

NEXT GENERATION MICROARRAYS TOUR

LIVE TOUR
2016

LABEL-FREE BIOMOLECULAR INTERACTION & KINETIC ANALYSIS MADE EASY

Berthold Technologies and Biametrics are inviting you to participate in and enjoy the Next Generation Microarrays Tour 2016

This free scientific seminar will focus on applications for next generation microarray technology enabling label-free biomolecular interaction & kinetic analysis and offers you opportunities for networking with your peers.

AGENDA

- 13.30 – 13.50 “No compromises. High throughput label-free molecular interaction”, Dr. F. Schleifenbaum, Berthold Technologies
- 13.50 – 14.20 bScreen System live demonstration
- 14.20 – 14.45 Coffee Break
- 14.45 – 15.15 Local Guest Speaker
- 15.15 – 15.35 “Application of label-free microarrays in drug development and clinical research”, Dr. G. Proll, Biametrics
- 15.35 – 16.30 Discussion & Snacks

Find out more at:

<https://www.berthold.com/bscreen>

Register now:

https://www.berthold.com/en/Microarrays_Tour_2016

October 11th

TÜBINGEN

Guest Speaker
Prof. Dr. G. Gauglitz
University of Tübingen

October 25th

BASEL

Guest Speaker
Dr. G. Roth
University of Freiburg

November 08th

MÜNCHEN-FREISING

Guest Speaker
t.b.a.

November 22nd

POTSDAM

Guest Speaker
Prof. Dr. M. Niedrig
Robert Koch Institute