

Tecan Synergence™

YOUR VISION EMPOWERED

COMPREHENSIVE OEM SERVICES FOR LAB AUTOMATION

Synergence is Tecan Partnering's unique approach to the development of OEM automated solutions for diagnostics and life sciences. Starting with your exact requirements, this service guides you seamlessly through your project, from product design and development to manufacturing, launch and post-launch activities.

Whatever your requirements of an OEM partner, we have the application knowledge and engineering expertise to turn concept into reality. With over 40 years of experience in laboratory automation and OEM instrument development, we have worked across the full spectrum of life sciences and diagnostics to take emerging technologies from research to routine testing. Our close relationship with our customers means that we understand their diverse workflows and applications, developing comprehensive, high quality solutions to help eliminate bottlenecks, improve productivity and ensure your product is a success in the market.

Excellence that comes from a truly synergistic partnership!

WHAT WE OFFER

Core offering

This flexible service includes all the 'elements' necessary to turn your vision for your OEM automation platform into a reality.

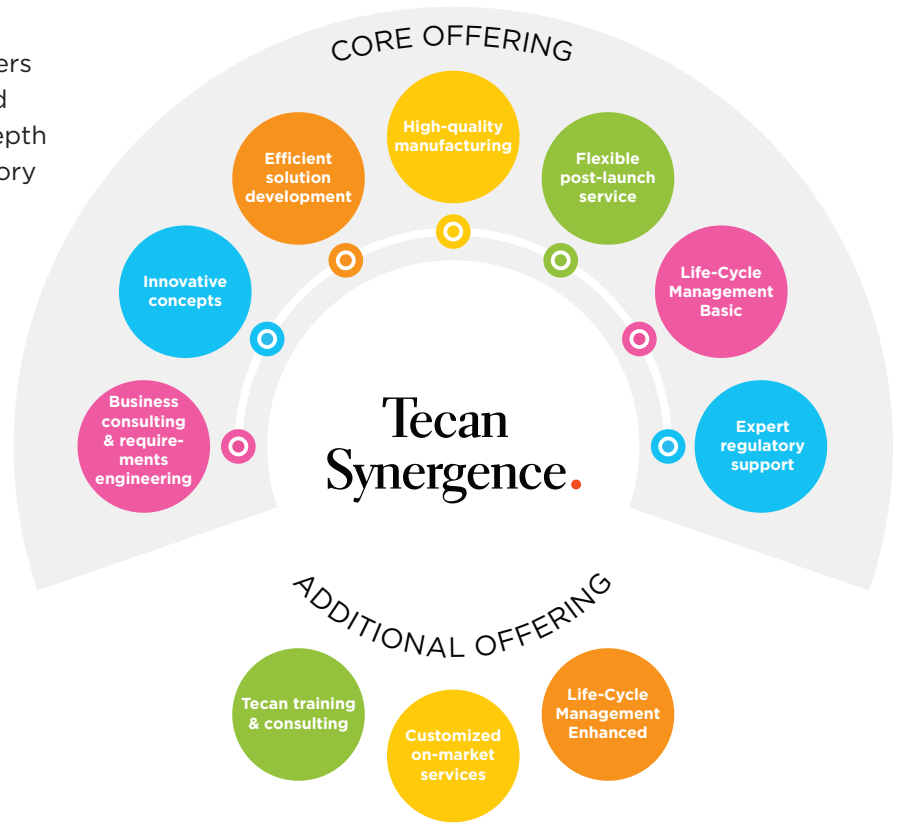
- Business consultancy and requirements determination
- Automation concept innovation
- Efficient solution development
- High quality manufacturing
- Flexible post-launch services
- Basic lifecycle management
- Expert regulatory support



Additional offerings

These supplementary services allow customers to complement their in-house know-how and capabilities by taking advantage of our in-depth market knowledge, global reach and regulatory expertise.

- Extended lifecycle management
- Instrument maintenance and repair
- Instrument refurbishment
- Product training
- Regulatory training and consultancy



DISCOVER THE TECAN PARTNERING ADVANTAGE

35+

OEM automated systems successfully launched to the IVD market worldwide

100+

audits completed by various major notified bodies worldwide

30+

IVD companies trust Synergence in bringing their vision to life

52

countries with Tecan Partnering representation

Our multidisciplinary team of scientists and engineers have extensive training experience across numerous life sciences and diagnostics market sectors, and can apply this expertise to applications in:

- Molecular diagnostics
- Immunoassays
- Mass spectrometry
- Histology and cytology
- Blood grouping/typing
- Biobanking
- Hematology
- Genomics
- Proteomics
- Drug discovery
- Cell biology
- Forensics
- Veterinary and agriculture
- Food sciences



LEARN MORE ABOUT OUR VALUED PARTNERSHIPS:

partnering.tecan.com/partners-and-case-studies

www.tecan.com/synergence

Australia +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170 **Italy** +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886 **Spain** +34 93 595 25 31 **Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +41 44 922 81 11

© 2019 Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit www.tecan.com