

## Member of the Supervisory Board of Sartorius AG

Prof. David Ebsworth, PhD, resident of Overath, Germany

Born in 1954

### Professional Profile:

- Since 2015 Provision of consultancy services to selected healthcare companies and financial investors, as well as angel investor in selected healthcare companies
- Since 2018: Pharma Investments SA, Luxembourg, Member of the Board "Conseil d'administration" (non-executive)
- Since 2018: Actimed Therapeutics Ltd, U.K., Chairman of the Board of Directors (non-executive)
- Since 2017: Kyowa Kirin International plc, U.K., Member of the Board of Directors (non-executive)
- Since 2016: Opterion Health AG, Switzerland, Chairman of the Board "Verwaltungsrat" (non-executive)
- Since 2016: Interpharma Investments Ltd, British Virgin Islands, Member of the Board of Directors (non-executive)
- Since 2016: Business School der University of Surrey, Guildford, U.K., Visiting Professor
- Since 2014 Verona Pharma plc, U.K. (stock market listed) - Chairman of the Board of Directors (non-executive)
- 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: President and General Manager, Pharmaceuticals
- 1995–1999: President, North American Pharmaceutical Division
- 1983–1995: Various positions (focal areas: Marketing, Sales, Business Operations)

#### **Educational Background:**

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

#### **Mandate i.S.d. § 125 Abs. 1 Satz 5 AktG**

- Verona Pharma plc, Großbritannien (börsennotiert) – Chairman of the Board of Directors (non-executive)
- Pharma Investments SA, Luxembourg – member of the board “Conseil d’administration” (non-executive)
- Actimed Therapeutics Ltd, U.K. – Chairman of the Board of Directors (non-executive)
- Kyowa Kirin International plc, U.K. – member of the Board of Directors (non-executive)
- Opterion Health AG, Switzerland – Chairman of the Board “Verwaltungsrat” (non-executive)
- Interpharma Investments Ltd, British Virgin Islands – member of the Board of Directors (non-executive)

**Major Activities Beyond Supervisory Board Position:**

- Consultant for various healthcare companies and financial investors
- Visiting Professor, Business School of the University of Surrey, Guildford, U.K.
- Positions held within the meaning of § 125, Subsection 1, sentence 5, of the German Stock Corporation Law (AktG)