



19 January 2022

## Updates in Goldcrest Patch I

### How to get the upgrade

Installation is easy enough for you to do yourself, but it is also perfectly okay to ask for our technical team to do next time they visit you. We can also offer to do it remotely if you allow remote access.

The software can also be accessed here:

- <https://ncx.sophion.net/index.php/s/0Ggfty6u1Gdw1Pn>
- pw: Qube4ever

This version is called Goldcrest Patch I and has the following version numbers:

- Viewpoint 2.6.70
- Viewpoint Maintenance 2.6.70
- Analyzer 6.6.70

### New features

- Default plans: Customer specific: Added block template for IC exchange with a single liquid exchange cycle.
- Genedata export can now also get baseline and full block periods from Rule Based Hill Fit methods
- Genedata export: A panel is shown where the user can select which liquid periods to use for baseline and full block
- Genedata exports can now only contain experiments that have been run with the same Assay (QPatch)/Experiment protocol (Qube)

### User interface changes

- Analyzer no longer prompts for matching compound plates. Instead, it relies on the automatic matching compound list matching
- Genedata export is now enabled as default

### Bug fixes

- Analyzer: Export to clipboard did not handle barcodes with line breaks very well
- Analyzer: Result tables did not always update when they only showed selected data
- Barcode reader: remove line feed from output format
- Data migration did not work when a protocol run had been data reduced shortly before
- SSD Switch: Dataserver could not always login after SSD Switch, because of an old database driver
- Analyzer: Could not export runs containing idle sweeps to CiPA or Genedata format
- Temperature: Combined measurement action could in some cases hang when requesting temperature
- Run down compensation: Fixed the most often occurring null pointer exception
- Calling listener error
- Transformed Boltzmann Fit method in I-V plot does not work when the voltage steps were given from higher to lower voltages
- Analyzer could not open some projects containing idle sweeps

- Analyzer: It was not always possible to select the correct liquid periods in data transfers
- Backup would hang if there was still a run running. It would not start again until the data PC software was restarted
- Data migration: Could not migrate projects that contained Boltzmann fits
- Data migration failed if the project had already been copied to the destination
- Boltzmann fit shows negative slope, no matter shape
- When changing colors in a sweep plot the changes are not implemented in the corresponding legend

### Performance and other changes

- Genedata & CiPA: Improved export speed on Qube by handling multiple experiments together

### Other

- Default plans: Customer specific: Added two worktable protocols for screening with special MTPs
- Default plans: Customer specific: Added IC exchange block template with different docking height
- Genedata & CiPA: Allows now exports containing compounds without concentration
- Genedata: Default is now to store in the new format (multi-xml, reduced data markup)
- Debug files now also contain S.M.A.R.T info from Qubes