



CellReporterXpress®

Image Acquisition and Analysis Software

Version 2.6.130

Known Issues (for MolDev Internal Use Only)



On-the-fly and post-acquisition analysis produce different results

The on-the-fly analysis of a multisite, non-stitched experiment will not perfectly match the analysis of the same experiment if the analysis is run post acquisition. The root cause of this is because each tile is analyzed individually in the on-the-fly analysis. Conversely, when analysis is run after acquisition, the entire image is analyzed at once. The workaround is to stitch multisite experiments or to always perform analysis the same way (that is, always on-the-fly or always post-acquisition).

Add Analysis to Experiment gets stuck trying to capture images when in Comparison mode

When a new analysis is added to an experiment, you can view it in single mode or comparison mode. Despite being in analysis rather than acquisition, if you change the locations, the software indicates it is “Capturing images”. If you have chosen Comparison mode, the software freezes with the message “Failure: Unknown server error”.

Existing IMEX versions are not removed during upgrade

When you install the latest version of the software, the CellReporterXpress Installation Utility should remove all IMEX icons, buttons, and functionality. However, if the user has an earlier version of IMEX on their system (either from 2.0 or 2.1 or from a pre-release version of 2.5), it is not removed as expected.

Live Preview stays on when the top door is opened

When live preview is on, and the top door is opened (either by pressing the button at the top right of the instrument or from the software), live preview does not turn off as expected and light can be seen inside the instrument.



WARNING! Do not look directly at the light inside the instrument.

Running two stitched protocols on the same host computer fails both

When the host computer is connected to more than one instrument, and you attempt to simultaneously run two acquisitions with stitched protocols, both acquisitions will fail. As a workaround, run stitched-protocol acquisitions one at a time.

Setting Z offset with transmitted light first sets the position for fluorescent channels

When an acquisition is configured to capture both a single fluorescent channel and transmitted light, and the transmitted light is set to be captured first, the focus offset value entered for the fluorescent channel will also be used for transmitted light. Further adjustments to the focus offset values yield unexpected results. As a workaround, set transmitted light to be captured last, set the focus offsets, and then set transmitted light to be captured first.

Added analysis for a single slide of a multislides experiment causes an error

When you add an analysis for a single slide with custom analysis regions to an existing multislides experiment, the software may stop responding or a timeout error may occur. As a workaround, include all analysis regions when you add an analysis.