



Automated

Chromatography.

DISCOVER OUR OPTIONS





Tecan's Freedom EVO® liquid handling platform and Fluent® Automation Workstation offer fully automated, parallel column chromatography using Repligen's OPUS® RoboColumn® system, completely eliminating manual processing steps. Miniaturization and parallel processing help to save precious biologics and reagents, reducing costs and shortening development times for new therapeutics.

Tecan's liquid displacement pipetting technology ensures successful parallel processing of up to eight RoboColumns held by the Te-Chrom™ module. The eluates are collected in microplates, and an optional shuttle system enables fraction collection. Large volumes of liquid can be drained directly to a waste container.

Explore our options for scalable, efficient and safe chromatography processing.

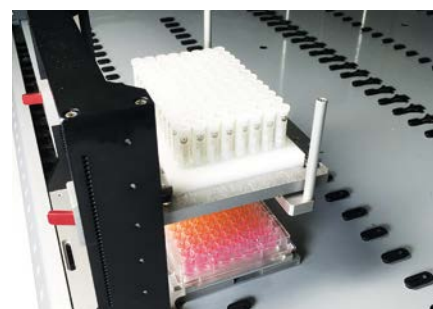
Te-Chrom on Freedom EVO.

INTEGRATION OPTIONS

Te-Chrom Stand alone option for batch elution

Applications:

Sample prep for mass spec, binding studies, sample concentration, PAT



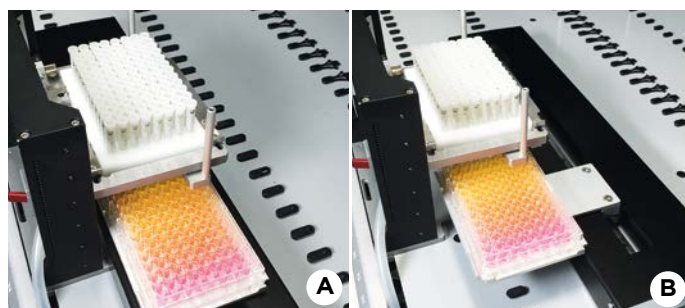
Te-Chrom with Te-Chrom Shuttle which holds the elution plate and allows to collect the RoboColumn flow through in fractions

Applications:

e.g. bioprocessing, method development, condition screening, resin screening, step gradient elution

Two mounting options for Te-Chrom shuttle:

- A) Below the RoboColumns, gives space for a 9 position microplate carrier in the back
- B) Behind the RoboColumns with plate shuttle extension



Te-Chrom with Te-Stack™ combines fraction collection with automated storage of up to 50 96-well elution plates.

Applications:

e.g. bioprocessing, method development, condition screening, resin screening, step gradient elution



Te-Chrom on Freedom EVO.



FREEDOM EVOWARE® SOFTWARE

Te-Chrom Wizard – offers easy chromatography process set-up, from equilibration and sample addition to washing and elution.

Buffer Creation Wizard – easy definition of buffer compositions, stock solutions and labware, allowing the preparation of up to 96 individual buffers in just 10 minutes.

EXPLORE OUR FREEDOM EVO-BASED SOLUTIONS

Antibody purification

Affinity chromatography-purified antibodies can be quickly prepared for downstream MS analysis or binding studies.

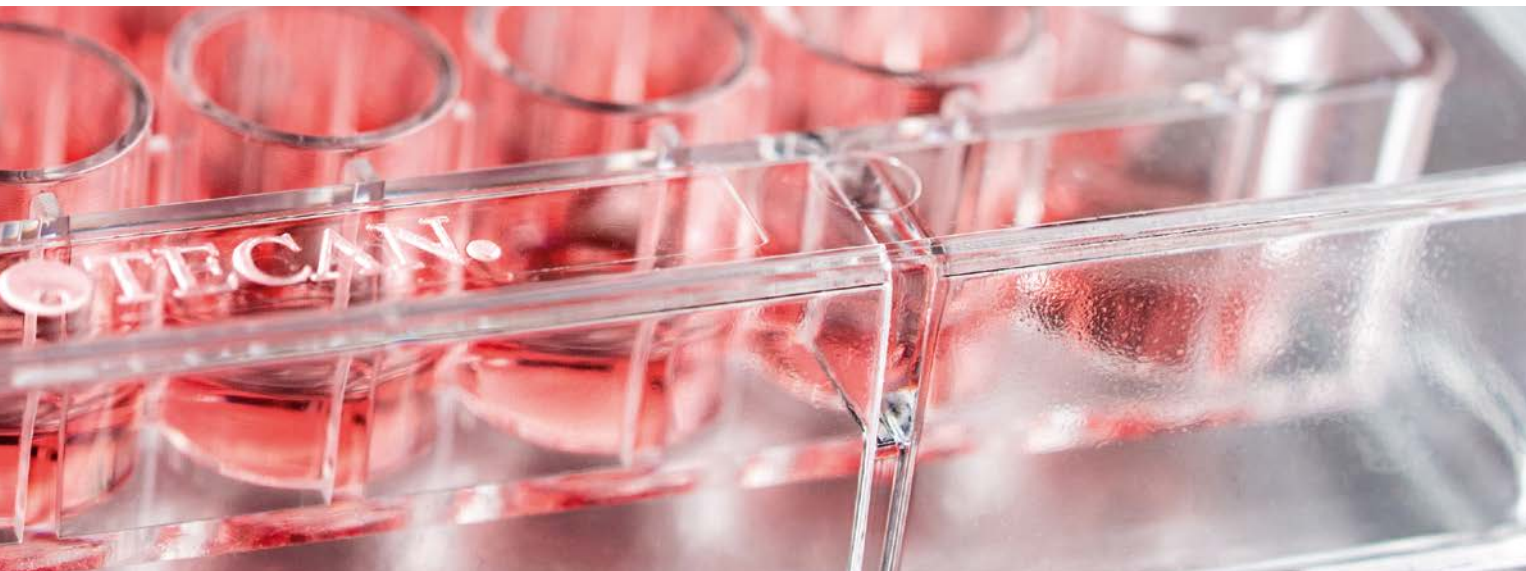
Step gradient elution

Fully automated plate handling allows simultaneous step-gradient elution, enabling you to collect individual fractions and find the optimal separation conditions for your proteins.

Condition screening

Increase walkaway time and add automated data collection and analysis to your workflow using an integrated microplate reader.





OPTIONAL DEVICES



Waste pump – to transfer large liquid volumes from the Te-Chrom waste tray to the liquid waste container.



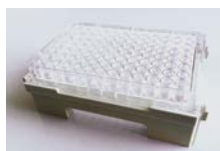
Infinite® 200 PRO reader integration – for absorbance measurements and protein quantification. Can be mounted on the left or right side, below the workdeck, or at the back of the Freedom EVO.



Nine-position microplate carrier – can be placed behind the Te-Chrom Shuttle.



Resting waste – holds the waste option in position with magnetic fixings.



Spacer for shallow-well microplates – adjusts to the same height as deep-well plates.



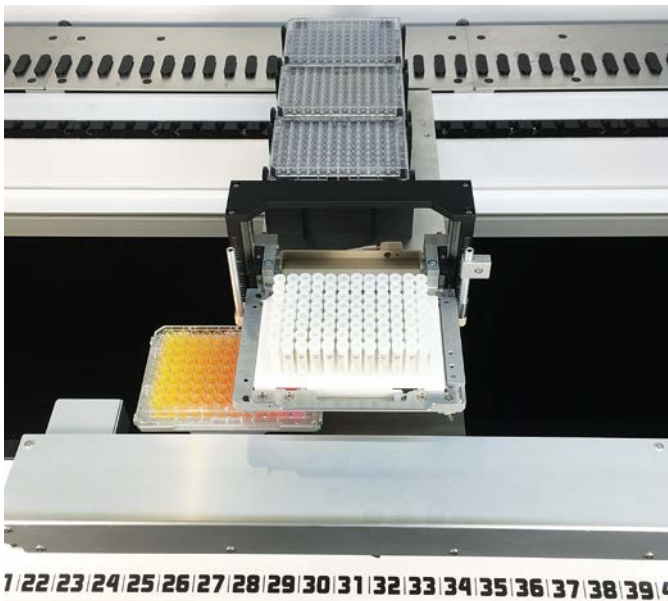
1 ml syringes – for pipetting volumes of up to 1 ml in a single stroke, with flow rates down to 16 cm/h.



Stainless steel fixed tips with conical ends – designed to ensure a good seal with the O-ring of the RoboColumns. Available in two sizes:

- **Standard** – for all RoboColumn/ elution plate combinations using Te-Chrom/Te-Chrom Shuttle.
- **Short** – for all RoboColumn/ elution plate combinations using Te-Chrom, Te-Chrom Shuttle and Te-Chrom with Te-Stack. (**NOTE:** Tips cannot reach the bottom of 50 ml centrifugation tubes.)

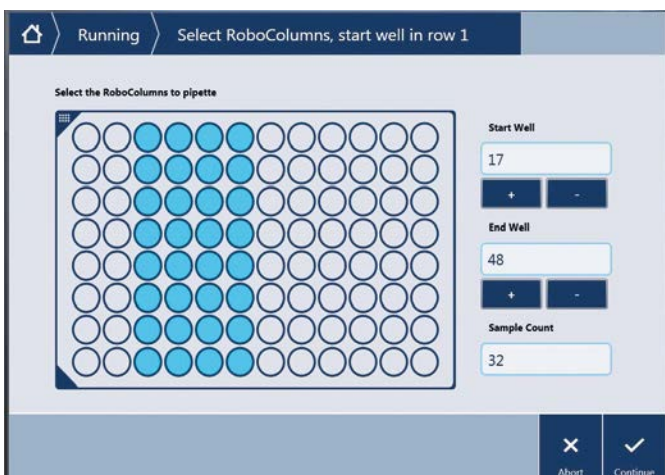
Te-Chrom on Fluent.



INTEGRATION

Both the Te-Chrom and Te-Chrom Shuttle are mounted on a seven-grid wide deck segment which additionally holds three nests for microplates. The Te-Chrom Shuttle offers a choice of batch elution or fraction collection.

Suitable for applications such as bioprocessing, method development, condition screening, resin screening, step gradient elution, sample prep for mass spec, binding studies, sample concentration and PAT.



FLUENTCONTROL® SOFTWARE

Chromatography process set-up - from equilibration and sample application to washing and elution - is straightforward using FluentControl's powerful and intuitive TouchTools™ touchscreen interface.



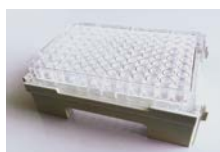
OPTIONAL DEVICES



Waste pump – to transfer large liquid volumes from the Te-Chrom waste tray to the liquid waste container.



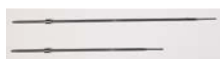
Infinite® 200 PRO reader integration – for absorbance measurements and protein quantification. Can be mounted on the left or right side, below the Dynamic Deck™, or at the back of the Fluent



Spacer for shallow-well microplates – adjusts to the same height as deep-well plates.



Resting waste – holds the waste option in position with magnetic fixings.



Stainless steel fixed tips with conical ends – designed to ensure a good seal with the O-ring of the RoboColumns. Available in two sizes:

- **Standard** – for all RoboColumn/ elution plate combinations using Te-Chrom/Te-Chrom Shuttle.
- **Short** – for all RoboColumn/ elution plate combinations using Te-Chrom, Te-Chrom Shuttle and Te-Chrom with Te-Stack. (**NOTE:** Tips cannot reach the bottom of 50 ml centrifugation tubes.)



1 or 5 ml syringes – for pipetting volumes of up to 4.8 ml in a single stroke, with flow rates down to 8 cm/h.

Automated chromatography. Miniaturized. Parallel.



Explore Repligen's **OPUS
RoboColumns**

www.repligen.com/robocolumn



Discover what Tecan offers
for your chromatography needs

www.tecan.com/parallelchromatography

Contact your local Tecan representative to define the chromatography workstation that fits your needs.

For research use only - not for use in diagnostic procedures.

.....
Australia +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170
Italy +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886 **Spain** +34 93 595 25 31
Switzerland +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +41 44 922 81 11
.....

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication, however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. All mentioned trademarks are protected by law. In general, the trademarks and designs referenced herein are trademarks, or registered trademarks, of Tecan Group Ltd., Männedorf, Switzerland. A complete list may be found at <http://www.tecan.com/trademarks>. Product names and company names that are not contained in the list but are noted herein may be the trademarks of their respective owners. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative.

Tecan is in major countries a registered trademark of Tecan Group Ltd., Männedorf, Switzerland.
© 2018 Tecan Trading AG, Switzerland, all rights reserved.

www.tecan.com

