

# TouchTools Suite™ 2.0

Since its launch last year, Tecan's user-friendly TouchTools Suite has made automated liquid handling more accessible to many scientists, by allowing direct real-time control of automated robotic pipetting systems through a touchscreen interface. It does not require any programming skills or pre-defined scripts to operate, and can virtually turn your robotic system into an automated hand pipette. The popular system has now been further enhanced by a set of added features to provide even more flexibility.

In TouchTools Suite 2.0, the unique, patent-pending Instant Pipetting™ module is now compatible with any Tecan pipetting arm, including all liquid handling (LiHa) arms, 8 Plus 1 Access™, MultiChannel Arm™ 384 and MultiChannel Arm 96. This makes straightforward procedures of a wider range of applications possible, including fast full plate replication. The new version is also able to control the tip spacing of 8-channel LiHa arms, which allows more complex pipetting patterns, even into 384-well plates.

Hit-Picking mode is another new feature that has been developed in response to many user requests. Hit-Picking mode displays the plate on the graphic user interface and allows the user to select individual wells through the touchscreen, commanding the automated platform to instantly access those positions.

Other improvements to the TouchTools Suite allow customized input screens to tailor the user interface to the laboratory's needs. A custom-loading guide is now able to display customized instructions saved as PowerPoint®, PDF or other formats. In the QuickStart module, the settings can be personalized, so that users can only access the methods that they are authorized to use.

Users of Tecan's robotic instruments can also look forward to a new version of the Common Notification System (CNS), due to launch in the near future, which will make the remote access and monitoring of the laboratory's automated systems from mobile devices even easier.

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

PowerPoint is a registered trademark of Microsoft Corporation in the United States and other countries. iPod touch is a trademark of Apple Inc., registered in the US and other countries.

App Store

