



Dr. Lin Zhang
Associate Research Fellow, Bioprocess R&D, Pfizer (USA)

Biography

Lin Zhang is an Associate Research Fellow in Pfizer's Global biologic group based in St. Louis, MO. Since joining Pfizer in 2002, Lin leads a lab responsible for the development, optimization and characterization of cell lines used for the production of biotechnology-derived therapeutic drug candidates. Lin's group also has responsibility for process development activities focusing on the evaluation and implementation of novel expression platform as well as enabling technologies for cell line development. Prior to Pfizer, Lin was a senior scientist at Sigma-Aldrich, where she led a lab that had a direct responsibility for the conception and development of products in the areas of protein expression and bioassay development from ground up. Lin received her Ph.D. in medicinal chemistry from Purdue University, West Lafayette, IN and a B.S degree in medicine from Dr. Bethune medical school, China.