

Curriculum Vitae | Susan Dexter

Member of the Board of Directors of Sartorius Stedim Biotech S.A.
Independent Director

Year of Birth: 1955

Nationality: US American

First appointment to the Board of Directors: April 2015

End of term: 2027 Shareholders' Meeting

Professional Profile

Since 2020	Sonnet Biotherapeutics, Inc., United States, Chief Technical Officer Non-clinical CMC Supply Chain. Responsible for product development for Sonnet's pipeline of biotherapeutic cytokine assets for treatment of solid tumor cancers
2008-2020	Latham Biopharm Group, United States, Managing Director; Due Diligence, Acting VP Business Development for multiple CMO's offering contract manufacturing services to the biotechnology life sciences industry, strategic consulting, single-use disposable technology implementation, project management and high-level business development and marketing, Advisor and speak for BioProcess International, Outsourced Pharma
2004-2008	Xcellerex, Inc. (now GE Healthcare), United States, Chief Business Officer; CMO services using fully integrated single-use bioprocessing technology, sales of single use bioprocessing technologies
1998-2004	Collaborative BioAlliance, Dow Chemical Company (Dow Biotechnology Contract Manufacturing Services), United States, Vice President, business development for microbial fermentation services, technologies and implementation of single use bioprocessing technologies
1986-1998	Celltech Biologics, Lonza Biologics, United States, business development, bioprocessing and manufacturing of biotechnology based biotherapeutics
1980-1986	Collaborative Research, United States, biotechnology sales in emerging markets for bioprocessing supplements and raw materials for biomanufacturing
1975-1980	University of Massachusetts Medical School, United States, research, mammalian cell culture, animal toxicology studies, basic research

Educational Background

- Degrees and Certifications: BS in Immunology and Marketing (double major, honors), American University, Washington, D.C., United States
- Harvard University Negotiation Course for Lawyers, Harvard University, Cambridge, Massachusetts, United States
- Finance for non-financial Managers, Harvard University through Dow Chemical Company internal training program

Other current directorships and positions outside the Group

- Member of the Board of Directors of ViroCell Biologics Ltd., United States
- Member of the Board of Directors of Virica Biotech Inc., Canada