



QXDx BCR-ABL %IS Reporter Tool

QXDx BCR-ABL %IS Reporter is a Microsoft Excel tool that will allow an end user to upload a file exported from QuantaSoft v 1.7 software and analyze the data to generate a summary report of patients tested for BCR-ABL.

QXDx BCR-ABL Reporter is available for download at: <http://strandls.com/bioinformatics/bcr-abl-reporter/>

The download will require the user to provide the following contact information: Name, institution, phone number, email (institution specific emails only, i.e. Yahoo/Google emails are not accepted) for traceability and notification in the event the tool is modified.

Key Features of the Tool

- Upload '.CSV' file from QuantaSoft v1.7
- Automated analysis
- Generate a Summary Report if the run passes
- Manual Result Generation – Analyze and pass a test to report result by overriding analysis rules (while not recommended, this will give the user additional flexibility)
- Samples can be distributed over any number of wells for sample-specific measurement with higher or lower desired sensitivity – the more wells used the higher the test sensitivity

Droplet Digital PCR

Key User Input

- Calibrators and controls minimum and maximum value card provided in the QXDx BCR-ABL %IS Kit
- CF Value – also part of QXDx BCR-ABL %IS Kit
- Minimum ABL copies per well: default value = 3000, gives user the flexibility to use higher/lower value based on their validation/verification

Key User Steps to Ensure Positive Experience

1. During plate setup
 - a. Name calibrators and controls wells appropriately (without '~'), example: "0.1% cal, 10% cal, Neg ctrl,..."
 - b. Name no template control appropriately – "NTC"
 - c. Name all sample wells appropriately (avoid keywords for controls/calibrators and NTCs)
2. Post droplet reader run and troubleshooting¹, export data to '.CSV' file and save file
3. Open QXDx BCR-ABL reporter file
 - a. Review/edit %IS values for calibrators and controls (in column C & D)
 - b. Modify CF value (as required by kit lot)
 - c. Click "Load Data File and Analyze"
 - d. Select '.CSV' file exported from QuantaSoft v1.7
4. Run summary generated w/BCR-ABL patients results, when quality rules passes
5. Print report

¹Refer to QXDx BCR-ABL IFU (P/N 12006672) for details on run and troubleshooting.



**Bio-Rad
Laboratories, Inc.**

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com

Life Science
Group

Website www.bio-rad.com **U.S.** 1-800-224-6783 **Australia** +61(2)9914-2800 **Austria** +43-1-877-8901 **Belgium** +32(3)710-53-00 **Brazil** +55(31)3689-6600 **Canada** 1-514-334-4372 **China** +86-21-61698500 **Czech Republic** +420-241-430-532 **Denmark** +45-4452-1000 **Finland** +358-9-804-22-00 **France** +33-1-47-95-60-00 **Germany** +49(0)89-318-840 **Greece** +30-210-7774396 **Hong Kong** +852-2789-3300 **Hungary** +36-1-459-6100 **India** +1800-180-1224 **Israel** +972-3-9636050 **Italy** +39-02-216091 **Japan** +81-3-6361-7070 **Korea** +82-2-3473-4460 **Mexico** +52(65)5488-7670 **The Netherlands** +31-318-540666 **New Zealand** +64-9-415-2280 **Norway** +47-23-38-41-30 **Poland** +48-22-3319999 **Portugal** +351-21-472-7700 **Russia** +7-495-721-1404 **Singapore** +65-6415-3170 **South Africa** +27-11-442-85-08 **Spain** +34-91-590-5200 **Sweden** +46-8-555-127-00 **Switzerland** +41(0)26-674-55-05/06 **Taiwan** +886-2-2578-7189 **Thailand** +662-651-8311 **United Kingdom** +44(0)20-8328-2000