

Spark[®] Cyto to empower CORE oncology research

Congratulations to Christophe Deben and the team at the University of Antwerp's Center for Oncological Research (CORE) on winning a fully-loaded Spark Cyto plate reader with live cell imaging and real-time cytometry! This competition, which ran from April to September 2019, required labs to submit a written proposal explaining how they would use the system to further their research and advance scientific understanding. A shortlist was selected from a strong field of over 100 entries by a panel of expert judges, and each finalist was invited to give a short webinar presentation providing further details of their planned research. Christophe's winning proposal described how he would use the Spark Cyto to investigate the influence of oxygen levels on chemotherapy-induced CD70 expression and antibody-dependent cellular cytotoxicity in non-small cell lung cancer cells. The judges were impressed by both the scope of this proposal and Christophe's vision of how

the unique features of the Spark Cyto could contribute to the growing field of immuno-oncology.

Christophe commented: "I'm very pleased to have been selected as the winner of this competition. When writing the project proposal, I let my imagination run wild in creating the experimental set-up, and it amazes me that nearly everything I thought of is possible with the Spark Cyto! The versatility of the instrument means that we will be able to use it for nearly every line of research running at CORE, and I can't wait to start working with it and discover all of its capabilities."

Look out for details of Christophe's work in a future issue of the Tecan Journal.

To find out more about the Spark Cyto, visit

www.tecan.com/sparkcyto



CORE received its Spark Cyto system in early December