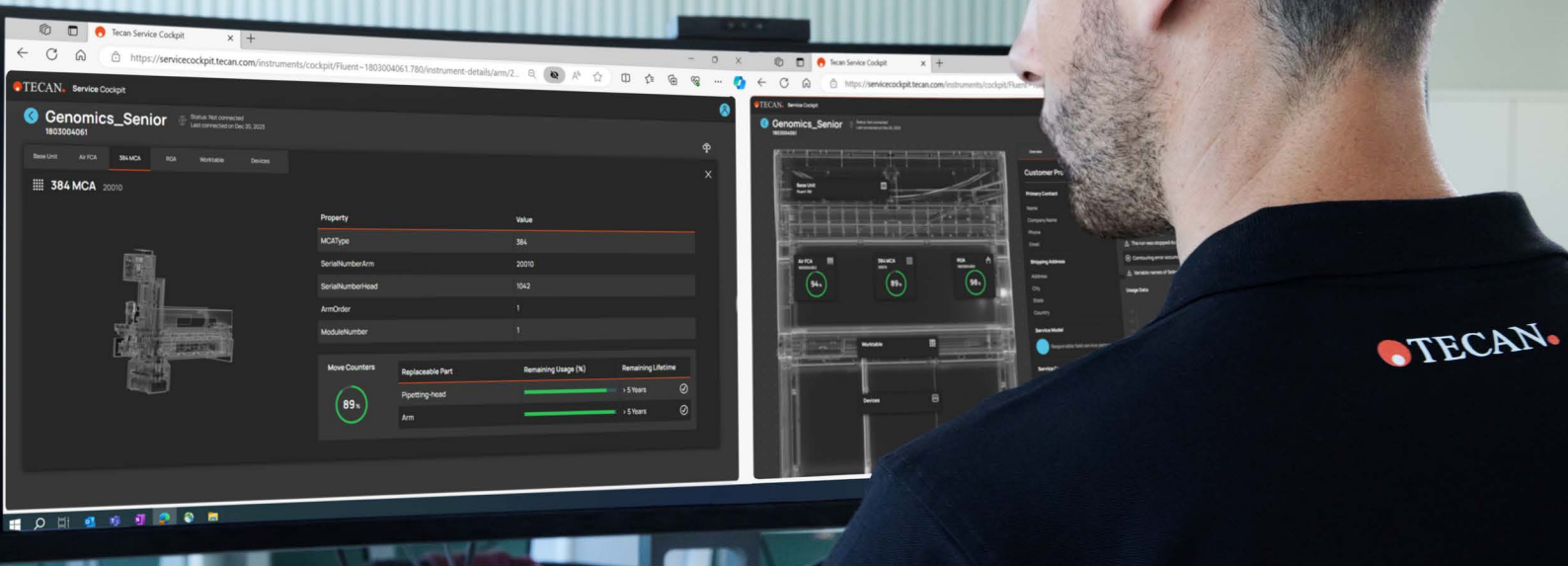


REMOTE DIAGNOSTICS FOR AN ENHANCED SERVICE EXPERIENCE.



Equipment downtime can have a major impact on your liquid handling instruments in terms of disruption and cost.

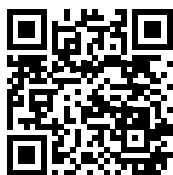
When connected to Introspect™, Tecan's digital productivity enhancement tool, you can unlock remote diagnosis and support with our behind-the-scenes hero Service Cockpit, Tecan's digital platform indicating the status of your instrument.

Service Cockpit improves your service experience and is your hidden ally. It offers insights into your liquid handling workstation, without compromising your data security.

Connect to Introspect, Tecan's insights platform, and automatically unlock service excellence with Service Cockpit.

Service Cockpit:

- Enables quicker responses and more effective solutions
- Minimizes questions and interactions
- Enhances overall productivity
- Empowers Tecan experts to troubleshoot issues using remote diagnostics
- Is ready to go when connected to Introspect



Learn more about remote diagnostics and how to sign up to Introspect:
www.tecan.com/remote-diagnostics

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

www.tecan.com

