

# Introspect™ software.


 CONNECT

## EFFICIENT, HOW SO?

Are you tracking productivity with manually maintained spreadsheets?

Introspect™ completely replaces these tools by automatically gathering information such as instrument runtime, consumables usage, error rates and more directly from your Freedom EVO® or Fluent® systems.

## INCREASED PRODUCTIVITY WITH UNPARALLELED INSIGHTS

Modern laboratory automation increases sample throughput and improves productivity by minimizing human error, enhancing precision, and delivering and documenting reproducible test results.

However, in today's busy labs, business-critical decisions on equipment purchases, service contract renewals, consumables spending, and staffing are often made based on incomplete information.

Why has a particular instrument run failed?

Why are some protocols less ef-

ficient, or generate more workflow errors?

Why do similar tasks take longer to complete dependent on the time of the day, or which day of the week they are run?

To optimize instrument uptime and increase throughput, labs need to get to the root of these problems and identify their underlying causes.

Assessing historical instrument usage patterns can highlight capacity gaps, helping labs to deploy resources more effectively and improving troubleshooting.

Tecan has developed Introspect, an easy and secure cloud-based reporting and analysis service, to help laboratories increase their productivity and get unparalleled insights from their instruments.

Data from connected automation workstations – such as Freedom EVO or Fluent systems – is displayed on intuitive dashboards, with real-time information on instrument uptime, consumables consumption and run success rates, completely replacing

time-consuming and less precise manually maintained spreadsheets.

In addition, both historical and archived instrument data can be uploaded to Introspect, allowing lab managers to study previously collected data and make informed decisions on productivity.

One or more instruments, across multiple locations, can be connected to Introspect, and dashboards viewed anywhere in the world through a web browser.

