

Infinite[®] 200 PRO offers new horizons in microplate detection

Building on the success of the Infinite 200 series, Tecan has developed the Infinite 200 PRO multimode microplate reader, with new enhancements that cater for the needs of today's scientists. The Infinite 200 PRO offers flexible, scalable detection solutions in an easy-to-use modular instrument, providing a full range of leading detection methods, plus innovations such as TR-FRET (HTRF[®]) functionality and improved cell-based applications. A variety of features designed to improve sensitivity and reproducibility are available – such as enhanced fluorescence intensity (FI) bottom reading with optimal read function (OR), linear/orbital shaking and temperature control – ensuring excellent performance and reliability for cell-based and biochemical measurements.

The Infinite 200 PRO uses advanced optics and high performance detectors, giving uncompromised performance in all detection modes, with automated z-focus adjustment for improved FI top measurements. Users can select from either Quad4 Monochromators™ or an intelligent filter slide system, adding a choice of modules to create the perfect reader for their needs, without losing the option to upgrade in the future.

A state-of-the-art injector module allows dispensing of up to two reagents per well, helping to replace a manual pipetting step or to trigger fast kinetic reactions. Its metal-free needles are ideal for ion studies, preventing interference of metal ions in reactions, and low dead volumes minimize wastage of substrates and buffers. The bulk reagent dispense function eliminates tedious pipetting steps for 6- to 384-well plates, with easily accessible prime/wash buttons to ensure straightforward maintenance.

The system is also compatible with Tecan's innovative NanoQuant Plate™, a quartz-based tool that can measure up to 16 samples simultaneously, from just 2 µl sample volumes. This unique tool extends the flexibility of the Infinite 200 PRO for micro-volume spectrophotometry applications, such as quantitation of nucleic acids for quality control and labeling efficiency, creating a truly multipurpose system.

To find out more about Tecan's Infinite 200 PRO, visit www.tecan.com/infinite
HTRF is a registered trademark of Cisbio Bioassays, France.

