

Haus Petersplatz / Department of Biomedicine; University Basel invites to

# Advances in Stool Diagnostics

Seminar, 15. August 2013 • 11:00 – 15:30  
Hotel Glockenhof • Zurich, Switzerland

Gastroenteritis is a serious medical and economic burden. Traditional serial selective testing practices are often slow and result in low diagnostic detection. This seminar discusses the potential of new techniques in stool diagnostics and provides the opportunity to share experiences with new applications in Gastro-intestinal pathogen detection.

Lunch and seminar are free but seats are limited. Please register by sending an email to:  
[jdaam@luminexcorp.com](mailto:jdaam@luminexcorp.com)

## Agenda

11:00 – 11:10

Welcome

11:10 – 11:50

**Implementation of a new Multiplex Test  
in a routine laboratory**

Prof. H. H. Hirsch, University Basel

11:50 – 12:30

**Syndromic Approach in the Diagnosis of  
Infectious Gastroenteritis**

Dr. J. Daam, Luminex

12:30 – 13:30

Lunch Buffet

13:30 – 14:10

***C. Difficile* - News on an old acquaintance**

Prof. L. von Müller, University of Homburg

14:10 – 14:50

**Gastroenteritic infections by Noro-  
and Rotavirus**

Dr. D. Hoffmann, Tech. University Munich

14:50 – 15:30

**New developments in Gastro-Diagnostics**

Chad Brueck, Luminex

### Seminar location:

Hotel Glockenhof

Sihlstrasse 31

CH-8001 Zurich

Supported by: **Luminex**<sup>®</sup>