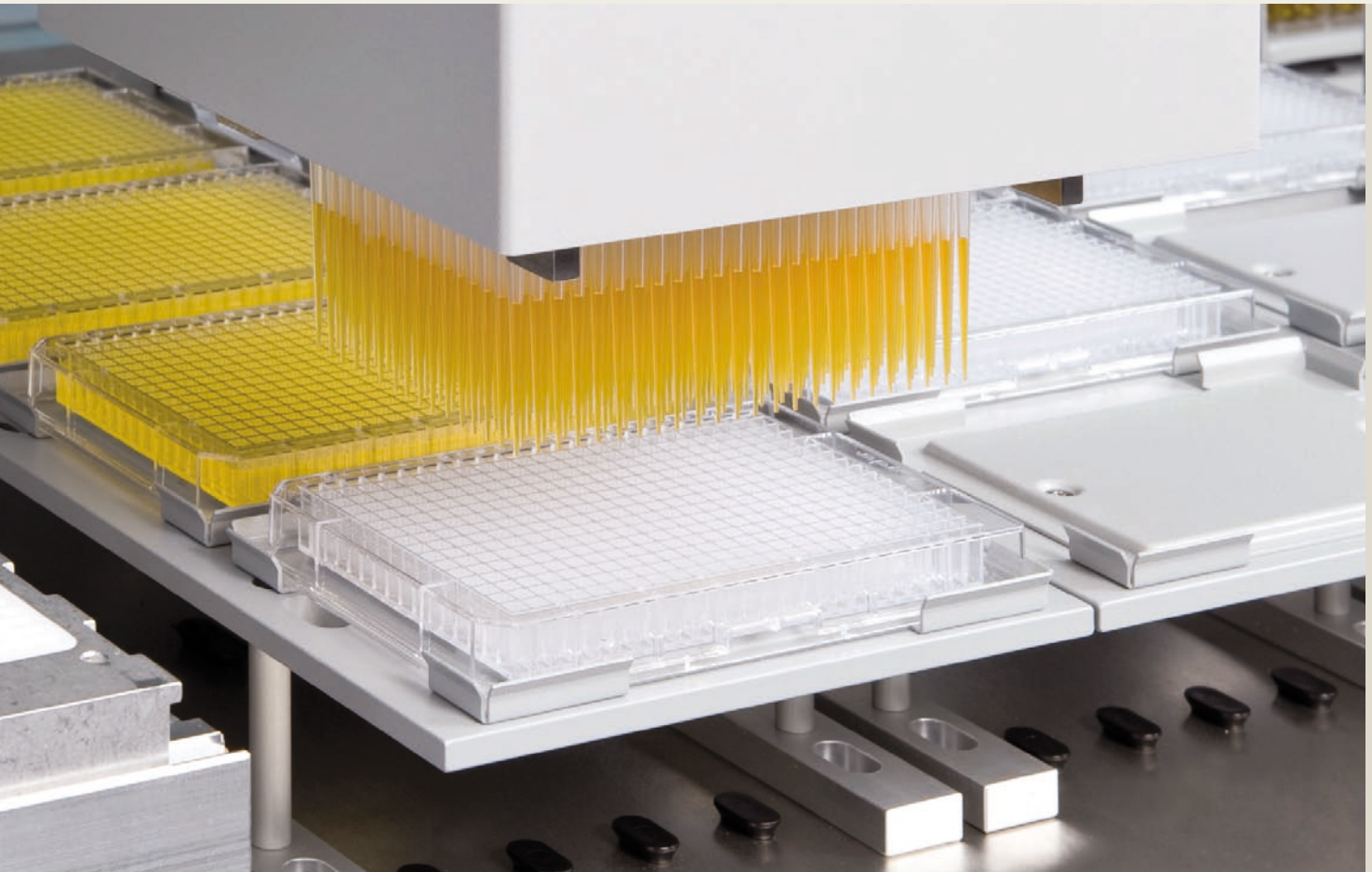


A new 384-channel head for the Freedom EVO[®]



Simultaneous pipetting of 384 wells

Tecan is launching a new 384-channel head for the Freedom EVO[®] liquid handling workstations, offering higher productivity to automated liquid handling processes in pharmaceutical and biotechnology research applications. The MultiChannel Arm™ 384 can be mounted onto Tecan's Freedom EVO 100, 150 or 200 platforms, increasing the efficiency and speed of pipetting processes for better throughput, and delivering a greater level of flexibility.

The head features a very broad volume range of 0.5 µl to 125 µl. It works with

both disposable tips and washable stainless steel tips, which can be automatically interchanged during a run. Alternatively, the head can rapidly mount 96 tips, or just a row of 24 or a column of 16 tips, for pipetting controls or conducting in-plate serial dilutions. An optional gripper can be mounted on the pipetting head to move plates or tip racks, for added functionality. This flexible tool is a great asset for all your pipetting applications.

In addition to the MultiChannel Arm, up to two other independent arms can be

fitted onto the Freedom EVO at the same time, offering multi-functionality and true parallel processing. The new 384-channel head is fully compatible with all Freedom EVO options, and many third party devices.

Not for clinical diagnostics use.