



Free for publication from 01.04.2021. If reprinted, please send us a copy.

New General Manager at Promega GmbH

Hans-Joachim Müller takes over business from April 1, 2021

Walldorf, Germany (01.04.2021). On April 01, 2021, Hans-Joachim Müller will take over as General Manager of Promega GmbH, replacing Peter Quick, who is retiring after more than 24 years with Promega GmbH and previous responsibilities at Johnson & Johnson Clinical Diagnostics, Kodak Diagnostics and Amersham. Peter Quick led the business of Promega GmbH since its foundation in 1997 and continuously expanded the company in the Life Science Research (LSR) market in Germany, Austria, Poland and Central Eastern Europe.

Peter Quick will continue to represent Promega's interests in the Life Science Research department of the Diagnostics Industry Association and the VDPG as a whole.

Peter Quick explains: "As a character with charisma, Hanseatic businessman in the best sense and molecular biologist with extensive experience in the life science research market. Hans-Joachim Müller will look after the interests of Promega GmbH, their customers, partners and employees with pleasure and commitment. Mr. Müller and the Promega teams look forward with great anticipation to the opportunity to fill the strong new location in Walldorf with life and, above all, to meet Promega's customers as soon as the pandemic allows."

Hans-Joachim Müller builds on 15 years of experience as a sales and marketing manager in the life science research market and another eight years of successful management of a radiopharmaceutical company.

In his new role, the Hamburg native, who has lived in Heidelberg since 1993, plans to expand Promega's market share in its key fields of LSR and in vitro diagnostics (IVD) by marketing its high-quality products. According to him, this is achieved primarily through an ambitious and motivated Promega team. "Employee satisfaction thus continues to be a high priority. Customers and partners also should have the desire and prefer to work with us."

Therefore, the focus will initially be on getting to know the team and looking at the market in more detail.

About Promega Corporation

Promega Corporation is a leader in providing innovative solutions and technical support to the life sciences industry. The company's product portfolio of over 4,000 products support a range of life science work across areas such as cell biology; DNA, RNA and protein analysis; drug development; human identification and molecular diagnostics. For over 40 years these tools and technologies have grown in their application and are used today by scientists and technicians in labs for academic and government research, forensics, pharmaceuticals, clinical diagnostics and agricultural and environmental testing. Promega is headquartered in Madison, WI, USA with branches in 16 countries and over 50 global distributors. For more information, visit www.promega.com.