

Tecan expands operations in China and Asia Pacific

- **Region headquarters opened in Shanghai**

Männedorf, Switzerland, October 29, 2008 – Tecan (SWX: TECN) 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, at a ceremony 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.

Under the new terms, the new Asia Pacific headquarters in Shanghai will now sell directly to many of the major multinational biotechnology, pharmaceutical and diagnostic companies flourishing in this rapidly developing area. Tecan's long-established Beijing operations will maintain close regional relationships with its customers in China. Similarly, a crucial well-established network of distributors will continue to cover the Asia Pacific region as before, including the important after-sales customer support. Tecan will also be able to support its distributors in these efforts by providing training, warehousing of spare parts and improved services out of Shanghai.

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."

About Tecan

Tecan (www.tecan.com) is a leading global provider of laboratory instruments and solutions in biopharmaceuticals, forensics, and clinical diagnostics. The company specializes in the development, production and distribution of automation solutions for laboratories in the life sciences sector. With its subsidiary REMP (www.remp.com), Tecan is the market leader in automated laboratory storage and logistics systems. Its clients include pharmaceutical and biotechnology companies, university research departments and diagnostic laboratories. Founded in Switzerland in



2

1980, the company has manufacturing, research and development sites in both North America and Europe and maintains a sales and service network in 52 countries. In 2007, Tecan generated sales of CHF 414.4 million (USD 345.3 million; EUR 252.7 million). Registered shares of Tecan Group are traded on the SWX Swiss Exchange (TK: TECN/Reuters: TECZn.S/Swiss security number: 1210019).

For further information:

Tecan Group

Dr. Rudolf Eugster
Chief Financial Officer

investor@tecan.com
www.tecan.com

Martin Brändle
Head of Corporate Communications &
Investor Relations
Tel. +41 (0) 44 922 84 30
Fax +41 (0) 44 922 88 89