

life
technologies

QUICK REFERENCE

Qubit[®] assays

Pub. no. MAN0010876

Rev. 1.0

Qubit 3.0
Fluorometer

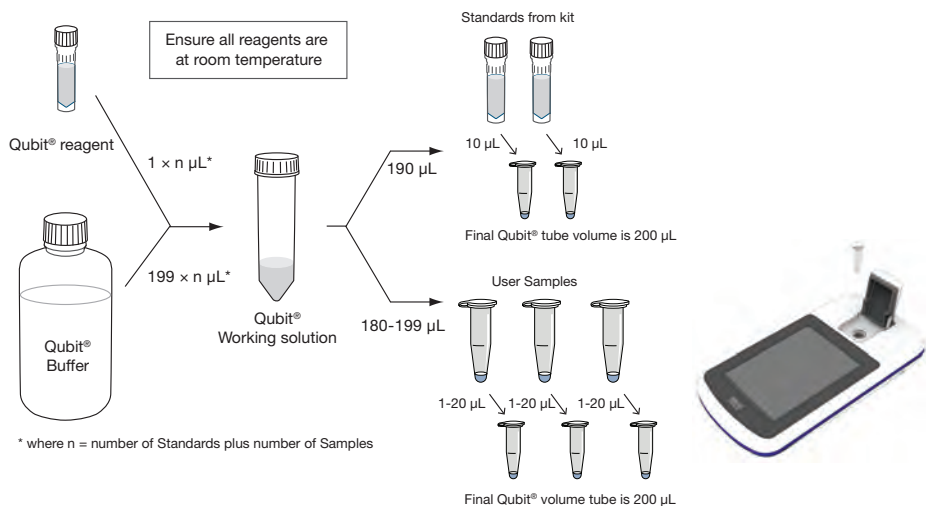
For Research Use Only. Not for use in diagnostic procedures.

NOTE: For best results, store the dye and the buffer at room temperature. Store the DNA, RNA, and protein standards at 4°C. Ensure that all assay reagents are at room temperature before you begin.

1. Set up two Assay Tubes for the standards (three for the protein assay) and one Assay Tube for each user sample.
2. Prepare the Qubit® Working Solution by diluting the Qubit® reagent 1:200 in Qubit® buffer. Prepare 200 µL of Working Solution for each standard and sample.
3. Prepare the Assay Tubes* according to the table below.
4. Vortex all tubes for 2–3 seconds.
5. Incubate the tubes for 2 minutes at room temperature (15 minutes for the Qubit® protein assay).
6. Insert the tubes in the Qubit® Fluorometer and take readings. For detailed instructions, refer to the Qubit® Fluorometer manual.

	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

* Use only thin-wall, clear 0.5 mL PCR tubes. Acceptable tubes include Qubit® assay tubes (set of 500, Cat. no. Q32856) or Oxygen PCR-05-C tubes (VWR, part no. 10011-830).



For more information, go to lifetechnologies.com/qubit

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 found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

DISCLAIMER - LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

© 2014 Thermo Fisher Scientific Inc. All rights reserved.

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.