



Dr. Gary Lye
Professor of Biochemical Engineering, University College of London (UK)

Biography

Gary Lye is Professor of Biochemical Engineering within the Advanced Centre for Biochemical Engineering at University College London (UCL). He received his PhD in Biotechnology from the University of Reading in 1992. Between 1993 and 1996 he was successively a Research Fellow and then Lecturer in Chemical Engineering at Imperial College London and the University of Edinburgh. He joined UCL in 1996. He is currently Chair of the Biochemical Engineering Subject Group of the UK Institution of Chemical Engineers (IChemE). He has broad research interests on the application of microscale and automation techniques to rapid bioprocess design, optimisation and scale-up. He currently leads the multidisciplinary Bioconversion-Chemistry-Engineering Interface (BiCE) programme at UCL which is collaborative with 13 leading international companies.