

30 years of Tecan

We are proud to have been true pioneers in the liquid handling market for three decades, and we really hope you value your Tecan as much as we love making them for you. As Tecan turns 30, here is how we celebrated across the world.



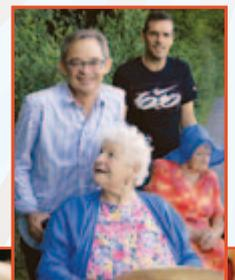
Happy Birthday Tecan! March 18th (1980)



30 years Tecan Party in Switzerland



30 years celebration in Japan with customers and Tecan employees



Tecan employees worldwide donate time to help charity organizations and assist on other occasions



Party on! 80's party with Tecan US



Salzburg 2010
Third Tecan
Symposium



For more information about 30 years of Tecan, visit www.tecan.com/history



Want to know more about our We love our Tecan campaign? Go to www.tecan.com/weloveourtecan



30 years celebrations in Shanghai with customers and Tecan employees