

A close-up, shallow depth-of-field photograph of several microcentrifuge tubes in a rack. The tubes are clear plastic with white caps and labels. The word "Cav" is visible on the labels. The background is dark and out of focus.

The OEM components of tomorrow. Today!

More reasons to Talk to Tecan

www.tecan.com/components

The new Cavro® generation from Tecan OEM Components



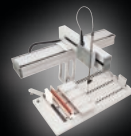
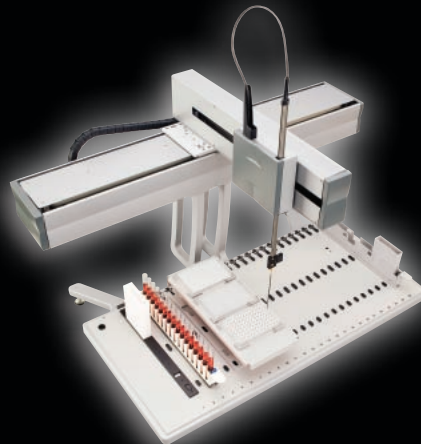
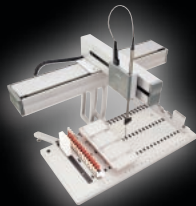
NEW Cavro® Centris Pump

Our new pump takes accuracy and precision to another level. Using long-life ceramics for the fluidic path outperforms conventional pumps in precision and longevity, even with aggressive solvents. Users can do more with one pump than ever before, offering a broader working volume range from a single syringe size.



NEW
Cavro® Omni Robot

Our new robot's innovative modularity offers superior reliability and flexibility, making it easily configurable for many different applications. It's also designed for easy integration, being both simple to install and to operate.



More reasons to Talk to Tecan

www.tecan.com/components

Tecan Systems, Inc.
2450 Zanker Road
San Jose, CA 95131
Tel +1 408.953.3100
Fax +1 408.953.3101
Toll Free +1 800.231.0711 (US)

www.tecan.com/components