



Dr. Wei-Shou Hu
Professor, University of Minnesota (USA)

Biography

Dr. Hu's research in cell culture technology deals with conventional and macroporous microcarriers, suspension culture, and cell-recycling systems. His recent work in cell culture involves the metabolic control of cell's physiological state and genome wide gene expression analysis of their regulation. He co-authored the textbook Bioseparations. He initiated the Engineering Foundation Conferences on Cell Culture Engineering more than a decade ago which has been among the most important forum of cell culture processing. Recent research activities of Dr. Hu also encompass metabolic engineering of antibiotic producing streptomycetes, and tissue engineering employing stem cells.