



CellReporterXpress®

Image Acquisition and Analysis Software

Version 2.5.769

Known Issues (for MolDev Internal Use Only)

Canceling an acquisition from the monitor page doesn't stop the instrument

When you cancel a running acquisition from the Monitor page, the software cancels the experiment and moves it from the IN PROGRESS tab to the SUCCEEDED tab. However, the experiment may not be canceled on the instrument. In this case, the instrument will finish the acquisition, and this can take a considerable amount of time. The only way to stop the acquisition on the instrument is to manually restart it by pressing the power button once.

Stitching timeouts during acquisitions with many sites

If you are running an acquisition with more than 900 sites, the acquisition will likely time out or you may have empty tiles in the data. Currently, the only workaround for this is to limit the number of wells in the acquisition protocol so that the total number of sites for the entire plate is less than 900.

Acquisition time outs with large stitched areas

If you acquire large stitched areas with multiple channels, the host computer saves the images two times: once in the experiment folder and again in the cache on the C drive of the host computer (which is used for stitching). Because the cache is not part of the storage calculation on the run protocol step, there is no indication of whether sufficient space is available to run the experiment. If the host computer runs out of memory, the acquisition will time out and fail, and no images will be available for the experiment. This issue is most likely to occur for slide acquisitions, but it can also occur with plates. As a workaround, confirm that there is enough space on the host computer for double the experiment size.

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

The on-the-fly analysis of a multi-site, 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 multi-site experiments or to always perform analysis the same way (that is, always on-the-fly or always post-acquisition).

Analysis Summary with multiple acquisition ROIs

If you analyze an experiment with multiple acquisition regions, the summary data only includes measurements for a single site. The workaround is to analyze one site at a time.

Segmentation overlays for Neurite Tracing and other analysis are missing in parts of the images

When analysis is tested, segmentation is drawn correctly. When analysis is run, the measurements values look correct, but the segmentation mask shows large gaps as if zoomed down in resolution and then zoomed back up.

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".

Exported colorimetric images > 2 GB are not down sampled correctly

When you export a stitched or tiled colorimetric image greater than 32K in the X or Y dimension or 2 GB in file size, the image is scaled down too much, resulting in an image with a lower resolution than expected.

Export Time Point(s) Images does not work

When you try to export images of individual time points from the time points tab on the Experiments page, a new browser window opens with the message "This site can't be reached, localhost refused to connect". This link was not updated during CellReporterXpress 2.5 development; it is attempting to export images through IMEX, which is no longer included with the software.

Existing IMEX versions are not removed during upgrade

When installing 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.