

# The power is in your hands with Freedom EVOware®

Tecan software continues to bring reliability and process security to users of automation. Freedom EVOware 2.2 software makes it easier than ever to control pipetting and scheduling operations of Freedom EVO® liquid handling platforms. Based on the same intuitive graphical user interface, version 2.2 of this software offers greater flexibility and ensures full compatibility with the latest hardware modules – all without compromising on operating speed and process security.

With Instant Pipetting™, Tecan introduces a completely new way to control pipetting robots, without the user needing any previous knowledge of programming. Instant Pipetting uses a touchscreen to enable users with any skill level to define applications and directly control pipetting with a liquid handling robot in real time. This innovative feature allows the user to perform common pipetting operations, like sample distribution, reagent addition and serial dilution, in a very easy and interactive way. All steps can be repeated to perform multiple iterations, or saved and recalled as a

### Software innovation

<b>Program</b>	Define your pipetting step easy and simple	<b>Instant Pipetting</b>
<b>RunControl</b>	Recall your pipetting easily via iPhone	<b>Phone RunControl</b>
<b>Monitor</b>	Monitor your instrument while performing your pipetting step	<b>CNS on iPhone or CNS gadget on desktop</b>
	Monitor your liquid channel and plates while performing your pipetting step	<b>iPhone monitoring applications</b>




script in Freedom EVOware or mobile phone applications, for increased ease and speed.

Once your assay is running, it is now easier than ever for you to stay connected to your application with the Common Notification System (CNS). This technology enables direct remote monitoring using your mobile phone device, your office desktop through a Windows

Vista® gadget, or via any networked computer with a web browser. The status of the system is immediately visible and if user interaction is required, it will be brought to your attention.

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

Vista is a registered trademark of Microsoft Corporation.

## Increased throughput and flexibility with Tecan quality MultiChannel Arm™ 96 nested disposable tips

Tecan's new disposable tip products for use with the MultiChannel Arm 96 option for Freedom EVO® automated liquid handling workstations offer higher throughput through improved logistics possibilities.

Tips without filters can be nested in passive stacks of up to 800 tips per SBS position on the deck, or externally in Te-Stack™ modules to provide a capacity of 3,840 tips per Te-Stack column, maximizing walk-away time and improving throughput. The nested design also offers environmental and logistical benefits, including reduced storage space requirements by 60 % compared to non-nested tip formats, lower transport costs and less wastage by using less packaging materials.

The fully SBS-compatible racks have been designed for increased throughput and reliability by optimizing the flexibility of the workstation deck space. Systems fitted with Te-Stack modules can now use the same SBS-format columns for tip racks and microplates, and platforms set up for non-SBS disposable tips can easily be converted.

All MCA 96™ disposable tips are produced in a cleanroom and are certified to be free of DNA, DNase/RNase, pyrogens and endotoxins. Non-nested disposable tips are also available in a sterile format.

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

Disposable tips stored externally in a Te-Stack module



Disposable tips nested in stacks on the workstation deck

