

Tecan US relocates to bigger and better premises

July this year saw Tecan US relocate from the Durham premises it had occupied for the past 21 years to a new, larger facility in the nearby Morrisville area of Raleigh. This new base offers scalable space and additional training rooms to meet the needs of continued growth over the coming years. With its high ceilings and clean lines, the new building has a very modern and sleek appearance, providing a pleasant environment for visitors and employees alike. Situated close to Raleigh-Durham International Airport, hotels, restaurants and shopping at Brier Creek, the new premises are convenient for both customers attending the Company's training facility and commuting employees.



Tecan US staff assembled outside the new premises



Cutting the ribbon at the opening ceremony

Careful planning enabled the relocation to be completed over the course of one weekend, as Peter Hornbach, Vice President of Operations, explained: "We closed slightly early on the Friday, and moved the IT portion of the business that evening. That was followed by 64 truckloads of equipment and, by mid afternoon on Saturday, we had actually moved everything in. We cleaned up on Sunday, and

were open for business as usual on Monday morning. Everyone was very motivated and worked together, and the move went really smoothly. Everyone is thrilled with the new base, and we are very proud of the premises."

To contact Tecan US, call +1 800 352 5128 or email sales-americas@tecan.com

Cornelias celebrates 25 years at Tecan

Celebrations were in order in June, as customer service representative Cornelias Rushing reached the landmark 25 years of service at Tecan US. Cornelias joined Tecan in 1987, and has seen many changes, as she explained: "When I joined Tecan, it had been working out of a garage at the home of a former President in Chapel Hill. I came on board shortly after the Company relocated and, since then, I have been involved in two more Company moves; it's been interesting, because each time Tecan goes up another

level. Before I joined the Company I was working in local government, so this was a really new area for me, seeing products such as the RSP 5000 and the MegaFlex. The product range has grown and improved over the 25 years I have been working here. When I started, the Company was just celebrating the acquisition of Cavro Scientific Instruments, and detection products were not even a part of Tecan. That's how far my days go back!"

"My actual anniversary date was Friday June 29th, and I had been told to make sure I was in the office on that day at a certain time, but nobody would tell me why. As it was my anniversary date, I just figured they would give me my long service award, I didn't expect anything extra. Instead, they surprised me with a beautiful cake and a




"...never in my wildest dreams did I imagine being here 25 years. It's been great, truly, a blessing."

reflections video, with pictures of me in different aspects over the years. Some of the photos were of events I had forgotten even happened, so it was really nice and I got a bit teary-eyed. The whole Company came together for the celebration and it was wonderful. I really, really appreciated it. I came here looking for a job that I could connect with, but never in my wildest dreams did I imagine being here 25 years. It's been great, truly, a blessing."