

# Around the clock monitoring with CNS

Tecan's Common Notification System (CNS) is now available for the Infinite® PRO series of microplate readers, allowing users to remotely monitor the status of any Infinite reader using Magellan™ software (v7.2). With CNS, users can now easily check if their Infinite reader is performing a measurement, awaiting a predefined user action, or ready for the next assay plate.

CNS allows readers to be remotely monitored through any networked computer or mobile device, offering increased walkaway times and giving laboratory scientists the freedom to perform other tasks while their assays are

running. This innovative feature is particularly useful for overnight operations, such as long-term microbial studies with the patent-pending Gas Control Module (GCM™), ensuring greater process security and alerting the user if there is a problem, for example running low on oxygen or carbon dioxide. Together with the existing Freedom EVO® Remote for monitoring of liquid handling instruments, this latest addition to the CNS portfolio provides complete peace of mind for walkaway laboratory automation.

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



CNS helps users to stay informed around the clock

# Tecan launches new options for Cavro® Omni Robot at AACCC's Clinical Lab Expo

Tecan has extended the flexibility of its popular Cavro Omni Robot with the introduction of embedded control functions and the option to select single axis configurations. Unveiled with the launch of the Cavro Omni Robot Version 4.0 at the AACCC's 2014 Clinical Lab Expo, these latest updates will make it even easier to configure the Cavro Omni Robot to suit specific applications or instrument designs.

The robot's new Embedded Command Processor Mode allows direct communication for precise control and coordination of axis movements and liquid handling operations. This OS-independent command schema allows line commands to be sent directly to the robot using virtually any computer or

custom control board, complementing the existing Windows®-based Command Processor Mode for greater integration flexibility.

The modular design of the Cavro Omni Robot allows users to choose from various lengths and orientations of all three axes, including a choice of single or dual arm X-axis configurations. To ensure there is a Cavro Omni for virtually every application, it is now also possible to choose any combination of X-Y or Y-Z axes, offering instrument designers the option to create their own axes without the time and expense of developing a complete liquid handling solution. These latest updates further extend the versatility of the Cavro Omni Robot and, together with the system's extensive choice

of liquid handling options and finishes, provide exceptional flexibility for OEM liquid handling applications.

To find out more about Tecan's Cavro Omni Robot, visit [www.tecan.com/omnirobot](http://www.tecan.com/omnirobot)



The Cavro Omni Robot offers exceptional integration flexibility

Perfection isn't instant.  
It evolves.

Tecan Cavro. 40 years of innovation.

