

Next-generation Infinite[®] F50

New state-of-the-art ELISA microplate reader with latest LED technology

The Infinite F50 is Tecan's next-generation, 8-channel absorbance reader that provides accurate, reproducible and fast measurements out of a compact, innovative design, and is ideal for a variety of ELISA applications. The Infinite F50 uses the latest LED technology to replace the traditional halogen lamp light source, providing longer service-free life, lower energy consumption and compact dimensions. LEDs typically last 10 times longer than halogen lamps, making inconvenient lamp changes a thing of the past.

The Infinite F50 comes complete with Magellan™ – Tecan's established reader control and data analysis software – to allow quick and easy measurements. The wizard-based software's interface in eight languages (Chinese, English, French, German, Italian, Portuguese, Russian and Spanish) provides full qualitative and quantitative EIA analysis, including endpoint, kinetic and multi-label measurements and all major curve fittings. A comprehensive QC function and a range of data import and export options are also available.

Both the Infinite F50 absorbance reader and Magellan software have been designed, tested, verified and validated according to ISO 13485:2003 quality system standards,



ensuring technical compliance and integrity of data. Furthermore, with the optional Magellan Tracker software, the Infinite F50 has been validated to meet the 98/79/EC IVD directive for *in vitro* diagnostic products and, by providing electronic signatures, records and multilevel user management, meets FDA's Title 21 CFR Part 11 to allow its use in diagnostic applications.

To find out more on Tecan's Infinite F50, visit www.tecan.com/infiniteF50

PowerScanner™ is the new flagship of Tecan's 'microarray suite' range

The PowerScanner is Tecan's open-format microarray scanner, designed for imaging of regular and high density microarrays of tissues, proteins and nucleic acids. This state-of-the-art system features:

- **High resolution** – from 2 to 40 µm, with two configurations available (4 µm system is upgradable to 2 µm)
- **Excellent sensitivity** – offering outstanding signal-to-noise ratio for reproducible detection
- **Large dynamic range** – plus multiple scans at varying gain
- **Unrivalled versatility** – allowing the widest possible range of applications
- **Flexible analysis** – including Array-Pro[®] Analyzer software

The highly customizable PowerScanner allows fully automated processing of up to 48 slides in one run, and features an open platform design to ensure compatibility with virtually all major brands of microarrays using transparent, opaque or reflective substrates. The system is configurable and upgradable according to individual resolution requirements, providing excellent resolution for imaging of both high- and low-density microarrays on a single platform.

To find out more on Tecan's PowerScanner, visit www.tecan.com/powerscanner



Array Pro Analyzer is a registered trademark of Media Cybernetics, Inc.