

## Infinite<sup>®</sup> M1000

### Top class performance with premium Quad4 Monochromators<sup>™</sup>

Reference 10001BO100395



Infinite M1000 with optional injector box



Infinite M1000 with integrated stacker option

The Infinite M1000 offers top-of-the-range quality and performance with robust and innovative new technology, and is the successor to Tecan's high performance Safire<sup>2™</sup> microplate reader. It has been specifically designed and built with the highest quality components and detection modules to ensure optimum performance, robustness and innovation for even the most demanding researchers in drug discovery, life science industry and advanced research institutions.

The high-end detection system gives excellent flexibility and performance through its next-generation premium Quad4 Monochromators, and the platform's modular concept allows upgrades to new detection modes at any time if further applications are required.

Apart from multi-channel absorbance measurements, the fully loaded platform offers you a wide range of detection modes such as fluorescence intensity top and bottom measurements, time-resolved fluorescence (TRF), fluorescence resonance energy transfer (FRET), time-resolved fluorescence resonance energy transfer (TR-FRET), fluorescence polarization (FP) and luminescence.

Tecan has introduced special on-board control functions for the Infinite M1000 that allow pre-programmed measurement workflows to be initiated at the touch of a button on the instrument itself, avoiding the need to go back to your PC in between workflows.

The Infinite M1000 has an optional state-of-the-art injector module, with variable volume and speed settings that allow the use of up to two injectors for dispensing reagents, to replace a manual pipetting step or trigger fast kinetic reactions in fluorescence, luminescence and absorbance modes.

The instrument can be easily combined with a stacker module for batch processing of up to 50 microplates. Alternatively, the Infinite M1000 can be fully integrated with an automated liquid handling workstation from Tecan's Freedom EVO<sup>®</sup> series.

Scientific instrumentation; not for use in human, clinical or diagnostic procedures. All systems or products are tailor-made one-offs that were exclusively developed on customer demand. All displayed pictures and images originate from development, test or installation phases and do not claim to represent the final installation at customer site.

Tecan, Infinite and Freedom EVO are in major countries registered trademarks and Safire<sup>2</sup> and Quad4 Monochromators are trademarks of Tecan Group Ltd., Männedorf, Switzerland.

© 2009, Tecan Trading AG, Switzerland, all rights reserved.