

Dako Omnis helping to accelerate cancer diagnostics



Dr Hermann Herbst

The Dako Omnis platform – developed and manufactured by Tecan – offers complete automation of both immunohistochemistry (IHC) and *in situ* hybridization (ISH) protocols. Following the instrument's launch in September 2013, this advanced staining platform is now helping clinical pathology

laboratories around the world to streamline their histology workflows, offering higher throughput and faster results for patients.

Featuring a wide range of functions designed to improve and simplify day-to-day operation – such as flexible loading, overnight running and fast turnaround times – the Dako Omnis is the result of combining Dako's expertise in tissue-based cancer diagnostics with Tecan's advanced technical knowledge. Britt Meelby Jensen, Dako General Manager and Agilent Vice President, commented: "Tecan's extensive experience in the development of liquid handling technologies and in-depth understanding of clinical laboratory

customer needs made the Company an obvious partner for the development of this advanced staining system."

To celebrate the success of Dako Omnis, Tecan invited Dr Hermann Herbst, Head of Pathology at Vivantes in Berlin, Germany – one of the first customers to receive their platform – to a special presentation on the shores of Lake Zurich. Dr Herbst described the changes to his laboratory's workflow since the system's introduction to an audience of Tecan employees and associates, highlighting the time savings and efficiency gains. Dr Herbst explained: "The Dako Omnis really is a state-of-the-art, fully automated IHC and ISH workstation, allowing us to run in either continuous or batch processing modes. It offers improved control over reagent use and staining and, by combining IHC and ISH on a single platform, it saves us considerable hands-on time in both processes."

To learn more about the Dako Omnis platform, visit www.dako.com/omnis

