



Dr. Michael West
CEO, BioTime (US)

Biography

Dr. West is the Chief Executive Officer of BioTime, Inc. and Embryome Sciences, Inc. of Emeryville, California and Adjunct Professor of Bioengineering at the University of California, Berkeley. BioTime, Inc. (OTCBB: BTIM) and its wholly-owned subsidiary Embryome Sciences are companies focused on human embryonic stem cell technologies. He received his Ph.D. from Baylor College of Medicine in 1989 concentrating on the biology of cellular aging. He has focused his academic and business career on the application of developmental biology to the age-related degenerative disease. He was the Founder of Geron Corporation of Menlo Park, California (Nasdaq: GERN) and from 1990 to 1998 he was a Director, and Vice President, where he initiated and managed programs in telomerase diagnostics, oligonucleotide-based telomerase inhibition as anti-tumor therapy, and the cloning and use of telomerase in telomerase-mediated therapy wherein telomerase is utilized to immortalize human cells. From 1995 to 1998 he organized and managed the research between Geron and its academic collaborators James Thomson and John Gearhart that led to the first isolation of human embryonic stem and human embryonic germ cells. From 1998 to 2007 he was President and Chief Scientific Officer at Advanced Cell Technology, Inc. (OTCBB: ACTC) where he managed programs in human somatic cell nuclear transfer, cell differentiation, and ACTCellerate, a technology for the multiplex derivation and characterization of human embryonic progenitor cell lines.