



**Dr. Lars J. Brandén**  
**Director Center for High Throughput Cell Biology, Yale University (USA)**

**Biography**

Lars J. Brandén, Ph.D., has recently moved to Yale University in New Haven, CT where he is establishing the Yale Laboratory for High Throughput Biology which he will be Director of. Prior to this move, Dr. Brandén was Associate Director for Chemical Genetics at the Judith P. Sulzberger Genome Center, Columbia University, in New York City, as well as Director of Automation for the Columbia Molecular Library Screening Center Network, in the Department of Biophysics and Cell Biology. Dr. Brandén received his undergraduate degree in molecular genetics from the University of Lund, Sweden, and completed his doctoral studies in cell and molecular biology at the Karolinska Institute, Stockholm, Sweden, under the guidance of Professor C.I. Edvard Smith. He pursued a postdoctoral research fellowship in the Laboratory of Cellular Biochemistry & Biophysics at the Sloan-Kettering Institute, Memorial Sloan-Kettering Cancer Center, New York City, under the direction of James E. Rothman, Ph.D.