

New partnership with FortéBio for bioprocessing and drug discovery assays

Tecan and FortéBio have worked together to integrate the new FortéBio Octet® 384 platform on the Freedom EVO® workstation for automated real-time, label-free assays to support bioprocessing and drug discovery workflows. Tecan has collaborated with FortéBio to deliver a robust, off-the-shelf integration that allows the user to get the most out of the Octet 384, allowing increased walkaway time as well as optional liquid handling for assay plate set-up.

Label-free, real-time methods avoid artifacts caused by label interference. FortéBio's Octet instruments provide reliable high throughput measurements for protein-protein, protein-peptide and protein-small molecule interactions using the innovative Dip and Read™ biosensor, that has been designed to analyze specific biomolecular interactions. The Octet RED384 performs label-free quantitation, kinetics and affinity measurements from protein-protein to protein-small molecule interaction analysis, including fragment or small molecule screening.

Tecan has designed a Freedom EVO deck extension plate to mount the Octet securely on the right hand side of the workstation. A robotic manipulator (RoMa) arm performs automated transport and handling of plates and sensor trays. An 8-channel liquid handling (LiHa) arm with fixed or disposable tips, or a MultiChannel Arm™ (MCA) with a 384- or 96-channel pipetting head can also be specified for liquid handling tasks.

Octet is a registered trademark and Dip and Read is a trademark of FortéBio Inc.

The logo for FortéBio, featuring the word "forté" in a blue, lowercase, serif font with a blue underline, and "BIO" in a blue, uppercase, sans-serif font to its right.

The integrated FortéBio Octet 384 platform on the Freedom EVO workstation for automated real-time, label-free assays

