



# Invitation.

## Capital Markets Day

Tecan Group AG  
Maennedorf, 6 June 2019

Tecan cordially invites analysts, institutional investors and media to its Capital Markets Day on 6 June 2019.

The event will be hosted by Dr. Achim von Leoprechting, CEO, along with other members of the management.

The 21st century has often been described as the Century of Biology. With an end-user as well as an OEM business, Tecan has a unique position to bridge from research to diagnostics settings and thereby help customers with the discovery of novel medications all the way to better diagnosis of diseases.

The day will provide an update on how Tecan technology is empowering key application areas, especially in Genomics, Cell Biology and Mass Spectrometry.

The focus in the afternoon will lie on product demonstrations and a site tour.

The morning presentations will also be webcast live and may be accessed here:  
[www.tecan.com/investor-overview](http://www.tecan.com/investor-overview)

### PROGRAM OUTLINE

**08:15 - 08:45**

Registration and Welcome Coffee

**08:45 - 13:00**

CEO, Management and  
Customer Presentations  
Q&A

**13:00 - 14:00**

Buffet Lunch

**14:00 - 15:30**

Product Demos  
Site Tour

All presentations and product demos will be in English, the site tour will be held in English as well as in German.

Please confirm your participation no later than 21 May 2019 by returning the reply form.

We look forward to welcoming you to this exciting event.

Kind regards,



Martin Braendle  
VP Communications & Investor Relations  
T +41 44 922 8430, F +41 44 922 8889  
[investor@tecan.com](mailto:investor@tecan.com)  
[www.tecan.com](http://www.tecan.com)



# Reply form.

## Capital Markets Day

Tecan Group AG  
Maennedorf, 6 June 2019

Date & time:

**6 June 2019, 08:45 - 15:30**

**Registration opens at 08:15**

Location:

**Tecan Headquarters**, Seestrasse 103, 8708 Maennedorf

- Yes, I will attend the Capital Markets Day
- Yes, I will attend lunch     No, I will not be able to attend lunch
  - Site Tour in English
  - Site Tour in German

- No, I will not be able to attend the Capital Markets Day

Please send your registration no later than 21 May 2019 to:  
[investor@tecan.com](mailto:investor@tecan.com)

Last name:

---

First name:

---

Job title:

---

Company:

---

Country:

---

Email:

---

Tecan Group Ltd.,  
Seestrasse 103,  
CH-8708 Maennedorf,  
Switzerland  
T +41 44 922 84 30  
F +41 44 922 88 89  
[info@tecan.com](mailto:info@tecan.com)  
[www.tecan.com](http://www.tecan.com)

 **TECAN.**

