

# Portelite™ Fluorimetric Total Nucleic Acid Quantitation Kit \*Optimized for CytoCite™ and Qubit™ Fluorometers\*

 Catalog number: 17631  
 Unit size: 200 Tests

| Component                             | Storage                                    | Amount                     |
|---------------------------------------|--------------------------------------------|----------------------------|
| Component A: Helixyte™ Green All      | Freeze (< -15 °C), Minimize light exposure | 1 vial (0.25 mL-200X DMSO) |
| Component B: Assay Buffer             | Refrigerated (2-8 °C)                      | 1 bottle (100 mL)          |
| Component C: Nucleic Acid Standard #1 | Refrigerated (2-8 °C)                      | 1 vial (1 mL, 0 ng/μL)     |
| Component D: Nucleic Acid Standard #2 | Refrigerated (2-8 °C)                      | 1 vial (1 mL, 10 ng/μL)    |

## OVERVIEW

Portelite™ Fluorimetric Total Nucleic Acid Quantitation Kit is designed to rapidly measure the total amounts of nucleic acids, including double-stranded DNA (dsDNA), single-stranded DNA (ssDNA) and RNA in a sample. The kit has all the essential reagents, including Helixyte™ Green ssDNA reagent, dilution buffer, and pre-diluted DNA standards. Helixyte™ Green All reagent is a sensitive fluorescent nucleic acid probe for measuring the total amounts of nucleic acids in a sample that may contain double-stranded DNA (dsDNA), single-stranded DNA (ssDNA), RNA and long oligonucleotides. Helixyte™ Green All reagent indiscriminately binds to dsDNA, ssDNA and RNA. Portelite™ Fluorimetric Total Nucleic Acid Quantitation Kit is optimized for measuring the total amounts of nucleic acids with CytoCite™ or Qubit® fluorometers.

## AT A GLANCE

### Protocol Summary

1. Prepare a Helixyte™ Green All working solution
2. Add 190 μL of 1X Helixyte™ Green All working solution into each 0.2 mL PCR tube
3. Add 10 μL of Nucleic Acid Standards or test samples into each tube
4. Incubate at room temperature for 2 minutes
5. Monitor the fluorescence intensity with CytoCite™ fluorometer or Qubit™ fluorometer

### Important

All kit components must be brought to room temperature before starting the experiment.

## KEY PARAMETERS

### Qubit Fluorometer

|                             |                 |
|-----------------------------|-----------------|
| Excitation                  | 480 nm          |
| Emission                    | 530 nm          |
| Instrument specification(s) | 0.2 mL PCR tube |

### CytoCite Fluorometer

|                             |                 |
|-----------------------------|-----------------|
| Excitation                  | 480 nm          |
| Emission                    | 530 nm          |
| Instrument specification(s) | 0.2 mL PCR tube |

## PREPARATION OF WORKING SOLUTION

### Helixyte™ Green All working solution

Make a 200-fold dilution of Helixyte™ Green All reagent (Component A) with Assay Buffer (Component B). For example, to prepare enough working solution for 5 samples, add 5 μL of Helixyte™ Green All (Component A) into 1 mL of Assay Buffer (Component B).

**Note** Protect the working solution from light by covering it with foil or placing it in the dark. It's recommended to prepare the solution in a plastic container rather

than a glass container, as the dye may adsorb to the glass surface. For best results, this solution should be used within a few hours after the dilution.

## SAMPLE EXPERIMENTAL PROTOCOL

The acceptable range for the sample volume could be 1~20 μL depending on the estimated concentration of the Nucleic Acid sample.

The following protocol is generated based on a sample volume of 10 μL.

1. Add 190 μL of 1X Helixyte™ Green All working solution into each CytoCite™ sample tube (#CCT100) or the equivalent 0.2 mL PCR tube.

**Note** Use thin-wall, polypropylene, clear 0.2 mL PCR tubes such as AAT Cat#CCT100.

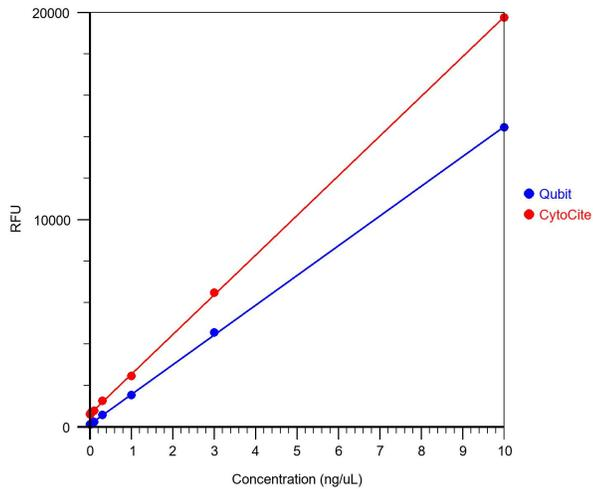
2. Add 10 μL of Nucleic Acid Standards or test samples into each tube, and then mix by vortexing for 2~3 seconds.
3. Incubate all tubes at room temperature for 2 minutes.
4. Insert the samples into CytoCite™ or Qubit™ and monitor the fluorescence intensity with the green fluorescence channel. Follow the appropriate procedures for CytoCite™ Fluorometer. See the link below for detailed instructions: <https://devices.aatbio.com/documentation/user-manual-for-cytocite-fluorometer>

### Preparation of Standard Calibration Curve

For Portelite™ assays, you have the choice to make a calibration curve with the Nucleic Acid Standards. Here is a brief protocol to generate a customized DNA standard curve.

1. Perform a 1:2 serial dilution: Add 10 ng/μL Nucleic Acid Standard #2 (Component D) into the Assay Buffer (Component B) to get 10, 5, 2.5, 1.25, 0.62, 0.31, 0.15 ng/μL DNA standard dilutions.
2. Add 190 μL of the Helixyte™ Green All working solution into each tube.
3. Add 10 μL of standards into a 0.2 mL PCR tube and then mix by vortexing for 2-3 seconds.
4. Incubate the reaction at room temperature for 2 minutes.
5. Insert the samples into CytoCite™ and monitor the fluorescence intensity with the green fluorescence channel.

## EXAMPLE DATA ANALYSIS AND FIGURES



Comparison of total nucleic acid dose response using the Qubit™ fluorometer (blue) or CytoCite™ fluorometer (red).

Image generated with Quesst™ Graph, ©2021 AAT Bioquest

**Figure 1.** Comparison of total nucleic acid dose response using the Qubit™ fluorometer (blue) or CytoCite™ fluorometer (red).

**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 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](mailto:info@aatbio.com) if you have any questions.