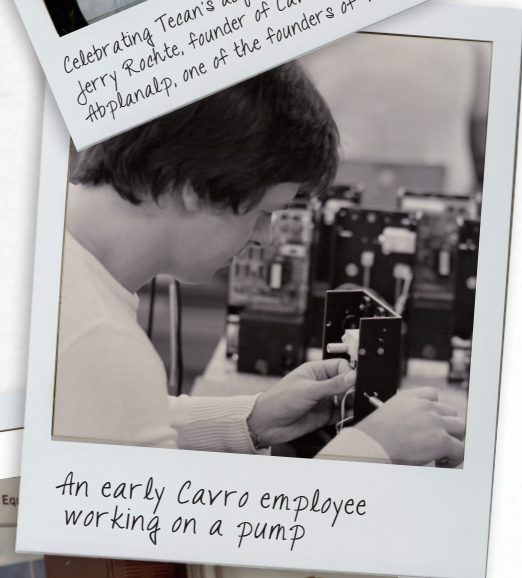


40 years of Cavro®

As we reach the milestone anniversary of 40 years of Cavro, Tecan Systems' Site Manager Nora Tryhorn takes stock and looks forward to a prosperous and successful future for Tecan's components business.



Celebrating Tecan's acquisition of Cavro in 1987. Jerry Roehle, founder of Cavro (left), and Heinz Abplanalp, one of the founders of Tecan



An early Cavro employee working on a pump



Jerry Roehle (left) and Bob Caveney at work in the early years of Cavro

Our business

Tecan's name is synonymous with excellence in the world of OEM components. For 40 years now Tecan Cavro brand components have provided innovative and reliable liquid handling for instrument designers for almost any application. Our outstanding Tecan Systems team of 104 employees is based in San Jose, in sunny California. Together, our vision is to be the best-in-class – a world class manufacturer and a design center of excellence for Tecan.

"I've worked in other companies and the people in Tecan are just different. Tecan Systems has really good employees who work hard, but also take the time to get to know each other."

Diana Morikawa, Director of Logistics and Production

Our products

Tecan Cavro components work robustly and efficiently in applications that are critical to society as a whole. We are very aware of the huge impact our products have on public health and safety.

"The products we make at Tecan touch so many lives. We are all very proud to work for a company that makes such great products."

Tawni Kestel, Purchasing Supervisor

Our customers

Talking to our customers during the entire development cycle of our products makes so much sense as we create exactly what is needed. We pride ourselves on our excellent communication channels and long-term relationships with our customers.

"I'm proud to be part of our service department, taking on all the responsibilities that are involved."

Bob Jackson, Service Technician



Early Cavro premises at 1234 Elko Drive, Sunnyvale, CA

Cavro milestone timeline

- 1972 Cavro Scientific Instruments founded
- 1980 Model 1500/2000 Pipettor/Dilutor
- 1985 IQ 190 Sample Processor
- 1986 SB 1200 Pump
- 1987 Tecan acquires Cavro
- 1989 RSP 9000
- 1990 XL 3000
- 1992 XL 300X
- 1994 XP 3000
- 1994 MSP 9000
- 1997 MiniWash Pump, Smart Valve, Smart Peristaltic Pump
- 1997 MSP 9500



Jerry Rochte (right) and a Cavro colleague working on a pump



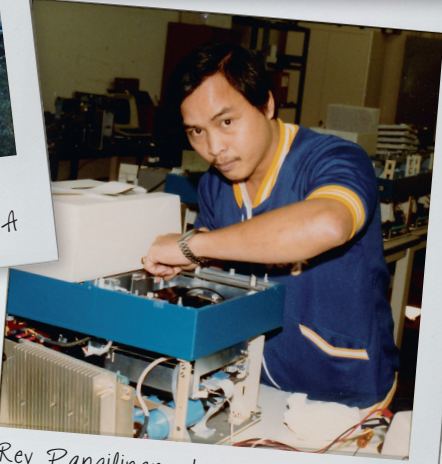
Nora Tryhorn in 1992 as Cavro's Customer Service Representative



The Tecan Systems' Sales Team. Left to right, Bruce Steffer (US), Jerry Maurice (US), Willi Krattenmacher (Europe), Chuck Fontana (US)



The manufacturing floor at Tecan Systems' previous Humboldt Court premises



Rey Pangilinan at work in 1984 - an early employee who is still working at the Company today!

Our quality

We've done a great job of improving our quality over the years and have seen significant improvements in our customer satisfaction levels since 2006. But there's

always room for improvement and we will continue to work hard as a team to make things better.

"My proudest achievement at Tecan Systems is the high level of productivity and quality accomplished by our relatively small R&D department."

Torleif Björnson, Director R&D

Our lean working principles

We started implementing lean manufacturing in 2009 and continue to strive for the utmost efficiency. But lean principles are based on a philosophy of constant improvement and we'll keep working to consistently find a better and faster way to do things.

"My proudest moment at Tecan was being the first person to build the new Cavro Omni Robot."

Kenny Tran, Track Assembler

Our people

For Tecan Systems, the secret of our success is undoubtedly our employees, several of whom have worked here for many years. We have a great team of extremely capable people who work hard and, most importantly, have a real passion for their jobs; everyone at Tecan Systems cares about the customer.

We asked our employees what's the best thing about working for Tecan Systems, and the answer was resounding....THE PEOPLE!!

"We are one big happy family, helping each other!"

Jose Rodriguez, Material Handler

To find out more about Tecan Cavro components, visit

tecان.com/components

