

Simplifying QC for Freedom EVO® workstations

Tecan has partnered with Artel – the worldwide leader in liquid handling quality assurance – to offer the QC Kit, a fully integrated pipetting performance verification solution for Freedom EVO liquid handling workstations.

The QC Kit will take advantage of Artel's proprietary Ratiometric Photometry™ technology to provide clinical, pharmaceutical and forensic laboratories with a simple and reliable way to meet increasing quality and regulatory demands for performance verification.

The QC Kit includes an Infinite® F50 or Infinite 200 PRO microplate reader, a calibrator plate, a microplate shaker, a barcode scanner and a Freedom EVOware® software add-on. Characterized verification microplates and color dye solutions will be available for ordering directly from Tecan. All color dye solutions are traceable to international reference standards, allowing direct comparison of precision and accuracy performance data between different sites and instruments. Liquid handling performance results can be obtained automatically in just 10 to 30 minutes – depending on the workstation configuration – without the need for sample preparation, standard curve generation or a carefully controlled testing environment.



QC Kit dye solution

To find out more about Tecan's QC Kit, visit www.tecan.com/qckit

Artel is a registered trademark and Ratiometric Photometry is a trademark of Artel.

The QC Kit is under development and will be available in the near future.

Tecan ends support for GENios™ plate reader

After its launch in 1999, the GENios plate reader became one of the most popular plate readers in the life science community. But, even the best things must come to an end, and Tecan has announced it is ending support for this well established, but past its prime, microplate reader.

Tecan is happy that thousands of satisfied customers have been using this robust instrument for decades, but it has now been replaced by a new generation of multimode microplate readers – the Infinite 200 PRO series – offering:

- Faster measurement of more samples in smaller volumes
- Measurement of small volume samples in 96- and 384-well microplates
- Predefined 'one-click' DNA measurements with user-friendly i-control™ software
- Automatic calculation of 260/280 purity ratios
- Measurement of up to 16 2 µl DNA drops in one run, with the patented NanoQuant Plate™
- Elimination of future compatibility problems with computers and printers

Talk to your Tecan sales representative to see how we can help you switch to a new generation of microplate readers, or visit www.tecan.com/dna-reader to learn more.

