

Tecan Appoints Global Head of Human Resources

- **Markus Schmid appointed Head of Corporate Human Resources & Internal Communications and member of the Management Board**

Männedorf, Switzerland, January 25, 2011 – The Board of Directors of Tecan Group (SIX Swiss Exchange: TECN) has appointed Markus Schmid as Head of Corporate Human Resources & Internal Communications and a member of the Management Board. He will assume the position at Tecan by September 1, 2011 at the latest. Schmid has been in charge of corporate human resources on an ad interim basis as an external consultant since October 2010.

Tecan CEO Thomas Bachmann commented: “We’re delighted to welcome Markus Schmid as our new Head of Corporate Human Resources & Internal Communications. With his extensive management and consulting experience in organizational and personnel development, he will make a positive and lasting impact on the implementation of the new organization introduced at the beginning of this year.”

Markus Schmid trained as a teacher and also studied psychology and journalism at the University of Freiburg, with a focus on occupational and organizational psychology. He has worked as a teacher and instructor at various educational levels and has held various consulting positions. In 1998 he became a partner and operations manager at MANRES AG, a consultancy that focuses on organizational development and transformation. There, he specialized in areas including executive and team coaching and organizational development consulting for prominent companies and public institutions. Markus Schmid is 42 years old and a Swiss national.

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 instruments and automated workflow solutions for laboratories in the life sciences sector. Its clients include pharmaceutical and biotechnology companies, university research departments, forensic and diagnostic laboratories. As an original equipment manufacturer, Tecan is also a leader in developing and manufacturing OEM instruments and components that are then distributed by partner companies. Founded in Switzerland in 1980, the company has manufacturing, research and development sites in both Europe and North America and maintains a sales and service network in 52 countries. In 2009, Tecan generated sales of CHF 392 million (USD 361.2 million; EUR 259.6 million). Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TK: TECN/Reuters: TECZn.S/ ISIN CH0012100191).

For further information:

Tecan Group

Martin Braendle

Head of Corporate Communications & Investor Relations

Tel. +41 (0) 44 922 84 30

Fax +41 (0) 44 922 88 89

investor@tecan.com

www.tecan.com