

Qubit® RNA Assay Kit

Online Specials

Cat. nos. Q32852, Q32855

Pub. no. MAN0006880

Revision 1.0

Detailed protocol is available online at www.lifetechnologies.com/manuals, and on the USB drive provided with the Qubit® 2.0 Fluorometer.

WARNING! For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from www.lifetechnologies.com/support.

Chemical	CAS	Warning
DMSO	67-68-5	Components of the product may be absorbed into the body through the skin.

1. Set up 2 Assay Tubes for the standards and 1 tube for each user sample.
2. Prepare Qubit® Working Solution by diluting the Qubit® RNA Reagent 1:200 in Qubit® RNA Buffer. Prepare 200 µL of Working Solution for each standard and sample.

For research use only. Not for use in diagnostic procedures.

3. Prepare the Assay Tubes (use 0.5-mL PCR tubes) according to the following table.

Volume	Standards	Samples
Working Solution (from step 2)	190 μ L	180–199 μ L
Standard (from kit)	10 μ L	—
User Sample	—	1–20 μ L
Total in each Assay Tube	200 μ L	200 μ L

4. Vortex standards and samples for 2–3 seconds and incubate at room temperature for 2 minutes.
5. Select RNA Assay on the Qubit® 2.0 Fluorometer to calibrate with standards and read the samples.

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.

© 2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate[s] or their respective owners.

For support visit lifetechnologies.com/support
lifetechnologies.com