

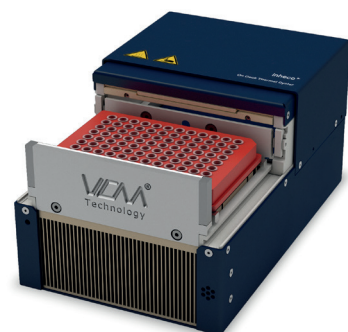
## Fully automate your genomics workflows

Integration of INHECO's On Deck Thermal Cycler (ODTC®) on the Freedom EVO®

Automate more of your genomics workflows – including reaction set-up for endpoint PCR or qPCR, hybridization, Sanger or next generation sequencing – with a dedicated on-deck cyclizer from INHECO. This cyclizer combines an ultra-compact design with exceptional thermal performance for even faster cycling times, and can be flexibly integrated into the Freedom EVO to suit your automation needs.

- Top thermal performance with patent-pending vapor chamber (VCM®) technology:
  - Faster heat transfer for faster cycling times
  - Excellent well-to-well uniformity
- Minimal footprint:
  - Most compact thermal cyclizer on the market
  - Flexible integration options (see Figure 1 overleaf)

PCR cycling methods can be set up with INHECO's script editor software and selected for a specific run using Freedom EVOware®. After the start of a run, Freedom EVOware drives the opening and closing of the ODTC and starts the PCR method for up to 10 integrated ODTC devices.



*ODTC with Bio-Rad low profile plate and auto sealing lid*



*ODTC closed*

Optimized for use with:



Bio-Rad Auto Sealing Lids (eg. Arched Auto-Sealing Lids #MSL-2022)

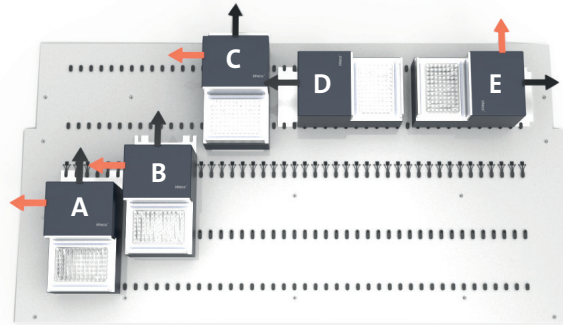


Bio-Rad low profile plates (eg. Hard-Shell® Low-Profile Thin-Wall 96-Well Skirted PCR Plates #HSP-9611)

### Integration options

The Inheco ODTC uses the SiLA standard for software control, and can be controlled by the Freedom EVOware SiLa driver V1.1 and higher.

During initial setup, the EVOware SiLa driver queries the device for its capabilities, and creates a new device driver in EVOware. The driver is capable of being replicated for controlling up to 10 devices, and all user interactions are shown in Tecan's Touch Tools user interface if installed.



The ODTC can be integrated in various positions on the worktable, depending on the workflow of your application. The direction of the ventilation exhaust has to be considered to ensure it blows away from the deck and not over reagents or samples.

### Suggested ventilation exhaust options

Position	ODTC SAP no. (ventilation direction)	Adapter plate SAP no.
<b>A</b> , front of the worktable	30101851 (left in red) or 30101852 (back)	30101077
<b>B</b> , back of the worktable (not under x-rail)	30101852 (back)	30101077
<b>C</b> , back of the worktable (under x-rail)	30101851 (left) or 30101852 (back)	30101078
<b>D</b> , back of the worktable, door right	30101852 (back)	30101078
<b>E</b> , back of the worktable, door left	30101851 (left) or 30101852 (back)	30101078

To find out more contact your local Tecan sales representative, visit [www.tecan.com](http://www.tecan.com), or contact [info@tecan.com](mailto:info@tecan.com).

**Australia** +61 3 9647 4100 **Austria** +43 62 46 89 33 **Belgium** +32 15 42 13 19 **China** +86 21 2206 3206 **Denmark** +45 70 23 44 50 **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 48 174 **Singapore** +65 644 41 886 **Spain** +34 935 95 25 31  
**Sweden** +46 31 75 44 000 **Switzerland** +41 44 922 81 11 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +43 62 46 89 33

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. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative. This brochure may contain reference to applications and products which are not available in all markets. Please check with your local sales representative. 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 [www.tecan.com/trademarks](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.

Tecan is a registered trademark of Tecan Group Ltd., Männedorf, Switzerland. Freedom EVO and Freedom EVOware are registered trademarks of Tecan Group Ltd., Männedorf, Switzerland. ODTC and VCM are registered trademarks of INHECO Industrial Heating & Cooling GmbH, Martinsried, Germany. Hard-Shell is a registered trademark of Eppendorf AG, Hamburg, Germany.

For research use only. Not for use in diagnostic procedures.

© 2015, Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit [www.tecan.com](http://www.tecan.com)

[www.tecan.com](http://www.tecan.com)