

Amplite™ Fluorimetric NADH Assay Kit

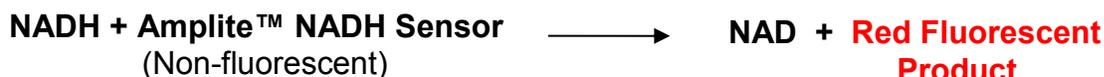
Red Fluorescence

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 15261 (400 assays)	Keep in freezer Avoid exposure to light	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⁺, the oxidized form of NADH. 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 done by monitoring the changes in NADH or NADPH absorption at 340 nm. The short UV wavelength of the traditional NAD/NADH and NADP/NADPH assays makes these methods to suffer low sensitivity and high interference. Due to the weak absorption of NAD and NADH, the UV absorption method requires large sample sizes, making the same NAD and NADH measurement unpractical if the availability of samples is limited.



This Amplite™ Fluorimetric NADH Assay Kit provides a convenient method for the detection of NADH. The enzymes in the system specifically recognize NADH in an enzyme recycling reaction. In addition, this assay has very low background since it is run in the red visible range that significantly reduces the interference resulted from biological samples. The Amplite™ Fluorimetric NADH Assay Kit provides a sensitive, one-step assay to detect as little as 100 pico-moles of NADH in a 100 μL assay volume (1 μM). 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 = 540/590 nm or an absorbance microplate reader at ~576 nm.

Kit Key Features

Broad Application:	Can be used for quantifying NADH in solutions and in cell extracts.
Sensitive:	Detect as low as 1 μM of NADH in solution.
Continuous:	Easily adapted to automation without a separation step.
Convenient:	Formulated to have minimal hands-on time. No wash is required.
Non-Radioactive:	No special requirements for waste treatment.

Kit Components

Components	Amount
Component A: NADH Recycling Enzyme Mixture	2 bottles (lyophilized powder)
Component B: NADH Assay Buffer	1 bottle (20 mL)
Component C: NADH Standard	1 vial (142 μg)

4.3 Monitor the fluorescence increase with a fluorescence plate reader at Ex/Em = 530 - 570/590 - 600 nm (optimal at Ex/Em = 540/590 nm).

Note1: 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.

*Note2: For cell based NADH measurements, ReadiUse™ mammalian cell lysis buffer *5X* (cat#20012) is recommended to use for lysing the cells.*

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. A NADH standard curve is shown in Figure 1.

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.

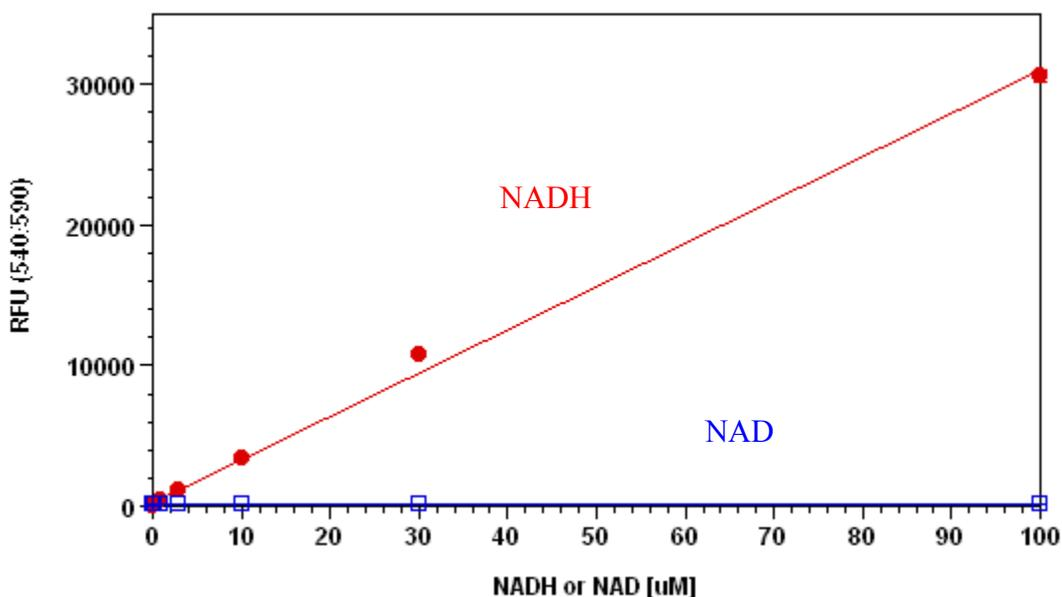


Figure 1. NADH dose response was measured with Amplite™ NADH Assay Kit in a 96-well black plate using a NOVStar microplate reader (BMG Labtech). As low as 1 µM (100 pmols/well) of NADH can be detected with 1 hour incubation (n=3) while there is no response from NAD.

References

1. Ziegenhorn J, Senn M, Bucher T. (1976) Molar absorptivities of beta-NADH and beta-NADPH. Clin Chem, 22, 151.
2. Ikegami T, Kameyama E, Yamamoto SY, Minami Y, Yubisui T. (2007) Structure and Properties of the Recombinant NADH-Cytochrome b(5) Reductase of Physarum polycephalum. BioSci Biotechnol Biochem.
3. Kimura N, Fukuwatari T, Sasaki R, Shibata K. (2006) Comparison of metabolic fates of nicotinamide, NAD⁺ and NADH administered orally and intraperitoneally; characterization of oral NADH. J Nutr Sci Vitaminol (Tokyo), 52, 142.
4. O'Donnell JM, Kudej RK, LaNoue KF, Vatner SF, Lewandowski ED. (2004) Limited transfer of cytosolic NADH into mitochondria at high cardiac workload. Am J Physiol Heart Circ Physiol, 286, H2237.

Warning: This kit is only sold to 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.