



**Dr. Guido Ströhlein, MBA**  
**CEO, ChromaCon AG (Switzerland)**

### **Biography**

Dr. Guido Ströhlein, MBA, holds the position of a chief executive officer at ChromaCon AG, a start-up company which brings new chromatographic processes (MCSGP-technology) for the purification of proteins and peptides to the market.

He received a MBA in Management, Technology and Economics from the ETH Zuerich, Switzerland. Guido earned his PhD under the guidance of Prof. Morbidelli at the Institute for Chemical- and Bioengineering, ETH Zurich, in the area of preparative bio-chromatography. He studied chemical engineering in Germany, England and the USA and gained industrial experience at DOW Chemical and Bayer in the area of bio-purifications and preparative chromatography.