

Amplite™ Luminometric Nitroreductase Assay Kit

 Catalog number: 12470
 Unit size: 100 Tests

Component	Storage	Amount
Component A: NTR Substrate	Desiccated, Freeze (< -15 °C), Minimize light exposure	1 vial
Component B: NTR Reaction Buffer	Freeze (< -15 °C)	25 mL
Component C: NTR Reaction Enzyme	Freeze (< -15 °C)	1 vial
Component D: NTR Detection Buffer	Freeze (< -15 °C)	5 mL
Component E: NTR Detection Mix 1	Freeze (< -15 °C)	1 vial
Component F: NTR Detection Mix 2	Freeze (< -15 °C)	1 vial
Component G: Nitroreductase Standard	Freeze (< -15 °C)	1 vial
Component H: DMSO	Freeze (< -15 °C)	100 μ L

OVERVIEW

Nitroreductases (NTR) are a family of evolutionarily related proteins involved in the reduction of nitrogen-containing compounds. They are absent in mammalian cells but widespread in bacteria. Nitroreductases play a crucial role in the reduction of nitroaromatic compounds via NAD(P)H-dependent reactions. NTR are also targets for developing novel antibiotics, degradation of pollutants, and cancer therapy. Although the quantification of NTR becomes extremely important for a number of biological applications, most of the current NTR assays are based on ELISA format with low sensitivity and low throughput. Amplite™ Luminometric Nitroreductase Assay Kit provides a highly sensitive and convenient method to quantify NTR activity. The kit uses a luciferin derivative that can be selectively reduced by NTR luciferin. The amount of generated luciferin is conveniently detected with the well-known luciferase detection system. The luminescent intensity generated is proportional to the NTR activity. This luminometric nitroreductase assay kit has been formulated to have minimal hands-on time and stable luminescence signal. It can detect as low as 20 ng/mL of NTR.

AT A GLANCE

Protocol summary

1. Prepare NTR Substrate stock solution
2. Add NTR standards or NTR test samples (50 μ L/well)
3. Add NTR Reaction working solution (50 μ L/well)
4. Incubate for 60 minutes at 37 °C
5. Prepare NTR Detection working solution
6. Add NTR Detection working solution (50 μ L/well)
7. Monitor luminescence intensity increase immediately

Important

Thaw all the kit components at room temperature before starting the experiment.

KEY PARAMETERS

Luminescence microplate reader

Recommended plate Solid white

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.

1. NTR substrate stock solution (100X)

Add 50 μ L of DMSO (Component H) into the vial of NTR Substrate (Component A) to make 100X NTR substrate stock solution.

Note Keep away from light.

2. NTR Reaction Enzyme stock solution (100X)

Add 50 μ L of NTR Reaction Buffer (Component B) into the vial of NTR Reaction Enzyme (Component C) to make 100X NTR Reaction Enzyme stock solution.

Note Keep away from light.

3. NTR Detection Mix 1 stock solution (100X)

Add 50 μ L of NTR Detection Buffer (Component D) into the vial of NTR Detection Mix 1 (Component E) to make 100X NTR Detection Mix 1 stock solution.

Note Keep away from light.

4. Nitroreductase Standard solution (250 μ g/mL)

Add 20 μ L of ddH₂O into the vial of Nitroreductase Standard (Component G) to make 250 μ g/mL Nitroreductase Standard solution.

PREPARATION OF STANDARD SOLUTION

For convenience, use the Serial Dilution Planner:

<https://www.aatbio.com/tools/serial-dilution/12470>

Nitroreductase Standard solution

Add 10 μ L of 250 μ g/mL Nitroreductase Standard solution to 240 μ L of NTR Reaction Buffer (Component B) to generate 10 μ g/mL nitroreductase standard solution (NS7). Then take 10 μ g/mL nitroreductase standard solution (NS7) and perform 1:3 serial dilutions in NTR Reaction Buffer (Component B) to get serially diluted Nitroreductase standards (NS2 - NS7). Note: Diluted NTR standard solution is unstable and should be used within 4 hours.

PREPARATION OF WORKING SOLUTION

1. NTR Reaction working solution

Add 50 μ L of 100X NTR substrate stock and 50 μ L of 100X NTR reaction enzyme stock solutions into 5 mL of NTR Reaction Buffer (Component B) to make a total volume of 5.1 mL NTR Reaction working solution.

Note Keep away from light.

2. NTR Detection working solution

Add 50 μ L of 100X Detection Mix 1 and 10 μ L of Detection Mix 2 (Component F) into 5 mL of NTR Detection Buffer (Component D) to make a total volume of 5.05 mL NTR Detection working solution.

Note Keep away from light.

SAMPLE EXPERIMENTAL PROTOCOL

Table 1. Layout of NTR standards and test samples in a solid white 96-well microplate. NS=NTR standards (NS7-NS1, 0.001 to 10 μ g/mL); BL=Blank Control; TS=Test Samples

BL	BL	TS	TS
NS1	NS1
NS2	NS2
NS3	NS3		
NS4	NS4		

NS5	NS5		
NS6	NS6		
NS7	NS7		

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.

Table 2. Reagent composition for each well

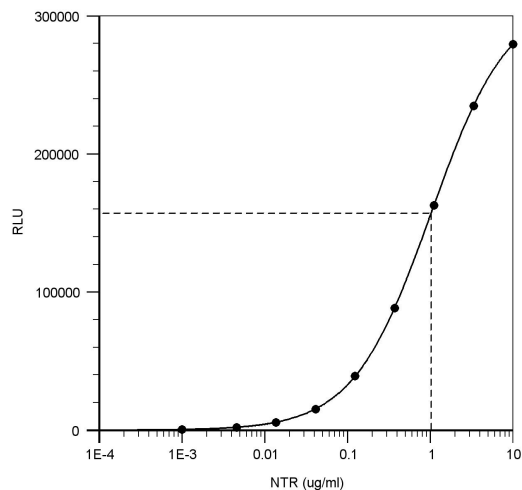
Well	Volume	Reagent
NS1-NS7	50 μ L	Serial Dilutions (10 to 0.001 μ g/mL)
BL	50 μ L	NTR Reaction Buffer (Component B)
TS	50 μ L	Test Sample

1. Prepare NTR standards (NS), blank controls (BL), and test samples (TS) according to the layout provided in Tables 1 and 2. For a 384-well plate, use 12.5 μ L of reagent per well instead of 50 μ L.

Note Treat cells or tissue samples as desired.

2. Add 50 μ L of NTR Reaction working solution to each well of NTR standard, blank control, and test samples. For a 384-well plate, add 12.5 μ L of NTR Reaction working solution into each well instead.
3. Incubate the reaction for 60 minutes at 37 °C, protected from light.
4. Add 50 μ L of NTR Detection working solution to each well to make the total assay volume 150 μ L/well. For a 384-well plate, add 12.5 μ L of Detection working solution into each well instead, for a total volume of 37.5 μ L/well.
5. Monitor the luminescence increase immediately with a luminescence microplate reader.

EXAMPLE DATA ANALYSIS AND FIGURES



Nitroreductase dose response was measured in a white solid 96-well plate with Amplitude™ Luminometric Nitroreductase Assay Kit using with a NOVO star plate reader (BMG Labtech). The kit can detect as low as 20 ng/mL of NTR (The data was obtained about 5 min after adding the NTR detection working solution).

Image generated with Quest™ Graph, ©2019 AAT Bioquest

Figure 1. Nitroreductase dose response was measured in a white solid 96-well plate with Amplitude™ Luminometric Nitroreductase Assay Kit using with a NOVO star plate reader (BMG Labtech). The kit can detect as low as 20 ng/mL of NTR (The data was obtained about 5 min after adding the NTR detection working solution).

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or