

Cavro® Ceramic Valves



- Compatible with a broad range of organic and inorganic solvents
- Resistant to abrasion
- Extended life
- Drop-in replacement for standard plastic valves


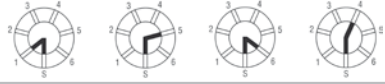
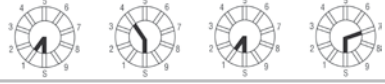
Tecan Systems now offers a line of Cavro® Ceramic Valves for the XCalibur™, XLP 6000™ and Smart Valve™ product lines. An innovative design (patent pending) eliminates the need for port seals, making these valves the first in the market with a 100% ceramic flow path. The ceramic material is 99.5% alumina which is 8 times harder than stainless steel and is compatible with a broad range of

organic and inorganic solvents. These unique characteristics make the new line of ceramic valves the ideal choice for applications requiring chemical inertness and/or resistance to abrasion.

The initial investment in ceramic valves is small given the significant long-term benefits they provide. The durability of the new valve reduces preventive maintenance requirements and instrument downtime due to planned or unplanned service.

Tecan offers an extra level of security by extending a 2-year warranty on ceramic valves against faulty material and/or workmanship.

Cavro® Ceramic Valve Specifications

Valve Configurations*	3-Port Distribution	
	6-Port Distribution	
	9-Port Distribution	
Port Fittings	1/4-28	
Nominal Port Diameter	1 mm (0.039 in)	
Port Depth	7.5 mm (0.295 in)	
Operating Pressure	4.14 bar (60 psi)	
Operating Temperature	15° to 40°C (59°F to 104°F)	
Storage Temperature	-20° to 65°C (-4°F to 149°F)	
Wetted Path	Ceramic (99.5% Alumina)	
Platform Compatibility	XCalibur, XLP 6000, Smart Valve Plus	

Ordering Information

Available Configurations*	Part Number
Ceramic Valves	
Valve 3+ 1/4-28 Ceramic	20737496
Valve 6+ 1/4-28 Ceramic	20737492
Valve 9+ 1/4-28 Ceramic	20737494
XCalibur	
XC 3+ 1/4-28 232/485 Ceramic	20740556
XC 3+ 1/4-28 CAN/485 Ceramic	20740558
XC 6+ 1/4-28 232/485 Ceramic	20740560
XC 6+ 1/4-28 CAN/485 Ceramic	20740562
XC 9+ 1/4-28 232/485 Ceramic	20740564
XC 9+ 1/4-28 CAN/485 Ceramic	20740566
XLP 6000 / XLP 3000	
XLP 3000 9+ 1/4-28 Ceramic	20740547
XLP 6000 3+ 1/4-28 Ceramic	20740549
XLP 6000 6+ 1/4-28 Ceramic	20740552
XLP 6000 9+ 1/4-28 Ceramic	20740554
Smart Valve Plus	
Smart Valve 3+ 1/4-28 Ceramic	20740570
Smart Valve 6+ 1/4-28 Ceramic	20740572
Smart Valve 9+ 1/4-28 Ceramic	20740574

* See your local Tecan representative for additional ceramic valve configurations

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication, 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 in this publication. Changes in this publication can be made at any time without notice. All mentioned tradeFor technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative.

394892 V1.0, 1-2007

© 2007, Tecan Trading AG, Switzerland, all rights reserved. Tecan® and Cavro® are in major countries registered trademarks of the Tecan Group Ltd., Männedorf, Switzerland. XCalibur™, XLP 6000™ and Smart Valve™ are trademarks of Tecan Group Ltd., Männedorf, Switzerland.

Austria +43 62 46 89 33 **Belgium** +32 15 42 13 19 **China** +86 10 5869 5936 **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 48 174 **Portugal** +351 21 000 82 16 **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 +41 44 922 8125

www.tecansystems.com