

Ignite the productivity of your lab with the Spark™ 10M reader

The new Spark 10M multimode microplate reader is designed to offer greater flexibility and increased productivity for cell biology and genomics customers. From microbiology research and cell-based assays to rapid DNA quantification, the all-new platform delivers a combination of advanced capabilities and exceptional ease of use to simplify your daily work.

At the heart of the Spark 10M are Tecan's unique Fusion Optics, which offer users the choice of filter- or Quad4 Monochromators™-based measurements – or even a combination of both – at the touch of a button. By allowing users to independently choose between filters and monochromators for both excitation and emission, this cutting-edge system means that laboratories no longer have to make a trade-off between flexibility and sensitivity.

This exceptional format flexibility is complemented by patent-pending High Speed Monochromators which allow complete absorbance scans – from 200 to 1,000 nm – in under five seconds. This option combines the speed of a diode array with the precision and flexibility of a monochromator, offering wide-ranging and detailed spectra with unsurpassed wavelength accuracy for rapid DNA and protein quantitation.

The Spark 10M reader has been developed from the outset with cell-based assays in mind, and includes a host of software and hardware features – including a built-in cell counter – designed to simplify cell biology protocols. Precise regulation of environmental conditions is crucial to any cell-based assay, and the system's fully integrated Gas Control Module (GCM™) offers simultaneous control of CO₂ and O₂ concentrations within the measurement



The Spark 10M reader is designed to improve productivity without compromising on flexibility

chamber. Combined with precise temperature regulation and plate shaking options, this provides predictable culture growth under physiological, hypoxic or anaerobic conditions, allowing long-term, walkaway studies. Productivity and reproducibility are further improved by the addition of a patent-pending humidity cassette to provide evaporation protection for any SLAS-format microplate, and automated lid handling to precisely balance gas exchange and minimize the risk of contamination for any measurement mode.

For even greater efficiency, the Spark 10M also features a bright-field cell counting module for automated, label-free cell counting and analysis. This innovative system offers accurate, reproducible cell counting and trypan blue-based cell viability analysis using disposable Cell Chip™ devices. To maximize sensitivity, even for low cell concentrations, the system allows flexible area selection and automated replicate processing, providing simple mean-value calculations, greater accuracy and more precise data.

All these measurement modes and options – along with an injector module with unique stirring and heating capabilities for rapid kinetic assays – are fully integrated and supported by the Spark 10M's user-oriented SparkControl™ software. Accessed via a PC or the instrument's intuitive touchscreen interface, SparkControl makes it easier than ever before to set up and run even complex assay protocols, with one-click operation and remote control for even greater efficiency and productivity.

To find out more about Tecan's Spark 10M multimode reader, visit www.tecan.com/spark



SparkControl makes it easy to run even complex assays

