

**EnzyFluo™ NAD<sup>+</sup>/NADH Assay Kit (EFND-100)**

**PROCEDURE FOR IN-PLATE NAD/NADH DETERMINATION IN 96-WELL PLATE**

Note: One EFND-100 kit is sufficient for 48 wells using this protocol. To measure a full 96-well plate, **two** EFND-100 kits are required.

- Cell preparation.** Plate (10 to 50 × 10<sup>3</sup>) and culture adherent cells in a clear-bottom black 96-well cell culture plate. Cells should be incubated for at least 6 hrs to ensure that cells are not labile to washes. Quantification of NAD and NADH requires cells from two separate wells. Replicates are recommended to account for cell growth variation. Eight wells should be left empty for the standard curves.
- Standard Curves.** Prepare 1 mL of 10 μM NAD Premix by mixing 10 μL 1 mM Standard and 990 μL distilled water. Dilute standard as follows.

NAD-quantification Standard wells:

No	Premix + H <sub>2</sub> O	[NAD] (pmol/well)
1	200 μL + 0 μL	100
2	120 μL + 80 μL	60
3	60 μL + 140 μL	30
4	0 μL + 200 μL	0

NADH-quantification Standard wells:

No	Premix + H <sub>2</sub> O	[Equivalent NADH] (pmol/well)
1	20 μL + 180 μL	10
2	12 μL + 188 μL	6
3	6 μL + 194 μL	3
4	0 μL + 200 μL	0

- Sample preparation.** Remove media and gently wash cells with room temperature PBS. Remove residual PBS.

Transfer 10 μL of each standard to the designated standard wells.

Add 50 μL of NAD extraction buffer to sample wells designated for NAD quantification plus both the NAD and NADH quantification standard wells. Add 50 μL of NADH extraction buffer to sample wells designated for NADH quantification. Heat plate at 37°C for 5 minutes. Add 50 μL of the opposite extraction buffer to neutralize the solution. Incubate plate at room temperature away from light for 10 minutes to equilibrate the temperature of the wells.

- Reagent Preparation.** Prepare sufficient Working Reagent by mixing for each reaction well, 14 μL Assay Buffer, 1 μL Enzyme A, 1 μL Enzyme B, 30 μL Lactate and 8 μL Probe. Fresh reconstitution is recommended.
- Reaction.** Add 50 μL Working Reagent per well quickly. Tap plate immediately to mix.
- Read fluorescence** at λ<sub>ex/em</sub> = 530/585 nm for time “zero” (F<sub>0</sub>). For NAD quantification, read F<sub>10</sub> after a 10-min incubation at room temperature. For NADH quantification, read F<sub>60</sub> after a 60-min incubation at room temperature. Protect plate from light during this incubation.

**CALCULATION**

For NAD quantification, calculate the ΔF's for standard and sample by subtracting F<sub>0</sub> from F<sub>10</sub>. Plot the ΔF's for NAD-quantification standards to determine the slope of the NAD standard curve. For NADH quantification, calculate the ΔF's for standard and sample by subtracting F<sub>0</sub> from F<sub>60</sub>. Plot the ΔF's for NADH-quantification standards to determine the slope of the NADH standard curve.

The NAD and NADH concentrations of the samples are computed as follows:

$$[\text{NAD(H)}] = \frac{\Delta F_{\text{SAMPLE}}}{\text{Slope (pmol/well)}^{-1}} \quad (\text{pmol/well})$$

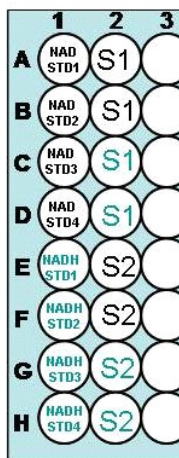
where ΔF<sub>SAMPLE</sub> is the change in fluorescence intensity values of the Sample (10 min – 0 min for NAD or 60 min – 0 min for NADH). Slope is the slope of the appropriate standard curve.

Note: If the sample ΔF values are higher than the ΔF value for the 100 pmol/well for NAD or 10 pmol/well NADH, reduce the cell count per well and repeat this assay. The detection limit of this assay is 0.05 pmol/well for either NAD or NADH.

**MATERIALS REQUIRED, BUT NOT PROVIDED**

Pipetting (multi-channel) devices, cell culture incubator, clear-bottom black 96-well cell culture plate and fluorescent plate reader capable of reading at λ<sub>ex/em</sub> = 530/585 nm.

**Example of Standards and Samples Layout in 96-Well Plate**



Add NAD Extraction Buffer to Wells for NAD/STD1, NAD/STD2, NAD/STD3, NAD/STD4, NADH/STD1, NADH/STD2, NADH/STD3, NADH/STD4, S1, S1, S2, S2,

Add NADH Extraction Buffer to Wells for S1, S1, S2, S2

