

Press Release

Tecan Schedules Meeting and Conference Call on March 4th to Discuss Full Year 2008 Financial Results

Männedorf, Switzerland, February 25, 2009 – Tecan Group AG (SIX Swiss Exchange: TECN) will hold an analyst and press meeting in Zurich, Switzerland, to discuss the 2008 annual results on Wednesday, March 4, 2009 at 10:00 AM CET. The results will be published on March 4 at 7:00 AM.

The meeting will also be accessible via telephone or webcast. The live webcast, including synchronized slides, will be available on the Tecan web site at www.tecan.com. A link to the webcast will be provided immediately prior to the event. A replay will be available via the web site following the live event.

The call will be conducted in English.

The dial-in numbers for the call are as follows:

Participants from Europe: +41 91 610 5600 or +44 207 107 0611 (UK)

Participants from the U.S.: +1 (1) 866 291 4166

Please dial in 15 minutes before the beginning of the meeting.

The replay of the conference call will be available for 24 hours, starting one hour after the live event.

The dial-in numbers for the replay are as follows:

For Europe: +41 (0)91 612 43 30 For the UK: +44 (0)207 108 62 33 For the US: +1 (1) 866 416 25 58

The code for the replay is 17826.

About Tecan

Tecan (www.tecan.com) is a leading global supplier of laboratory instruments and solutions for the biopharma, forensic and diagnostic industries. The company specializes in the development, production and distribution of automation solutions for life science laboratories. Through its REMP subsidiary (www.remp.com), Tecan is the premier supplier of automated laboratory storage and logistics systems. Tecan's clients include pharmaceutical and biotech companies, university research departments, diagnostic laboratories and companies as well as forensics laboratories. Founded in Switzerland in 1980, the company has production, research and development sites in both North America and Europe and maintains a sales and service network in 52 countries. In 2007, Tecan achieved sales of CHF 414.4 million (USD 345.3 million; EUR 252.7 million). Registered shares of the Tecan Group are traded on the SIX Swiss Exchange (TK: TECN/Reuters: TECZn.S/Valor: 1210019).



2

For further information, please contact:

Tecan Group

Dr. Rudolf Eugster Chief Financial Officer

investor@tecan.com www.tecan.com Martin Braendle Head of Corporate Communications & Investor Relations Tel. +41 (0) 44 922 84 30 Fax +41 (0) 44 922 88 89