

Curriculum Vitae | Prof. David Raymond Ebsworth, Ph.D.

Member of the Supervisory Board of Sartorius AG

Year of birth: 1954

Nationality: British

Place of residence: Overath, Germany

Member since: January 2020

Professional Profile

- | | |
|------------|--|
| Since 2015 | Provision of independent consultancy services to selected healthcare companies and financial investors, as well as angel investor in selected healthcare companies |
| 2009–2015 | Galenica AG, Switzerland

2014–2015: Project support for CEO of Vifor Pharma and CEO of Galenica Santé until retirement

2012–2014: CEO and Chairman of the Corporate Executive Committee (Zurich and Bern, Switzerland)

2009–2014: CEO of Vifor Pharma AG and member of the Corporate Executive Committee of Galenica AG |
| 2003–2009 | Various positions as non-executive director and consultant to the industry and financial investors |
| 2002–2003 | Oxford Glycosciences Ltd, U.K., CEO |
| 1983–2002 | Bayer AG, Germany

2000–2002: Business Unit Manager, Pharmaceuticals

1995–1999: President, North American Pharmaceutical Division

1983–1995: Various positions (focal areas: Marketing, Sales, Business Operations) |

Educational Background

- Doctor of Humane Letters (2000), Honorary Degree, University of New Haven, Connecticut, USA
- Ph.D. in Comparative Industrial Relations (1980), University of Surrey, Guildford, U.K.
- B.Sc. in Chemistry and German (1976), University of Surrey, Guildford, U.K.

Membership in Other Supervisory Boards Mandated by German Law

- Synlab AG, Munich, Germany (stock exchange listed) – Chairman of the Supervisory Board

Membership in Comparable Supervisory Committees of Business Enterprises Within Germany and Abroad

- Verona Pharma plc, Großbritannien (stock exchange listed) – Chairman of the Board of Directors (non-executive)
- Actimed Therapeutics Ltd, U.K. – Chairman of the Board of Directors (non-executive)
- Kyowa Kirin International plc, U.K. – Chairman of the Board of Directors (non-executive)
- Opterion Health AG, Switzerland – Chairman of the Board “Verwaltungsrat” (non-executive)

Other Major Activities Beyond Supervisory Board Position

- Visiting Professor at the Business School of the University of Surrey, Guildford, U.K.