

Portelite™ Fluorimetric RNA Quantification Kit *5-100 ng High Sensitivity*

 Catalog number: 17695
 Unit size: 100 Tests

Component	Storage	Amount (Cat No. 17695)
Component A: StrandBrite™ Red RNA HS	Freeze (< -15 °C), Minimize light exposure	1 vial (50 µL, 100X in DMSO)
Component B: RNA Assay Buffer	Refrigerated (2-8 °C)	1 bottle (25 mL)
Component C: RNA Standard #1	Freeze (< -15 °C)	1 vial (RNA: 0 ng/µL)
Component D: RNA Standard #2	Freeze (< -15 °C)	4 vials (4 × 250 µL) (RNA: 10 ng / µL)

OVERVIEW

The Portelite™ Fluorimetric RNA Quantification Assay Kit offers a highly sensitive method for quantifying RNA in the range of 5 to 100 ng. Utilizing a proprietary RNA-selective fluorescent dye, this assay provides superior accuracy and specificity over traditional absorbance-based methods by minimizing interference from DNA, proteins, and free nucleotides.

Designed for researchers working in molecular biology, transcriptomics, and RNA-based workflows, the Portelite™ Fluorimetric RNA Quantification Assay Kit streamlines RNA quantification with minimal sample input and user-friendly steps. Its ready-to-use format minimizes preparation time while ensuring consistent, reproducible results across multiple sample types. This kit is compatible with multiple downstream applications of RNA such as RT-qPCR, RNA-seq, or cDNA synthesis.

AT A GLANCE
Protocol summary

1. Add StrandBrite™ Red DNA HS reagent working solution (200 µL).
2. Add test samples (10 µL).
3. Incubate at room temperature for 5 minutes.
4. Monitor fluorescence intensity in Qubit™ fluorometer using Red filter.

Important: The following protocol is provided as an example for quantifying total RNA with StrandBrite™ Red RNA HS. Warm all the components to room temperature before opening. No data are available for the mutagenicity or toxicity of StrandBrite™ Red RNA HS stain. Because this reagent binds to nucleic acids, it should be treated as a potential mutagen and handled with appropriate care. The DMSO stock solution should be handled with particular caution as DMSO is known to facilitate the entry of organic molecules into tissues.

KEY PARAMETERS
Qubit Fluorometer

Emission	665-720 nm
Excitation	635 nm
Instrument specification(s)	0.2 mL PCR vial

PREPARATION OF WORKING SOLUTION

Add 2.5 µL StrandBrite™ Red RNA HS (Component A) into 1 mL of RNA Assay Buffer (Component B) and mix well. Protect the working solution from light by covering it with foil or placing it in the dark.

Note: 1 mL of working solution is enough for 5 tests.

Note: We recommend preparing this solution in a plastic container rather than glass, as the dye may adsorb to glass surfaces. For best

results, this solution should be used within a few hours of its preparation.

SAMPLE EXPERIMENTAL PROTOCOL

The acceptable sample volume could range from 1–20 µL depending on the estimated concentration of the RNA sample. The recommended sample volume is 10 µL with the RNA concentration in 0.5–10 ng/µL range. If other sample volume is being used, please adjust the dilution factor in the concentration calculations.

The following protocol is generated based on 10 µL sample volume with the RNA concentration in 0.5-10 ng/µL range.

1. Add 200 µL StrandBrite™ Red RNA HS working solution into each CytoCite™ sample tube (AAT cat# CCT100) or 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 RNA standard #1 and #2 or test samples into each tube, and then mix by vortexing 2-3 seconds.
3. Incubate all tubes at room temperature for 2 minutes.
4. Insert the samples into Qubit™ and monitor the fluorescence with red fluorescence channel.

Preparation of standard calibration curve (Optional)

For StrandBrite™ assays, you have the choice to make a calibration curve with the RNA standards. Here is a brief protocol to generate a customized RNA standard curve:

1. Perform dilution of RNA standard (Component D) to 10, 8, 6, 4, 2, 1, 0.5 ng/µL in RNA Assay Buffer (Component B).
2. Add 200 µL StrandBrite™ Red DNA BR working solution into each tube.
3. Add 10 µL RNA standards or test samples into each tube, and then mix by vortexing 2–3 seconds.
4. Incubate the reaction at room temperature for 2 minutes.
5. Insert the samples into Qubit™ and monitor the fluorescence with red fluorescence channel.

EXAMPLE DATA ANALYSIS AND FIGURES

We recommend using the Online Linear Regression Calculator which can be found at:

<https://www.aatbio.com/tools/linear-logarithmic-semi-log-regression-onlinecalculator/>

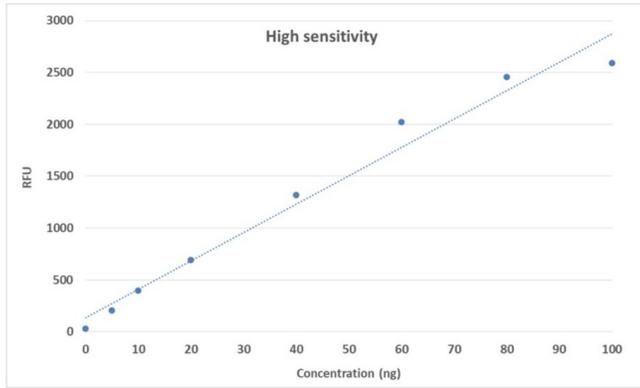


Figure 1. RNA dose response generated with StrandBrite™ Fluorimetric RNA Quantitation Kit *High Sensitivity*.

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 if you have any questions.