



A laptop is not included with the Cavro Integration Kit.

## Cavro® Integration Kit

Have your components running in minutes

In today's fast-paced environment, time-to-market is a major consideration for any instrument designer. Selecting proven, off-the-shelf components is a sure way to accelerate your instrument development.

You can also speed up and simplify validation of your instruments by choosing OEM components that have already been extensively tested and validated. Tecan Cavro brand components are all constructed to meet ISO 13485 and ISO 9001 requirements, and also comply with national and regional standards, including Europe's 2006 RoHS Directive and the US FDA's QS regulations.

The Cavro Integration Kit provides all the necessary items for you to start evaluating your Tecan Cavro components straight out of the box, so you can focus on your application and OEM development. The kit is designed to be easy to use, so that you can be up and running in a matter of minutes. The Cavro Integration Kit operates and supports all Tecan Cavro components – you can even run pumps and robots simultaneously.

The kit includes a power supply, cables, a communications interface and software; you can command and operate all connected devices via a PC through the intuitive graphical user interface.



The Cavro Fusion Software included in the Cavro Integration Kit automatically identifies which communications interface port your Tecan Cavro component is connected to. This lets you focus on operating the component, rather than connecting it.

## Easy evaluation of your components

The Cavro Integration Kit is available in three configurations:

- The **Basic** kit contains a 3 A power supply and operates up to three pumps via RS-485.
- The **Power** kit comes with a 5 A power supply and is able to operate up to three pumps via RS-232, or one RSP plus one to two pumps. A Cavro RSP interface cable is also included in the **Power** kit.

In the **Basic** and **Power** kits, the communications interface and the power supply for your Tecan Cavro component are combined in the Cavro Hub, allowing easy connection of your component through a USB port.

- The **Ethernet** kit also contains a 5 A power supply and supports a single-arm Cavro Omni Robot with up to two pumps.

### Specifications

	Cavro Integration Kit – Basic	Cavro Integration Kit – Power	Cavro Integration Kit – Ethernet
<b>Operation</b>	1-3 pumps	1-3 pumps Or 1 single-arm Cavro RSP and 1-2 pumps Or 1 dual-arm Cavro RSP and 1 pump	1 single-arm Cavro Omni Robot and 1-2 pumps
<b>Power supply</b>	24 V/3 A	24 V/5 A	24 V/5 A
<b>Cavro Hub with USB cable</b>	●	●	
<b>Three-pump cable (to connect with Cavro Hub)</b>	●	●	
<b>Cavro RSP interface cable</b>		●	
<b>Ethernet cable</b>			●
<b>Two-pump cable (to connect with Cavro Omni Robot)</b>			●
<b>Cavro Fusion Software</b>	●	●	●
<b>Quick start guide</b>	●	●	●

### Minimum PC requirements

Operating system Windows® XP, Windows 7

### Spares

Power supply (3 A)

Power supply (5 A)

Cavro Hub (USB/ RS-232)

Three-pump cable (to connect with Cavro Hub)

Cavro RSP interface cable

Ethernet cable

Two-pump cable (to connect with Cavro Omni Robot)

USB A-B 0.9 m cable

Cavro Fusion Software

### Tecan – Who we are

Tecan ([www.tecan.com](http://www.tecan.com)) is a leading global provider of laboratory instruments and solutions in biopharmaceuticals, forensics and clinical diagnostics. The company specializes in the development, production and distribution of automated workflow solutions for laboratories in the life sciences sector. Its clients include pharmaceutical and biotechnology companies, university research departments, forensic and diagnostic laboratories. As an original equipment manufacturer (OEM), Tecan is also a leader in developing and manufacturing OEM instruments and components that are then distributed by partner companies. Founded in Switzerland in 1980, the company has manufacturing, research and development sites in both Europe and North America and maintains a sales and service network in 52 countries.

**Australia** +61 3 9647 4100 **Austria** +43 62 46 89 33 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **Denmark** +45 70 23 44 50 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170  
**Italy** +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Singapore** +65 644 41 886 **Spain** +34 93 490 01 74 **Sweden** +46 31 75 44 000  
**Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +43 62 46 89 33

### Safety and regulatory compliance

Tecan Cavro OEM components are designed and manufactured to the highest quality, meeting EU RoHS directives. Many Tecan Cavro components are UL-recognized components. Tecan Systems' Quality Management System (QMS) is compliant with ISO 9001, ISO 13485, and the Quality System Regulations (QSR).

Tecan Group Ltd. makes every effort to include accurate and up-to-date information, however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided. Changes can be made at any time without notice. All mentioned trademarks are protected by law. For technical details and detailed procedures of the specifications provided please contact your Tecan representative. This may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

Windows is a registered trademark of Microsoft Corporation in the United States and/or other countries.

© 2014, Tecan Trading AG, Switzerland, all rights reserved.

[www.tecan.com/components](http://www.tecan.com/components)