

NADP/NADPH Quantification Colorimetric Kit

(Catalog #K347-100; 100 assays; Store kit at -20°C)

I. Introduction:

Assays of nicotinamide nucleotides are of continual interest in the studies of energy transforming and redox state of cells or tissue. The **NADP/NADPH Quantification Kit** provides a convenient tool for sensitive detection of the intracellular nucleotides: NADP, NADPH and their ratio. The enzymes in the system specifically recognize NADP/NADPH in an enzyme cycling reaction (It does not recognize NAD⁺/NADH). There is no need to purify NADP/NADPH from sample mix. The enzyme cycling reaction significantly increases detection sensitivity. Results can be quantified using plate reader at OD_{450 nm}.

II. Kit Contents:

Components	K347-100	Cap Code	Part No.
NADP/NADPH Extraction Buffer	50 ml	NM	K347-100-1
NADP Cycling Buffer	15 ml	WM	K347-100-2
NADP Cycling Enzyme Mix	0.2 ml	Green	K347-100-3
NADPH Developer	1 vial	Purple	K347-100-4
Stop Solution	1.2 ml	Red	K347-100-5
NADPH Standard (MW: 833.36)	166.7 µg	Yellow	K347-100-6

III. NADP/NADPH Assay Protocol:

A. Reagent Reconstitution and General Consideration:

- Reconstitute NADPH developer with 1.2 ml of ddH₂O. Pipet up and down several times to completely dissolve the pellet into solution. Aliquot enough NADP Cycling Enzyme mix (2 µl per assay) for the number of assays to be performed in each experiment and aliquot and freeze the stock solution immediately at -20°C for future use. The reconstituted enzymes are stable for up to 2 months at -20°C.
- Reconstitute NADPH standard with 200 µl pure DMSO to generate 1 nmol/µl NADPH standard stock solution.
- Ensure that the NADP Cycling Buffer is at room temperature before use. The optimal temperature is 22°C. Keep other enzymes on ice during the assay and protect from light as much as possible.

B. Sample Preparation:

- For cell samples*, wash cells with cold PBS. Pellet 4 x 10⁶ cells for each assay in a microcentrifuge tube (2000 rpm for 5 min). Lyse the cells with 800 µl of NADP/NADPH Extraction Buffer in a microfuge tube and keep on ice for 10 min. Spin down at 10,000 x g for 10 min, and collect the supernatant. Transfer the extracted NADP/NADPH solution into a new labeled tube.
- For tissue samples*, weight ~50 mg tissue for each assay, wash with cold PBS, homogenize with 500 µl of NADP/NADPH Extraction Buffer in a microcentrifuge tube. Keep on ice for 10 min. Spin the sample at 10000 x g for 10 min. Transfer the extracted NADP/NADPH solution into a new labeled tube.

***Note:** Cell or tissue lysates may contain enzymes that consume NADPH rapidly. We suggest removing these enzymes from the sample either by filtering the samples through 10 kDa molecular weight cut off filters (BioVision Cat # 1997-25) or deproteinizing the sample using Deproteinizing Sample preparation Kit (BioVision Cat# K808) before performing the assays.

C. NADP/NADPH Assay Protocol:

- Standard Curve:** Dilute 10 µl of the 1 nmol/µl NADPH Standard with 990 µl NADP/NADPH Extraction Buffer to generate 10 pmol/µl Standard NADPH (**Note:** Diluted NADPH solution is unstable, must be used within 4 hr). Add 0, 2, 4, 6, 8, 10 µl of the diluted NADPH Standard into labeled 96-well plate in duplicate to generate 0, 20, 40, 60, 80, 100 pmol/well Standard. Make the final volume to 50 µl with NADP/NADPH Extraction Buffer.

Samples: To detect total NADP/NADPH (NADPt), transfer 50 µl of extracted samples into labeled 96-well plate in duplicates. **Note:** Several sample dilutions should be performed to ensure the reading can be within the Standard Curve range.

Decompose of NADP from extraction: To detect NADPH only, aliquot 200 µl samples into eppendorf tubes. Heat samples to 60°C for 30 min in a water bath or a heating block. Under the conditions, all NADP will be decomposed while NADPH will still be intact. Cool samples on ice. Quick spin samples if precipitates occur.

Transfer 50 µl of NADPH samples into labeled 96-well plate in duplicates (**Note:** several sample dilutions should be performed to ensure the reading can be within the standard curve range).

- Prepare a NADP Cycling Mix for each reaction:

NADP Cycling Buffer Mix: 98 µl
NADP Cycling Enzyme Mix: 2 µl

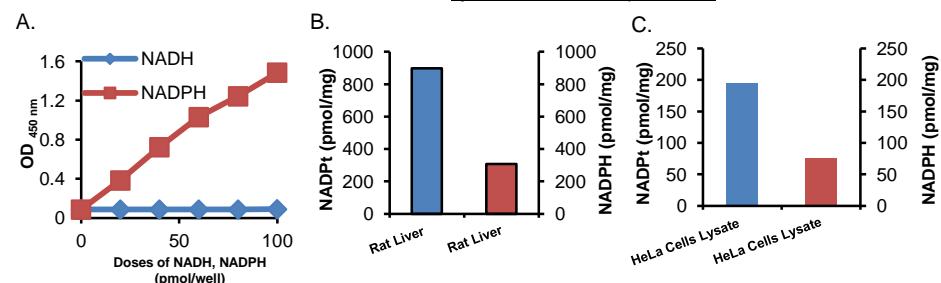
Mix well and add 100 µl of the mix into each well, mix well.

- Incubate the plate at RT for 5 min to convert NADP to NADPH.
- Add 10 µl NADPH developers into each well. Let the reaction develop for 1 to 4 hs at RT. Read the plate at OD_{450 nm}.

Note: The signal increases as the reaction time. The plate can be read multiple times while the color is in developing. The reaction can be stopped by addition of 10 µl Stop Solution each well and mix well. The color should be stable within 48 hr in a sealed plate, after the reactions are stopped.

- Calculation:** Subtract 0 Standard reading from all readings. Apply the sample OD_{450 nm} reading to standard curve. The amount of NADPt or NADPH can be expressed in pmol/10⁶ cells or ng/mg protein (NADPH molecular weight 745.4).

NADP/NADPH Ratio is calculated as: $\frac{(NADPt - NADPH)}{NADPH}$



Figures: A) NADH Standard Curve. Measurement of NADPt and NADPH in rat liver lysate (20 µg) (B) HeLa cell lysate (80 µg) (C). Assays were performed following the kit protocol.

FOR RESEARCH USE ONLY! Not to be used on humans.

Problems	Cause	Solution
Assay not working	<ul style="list-style-type: none"> • Use of ice-cold buffer • Omission of a step in the protocol • Plate read at incorrect wavelength • Use of a different 96-well plate 	<ul style="list-style-type: none"> • Buffer must be at room temperature • Refer and follow the data sheet precisely • Check the wavelength in the data sheet and the filter settings of the instrument • Fluorescence: Black plates ; Luminescence: White plates; Colorimeters: Clear plates
Samples with erratic readings	<ul style="list-style-type: none"> • Use of an incompatible sample type • Samples prepared in a different buffer • Samples were not deproteinized (if indicated in datasheet) • Cell/ tissue samples were not completely homogenized • Samples used after multiple free-thaw cycles • Presence of interfering substance in the sample • Use of old or inappropriately stored samples 	<ul style="list-style-type: none"> • Refer data sheet for details about incompatible samples • Use the buffer provided in the kit or refer data sheet for instructions • Use the 10 kDa spin cut-off filter or PCA precipitation as indicated • Use Dounce homogenizer (increase the number of strokes); observe for lysis under microscope • Aliquot and freeze samples if needed to use multiple times • Troubleshoot if needed, deproteinize samples • Use fresh samples or store at correct temperatures till use
Lower/ Higher readings in Samples and Standards	<ul style="list-style-type: none"> • Improperly thawed components • Use of expired kit or improperly stored reagents • Allowing the reagents to sit for extended times on ice • Incorrect incubation times or temperatures • Incorrect volumes used 	<ul style="list-style-type: none"> • Thaw all components completely and mix gently before use • Always check the expiry date and store the components appropriately • Always thaw and prepare fresh reaction mix before use • Refer datasheet & verify correct incubation times and temperatures • Use calibrated pipettes and aliquot correctly
Readings do not follow a linear pattern for Standard curve	<ul style="list-style-type: none"> • Use of partially thawed components • Pipetting errors in the standard • Pipetting errors in the reaction mix • Air bubbles formed in well • Standard stock is at an incorrect concentration • Calculation errors • Substituting reagents from older kits/ lots 	<ul style="list-style-type: none"> • Thaw and resuspend all components before preparing the reaction mix • Avoid pipetting small volumes • Prepare a master reaction mix whenever possible • Pipette gently against the wall of the tubes • Always refer the dilutions in the data sheet • Recheck calculations after referring the data sheet • Use fresh components from the same kit
Unanticipated results	<ul style="list-style-type: none"> • Measured at incorrect wavelength • Samples contain interfering substances • Use of incompatible sample type • Sample readings above/below the linear range 	<ul style="list-style-type: none"> • Check the equipment and the filter setting • Troubleshoot if it interferes with the kit • Refer data sheet to check if sample is compatible with the kit or optimization is needed • Concentrate/ Dilute sample so as to be in the linear range
<p>Note: The most probable list of causes is under each problem section. Causes/ Solutions may overlap with other problems.</p>		