

Quant-iT™ dsDNA Broad-Range Assay Kit

Catalog Number Q33130

Pub. No. MAN0002341 Rev. B.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](https://www.thermofisher.com/support).

Product description

The Quant-iT™ dsDNA Broad-Range Assay Kit makes DNA quantitation easy and accurate. The assay is highly selective for double-stranded DNA over RNA, and in the range of 4–1,000 ng, the fluorescence signal is linear with DNA (see “Assay selectivity” on page 1). Common contaminants, such as salts, solvents, detergents, or protein are well tolerated in the assay (Table 1).

In addition to the Quant-iT™ dsDNA Broad-Range Assay Kit described here, we also offer the Quant-iT™ dsDNA High-Sensitivity Assay Kit (Cat. No. [Q33120](#)). The Quant-iT™ dsDNA High-Sensitivity Kit is designed for assaying samples containing 0.2–100 ng of DNA.

Contents and storage

Component	Amount	Concentration	Storage ^[1]
Quant-iT™ dsDNA BR reagent (Component A)	1.0 mL	200X in DMSO	2°C to 8°C Desiccate Protect from light
Quant-iT™ dsDNA BR buffer (Component B)	250 mL	Not applicable	≤30°C
λ dsDNA BR standards (Component C)	set of 8 (500 µL each)	0, 5, 10, 20, 40, 60, 80, and 100 ng/µL	2°C to 8°C ^[2]

Number of labelings: 1,000, with a 200 µL assay volume in a 96-well microplate format. The Quant-iT™ dsDNA BR assay can be adapted for use in cuvettes or 384-well microplates.

Approximate fluorescence excitation/emission maxima: 510/527 nm (see “Spectral data” on page 2)

^[1] When stored as directed, kit contents are stable for at least 6 months.

^[2] For long-term storage, the dsDNA standards can be stored at ≤–20°C.

Required materials not supplied

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

Critical assay parameters

Assay temperature

Quant-iT™ assays deliver optimal performance when all solutions are at room temperature; temperature fluctuations can influence the accuracy of the assay.

Assay selectivity

The Quant-iT™ dsDNA BR assay is highly selective for double-stranded DNA over RNA (Figure 1).

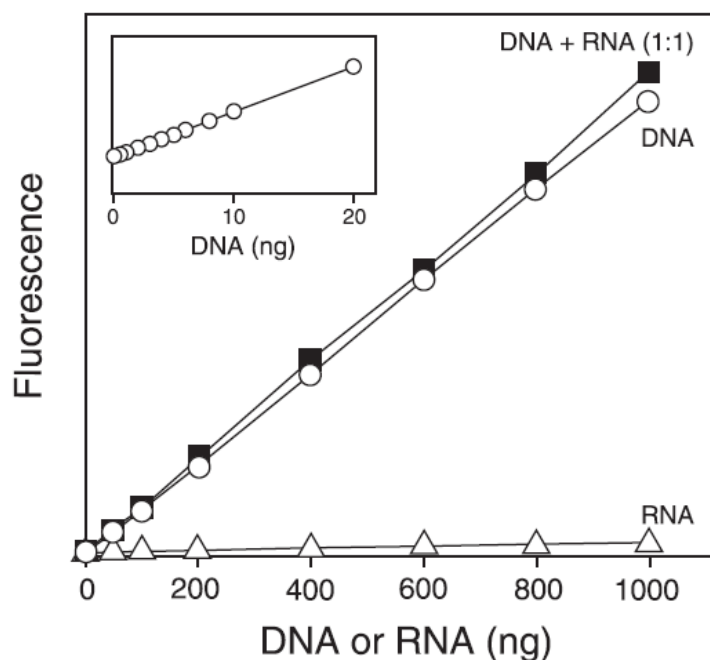


Figure 1 DNA selectivity and sensitivity of the Quant-iT™ dsDNA BR assay.

Triplicate 10 μ L samples of λ DNA (O), *E. coli* rRNA (Δ), or a 1:1 mixture of DNA and RNA (■) were assayed in the Quant-iT™ dsDNA BR assay. Fluorescence was measured at 485/530 nm and plotted versus the mass of nucleic acid for the DNA alone or RNA alone, or versus the mass of the DNA component in the 1:1 mixture. The variation (CV) of replicate DNA determinations was $\leq 3\%$. The inset, a separate experiment with octuplicate determinations, shows the sensitivity of the assay for DNA. Background fluorescence has not been subtracted.

Incubation time

To allow the Quant-iT™ dsDNA BR assay to reach maximum fluorescence, incubate for 2 minutes after mixing the sample or standard with the working solution. After this incubation period, the fluorescence signal is stable for 3 hours at room temperature when the samples and standards are protected from light.

Spectral data

The Quant-iT™ dsDNA BR reagent has excitation and emission maxima of 510/527 nm when bound to DNA.

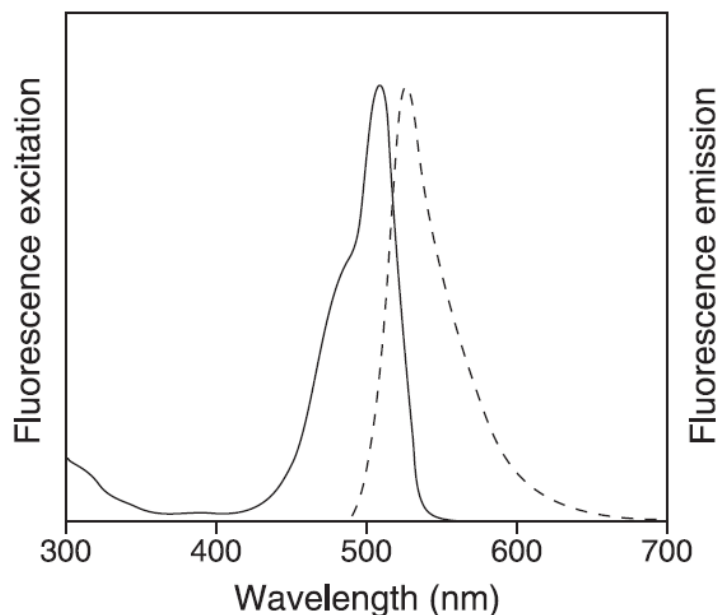


Figure 2 Excitation and emission maxima for the Quant-iT™ dsDNA BR reagent bound to DNA.

Photostability of the Quant-iT™ reagent

Avoid multiple sequential fluorescence reads to minimize reagent photobleaching.

Effects of contaminating substances

A number of common contaminants have been tested in the Quant-iT™ dsDNA BR assay, and most are well tolerated. For untested contaminating substances and in general, the standards should be assayed under the same conditions as the unknowns for highest accuracy. For example, if the experimental samples are in an unusual buffer and if 10 µL volumes of these samples are used, then add 10 µL volumes of the unusual buffer (lacking DNA) to the assays of the standards.

Table 1 Effects of contaminants in the Quant-iT™ dsDNA Broad-Range Assay.

Contaminant	Final concentration in the assay	Concentration in 20 µL sample	Concentration in 10 µL sample	Result ^[1]
Sodium chloride	10 mM	100 mM	200 mM	OK ^[2]
Magnesium chloride	2 mM	20 mM	40 mM	OK ^[2]
Sodium acetate	10 mM	100 mM	200 mM	OK
Ammonium acetate	10 mM	100 mM	200 mM	OK ^[2]
Potassium phosphate, pH 7.4	5 mM	50 mM	100 mM	OK ^[2]
Ethanol	1%	10%	20%	OK
Phenol	0.1%	1%	2%	OK
Chloroform ^[3]	0.2%	2%	4%	OK
SDS	0.01%	0.1%	0.2%	OK
Triton™ X-100	0.001%	0.01%	0.02%	OK ^[2]
dNTPs ^[4]	100 µM	1 mM	2 mM	OK
BSA	20 µg/mL	200 µg/mL	400 µg/mL	OK ^[2]
IgG	10 µg/mL	100 µg/mL	200 µg/mL	OK

^[1] Results are given as OK, usually less than 10% perturbation, or as NR (not recommended).

^[2] An acceptable result, but with some distortion of the standard curve; for best results, add the same amount of contaminant to the standard samples.

^[3] Immiscible.

^[4] A mixture of dATP, dCTP, dGTP, and dTTP.

Prepare and read samples and standards using the Quant-iT™ dsDNA Broad-Range Assay Kit with a fluorescence microplate reader

This protocol describes the use of the Quant-iT™ dsDNA Broad-Range Assay Kit with a fluorescence microplate reader equipped with excitation and emission filters appropriate for fluorescein or Alexa Fluor™ 488 dye. For an overview of this procedure, see Figure 3.

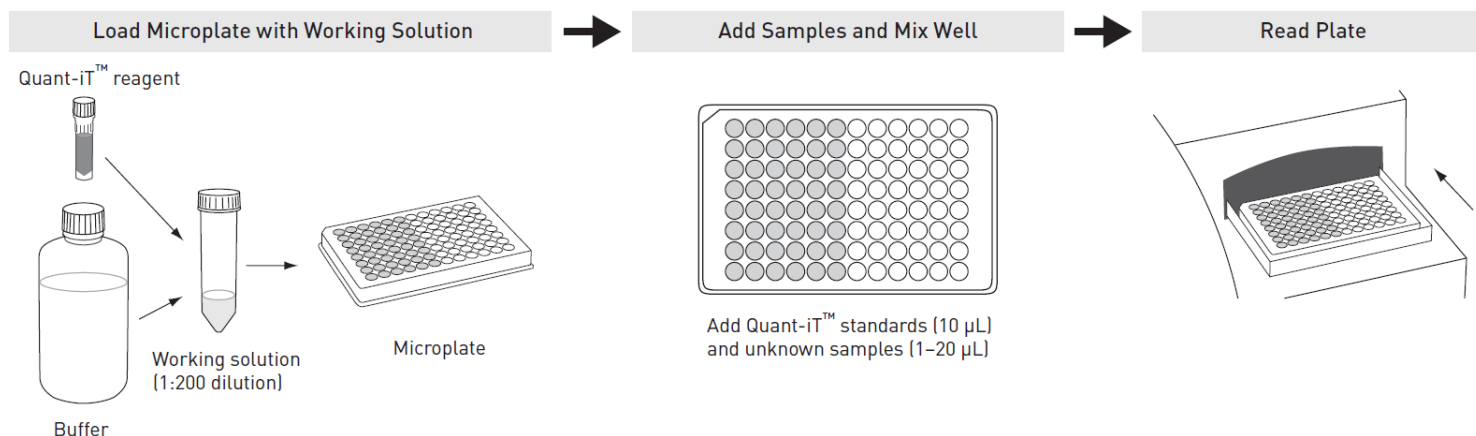


Figure 3 Overview of the Quant-iT™ dsDNA Broad-Range assay.

IMPORTANT! For best results, ensure that all materials and reagents are at room temperature.

1. Make a working solution by diluting Quant-iT™ dsDNA BR reagent 1:200 in Quant-iT™ dsDNA BR buffer. For example, for ~100 assays combine 100 µL of Quant-iT™ dsDNA BR reagent (Component A) and 20 mL of Quant-iT™ dsDNA BR buffer (Component B) in a disposable plastic container and mix well. The Quant-iT™ working solution is stable for at least 3 hours at room temperature, protected from light.

IMPORTANT! Do not use glass containers. Do not use buffers other than the Quant-iT™ dsDNA BR buffer to make the working solution.

2. Load 200 µL of the working solution into each microplate well.
3. Add 10 µL of each λ DNA standard (Component C) to separate wells and mix well. Take care not to introduce nucleases into the tubes of DNA standard as you remove aliquots for the assay. Duplicates or triplicates of the standards are recommended.
4. Add 1–20 µL of each unknown DNA sample to separate wells and mix well. Duplicates or triplicates of the unknown samples are recommended.
5. If possible, mix the 96-well plate using a plate mixer or using the plate reader for about 3–10 seconds. After mixing, allow the plate to incubate at room temperature for 2 minutes.
6. Measure the fluorescence using a microplate reader (excitation/emission maxima are 510/527 nm; see “Spectral data” on page 2). Standard fluorescein wavelengths (excitation/emission at ~480/530 nm) are appropriate for this dye. The fluorescence signal is stable for 3 hours at room temperature, when protected from light.
7. Use a standard curve to determine the DNA amounts. For the λ DNA standards, plot amount vs. fluorescence, and fit a straight line to the data points.

Data analysis considerations – standard curves and extended ranges

The fluorescence of the Quant-iT™ dsDNA BR reagent bound to dsDNA is extremely linear up to 1,000 ng. For best results at the low end of the standard curve, the line should be forced through the background point (or through zero, if background has been subtracted). When 10 µL volumes of the standards are used, the lowest DNA-containing standard represents 50 ng of DNA; nevertheless, highly accurate determinations of DNA down to 2 ng are attained using the standard curve as described above.

To assess the reliability of the assay in the low range, use smaller volumes of the standards; for example, 2 µL volumes for a standard curve ranging from 0–200 ng (Figure 4A). Alternatively, dilute the standards in buffer for an even tighter range (Figure 4A, inset).

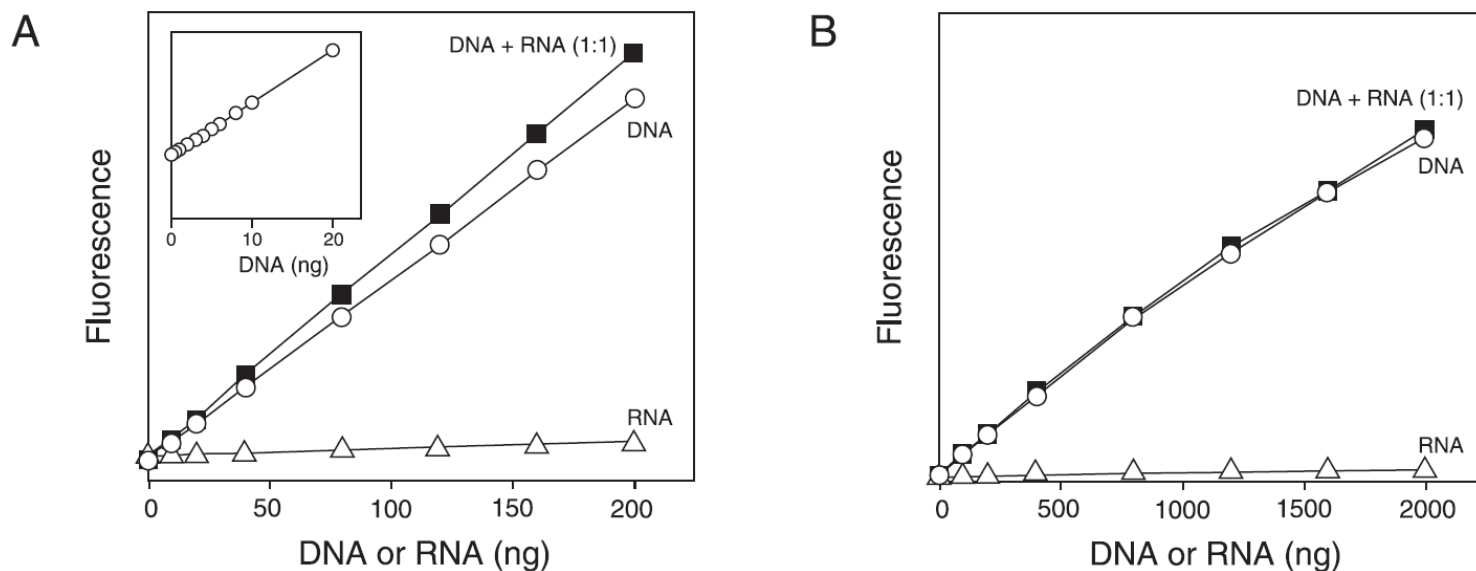


Figure 4 Extended ranges for the Quant-iT™ dsDNA BR assay.

Triplicate 2 µL (Panel A) or 20 µL samples (Panel B) of λ DNA (○), *E. coli* rRNA (Δ), or a 1:1 mixture of DNA and RNA (■) were assayed in the Quant-iT™ dsDNA BR assay. Fluorescence was measured at 485/530 nm and plotted versus the mass of nucleic acid for the DNA alone or RNA alone, or versus the mass of the DNA component in the 1:1 mixture. The inset (Panel A), a separate experiment with octuplicate determinations, shows the sensitivity of the assay for DNA. Background fluorescence has not been subtracted.

If desired, the utility of the Quant-iT™ dsDNA BR assay can be extended beyond 1,000 ng, up to 2,000 ng (Figure 4B). For standards in this range, use 20 µL volumes of the provided standards. Note that the standard curve may not be linear in the range 1,600–2,000 ng.

Related products

Table 2 Bulk Reagents and Kits

Product	Quantity	Cat. No.
Quant-iT™ PicoGreen™ dsDNA Assay Kit	1 mL assay kit 10 x 100 µL	P7589 P11496
Quant-iT™ PicoGreen™ dsDNA Reagent	1 mL reagent 10 x 100 µL	P7581 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 3 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.



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. MAN0002341

Revision	Date	Description
B.0	8 March 2022	The format and content were updated.
A.0	16 February 2015	New document for the Quant-iT™ dsDNA Broad-Range Assay Kit.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, 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.