

Tecan announces changes in management team

Männedorf (Switzerland), 31 March 2008 — Bernhard Iseli, a Member of the Executive Committee and head of the Business Segment Sample Management has decided to resign from his position effective June 30th 2008 in order to focus on new challenges outside of Tecan.

Bernhard Iseli has been with REMP since 2002 which was acquired by Tecan in the summer of 2005. REMP today encompasses the Business Segment Sample Management within the Tecan Group. Tecan's Board of Directors and the Management regret Mr. Iseli's decision and thank him for his significant contributions to both, the Business Segment and the Group.

A search for a successor is currently underway.

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 SWX Swiss Exchange (TK: TECN/Reuters: TECZn.S/Valor: 1210019).

For further information, please contact:

Tecan Group

Christopher Hanan
Head Business Development & Corporate Communications

Tel. +41 (0) 44 922 84 30

Fax +41 (0) 44 922 88 89

www.tecan.com