



Dr. Jürgen Hubbuch
Professor, Forschungszentrum Jülich (Germany)

Biography

- 10/90 – 01/97 M.Sc. / Diplom Ingenieur in Chemical Engineering, Faculty of Chemical Engineering University of Karlsruhe (TH), Karlsruhe, Germany
- 10/93 – 10/95 Master of Philosophy (M.Phil.), Heriot-Watt University, Edinburg, Scotland
- 03/97 – 11/00 Doctor of Philosophy (Ph.D.), BioCentrum, Center for Process Biotechnology, Prof. O.R.T. Thomas Technical University of Denmark, Lyngby, Denmark
- 01/01 – 12/02 Group leader - Downstream Processing, Institute of Enzyme Technology, Prof. M.-R. Kula, Heinrich-Heine University Düsseldorf, Düsseldorf, Deutschland
- 01/03 – 01/08 Group leader – Bioseparation, Institute of Biotechnology 2, Prof. C. Wandrey Forschungszentrum Jülich GmbH, Jülich, Deutschland
- 02/08 – Professor, Institute of Process Engineering in Life Sciences Section IV: Biomolecular Separation Engineering, University of Karlsruhe