

NGS library preparation at the touch of a button

Tecan has launched the Freedom EVO® NGS workstation to simplify NGS sample preparation and PCR set-up. Offering user-friendly, walkaway automation of library preparation at the touch of a button, it eliminates the need for extensive manual processing.



The Freedom EVO NGS workstation guides the user through instrument set-up

NGS is the fastest growing application in genomics, yet many sample preparation protocols can be labor intensive and time consuming. The Freedom EVO NGS workstation has been designed to offer robust and reliable automation of library preparation, quantification, qPCR set-up, normalization, pooling and capture. It includes all the modules required for precise set-up of NGS sequencing libraries and, with one-touch method selection and step-by-step user instructions via Tecan's intuitive TouchTools™ graphical interface, it reduces hands-on time and increases productivity – even for inexperienced users.

To further streamline processing, Tecan has worked with leading industry partners to supply ready-to-run protocols for a number of popular NGS technologies. Protocols for Illumina's TruSeq® Nano DNA, TruSeq Stranded mRNA* and Nextera® Rapid Capture Exome* chemistries have been developed in collaboration with Illumina and the High-Throughput Sequencing Facility, University of North Carolina at Chapel Hill, with thorough testing and sequencing data analysis to ensure high quality results.

In addition, the system is supplied with protocols covering the Ion Torrent™ library

preparation workflow for Ion AmpliSeq™ panels, developed and supported by Thermo Fisher Scientific, helping you to hit the ground running and simplifying day-to-day activities.

To find out more about Tecan's Freedom EVO NGS workstation, visit www.tecan.com/ngs

*Illumina-qualified protocol.