

BIOPHARMA SUMMIT

Shaping the
future together



14th November

IZB Martinsried | IZB Martinsried, 82152
Planegg/Martinsried, Germany

to register please follow:

www.thermofisher.com/biopharmasummit

- **Summit Agenda**

- 8:45am Arrival and Registration (Breakfast available)
- 9:15 Meeting welcome and overview
- 09:30 A multi-attribute method (MAM) for advanced QA/QC characterization
- 10:00 Automation of differential protease digestions with Trypsin and Chymotrypsin using SMART resins
- 10:30 Networking break
- 11:00 Taking charge variant & aggregate analysis from research into the routine environment
- 11:30 Structural Proteomics in BioPharma
- 12:00 ADC Characterzation on a single high-resolution accurate-mass platform
- 12:30 [Native MS for highly complex glycosylationanalysis of biopharmaceuticals](#)
[Professor Christian Huber, Christian Doppler Laboratory for Biosimilar Characterization](#)
- 12:30 Networking lunch