

**Analyst and Media Conference  
Full-Year Financial Results 2013  
Tecan Group AG  
Zurich, 11 March 2014**

SIX Swiss Exchange  
Convention Point, Auditorium  
Selnaustrasse 30  
8021 Zurich

- Yes, I will attend the conference in Zurich
- I will stay for the Aperitif
- I will participate in the webcast / conference call
- No, I am not able to participate

Name: \_\_\_\_\_

Surname: \_\_\_\_\_

Company: \_\_\_\_\_

Address: \_\_\_\_\_

Phone: \_\_\_\_\_

Fax: \_\_\_\_\_

Email: \_\_\_\_\_

**Please send your response by email or fax no later than February 27, 2014 to:**

Martin Braendle, [investor@tecan.com](mailto:investor@tecan.com) or Fax +41 (0)44 922 88 89