

Qubit Flex
Fluorometer

Qubit™ Flex Fluorometer

Quick Reference

Cat. No. Q33327 | Pub. Part No. 100086216 | Pub. No. MAN0018187 | Rev. A.0

Gather Your Materials



**Single Channel
Pipette (1-20 µL)**



**Multichannel
Pipette (200 µL)**



**Qubit
Assay**



**Qubit Flex
Tube Strips**



**Qubit Flex
Reservoir (10 mL)**

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.

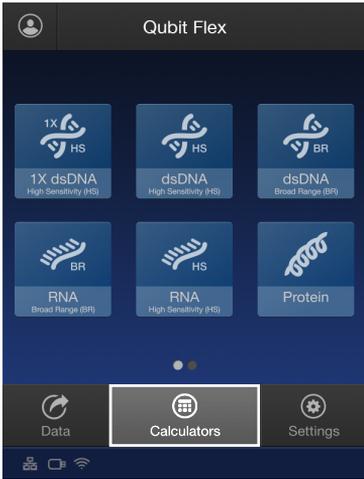
Manufacturer:
Life Technologies Holdings Pte Ltd | Block 33 | Marsiling Industrial Estate Road 3 | #07-06, Singapore 739256

Find out more at thermofisher.com/qubit

ThermoFisher
SCIENTIFIC

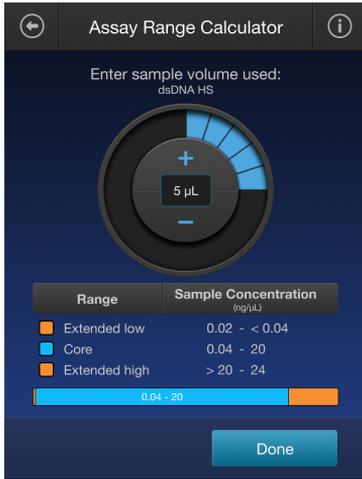
For Research Use Only. Not for use in diagnostic procedures. © 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

1 Before You Begin



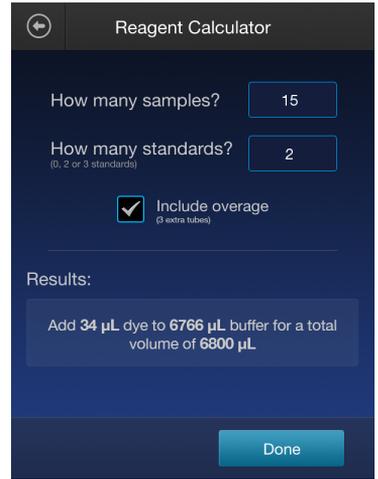
Select **Calculators** from the Qubit Flex home screen.

These two calculators will help you prepare your assays.



Use the **Assay Range Calculator** to show how the assay range changes based on the sample volume used (1-20 µL). Tap **i** to view range details.

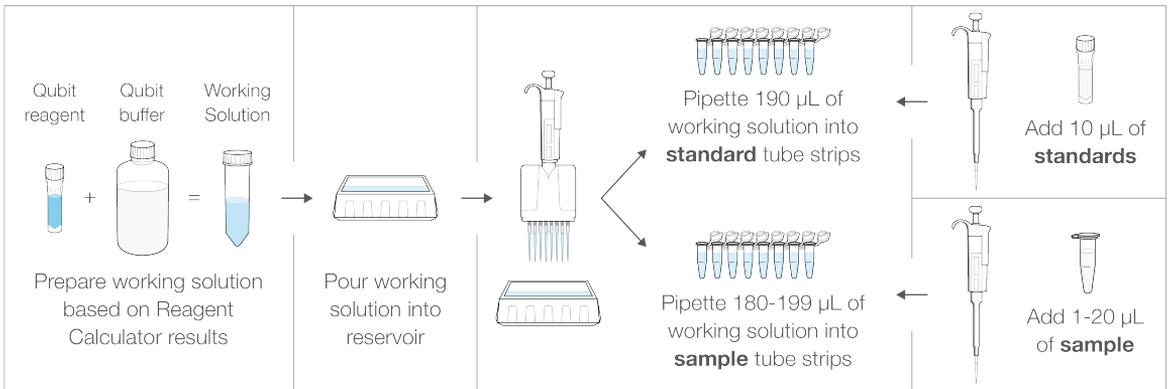
Core range = High accuracy
Extended range = Moderate accuracy



Use the **Reagent Calculator** to determine the amount of working solution to prepare.

One sample = One tube of the tube strip
One standard = All 8 tubes of the tube strip
Overage = 3 extra tubes

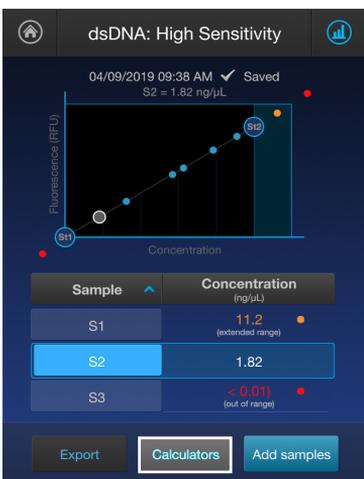
2 Prepare the Assay



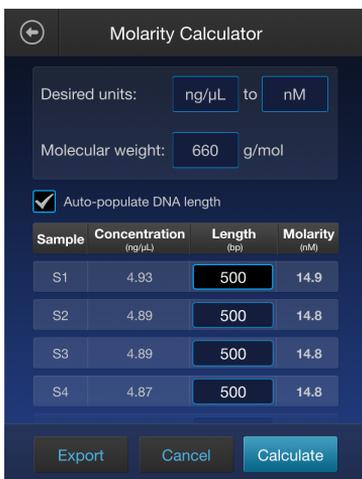
Vortex all tubes for 2-3 seconds, check for bubbles, then incubate for 2 minutes (15 minutes for protein) at room temperature

3 Read Standards & Run Samples

4 Analyze the Results



Open circles on graph = Standards
Blue = Samples in core range
Orange = Samples in extended range
Red = Samples outside of range



Use the **Molarity Calculator** to convert values to molarity based on nucleic acid length.

Sample	Add sample (µL)	Add buffer (µL)
S1	2.3	7.7
S2	2.2	7.8
S3	2.3	7.7
S4	2.2	7.8
S5	2.2	7.8
S6	2.3	7.7
S7	2.2	7.8
S8	2.3	7.7

Use the **Normalization Calculator** to determine how to dilute the samples to the same concentration.