

BioQule V1.0.0 Release Notes

Bug Fixes

1. Pressing the Return key on windows other than the Barcode Entry window no longer transports users to different windows or produces popups
2. A script which BioQule cannot run no longer crashes the software. It lets the user know that the script can't be run and transports the user back to the Main Window
3. Accidentally disconnecting the USB connection from BioQule to computer no longer crashes the Software. Instead it lets the user know that the USB has disconnected and returns the user to the Device Select Page.
4. The Abort button no longer causes the software to crash if the user runs and aborts a run right after aborting a run instead of restarting. Instead the software automatically restarts upon abort.
5. If BioQule results folder does not exist or has been accidentally deleted the software can create it.
6. The Information put in the table does not get preserved when you cancel from the final information page and rerun using the same instance of the software
7. Popups are no longer scalable
8. Can no longer select multiple devices in the Select Device page

Added Features

1. Can now change the directory that the results get saved in on the Enter Information Page
2. The Results CSV also contains information about the run: Computer ID, User, Time Start, Time Stop, Assay Run, Box ID ...