

# Customers benefit as Tecan Australia becomes a subsidiary

Customers in Australia, New Zealand and the Pacific Islands have welcomed the news that in early 2012 Tecan Australia became a fully owned Tecan subsidiary. A new facility in Melbourne and local presence in several major cities reflects Tecan's uniquely strong commitment and continuing investment in the Australasian market. The new Melbourne base houses a comprehensive sales and support organization, complete with a workshop and additional facilities for customer training and demonstrations. Especially important for customers spread over such a wide geographical territory, the new headquarters includes additional warehouse facilities and stock levels to quickly support customers in need.



Tecan Australia also provides a local presence at scientific meetings and exhibitions

Roland Durner, General Manager for Tecan Australia, said: "This is great news for our customers and employees alike. Our new facility, backed by strong operating procedures and with Tecan's renowned emphasis on regulatory and quality requirements, will help us to improve every element of our business. Our existing team of highly trained sales and support staff will now be even better placed to provide a dedicated service to our customers in research, clinical diagnostics, forensics, and food and agriculture in the region."

To find contact information on Tecan Australia, visit [www.tecan.com.au](http://www.tecan.com.au)

To hear more about Roland's views on the challenges of the Australasian market, read this month's Leading the debate on page 27.

## Meet the team



Roland Durner



Luke Danielewski



Adam Gangur



James Hunter



Biju Kishor



John Ince



Clair Joyce



Josh Ardley



Michael Zhen

