

China and Asia Pacific operations expanded

Tecan has opened its new Asia Pacific region headquarters, based in the heart of Shanghai's pharmaceutical and life science center. The launch of the regional headquarters reflects the Company's significant growth and strong commitment to the market in China, as well as the whole Asia Pacific region, and will improve service and support for customers and distribution partners alike.

The new office was opened by Thomas Bachmann, CEO of the Tecan Group, on the 28th October, 2008. The ceremony was attended by key customers from around the Asia Pacific region, major distributors, officials and representatives from key Chinese Provinces, the Swiss and Austrian Embassies in China and the Tecan Group. Guests were welcomed to the Shanghai office by the CEO and were treated to a traditional Chinese ceremonial inauguration of the facility, followed by a celebratory dinner.

Mr Bachmann said: "We are delighted to be making this positive move into what is almost certainly one of the fastest growing regions in life sciences today. With a stronger presence and new facilities, we will be able to improve sales coverage, service and support to all of our customers and partners in the entire Asia Pacific region."



Thomas Bachmann thanks Peter Grünig for building up Tecan (Shanghai) Trading Co., Ltd.



Dinner with our Chinese guests



Kirsti Tavast (Head of Market Unit Asia Pacific) and Shinya Dobashi (new General Manager Tecan Japan Co.) in front of the Asia Pacific map

