



## Release Notes for the QX200 Droplet Reader and QX Manager Software

### Version 1.2

March 2021 (*release notes only*)

## Contents

|                                   |   |
|-----------------------------------|---|
| Introduction .....                | 1 |
| Supported Operating System.....   | 1 |
| Upgrading to New Versions .....   | 2 |
| New Features.....                 | 2 |
| Fixed Issues.....                 | 3 |
| Known Issue .....                 | 3 |
| Documentation.....                | 3 |
| Contacting Technical Support..... | 4 |
| Legal Notices .....               | 4 |

## Introduction

QX Manager Software, when connected to your QX200 Droplet Reader, provides the necessary functionality to create, run, and analyze experiments on your samples.

## Supported Operating System

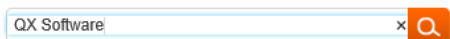
QX Manager Software is supported on a 64-bit Windows 10 operating system.

## Upgrading to New Versions

This section explains how to download and install Standard Edition. If you are using Regulatory Edition, contact Bio-Rad Technical Support.

### To upgrade the software to v1.2

1. Enter the following URL in your browser: <http://www.bio-rad.com>
2. Log into the website with your Bio-Rad user name and password.
3. Enter QX Software into the Search field and click the search icon.



4. Click QX Software to open the corresponding page.
5. Scroll to the Download section and for QX Manager Software, Standard Edition, version 1.2, click Download Software.
6. In the Software Download window, select the *I agree with the compliance requirements* checkbox, and then click Download.
7. A zip file is downloaded and appears in your Downloads folder. The download process can take a few minutes.
8. Open the Downloads folder and select the QXManagerStandard\_v1.2 zip file.
9. Right-click and select Extract All.
10. Select a destination folder and then select the *Show extracted files when complete* checkbox, and then click Extract.
11. Open the folder containing the extracted files, and locate the .exe file.
12. Double-click the file to begin the installation.
13. Follow the prompts to install QX Manager Software, Standard Edition, v1.2.

## New Features

### Automatic Thresholding for Positive Control Wells

QX Manager automatically generates thresholds for each channel, based on positive controls currently selected. QX Manager then applies those thresholds to all currently selected wells.

## Fixed Issues

- Entering commas in the Well Notes field, or entering a note on the next line, no longer pushes information entered for a well to the next well row/column in the .csv file.
- QX Manager Software no longer displays an error message that cannot be cleared when you open a data file in which the Drop-Off (DOF) experiment type is selected.
- QX Manager Software no longer displays NO CALL instead of 0 (zero) for experiments in which the Advanced Classification Mode experiment type is selected. The calculated concentration is now correct.
- Users can now change well attributes in the plate editor in a .qlp file.
- QX Manager Software no longer crashes when users select multiple wells for thresholding.
- Data now appears when users export the plate setup or analysis data to reports.

## Known Issue

- QX Manager Software incorrectly uses unselected positive control wells to calculate positive control-based automatic thresholding.

**Workaround:** For any positive control well that QX Manager should exclude from the automatic thresholding calculation, change the sample type from Pos-Ctrl to Unknown.

## Documentation

Information about the instrument and software is available in the following resources:

- QX200 Droplet Reader and QX Manager Software, Standard Edition, User Guide
- QX200 Droplet Reader and QX Manager Software, Regulatory Edition, User Guide

## Contacting Technical Support

The Bio-Rad Technical Support department in the U.S. is open Monday through Friday, 5:00 AM to 5:00 PM, Pacific Time.

**Phone:** 1-800-424-6723, option 2

**Email:** [Support@bio-rad.com](mailto:Support@bio-rad.com) (U.S./Canada Only)

For technical assistance outside the U.S. and Canada, contact your local technical support office or click the Contact Us link at [www.bio-rad.com](http://www.bio-rad.com).

## Legal Notices

No part of this publication may be reproduced or transmitted in any form or by any means, electronic or mechanical, including photocopy, recording, or any information storage or retrieval system, without permission in writing from Bio-Rad Laboratories, Inc.

Bio-Rad reserves the right to modify its products and services at any time. This guide is subject to change without notice. Although prepared to ensure accuracy, Bio-Rad assumes no liability for errors or omissions, or for any damage resulting from the application or use of this information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.

Purchase of Digital PCR products from Bio-Rad Laboratories includes a limited, non-transferable right under our intellectual property for you to use the products for your internal research purposes only. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for you to distribute or resell the products or to use the products for commercial applications, such as contract services or fee for services. No rights are granted for diagnostic uses. Information concerning a license for such uses can be obtained from Bio-Rad Laboratories. It is the responsibility of the purchaser/end user to acquire any additional intellectual property rights that may be required. The QX200 Droplet Digital PCR System and/or its use is covered by claims of U.S. patents, and/or pending U.S. and non-U.S. patent applications owned by or under license to Bio-Rad Laboratories, Inc. See [www.bio-rad.com/en-us/trademarks](http://www.bio-rad.com/en-us/trademarks) for details.

All trademarks used herein are the property of their respective owner.

Copyright ©2021 Bio-Rad Laboratories, Inc.