



Dr. Jean-Philippe Stephan
Research Scientist Advanced Automation Technologies, Genentech (USA)

Biography

Jean-Philippe Stephan, originally from Rennes, Brittany, France, received his Ph.D. in Developmental Biology from Pierre and Marie Curie University (Paris VI, France). After a postdoctoral fellowship in Genentech Protein Chemistry Department with Dr. Jennie Mather, he was hired as a Research Scientist in the Assay and Automation Technology department (AAT) at Genentech, Inc. Since then Dr. Stephan has had a long-term interest in the development, evaluation and implementation of new technologies that could be applied to the discovery and characterization of therapeutic antibodies. As Research Senior Scientist in AAT, Dr. Stephan supervise a group that support multiple projects from early Research to early Development and manage an automated cell culture lab using the Tecan Cellerity platform.