

Qubit 4
Fluorometer

QUICK REFERENCE

Qubit™ Assays

Pub. No. MAN0017210

Rev. B.0

invitrogen
by Thermo Fisher Scientific

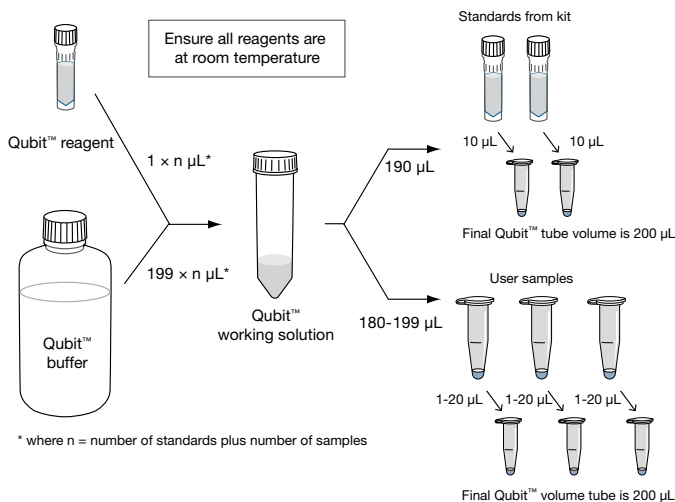
For Research Use Only. Not for use in diagnostic procedures.

- Set up two assay tubes for the standards (three for the protein or RNA IQ assay) and one assay tube for each sample.
- 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.†
- Prepare the assay tubes* according to the table below.
- Vortex all tubes for 2–3 seconds.
- Incubate the tubes for 2 minutes at room temperature (15 minutes for the Qubit™ protein assay).
- 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

† Qubit 1X dsDNA HS Working Solution (Cat. Nos. Q33230 and Q33231) needs to be prepared only once, and subsequently can be stored as a ready-to-use working solution.

* Use only thin-wall, clear 0.5 mL PCR tubes. Acceptable tubes include Qubit™ assay tubes (set of 500, Cat. No. Q32856).



Find out more at thermofisher.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.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 document is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC 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.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.



Manufacturer: Multiple Life Technologies Corporation manufacturing sites are responsible for manufacturing the products associated with the workflow covered in this guide.

ThermoFisher
SCIENTIFIC