

Materials Required

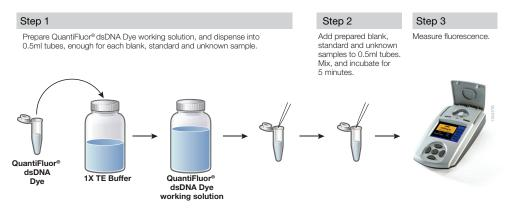
- QuantiFluor® dsDNA System (Cat.# E2670)
- Quantus[™] Fluorometer (Cat.# E6150)
- thin-walled 0.5ml PCR tubes (Cat.# E4941 or Axygen Cat.# PCR-05-C)
- nuclease-free water

Warm all assay components to room temperature before use.

Caution: We recommend the use of gloves, lab coats and eye protection when working with these or any chemical reagents.

The *Quantus*[™] *Fluorometer Operating Manual* #TM396 and *QuantiFluor*[®] *dsDNA System Technical Manual* #TM346 are available at: **www.promega.com/protocols**

Single-Tube Format Protocol



Note: If the Quantus[™] Fluorometer was previously calibrated, you may not need to calibrate it again. Therefore, do not prepare blank and standard samples, and skip Steps 2, 3 and 7.

- 1. Prepare 1X TE Buffer: Dilute the 20X TE Buffer 20-fold with nuclease-free water.
- 2. Prepare Working Solution: Dilute the QuantiFluor® dsDNA Dye 1:400 in 1X TE buffer, and mix.
- 3. **Prepare Blank:** Mix 2µl of 1X TE buffer with 200µl of QuantiFluor[®] dsDNA Dye working solution in an empty 0.5ml PCR tube. Vortex well and protect tube from light.
- 4. **Prepare Standard:** Add 2µl of the provided DNA Standard (100ng/µl) to 200µl of QuantiFluor[®] dsDNA Dye working solution in an empty 0.5ml PCR tube. Vortex well and protect tube from light.
- 5. **Prepare Unknown(s):** Add 1–20µl of unknown samples to 200µl of QuantiFluor[®] dsDNA Dye working solution in 0.5ml PCR tubes. Vortex well, and protect tube from light.
- 6. Incubate the prepared samples at room temperature for 5 minutes, protected from light.
- 7. Select the dsDNA protocol on the Quantus[™] Fluorometer.
- 8. If needed, calibrate the Quantus[™] Fluorometer by reading the blank (Step 3) and standard (Step 4) samples in the Calibration screen, then select "Save".
- 9. Enter the volume of the unknown sample (1–20µl used in Step 5) and desired concentration units.
- 10.Measure fluorescence of the unknown sample and record the final sample concentration results.

Instructions for Use of Product E2670.



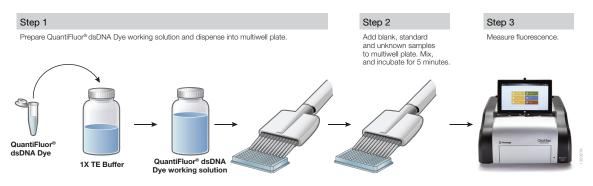
Materials Required

- multiwell detection instrument capable of measuring fluorescence (e.g., GloMax® Discover System [Cat.# GM3000])
- Nuclease-Free Water (Cat.# P1195)
- black, flat-bottom 96-well plates
- 1.5ml tubes

Warm all assay components to room temperature before use.

The *QuantiFluor® dsDNA System Technical Manual* #TM346 and *GloMax® Discover System Operating Manual* #TM397 are available at: **www.promega.com/protocols**

Multiwell Plate Protocol



We recommend preparing a standard curve that extends above and below the likely concentration range for your unknown samples.

- 1. Prepare 1X TE Buffer: Dilute the 20X TE Buffer 20-fold with nuclease-free water.
- 2. Prepare Working Solution: Dilute the QuantiFluor® dsDNA Dye 1:400 in 1X TE buffer, and mix.
- 3. **Prepare a Standard Curve:** Using dsDNA standards, prepare samples that result in 0.05–200ng/well when dispensing 10µl of standard to each well.
- 4. Pipet 200µl of QuantiFluor® dsDNA Dye working solution into each well.
- 5. Dispense 10µl of the prepared dsDNA standards prepared as shown in Figure 1.
- 6. For the blank, pipet 10µl of 1X TE Buffer into row H.
- 7. Add 1–20µl of unknown sample to the remaining wells, recording the dilution factor.
- 8. Mix the plate thoroughly.
- 9. Incubate assays for 5 minutes at room temperature, protected from light.
- 10.Measure fluorescence (504nm_{Ex}/531nm_{Em}) using a plate reader. For the GloMax[®] Discover System, select "QuantiFluor dsDNA System."
- 11.Calculate the dsDNA concentration by copying and pasting your raw fluorescence data into our online tool: www.promega.com/resources/tools/ quantifluor-dye-systems-data-analysis-workbook/

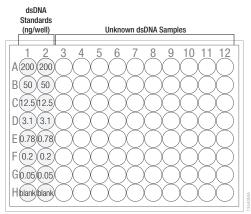


Figure 1. Dispense standard dilutions and blank samples in duplicate into Columns 1 and 2 of a multiwell plate.

PROMEGA CORPORATION • 2800 WOODS HOLLOW ROAD • MADISON, WI 53711-5399 USA • TELEPHONE 608-274-4330