

Process Intensification in the World of Chromatography – Training Content

- Introduction to MCC theory & benefits
- Transferring a batch process to MCC: steps involved and the use of models
- Key points to consider for transfer to MCC - avoiding issues: common mistakes and experimental set-up (number of cycles, how to start and end the process, sampling strategy, scale up considerations...)
- Recipe writing on the BioSMB PD system: demonstration and hands-on
- Lab session: execution of a real MCC experiment on the BioSMB PD system