invitrogen USER GUIDE

Quant-iT™ RiboGreen™ RNA Reagent and Kit

Catalog Numbers R11490, R11491, T11493

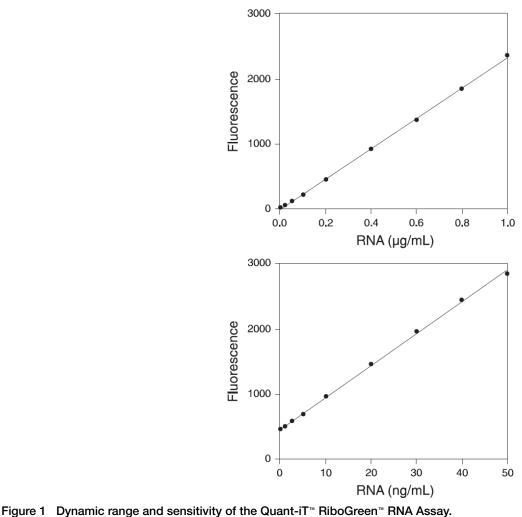
Pub. No. MAN0002073 Rev. A.0



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from **thermofisher.com/support**.

Product description

The Quant-iT[™] RiboGreen[™] RNA Reagent is an ultrasensitive fluorescent nucleic acid stain for quantitating RNA in solution. Detecting and quantitating small amounts of RNA is important in many applications including measuring yields of *in vitro* transcribed RNA and measuring RNA concentrations before performing Northern blot analysis, S1 nuclease assays, RNase protection assays, cDNA library preparation, reverse transcription PCR, and differential display PCR.



For the high-range assay (top panel), the Quant-iT RiboGreen RNA Reagent was diluted 200-fold into 10 mM Tris-HCl, 1 mM EDTA, pH 7.5 (TE) and 100 µL of the reagent solution was added to microplate wells containing 100 µL of ribosomal RNA in TE. For the low-range assay (bottom panel), the Quant-iT RiboGreen RNA Reagent was diluted 2,000-fold into TE and 100 µL of the reagent solution was added to microplate wells containing 100 µL of ribosomal RNA in TE. Samples were excited at 485 ± 10 nm and the fluorescence emission intensity was measured at 530 ± 12.5 nm using a fluorescence microplate reader. Fluorescence emission intensity was then plotted versus RNA concentration.

The Quant-iT[™] RiboGreen[™] RNA Reagent enables quantitation of as little as 1 ng/mL RNA (200 pg RNA in a 200 µL assay volume) with a fluorescence microplate reader using fluorescein excitation and emission wavelengths. The linear range of the Quant-iT[™] RiboGreen[™] RNA

Reagent extends over three orders of magnitude in RNA concentration (1 ng/mL to 1 μg/mL) using two dye concentrations (Figure 1). The high-range assay allows quantitation of 20 ng/mL to 1 μg/mL RNA, and the low-range assay allows quantitation of 1 ng/mL to 50 ng/mL RNA. This linearity is maintained in the presence of several compounds commonly found to contaminate nucleic acid preparations, including nucleotides, salts, urea, ethanol, chloroform, detergents, proteins, and agarose. Although the Quant-iT[™] RiboGreen RNA Reagent also binds to DNA, pretreatment of mixed samples with DNase can be used to generate an RNA-selective assay (see "Eliminate DNA from samples" on page 5).

Contents and storage

Component	Quant-iT [™] RiboGreen [™] RNA Reagent ^[1]	Quant-iT™ RiboGreen™ RNA Assay Kit	Concentration	Storage ^[2]
	Cat. No. R11491	Cat. No. R11490		
Quant-iT™ RiboGreen™ RNA	1 mL	1 mL	Solution in DMSO	2°C to 8°C ^[3]
Reagent (Component A)				Desiccate
				Protect from light
20X TE Buffer, RNase-free (Component B)	Not applicable	25 mL	200 mM Tris-HCl, 20 mM EDTA, pH 7.5 in DEPC- treated water	≤30°C
Ribosomal RNA standard, 16S and 23S rRNA from <i>E. coli</i> (Component C)	Not applicable	5 × 200 μL	100 μg/mL in TE buffer	2°C to 8°C ^[4] Avoid freeze-thaw cycles

Number of labelings: 2,000 to 20,000 with an assay volume of 200 μL in a 96-well microplate format. The Quant-iT™ RiboGreen™ RNA Assay can be adapted for use in cuvettes or 384-well microplates.

Approximate fluorescence excitation/emission maxima: 500/525 nm, bound to nucleic acid.

- [1] Stand-alone reagent does not include Components B and C.
- [2] When stored as directed, products are stable for at least 6 months.
- [3] For long-term storage, the Quant-iT™ RiboGreen™ RNA Reagent can be stored at ≤-20°C
- [4] For long-term storage, store the rRNA standards at ≤-20°C or -70°C.

Required materials not supplied

- Nuclease-free pipettors and tips
- Nuclease-free water
- Microplates for Fluorescence-based Assays, 96-well (Cat. No. M33089)

Prepare the assay buffer

Prepare the 1X TE working solution by diluting the concentrated buffer (Component B) 20-fold with nuclease-free water. Prepare nuclease-free water by treating distilled, deionized water with 0.1% diethyl pyrocarbonate (DEPC), incubating for several hours at 37°C, and autoclaving for at least 15 minutes at 15 lbs/sg. inch to sterilize the water and eliminate DEPC.

IMPORTANT! TE buffer (10 mM Tris-HCl, 1 mM EDTA, pH 7.5) is used to prepare the Quant-iT[™] RiboGreen[™] RNA Reagent and for diluting RNA standards and samples. Make sure the TE solution is free of contaminating nucleases and nucleic acids. The 20X TE buffer included in the Quant-iT[™] RiboGreen[™] RNA Assay Kit is nuclease-free and nucleic acid–free.

Prepare the reagent

Two different dye concentrations are required to achieve the full linear dynamic range of the Quant-iT[™] RiboGreen[™] RNA Assay. Before preparing the working solution of the Quant-iT[™] RiboGreen[™] RNA Reagent, decide whether you wish to perform the **high-range** assay (20 ng/mL to 1 μ g/mL RNA), **low-range** assay (1 ng/mL to 50 ng/mL RNA), or both.

On the day of the experiment, allow the Quant-iT[™] RiboGreen[™] RNA Reagent to warm to room temperature before opening the vial, then prepare an aqueous working solution of the Quant-iT[™] RiboGreen[™] RNA Reagent by diluting the concentrated DMSO stock solution (Component A) into TE, 200-fold for the **high-range** assay or 2,000-fold for the **low-range** assay. For microplate assays of a total 200 µL assay volume, you need 100 µL of the Quant-iT[™] RiboGreen[™] RNA Reagent working solution per sample.

For example, to prepare enough working solution to assay 100 samples in 200 μL volumes, add 50 μL Quant-iT[™] RiboGreen[™] RNA Reagent to 9.95 mL TE for the **high-range** assay or add 5 μL Quant-iT[™] RiboGreen[™] RNA Reagent to 9.995 mL TE for the **low-range** assay.

Note: Allow the Quant-iT[™] RiboGreen RNA Reagent to warm to room temperature before opening the vial. Cold DMSO solutions absorb moisture from warmer, room temperature air, resulting in loss of efficacy for the reagent. Always store the DMSO stock solution in the

presence of desiccant when not in use. We recommend preparing the working solution in sterile, disposable polypropylene plasticware rather than glassware, as the reagent may adsorb to glass surfaces. Protect the working solution from light, as the Quant-iT[™] RiboGreen RNA Reagent is susceptible to photodegradation. **For best results, use the working solution within a few hours of preparation.**

Prepare the RNA standard curve

 Prepare a 2 μg/mL solution of RNA in TE using nuclease-free plasticware. Determine the RNA concentration on the basis of absorbance at 260 nm (A₂₆₀) in a cuvette with a 1 cm pathlength; an A₂₆₀ of 0.05 corresponds to 2 μg/mL RNA.

The ribosomal RNA standard (Component C), provided at 100 μ g/mL in the Quant-iT[™] RiboGreen[™] RNA Assay Kit, is diluted 50-fold in TE to make the 2 μ g/mL working solution. For example, 4 μ L of the RNA standard mixed with 196 μ L of TE is sufficient for the standard curve described in step 2.

Note: For a standard curve, we commonly use 16S and 23S ribosomal RNA, although any purified RNA preparation may be used. It is sometimes preferable to prepare the standard curve with RNA similar to the type being assayed. However, most single-stranded RNA molecules yield approximately equivalent signals.

Note: The RNA solution used to prepare the standard curve should be treated the same way as the experimental samples and should contain similar levels of contaminants. See "Effects of common contaminants" on page 4 for a list of contaminants tested in the Quant-iT™ RiboGreen™ assay.

2. For the **high-range** standard curve, dilute the 2 μg/mL RNA solution into microplate wells as shown in Table 1. For the **low-range** standard curve, dilute the 2 μg/mL RNA solution 20-fold into TE to make a 100 ng/mL RNA stock solution, then prepare the dilution series shown in Table 2.

Table 1 Protocol for preparing a high-range standard curve.

Volume of TE buffer	Volume of 2 μg/mL RNA stock	Volume of 200-fold diluted Quant-iT [™] RiboGreen [™] RNA Reagent	Final RNA concentration in the assay
0 μL	100 µL	100 μL	1 μg/mL
50 μL	50 μL	100 μL	500 ng/mL
90 µL	10 μL	100 μL	100 ng/mL
98 µL	2 μL	100 μL	20 ng/mL
100 μL	OμL	100 μL	blank

Table 2 Protocol for preparing a low-range standard curve.

Volume of TE buffer	Volume of 100 ng/mL RNA stock	Volume of 2,000-fold diluted Quant-iT™ RiboGreen™ RNA Reagent	Final RNA concentration in the assay
0 μL	100 µL	100 µL	50 ng/mL
50 μL	50 μL	100 μL	25 ng/mL
90 μL	10 μL	100 µL	5 ng/mL
98 µL	2 μL	100 μL	1 ng/mL
100 μL	OμL	100 μL	blank

- 3. Add 100 µL of the appropriate aqueous working solution of Quant-iT[™] RiboGreen[™] RNA Reagent (prepared in "Prepare the reagent" on page 2) to each microplate well. Use the high-range working solution for performing the high-range assay, and use the low-range working solution for performing the low-range assay. Mix well and incubate for 2–5 minutes at room temperature, protected from light.
- 4. Measure the sample fluorescence using a fluorescence microplate reader and standard fluorescein wavelengths (excitation ~480 nm, emission ~520 nm).

Note: To ensure that the sample readings remain in the detection range, set the instrument's gain so that the sample containing the highest RNA concentration yields a fluorescence intensity near the microplate reader's maximum. For optimal detection sensitivity, the instrument gain can be increased for the low-range assay relative to the high-range assay. To minimize photobleaching effects, keep the time for fluorescence measurement constant for all samples.

5. Subtract the fluorescence value of the reagent blank from that of each of the samples. Use corrected data to generate a standard curve of fluorescence versus RNA concentration (see Figure 1).

Analyze samples

- 1. Dilute the experimental RNA solution in TE to a final volume of 100 µL in microplate wells.
 - Note: You can alter the amount of sample diluted, provided that the final volume remains $100 \, \mu L$. High dilutions of the experimental sample may serve to diminish the interfering effect of certain contaminants. However, extremely small sample volumes should be avoided because they are difficult to pipet accurately. In addition, the level of assay contaminants should be kept as uniform as possible throughout an experiment, to minimize sample-to-sample signal variation. For example, if a series of RNA samples contain widely differing salt concentrations, then they cannot be compared to a single standard curve. To avoid this problem, simply adjust the concentration of contaminants to be the same in all samples, if possible (see "Effects of common contaminants" on page 4).
- 2. Add 100 µL of the aqueous working solution of the Quant-iT[™] RiboGreen RNA Reagent (prepared in "Prepare the reagent" on page 2) to each sample. Incubate for 2–5 minutes at room temperature, protected from light.
- 3. Measure the fluorescence of the samples using the same instrument parameters used to generate the standard curve (see step 4). To minimize photobleaching effects, keep the time for fluorescence measurement constant for all samples.
- 4. Subtract the fluorescence value of the reagent blank from that of each of the samples. Determine the RNA concentration of the sample from the standard curve generated in "Prepare the RNA standard curve" on page 3.
- 5. The assay can be repeated using a different dilution of the sample to confirm the quantitation results.

Effects of common contaminants

The Quant-iT RiboGreen Assay remains linear in the presence of several compounds that commonly contaminate nucleic acid preparations, although the signal intensity may be affected (Table 3). For the highest accuracy, the standards should be prepared under the same conditions as the experimental samples and contain similar levels of contaminants.

Table 3 Effects of common contaminants on the signal intensity of the assay.

Compound	Maximum acceptable concentration	% Signal change ^[1]
Salts		
Ammonium acetate	20 mM	4% decrease
Sodium acetate	20 mM	11% decrease
Sodium chloride	20 mM	15% decrease
Zinc chloride	1 mM	9% decrease
Magnesium chloride	0.5 mM	9% decrease
Calcium chloride	0.1 mM	2% increase
Cesium chloride	10 mM	8% decrease
Guanidinium thiocyanate	10 mM	9% decrease
Urea	3 M	13% decrease
Organic solvents		
Phenol	0.5%	5% decrease
Ethanol	20%	10% decrease
Chloroform	2%	15% increase
Detergents		
Sodium dodecyl sulfate	0.05%	10% decrease
Triton™ X-100	0.5%	8% decrease
Proteins		
Bovine serum albumin	0.2%	11% decrease
lgG	0.02%	4% decrease

Compound	Maximum acceptable concentration	% Signal change ^[1]
Other compounds		
Formamide	10%	12% decrease
Sucrose	>500 mM	4% decrease
Boric acid	100 mM	15% decrease
Polyethylene glycol	10%	10% decrease
Agarose	0.2%	3% increase

^[1] The compounds were incubated at the indicated concentrations with the Quant-IT™ RiboGreen™ RNA Reagent in the presence of 1.0 mg/mL ribosomal RNA. All samples were assayed in a final volume of 200 µL in 96-well microplates using a fluorescence microplate reader. Samples were excited at 485 ± 10 nm and fluorescence intensity was measured at 530 + 12.5 nm

Eliminate DNA from samples

The Quant-iT[™] RiboGreen RNA Reagent also binds to DNA. Fluorescence in samples that is due to the Quant-iT RiboGreen RNA Reagent binding to DNA can be eliminated by pre-treating the sample with RNase-free DNase, ensuring that the entire sample fluorescence is due to dye bound to RNA.

- 1. Prepare 10X DNase digestion buffer: nuclease-free 200 mM Tris-HCl, pH 7.5, containing 100 mM MgCl₂ and 20 mM CaCl₂.
- 2. Add 0.11 volume of 10X DNase digestion buffer to each DNA-containing sample (for example, to a 9 μL sample, add 1 μL 10X buffer).
- 3. Add \sim 5 units of RNase-free DNase I per μg of DNA estimated to be in the sample.
- 4. Incubate the sample at 37°C for 90 minutes.
- 5. Dilute the sample at least 10-fold into TE to diminish effects of the digestion buffer salts on the Quant-iT[™] RiboGreen[™] assay procedure.
- 6. Perform the Quant-iT[™] RiboGreen[™] assay as described (see "Analyze samples" on page 4).

Related products

Table 4 Bulk Reagents and Kits

Product	Quantity	Cat. No.
Quant-iT™ PicoGreen™ dsDNA Assay Kit	1 mL assay kit	P7589
	10 x 100 μL	P11496
Quant-iT™ PicoGreen™ dsDNA Reagent	1 mL reagent	P7581
	10 x 100 μL	P11495
TE Buffer (20X), RNase-free	100 mL	T11493
Quant-iT™ RiboGreen™ RNA Assay Kit	1 mL assay kit	R11490
Quant-iT™ RiboGreen™ RNA Reagent	1 mL reagent	R11491
Quant-iT™ RediPlate™ 96 RiboGreen™ RNA Quantitation Kit	1 plate	R32700
Quant-iT™ OliGreen™ ssDNA Assay Kit	1 mL assay kit	O11492
Quant-iT™ OliGreen™ ssDNA Assay Reagent	1 mL reagent	O7582

Table 5 Microplate Reader Assays

Product	Dynamic Range	Quantity	Cat. No.
Quant-iT™ 1X dsDNA Assay Kit, High Sensitivity	200 pg-100 ng	1,000 reactions	Q33232
Quant-iT™ 1X dsDNA Assay Kit, Broad-Range	4 ng–2 μg	1,000 reactions	Q33267
Quant-iT™ DNA Assay Kit, High Sensitivity	200 pg-100 ng	1,000 reactions	Q33120
Quant-iT™ DNA Assay Kit, Broad- Range	4 ng−1 μg	1,000 reactions	Q33130
Quant-iT™ RNA Assay Kit	5-100 ng	1,000 reactions	Q33140
Quant-iT™ RNA Reagent	5-100 ng	1,000 reactions	Q32884
Quant-iT™ RNA Assay Kit, Broad Range	20 ng−1 μg	1,000 reactions	Q10213
Quant-iT™ RNA XR Assay Kit	200 ng-10 μg	1,000 reactions	Q33225
Quant-iT™ microRNA Assay Kit	1-100 ng	1,000 reactions	Q32882
Quant-iT™ Protein Assay Kit	250 ng-5 μg	1,000 reactions	Q33210
Microplates for Fluorescence-based Assays, 96-well	-	10 plates	M33089

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Life Technologies Corporation | 29851 Willow Creek Road | Eugene, Oregon 97402 USA For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0002073

Revision	Date	Description
A.0		The format and content were updated. The version numbering was reset to A.0 in conformance with internal document control.
1.00	10.line 2008	New document for the Quant-iT™ RiboGreen™ RNA Assay Kit and Quant-iT™ RiboGreen™ RNA Reagent.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific, Inc. and its subsidiaries unless otherwise specified.



15 March 2022