

Sias joins Tecan Partnering

Sias – a leading OEM supplier of laboratory automation solutions – has joined the Tecan Group, adding complementary new platforms, capabilities and expertise to customers of Tecan Partnering. A dedicated OEM supplier, Sias is committed to the highest quality standards and building close relationships with its customers to ensure projects are delivered in the shortest possible time.

sias
Empowered with Tecan ●

Complete automation of laboratory workflows has become increasingly important over the last decade, particularly in the *in vitro* diagnostics and contract research sectors. The accuracy, reproducibility and comprehensive audit trails offered by modern laboratory automation solutions provide the high level of process security necessary to meet stringent quality control and regulatory standards.

Developing automation systems capable of delivering these features is both time consuming and complex, requiring extensive liquid handling expertise to ensure reliable, error-free processing. To overcome this issue, many leading providers choose to work with a knowledgeable OEM laboratory automation partner – such as Tecan Partnering – to reduce both the time and cost of developing new products. Founded in 2000, Sias has grown into a leading supplier in the sector, combining a flexible approach with open and efficient communication to establish itself a reputation for exceptional customer satisfaction. The common history of a customer-focused approach, together with its close proximity to Tecan's Swiss facilities, make Sias an excellent fit with Tecan's existing Partnering portfolio, allowing close cooperation

and providing a healthy funnel for new development projects. Dr Achim von Leoprechting, Head of Tecan Partnering, explained: "Sias' core expertise in automating immunoassays and molecular diagnostics workflows – as well as its strong client base in China – perfectly match our corporate strategy.

“ We are pleased to have reached this agreement and look forward to welcoming the Sias team into Tecan. Sias will further extend our leading Partnering portfolio, providing access to complementary automation platforms which are particularly well suited to small to medium-sized OEM systems for diagnostic and research applications. ”

In addition, the close proximity of the two Swiss facilities will allow synergies to be implemented.”

The company's modular platforms and robotic components are ideally suited for low- to medium-throughput applications, further extending Tecan's capabilities into new markets and applications. Sias' existing customers will also be able to benefit from Tecan's global sales infrastructure and after-sales support network. Andre Guenter, CEO of Sias, commented: "This agreement will be mutually beneficial for both companies and for our customers. Sias has a

dedicated, highly trained team, innovative technologies and leading products, as well as a strong pipeline of new development projects and dedicated business partners all over the world. Tecan's global presence will allow us to grow even faster, leveraging local sales and service organizations to the advantage of our customers. We will also greatly benefit from the combined capabilities of a larger development organization.”

Tecan CEO Dr David Martyr added: "We are pleased to have reached this agreement and look forward to welcoming the Sias team into Tecan. Sias will further extend our leading Partnering portfolio, providing access to complementary automation

platforms which are particularly well suited to small to medium-sized OEM systems for diagnostic and research applications.”

To find out more about Sias, visit www.sias.biz

To find out more about Tecan Partnering, visit partnering.tecan.com

