

Infinite[®] 200 PRO and Gas Control Module used to investigate cancer therapies

Researchers in Salzburg, Austria, have been using an Infinite 200 PRO multimode reader and Gas Control Module (GCM™) to study the efficacy of novel photodynamic therapies for skin cancers.

Photodynamic therapies based on δ -aminolevulinic acid (ALA) potentially offer a highly selective approach for treatment of dermatological pre-cancers and early malignancies by causing a build up of the photosensitizing agent protoporphyrin IX in neoplastic cells. A collaborative team from

Paracelsus Medical University, the University of Salzburg and the Norwegian University of Science and Technology – along with representatives from Tecan – has recently published a study entitled *Real-time analysis of endogenous protoporphyrin IX fluorescence from δ -aminolevulinic acid and its derivatives reveals distinct time- and dose-dependent characteristics in vitro*¹. This study takes advantage of the in-reader incubation capabilities of the Infinite 200 PRO and GCM to investigate the characteristics of novel ALA derivatives which might improve the pro-drug's cellular uptake.

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



The GCM enables incubation of cell-based assays within the Infinite 200 PRO's measurement chamber

¹ Kiesslich et al. *J Biomed Opt*, 2014, 19(8), 85007

Fluent™ Laboratory Automation Solution tours European exhibitions

Tecan has been showcasing the recently launched Fluent Laboratory Automation Solution at exhibitions across Europe, giving scientists the opportunity to see firsthand how this innovative system could help to streamline their workflows and take the complexity out of cell biology research. The latest addition to the Company's world-class laboratory automation portfolio, the Fluent workstation brings together all the modules and devices required for automation of a wide range of cell-based assays in a single efficient and easy-to-use system, and has already been on display at ELRIG Drug Discovery in the UK, MipTec in Switzerland, and World of Technology & Science in the Netherlands. Visitors to these shows had the chance to discuss the potential benefits of this powerful solution with Tecan's applications specialists, as well as to learn

about the Company's wide range of advanced automation solutions for cell-based and biochemical assays.

To learn more about Tecan's Fluent Laboratory Automation Solution, go to www.tecan.com/fluent

