

# CUSTOMER TRAINING COURSES AND SEMINARS

Process Control

Tower-Sens LB 490

3<sup>rd</sup> Extension Module

2<sup>nd</sup> Extension Module

length



 **BERTHOLD**





## Training and seminars for process measurement technology

With the purchase of your Berthold measurement application you have acquired a high-tech product. Regardless of whether you are measuring dose rates, contamination or activities: The particular measuring methods always require exact adherence to all parameters involved in the process. This applies to assembly at the measuring point as well as to configuration and calibration.

## Our service for you

We are happy to conduct our product training and radiation protection courses at your premises. Feel free to contact us so that we can coordinate location, whether at ours or yours, topics for training and the date together. On request, we can help you find accommodation. Dates and more information about our courses and training can be found on our website.

## Practical product training

Our product trainings give you safety and certainty in your daily work and ensure that you can fully use the potential of your Berthold product. Experienced service technicians, as well as specialists from development and product management convey in-depth practical knowledge that is tailored to the applications and needs of the participants.

### Our course offers

1.5-day product training in small groups with 4 to max. 10 participants. For example, for DuoSeries LB 47x, UniProbe LB 490 or Micro-Polar LB 56x, for any application (level, density, concentration). More offers on request.

### Product training process (example)

- Radiation protection basics
- Measurement basics
- Handling / operation / calibration
- Overview of the various measurement functions
- Intensifying customer-specific applications
- Practical exercises on the device
- Discussion







## Contact us

Have we sparked your interest? We look forward to hearing from you. We will be happy to advise you noncommittally on our training courses and will gladly accept your registration for any of our courses

### Contact product training:

Gerlinde Thal  
gerlinde.thal@berthold.com  
Phone: +49 (0) 7081 177 231  
Fax: +49 (0) 7081 177 339

### Berthold Technologies GmbH & Co. KG

Calmbacher Straße 22 · 75323 Bad Wildbad · Germany  
+49 7081 1770 · [industry@berthold.com](mailto:industry@berthold.com) · [www.berthold.com](http://www.berthold.com)

