

# New Active Worktable



Active Worktable has been designed for simplicity in operation, bringing extra process security, control and monitoring to research scientists.

The new Active Worktable option for the Freedom EVO® 100, 150 and 200 platforms, launching at Lab Automation 2010, offers benefits of extra process security, control and monitoring to research scientists. It brings the functionality available in Freedom EVOLyzer® and Freedom EVO Clinical workstations into the researchers' domain, with added flexibility to run their own scripts for various applications.

Active Worktable has been designed for simplicity in operation, keeping the user informed of what is on the deck through a graphical user interface and, with red and green lights to signal carrier status. All automated applications can benefit from the quick operational status update that is available at a glance.

## Worktable monitoring ensures safety

The Active Worktable closely monitors all user interactions with the worktable, and confirms all loading and unloading actions using a magnetic sensor on each individual grid on the worktable. Active Worktable notifies the user what needs to be done, and where carriers need to be positioned with multicolored LEDs on each grid. It also monitors the worktable during a run and, if it detects that a required carrier has been removed, it will stop the system to avoid false results. Optional PosID™ adds automatic barcode scanning and evaluation of the samples, loaded plates and reagent containers to ensure further process security.

## Simple interaction with the operator

The Active Worktable's integrated system of green and red lights simplifies the placement of sample and reagent racks on the instrument. The lights, together with an acoustic signal, alert the user whenever interaction with the system is needed, making it easy for the operator to know the status of the workstation.

Active Worktable has been tightly integrated into Freedom EVOware® Sample Oriented, so customers are offered the flexibility in the control of added modules as well as wizard-based sample pipetting.

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