

The Infinite[®] M1000

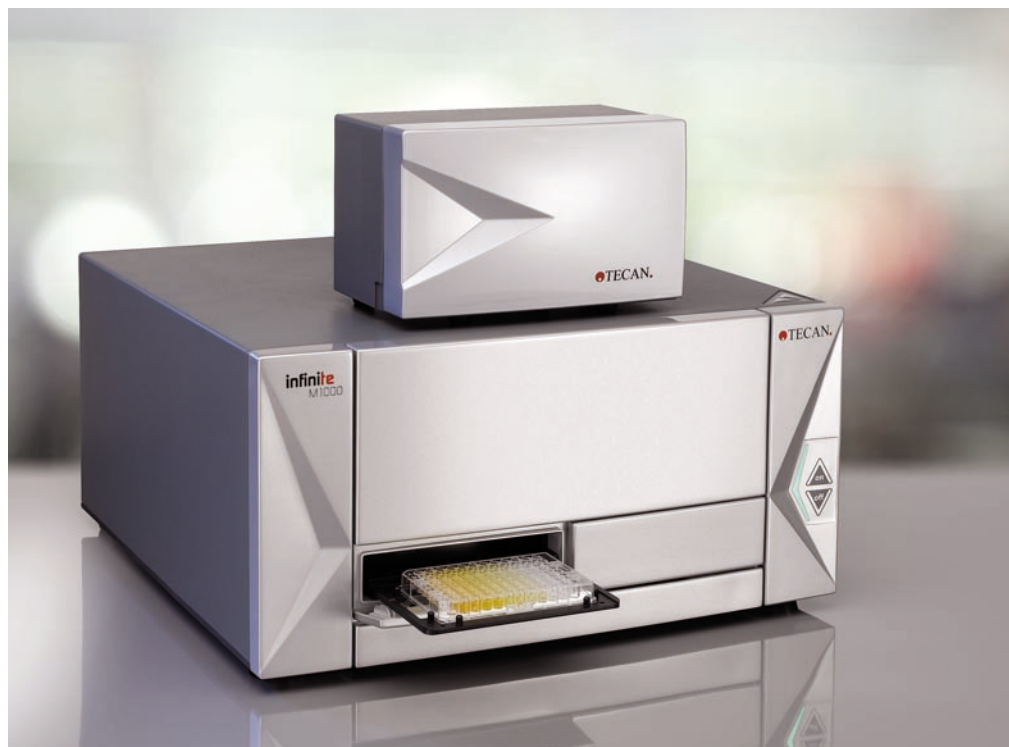
Tecan's multimode flagship microplate reader, is launched at LabAutomation 2008

The Infinite[®] M1000 brings flexibility, sensitivity and speed in one high-end, multimode reader, bridging the gap between research, assay development and screening.

Tecan is launching the Infinite M1000, its new flagship microplate reader, at LabAutomation 2008, booth 305, 27-30 January in Palm Springs, USA. The high-end detection system offers top-of-the-range quality and performance with robust and innovative new technology.

The Infinite M1000 joins the Infinite 200 and Infinite 500 series of detection systems, and is the successor to Tecan's high performance Safire²™ microplate reader. The new platform gives excellent flexibility through its next-generation premium quad4 monochromators™, but it also provides the uncompromising sensitivity and speed typically associated with filter-based systems.

The instrument's special combination of flexibility, sensitivity and speed is ideal for bridging the drug discovery gap between research, assay development and screening in the biopharmaceutical industry, as well as for advanced research laboratories with multiple users and ever-changing applications. The platform's modular concept allows upgrades to new detection modes at any time if further applications are required.



The Infinite M1000 shown with optional injector box

The Infinite M1000 has an optional state-of-the-art injector module that allows 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 injectors have variable volume and speed settings and can be used in combination with the ratio mode to allow fast switching of wavelengths for a wide range of applications. The instrument can be easily combined with a stacker module for batch processing of up to 50 microplates.

Tecan has introduced special on-board control functions for the Infinite M1000 that allow, for example, 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.

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

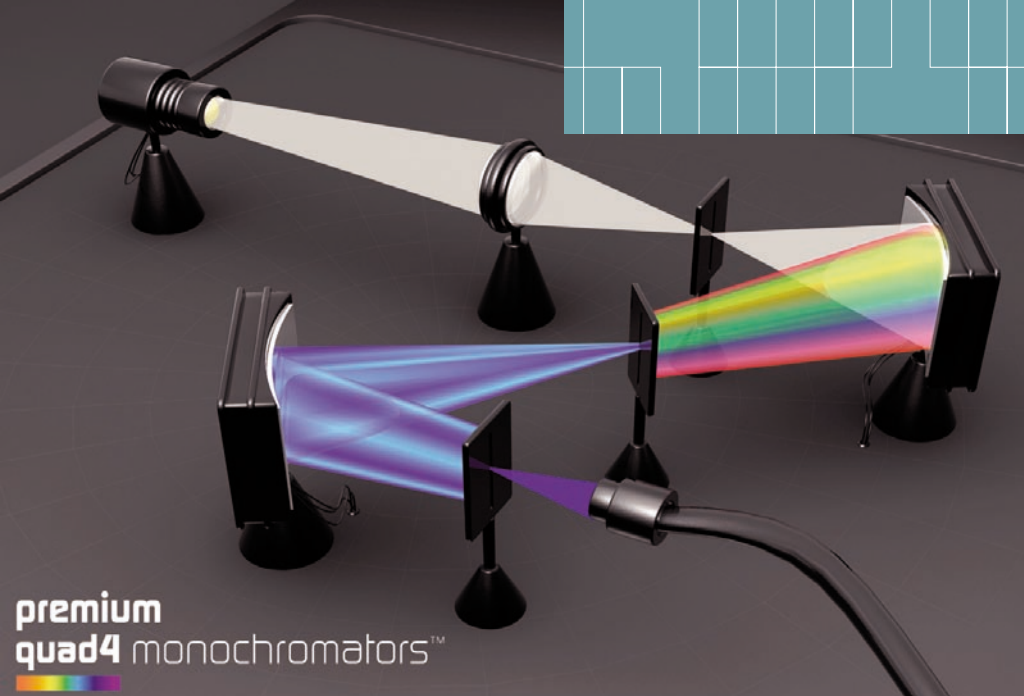


Image of the quad4
monochromators



A stacker module can be used for batch processing

Infinite M1000 – features at a glance

- High-end multimode microplate reader
- Equipped with premium quad4 monochromators
- Provides a wide range of detection modes and plate formats
- Optimized for luminescence and TR-FRET-based assays
- Modular and upgradeable
- Offers injector, stacker and barcode options
- Includes special on-board control functions for plate in/out and starting measurements
- Controlled by easy to use, workflow-oriented i-Control™ software
- Compatible with Magellan™ software for comprehensive data evaluation and processing

resonance energy transfer (FRET) and fluorescence polarization (FP). Furthermore, the Infinite M1000 has been especially optimized for TR-FRET based assays as well as glow luminescence, fast luminescence and dual color luminescence assays.

The Infinite M1000 supports a broad spectrum of applications including the latest biomolecular assays for primary and secondary screening; receptor-ligand binding studies and other molecular interaction assays; kinase assays; protease

assays; G protein-coupled receptor assays; cell-based assays; DNA/RNA quantification; and applications based on UV fluorometry.

If you would like to learn more about our new flagship microplate reader, see www.tecan.com/highend or please visit us at the Tecan booth during LabAutomation 2008 – we look forward to seeing you!

For availability of products in your country for the identified application areas, please contact your local Tecan sales organization. For further reference, please refer to www.tecan.com

Product names and configurations as well as technical details are subject to change and the current product offerings may differ from those outlined in the text.