

Customers celebrate the launch of Tecan Australia



Tecan recently invited customers from around Australasia to mark the inauguration of Tecan Australia as a fully owned Tecan subsidiary, with a series of launch events in Sydney, Brisbane and Melbourne.

The creation of Tecan Australia reflects Tecan's uniquely strong commitment to the Australasian market, and will provide customers in Australia, New Zealand and surrounding territories with an even better service, thanks to a new facility in Melbourne and local presence in several major cities.

The launch events gave customers a rare opportunity to discuss their applications and automation issues with Tecan staff and their fellow scientists in a relaxed environment, and proved so successful that Tecan Australia will soon be establishing regular user group meetings to keep these valuable communication channels open.



The new Melbourne facility offers customers in Australia a local service



The launch parties brought together Tecan users from around Australasia



The 35th floor dining experience offered a good view of Melbourne



The Sydney Harbour Bridge offered a spectacular backdrop for the first event