

# Higher throughput with CrysScreen™

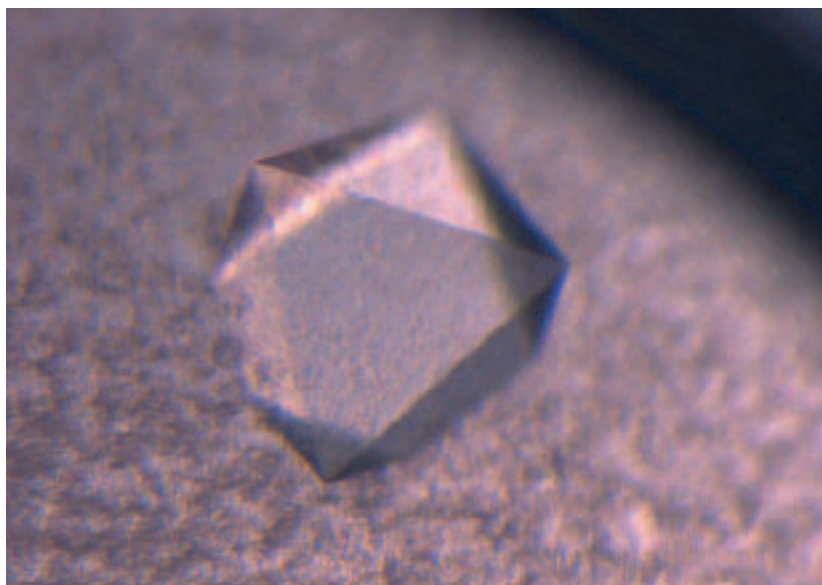


CrysScreen software provides crystallographers with an intuitive user interface to help manage complex protein crystallography protocols on Freedom EVO® workstations. The software guides users through automation of protein crystallography trials, including buffer creation, set-up of experimental conditions, running experiments, data management and results analysis.

The latest release of the software offers higher throughput than ever before, with an enhanced worklist generator to allow parallel pipetting tasks on a single instrument, and straightforward control and review of project management using the powerful Project Tree. To ensure compatibility with your laboratory workflow, CrysScreen is closely integrated with Tecan's Freedom EVOware® instrument control software, and has now been updated to guarantee full compatibility with Windows Vista® for simple control of hanging drop, sitting drop and microbatch under oil crystallization experiments.

To find out more on Tecan's CrysScreen, visit [www.tecan.com/cryscreen](http://www.tecan.com/cryscreen)

Windows Vista is a registered trademark of Microsoft Corporation.



## FE500pro™ – pre-analytics made easy

Automation is fast becoming the method of choice for clinical laboratories facing demands for increased throughput and traceability. The FE500pro is Tecan's front-end solution for automation of pre-analytical processing, improving process security and safety, while reducing costs and allowing staff to concentrate on other tasks.

The FE500pro offers walkaway processing, incorporating a wide range of functions onto a compact system, including:

- sample registration and pre-sorting
- centrifugation
- liquid level verification
- clot detection

- de-capping
- secondary tube labeling
- aliquoting
- sorting of tubes into analyzer racks

The FE500pro has an enclosed worktable to ensure biosafety and minimize risk of sample contamination, and the platform's twin lane conveyor maximizes processing flexibility. The system uses automatic barcode and liquid level verification to guarantee security of patients' samples during processing, and communicates directly with your LIMS to generate the correct number of secondary tubes for your workflow.

To find out more on Tecan's FE500pro, visit [www.tecan.com/fe500](http://www.tecan.com/fe500)

The FE500pro workstation.

