

Infinite[®] 200 PRO provides a comprehensive solution for cell-based measurements

The new Gas Control Module (GCM™) for the Infinite 200 PRO is a breakthrough for cell-based assays, offering precise regulation of oxygen or carbon dioxide levels within the reader chamber for the first time. Variations in environmental conditions can lead to inconsistent and unreliable data for cell-based optical studies, due to changes in the pH and color of the media during incubation. The GCM provides a more stable culture environment over time, making it ideally suited to *in vitro* investigation of eukaryotic cell lines and allowing the duration of cell-based experiments to be extended without adversely affecting results.

The GCM also ensures greater biological relevance for a wide range of studies, by allowing assays to be performed under hypoxic or physiological conditions. This will extend the experimental window of microplate-based investigations using anaerobic or facultative anaerobic bacteria, allowing incubation within the reader chamber. It also allows the Infinite 200 PRO to be used for a range of biomedical studies, providing researchers with a simple mechanism for studying cell behavior under changing environmental conditions, mimicking events such as ischemia or oxidative shock for *in vitro* investigation.

To find out more on Tecan's Infinite 200 PRO and Gas Control Module, visit www.tecan.com/gcm



Infinite 200 PRO with Gas Control Module

Integrated Adaptive Focused Acoustics™ for the Freedom EVO

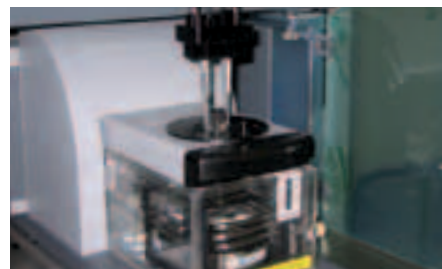
Tecan is collaborating with Covaris, Inc. to integrate Adaptive Focused Acoustics (AFA) non-contact acoustic sample preparation technology into Freedom EVO workstations. High quality, reproducible sample preparation is the critical first step in improving analytical results, and Covaris' highly controlled AFA system sets the standard for today's rapidly advancing analytical technologies and sophisticated assays. This joint offering will be available to customers worldwide, providing a scalable, turnkey solution for automated sample preparation, compound management and next generation sequencing applications.

Kevin Moore, Tecan's Director of Market and Application Management for BioPharma, commented: "AFA is gaining widespread acceptance in the compound management

community for rapid thawing, mixing and dissolution of samples, as well as for DNA shearing for next generation sequencing. We are extremely pleased to be working together with Covaris to provide valuable automated solutions to our customers working in these areas."

To find out more about this application, visit www.tecan.com/covaris

Adaptive Focused Acoustics is a trademark of Covaris, Inc.



Adaptive Focused Acoustics (AFA) sample preparation technology on the Freedom EVO workstation

