		
NS7	NS7		

Table 2. Reagent composition for each well

Well	Volume	Reagent
NS1 - NS7	50 μ L	serial dilution (3.13 to 200 μ M)
BL	50 μ L	PBS
TS	50 μ L	sample

chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.

1. Prepare NADPH standards (NS), blank controls (BL), and test samples (TS) into a white wall clear bottom 96-well microplate according to Table 1 and Table 2. For a 384-well plate, add 25 μ L of reagent per well instead of 50 μ L.

Note: Prepare cells or tissue samples as desired. Lysis Buffer (Component D) can be used for lysing the cells for convenience.

2. Add 50 μ L of the NADH/NADPH probe working solution into each well of NADPH standard, blank control, and test samples to make the total NADPH assay volume of 100 μ L/well. For a 384-well plate, add 25 μ L of the NADH/NADPH probe working solution into each well instead, for a total volume of 50 μ L/well.
3. Incubate the reaction at room temperature for 15 minutes to 2 hours; protect from light.
4. Monitor the absorbance increase with an absorbance plate reader at 460 nm.

EXAMPLE DATA ANALYSIS AND FIGURES

The reading (Absorbance) obtained from the blank standard well is used as a negative control. Subtract this value from the other standards' readings to obtain the base-line corrected values. Then, plot the standards' readings to obtain a standard curve and equation. This equation can be used to calculate NADPH samples. We recommend using the Online Linear Regression Calculator which can be found at:

<https://www.aatbio.com/tools/linear-logarithmic-semi-log-regression-online-calculator>

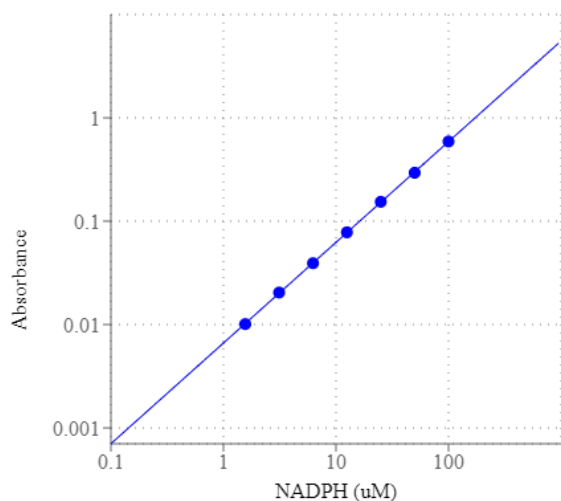


Figure 1. NADPH dose response was measured with Amplite® Colorimetric NADH and NADPH Assay Kit in a 96-well white/clear bottom plate using a SpectraMax microplate reader (Molecular devices).

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous