

Release Notes

v2.1.0 (19 March 2024)

New Features

- Volume pairs can be manually registered to support overlaying volumes that were not originally acquired in the same location
- 2Dt volumes support the creation of markups for quantifying data
- A Keyboard Shortcuts reference has been added to the Help menu
- IVIS 3D FLIT volumes can be loaded for basic visualizing

Improvements

- User interface updated to adopt the Revvity design system to have a consistent user experience across various Revvity software applications
- An all new application icon corresponding to the Revvity Life Sciences software portfolio
- Shear Wave Elastography scan plan settings includes a "Total Number of Steps" option to maintain the same number of acquisition locations while adjusting the ROI size
- Optimized Linear Array TGC values to help avoid datasets appearing too bright under default conditions
- Caliper properties can be controlled in a new section in the "Utilities" area of analysis allowing for customization of color and glyph type
- Study report can now be generated to include information for secondary volumes
- Volume rendering is supported for 3D volumes that are secondary volumes
- Shear Wave Elastography processing dialog updates only when showing one type of scan
- Easier viewing of log file information inside the Report a Problem dialog
- Markup statistics for microCT volumes can be exported using the "Export Stats From Folder" functionality
- The Wobbler transducer driver was updated to a Microsoft generic driver to avoid CVEs (CVE-2017-14075, CVE-2017-14153, CVE-2017-14344) impacting the old version
- Support for reverting Linear Array firmware to a version compatible with SonoEQ 1.12.x and earlier has been removed
- Increased tooltip usage throughout application for improved learnability

Fixes

- Acquisition 3D Scan ROI size and position were not being remembered properly when changing imaging modes
- M-Mode artifacts were being generated when using Linear Array outside of M-Mode
- Shear Wave Elastography lower threshold value was not scaled appropriately when changing representation
- Secondary statistics were sometimes not saved for 3D Shear Wave Elastography and B-Mode volume pairs
- Certain keyboard shortcuts could be executed resulting in a running acquisition not processing new instructions

- Stats CSV files could have content shifted under the incorrect header if a comma was present in the Volume Filepath field
- Heart Finder scans could be started in the 3D Scan tab with the ROI hidden
- Frame spacing was not restored after running Heart Finder scans
- Slice offsets were not being centered on a loaded segmentation
- Irrelevant metadata was being saved with certain acquisitions
- Fix the radiance scalar bar being shown in acquisition when swapping from analysis
- Entering Research Imaging Mode would launch the Quality Control window
- Servicing Dashboard Quality Control individual test buttons could be pressed while in the middle of a QC run
- Servicing Dashboard Quality Control was not applying appropriate settings for the brightness tests