

Quant-iT™ Assays

Abbreviated Protocol

IMPORTANT: Ensure all assay reagents are at **room temperature** before you begin.

1. Set up your tubes: you'll need two tubes for the standards (three for the protein assay) and one tube for each of your samples.
2. Make the Quant-iT™ **Working Solution** by diluting the Quant-iT™ reagent 1:200 in Quant-iT™ buffer. 200 μL of **Working Solution** are required for each sample and standard.
3. Prepare Assay Tubes according to the table below.

	Standard Assay Tubes	User Sample Assay Tubes
Volume of Working Solution (from step 2) to add	190 μL	180-199 μL
Volume of Standard (from kit) to add	10 μL	—
Volume of User Sample to add	—	1–20 μL
Total Volume in each Assay Tube	200 μL	200 μL

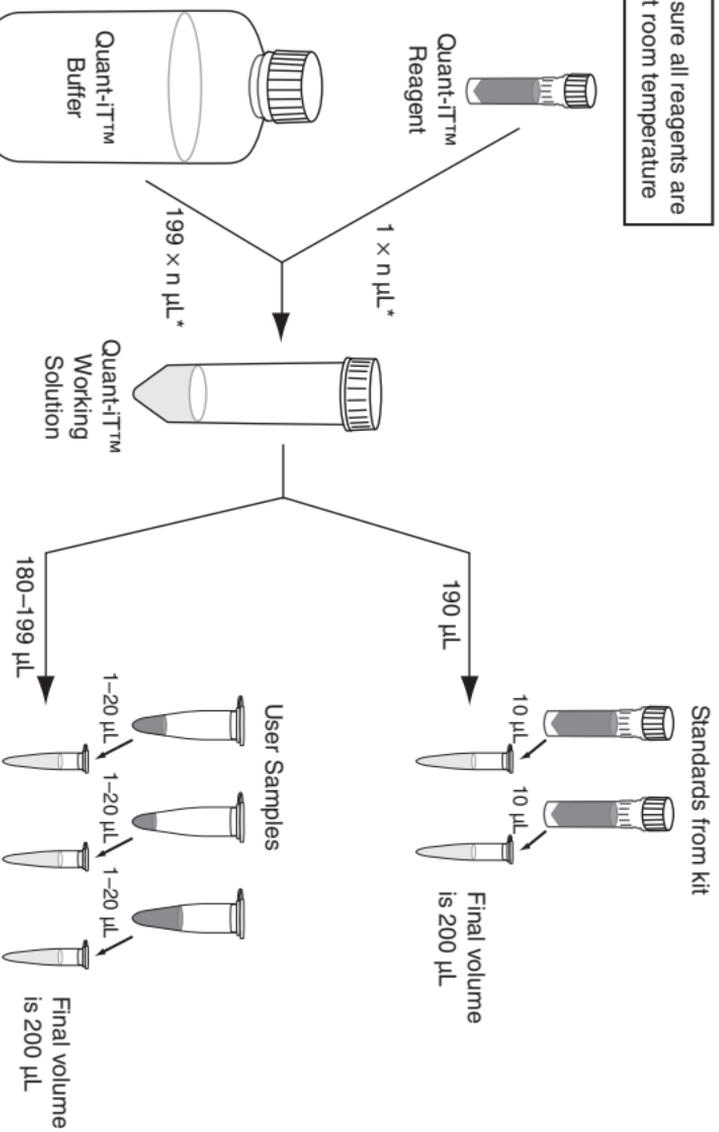
4. Vortex all tubes for 2–3 seconds.
5. Incubate the tubes for 2 minutes at room temperature (15 minutes for the Quant-iT™ protein assay).
6. Read tubes in Qubit™ fluorometer.
7. Multiply by the dilution factor (see Manual) to determine concentration of your original sample. Alternatively, choose **Calculate sample concentration** to have the Qubit™ fluorometer perform this multiplication for you.

* Use only thin-wall, clear 0.5 mL PCR tubes. Acceptable tubes include Qubit™ assay tubes (500, Invitrogen Cat. no. Q32856) or Axygen PCR-05-C tubes (VWR, part number 10011-830).

probes.invitrogen.com/qubit



Ensure all reagents are at room temperature



* where n = number of Standards plus number of Samples

 **invitrogen**[™]

Vortex all assay tubes for 2-3 seconds

Incubate at room temperature for 2 minutes (15 minutes for Quant-iTTM protein assay)

Read tubes in Qubit[™] fluorometer

