

Table 2. Reagent composition for each well*

RNA Standard	Blank Control	Test Sample
Serial dilutions* (100 μ L)	TE: 100 μ L	100 μ L

*Note: Add the serially dilutions of RNA standards from 15.6 to 1000 ng/mL into wells from DS1 to DS7 in duplicate.

3. Run dsDNA assay:

3.1 Add 100 μ L of StrandBrite™ Green working solution (from Step 2) to each well of the RNA standard, blank control, and test samples (see Step 3) to make the total RNA assay volume of 200 μ L/well.

Note: For a 384-well plate, add 25 μ L sample and 25 μ L of StrandBrite™ Green working solution per well.

3.2 Incubate the reaction at room temperature for 5 to 10 minutes, protected from light.

3.3 Monitor the fluorescence increase with a spectrofluorometer at Ex/Em = 490/525 nm (cutoff at 515 nm).

Note: To minimize photobleaching effects, keep the time for fluorescence measurement constant for all samples.

3.4 The fluorescence in blank wells (with the assay buffer only) is used as a control, and is subtracted from the values for those cuvettes with RNA standard or test samples. The RNA concentration of the sample is determined according to the RNA standard curve.

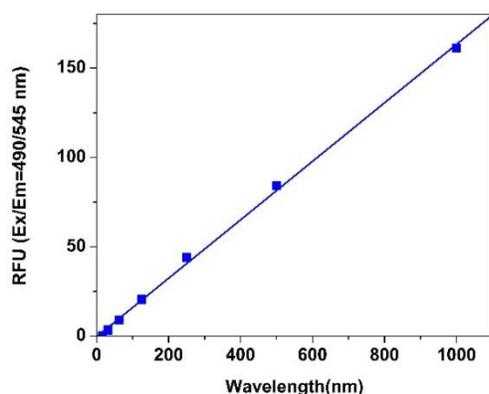


Figure 2. RNA dose response with StrandBrite™ Green in a solid black 96-well microplate and measured using a Gemini microplate reader (Molecular Devices).

References

1. Jones LJ, Yue ST, Cheung CY, Singer VL. (1996) RNA quantitation by fluorescence-based solution assay: RiboGreen reagent characterization. *Anal Biochem* 15, 265(2):368-74.
2. Le Pecq JB, Paoletti C. (1993) A new fluorometric method for RNA and DNA determination. *Anal Biochem* 17(1):100-7