

# The future of microplate detection has arrived

Tecan has recently updated its Infinite® M1000 high-end multimode microplate reader with new features designed to offer the same top class performance for an even wider range of applications. First launched in 2008, the Infinite M1000 uses Tecan's premium Quad4 Monochromators™ technology to offer exceptional speed, sensitivity and flexibility for academic, biotechnology and pharmaceutical research applications.

Already highly acclaimed for its scanning capabilities, the Infinite M1000 now offers luminescence scanning to complement its existing 3D and ultra-fast absorbance and fluorescence scanning functions. Luciferase reporter mechanisms have become commonplace across a broad range

of disciplines, making development of recombinant luciferases and novel substrates an important area of research. This new feature allows the emission spectra of stable luminescent signals to be recorded at 1 nm resolution using the instrument's emission monochromators, making the Infinite M1000 an ideal platform for luminescence-based assay development.

The format flexibility of the Infinite M1000 has also been expanded, allowing Tecan's patent pending NanoQuant Plate™ for very low volume absorbance measurements to be used alongside virtually all 6- to 1,536-well plate formats. A new double orbital shaking option – moving the plate transporter in a figure-of-eight motion – offers enhanced mixing of reagents and biomolecules, as well as improved suspension of non-adherent cell cultures. For adherent cell cultures, the introduction of an Optimal Read (OR) function has further enhanced measurement reliability in bottom reading fluorescence intensity mode, by performing multiple, spatially separated measurements to achieve maximum illumination and sensitivity.

The new Infinite M1000 also features a redesigned injector module, offering up

to two injectors for precisely controlled pipetting for triggering fast kinetic reactions and bulk dispensing. This state-of-the-art option is available with various syringe sizes and features a tilting reagent carrier to ensure low dead volumes and minimize wastage of substrates and buffers, as well as a bulk dispense function to eliminate tedious pipetting steps for 6- to 384-well plates, saving you both time and reagents for plate-based kinetic studies. The modular, upgradeable nature of Infinite series readers means that customers who purchase the new Infinite M1000 will also be able to benefit from the soon to be launched AlphaScreen® option, making the Infinite M1000 a truly flexible solution to meet your changing laboratory needs.

To find out more about Tecan's Infinite M1000, visit [www.tecan.com/infiniteM1000](http://www.tecan.com/infiniteM1000)

NOTE: New features are only available for the updated Infinite M1000 (REF 30061442) with V2.0 or higher firmware and V1.8 or higher i-control™ software. For further information please contact your local sales representative.

AlphaScreen is a registered trademark of Perkin Elmer, USA.

