



FLUENT®

Product Catalog



This Tecan product catalog contains a wide range of Fluent Options, but is not comprehensive. If something additional is needed, please contact us.

Table of contents

 Platform and arms	3
Fluent series	4
Liquid handling arms	5
Liquid Flexible Channel Arm	5
Air Flexible Channel Arm	6
Multiple Channel Arm 96	8
Multiple Channel Arm 384	8-11
Grippers	12
Robotic Gripper Arm	12-13
Finger Exchange System	14
 Deck segments	15
Wash and waste segments.....	17
Trough segments	18
Microplate segments	19-20
Plates and tubes adapters	21
Fluent tube runners	22
Fluent ID tube runners	23
Trough runners	24
Deck segments with grid pins	25

 Modules	26
Containment	27-28
Barcode scanning	29-30
Logistics and storage	31-33
Heating, cooling, shaking	34-36
Separation and purification	37-39
Frida Reader	40
Mix and Pierce Workstation (tube rotator)	41
Incubators	42
Microbiology modules	43-46
Microplate sealer and seal remover	47
 Integrated and stand-alone devices	48
Microplate washers	49-50
Microplate readers	51-55
Positive pressure workstation	56
Digital dispensers	57
 Plastic consumables	59
Disposable tips	60
LiHa tips	60-62
MCA384 tips	63
MCA96 tips	63
Labware	64
Troughs	64
 Order information	65

Platform and arms

Fluent series	4
Liquid handling arms	5
Liquid Flexible Channel Arm	5
Air Flexible Channel Arm	6
Multiple Channel Arm 96	8
Multiple Channel Arm 384	8-11
Grippers	12
Robotic Gripper Arm	12-13
Finger Exchange System	14



Fluent series

Fluent 480



- Bench-top or cabinet mounted
- Up to two independent robotic arms in any order
- Compact size / medium capacity
- 32 deck grids accommodate up to 30 microplates (ANSI/SBS/SLAS format)

Fluent 780



- Bench-top or cabinet mounted
- Up to three independent robotic arms in any order
- Medium size / high deck capacity
- 52 deck grids accommodate up to 48 microplates (ANSI/SBS/SLAS format)

Fluent 1080



- Bench-top or cabinet mounted
- Up to three independent robotic arms in any order
- Largest system / highest deck capacity
- 72 deck grids accommodate up to 72 microplates (ANSI/SBS/SLAS format)



Liquid Flexible Channel Arm™ (L-FCA)

The L-FCA is an optional arm for the Fluent, which is characterized by a liquid-filled system, controlled by syringe dilutors. The L-FCA has 8 channels with independent vertical movement (Z-axis) for liquid detection, and automatic spacing between channels of 9-38 mm – allowing different labware formats to be addressed. Each channel can use fixed tips or disposable tips (DiTi). Fixed tips can be used for piercing and are washable. Disposable tips help pipetting of carryover-critical steps. Fixed and disposable tip channels can be combined on one arm.

The L-FCA can pipette liquids within different volume ranges, depending on the tips used and the syringe size(s). The volume range covers below 1 µL to over 5 mL, depending on configuration*.



Check out
fixed tips

Part number	Part name	Description
30042145	ARM Fluent FLEXIBLE CHANNEL 8 CHA.LIQUID	L-FCA ordered for new instrument.
30042102	UPGRADE ARM FLUENT FCA 8 CHA.LIQUID	L-FCA for upgrade in field. For fixed tips. Does not include tip ejection system. For 2014 Fluent design.
30042108	UPGRADE ARM FLUENT L-FCA 8 CHA.EJECT SYS	L-FCA for upgrade in field. For disposable tips. Includes tip ejection system. For 2014 Fluent design.
30042146	UPGRADE ARM FLUENT FCA 8 CHA.LIQUID	L-FCA for upgrade in field. For fixed tips. Does not include tip ejection system. For 2018 Fluent design.
30042147	UPGRADE ARM FLUENT L-FCA 8 CHA.EJECT SYS	L-FCA for upgrade in field. For disposable tips. Include tip ejection system. For 2018 Fluent design.



* Find more information in the [specification sheet](#).

Air Flexible Channel Arm™ (A-FCA)

The A-FCA is an optional arm for the Fluent, characterized by an air-filled system, powered by air displacement technology. The A-FCA has 8 channels with independent vertical movement (Z-axis) for liquid detection, and automatic spacing between channels of 9 mm - 38 mm – allowing different labware formats to be addressed. The channels use disposable tips (DiTi) only.

The A-FCA pipets liquids from 0.5 µL – 1,000 µL using different disposable tip sizes to optimize accuracy and precision*.



Check out
disposable tips

* Find more information in the [specification sheet](#).

Part number	Part name	Description
30042201	ARM FLUENT FLEXIBLE CHANNEL 8 CHA AIR MS	A-FCA ordered for new instrument.
30042203	UPGRADE ARM FLUENT FCA 8 CHA. AIR MS	A-FCA for upgrade in field. For disposable tips. Includes tip ejection system. For 2014 & 2018 Fluent design.



Accessories for L-FCA

Fixed tips developed for Fluent

Fixed tips are intended for multiple transfer cycles. The tips are made of stainless steel with a polymer coating. Fixed tips are washable in a Fluent wash station and use thereof helps to reduce consumables costs and logistics. Fixed tips allow piercing through foils and seals and access deep tubes or bottles.



30042195



30042193

Part number	Part name
30042195	TIP STD. PFA COATED SILVER WITH LOCK NUT
30042193	30042196 WITH LOCKNUT
30042196	TIP LOW VOLUME WITH ADJUSTABLE LOCK NUT

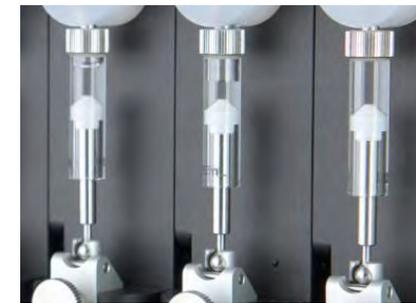
Syringes

The size of the syringe dictates the largest volume that can be handled in a single aspiration or dispense action on any given channel. Syringe sizes can be mixed on an instrument to offer a wide optimized liquid handling range across the pipetting channels.

The syringes are available in three sizes: 5,000 μL , 1,250 μL (standard) and 250 μL .



30042155



30042158

Part number	Part name
30042155	SYRINGE CXP 250 μl
30042158	SYRINGE CXP 1250 μl
30042160	SYRINGE CXP 5000 μl



Multiple Channel Arm™ (MCA) 96

The MCA 96 is an optional arm for the Fluent. The channels of the MCA 96 head are compatible with disposable tips for a volume range from 1 µL to 1,000 µL*. Capacitive liquid level detection is possible at any tip position, even if only one tip is picked up.

The MCA 96 can be upgraded with an optional gripper with the Finger Exchange System™ (FES) for simple labware (ANSI/SBS/SLAS format) handling tasks.



Part number	Part name
30042300	ARM Fluent MULTIPLE CHANNEL 96
30042306	GRIPPER MULTIPLE CHANNEL 96
30042314	SET FINGER GRIPPER MULTIPLE CHANNEL 96
30042301	UPGRADE ARM FLUENT MULTIPLE CHANNEL 96**
30042307	UPGRADE GRIPPER MULTIPLE CHANNEL 96**

* Find more information in the [MCA 96 brochure](#).

Multiple Channel Arm™ (MCA) 384

The MCA 384 is an optional arm for the Fluent. The channels of the MCA 384 head are compatible with disposable and fixed tips. The pipetting head features a very broad volume range from 0.25 µL to 125 µL in the 384-well format and from 0.5 µL to 500 µL in the 96-well format*.



Part number	Part name
30042350	ARM Fluent MULTIPLE CHANNEL 384
30042351	UPGRADE ARM FLUENT MULTIPLE CHANNEL 384**

* Find more information in the [MCA 384 brochure](#).

** For upgrade of Fluent in field.

Check out
disposable tips



Flexible Channel Arm™ (FCA) - Gripper

The FCA gripper is a tool for simple plate movements (ANSI/SBS/SLAS format). It is recommended for genomic workflows such as nucleic acid extraction and clean up. The FCA gripper is ideally suited for instruments with only a Flexible Channel Arm. Centric gripper fingers are mounted automatically between pipetting steps for plate transfers to and from the following locations:

- Pipetting positions*
- Magnetic separation tool*
- Plate shaker*

The FCA gripper is only available for disposable tip systems.

Part number	Part name
30042450	SET FINGERS FCA GRIPPER CENTRIC
30042451	STATION DOCKING FCA GRIPPER

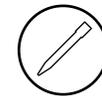


30042450



30042451

* Standard Fluent® microplate nests; Magnum EX/FLX® (Alpaqua® LLC, MA 01915, USA); Bioshake ELM series (Q Instruments, 07749 Jena, Germany).



Accessories for MCA 384

Segments

In addition to the standard microplate nest segments, runners and the disposable tips (DiTi) nest segments, the MCA 384 makes use of the system segment and an active DiTi nest segment, specifically designed for the MCA 384.



30042742



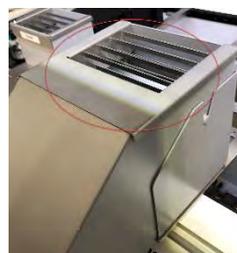
30042743



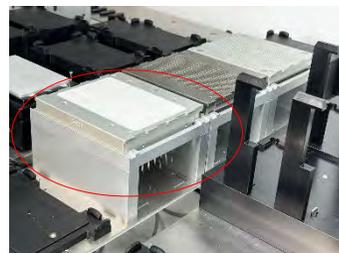
30032026



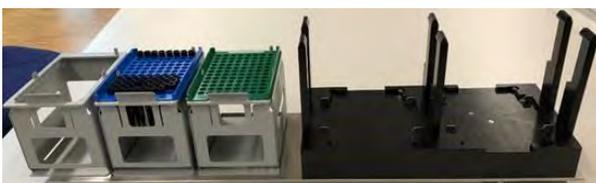
30042825



30042826



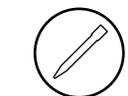
30032027



30042305 (comes without the DiTi nests)

Part number	Part name	Description
30042742	SEGMENT MCA 384 DITI ACTIVE LAND.2	Picking up 384 DiTis with the MCA 384 pipetting head.
30042743	SEGMENT SYSTEM BASE MCA Fluent	Special segment for the MCA 384 to access a wash station, an adapter plate or DiTi boxes.
30032026	NEST SBS MCA384 SYSTEM CARRIER	Nest to place DiTi boxes at the correct height.
30042825	GUIDE DROP TIP 384 WASTE THRU	The waste guide is needed to correctly dispose MCA 384 tips directly to waste.
30042826	GUIDE DROP TIP 384 WASTE FRONT	The waste guide is needed to correctly dispose MCA 384 tips directly to waste.
30032027	RACK ADAPTER MCA 384	Rack to place MCA 384 head adapters.
30042305	SEGMENT MCA 384 DITI ACT.FRONT RACK2 REAR	Picking up 384 DiTis with the MCA 384 pipetting head.

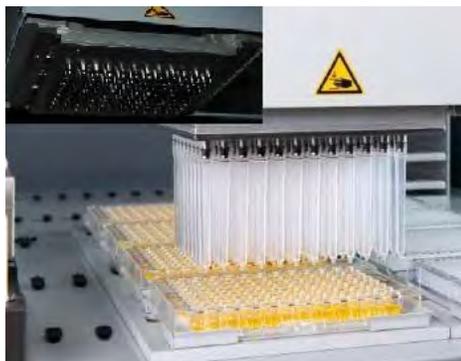
Check out more deck segments



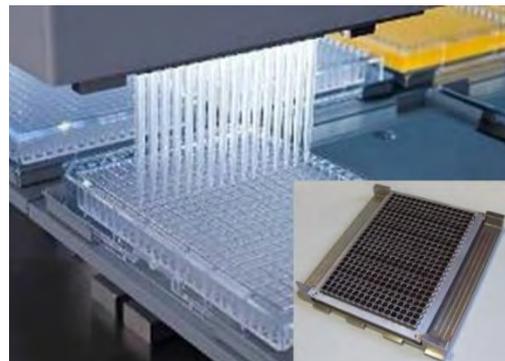
Accessories for MCA 384

Adapter plates

The MCA 384 head can be equipped with a variety of MCA 384 adapters. If required during a pipetting run, the MCA 384 head can automatically mount the appropriate MCA 384 adapter from one or more adapter racks that are situated on the MCA system segment, and return the adapter to the adapter rack at the completion of the step/run.



30032062



30032060

Pictures are a selection of MCA 384 adapter plates. For more information talk to your Tecan sales representative.

Part number	Part name	Description
30032060	ADAPTER 1:1 COMBO 384 FOR MCA 384	Head adapter for 384 tips for the multiple channel arm 384. Row- or column-wise DiTi picking possible.
30032061	ADAPTER 1:1 384 FOR MCA 384	Head adapter for 384 tips for the multiple channel arm 384. No row- or column-wise DiTi picking possible.
30032063	ADAPTER COMBO 96 DiTi MCA 384	Head adapter for 96 tips for the MCA 384.
30032048	ADAPTER 96 MCA 96 DITI MCA 384	Head adapter for usage of MCA 96 tips for the MCA 384. Only use the ANSI/SBS/SLAS box for 96 tips.
30032062	ADAPTER EXT VOL 96 MCA 96 DITI MCA 384	Head adapter for usage of MCA 96 tips for the MCA 384. Only use the ANSI/SBS/SLAS box for 96 tips. Max. pipetting capacity of 125 µL each.
30051709	SET ADAPTER COMBO 384 AND EVA	Head adapter for usage of MCA 96 tips for the MCA 384. Only use the ANSI/SBS/SLAS box for 96 tips. Increased pipetting capacity of 500 µL each.
30032064	ADAPTER+384 FIXED TIPS LC FOR MCA 384	Head adapter with 384 fixed tips for the multiple channel arm 384. Volume range: 3 µL – 125 µL (water), 2 µL – 125 µL (DMSO).
30032065	ADAPTER+384 FIXED TIPS SC FOR MCA 384	Head adapter with 384 fixed tips for the multiple channel arm 384. Volume range: 1 µL – 15 µL (water), 0.5 µL – 15 µL (DMSO).
30032066	ADAPTER+96 FIXED TIPS LC FOR MCA 384	Head adapter with 96 fixed tips for the multiple channel arm 384. Volume range: 5 µL – 125 µL (water and DMSO).
30032067	ADAPTER+96 FIXED TIPS SC FOR MCA 384	Head adapter with 96 fixed tips for the multiple channel arm 384. Volume range: 0.5 µL – 15 µL (water and DMSO).

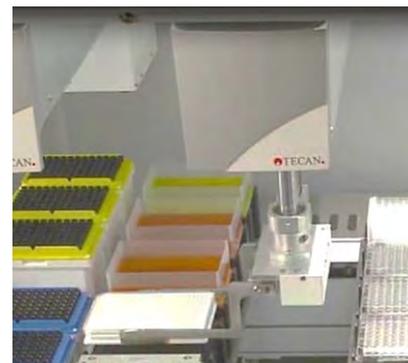


Robotic Gripper Arm™ (RGA)

The Fluent instrument can be equipped with one Robotic Gripper Arm. This robotic arm can transfer a variety of labware to and from the deck, storage positions, and devices such as plate shakers or microplate readers.

An RGA can be fitted with either a regular gripper head for fixed gripper fingers or with a Finger Exchange System™ (FES) head for exchangeable fingers. The RGA's technical data and work ranges are identical for both types of gripper heads. The RGA can also be 'standard' length or have an extended reach (long Z) version that can reach beneath the deck, for access to cabinet-integrated devices.

Part number	Part name
30042400	Robotic Gripper Arm Standard Z axis with Regular Gripper Head
30042405	Robotic Gripper Arm Long-Z** axis with Regular Gripper Head
30042401	Robotic Gripper Arm Standard Z axis with Finger Exchange System Gripper Head
30042406	Robotic Gripper Arm Long-Z** axis with Finger Exchange System Gripper Head



30042400



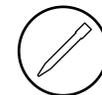
30042405



30042401



30042406

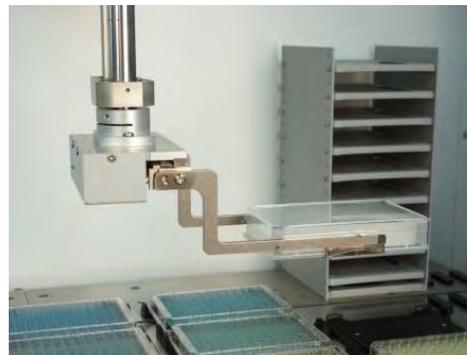
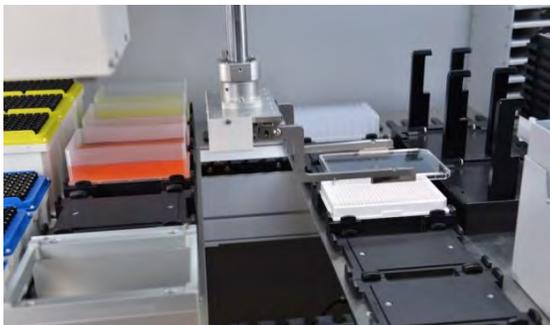


Check out
Finger Exchange System

Robotic Gripper Arm (RGA) Fingers

Various gripper fingers are available for the RGA, depending on the application. These include:

- Eccentric, eccentric long fingers or centric fingers for plates, deep-well plates and lids
- Tube fingers for tubes or vials



30042430



30042431



30042432



30042433



30042425

Part number	Part name	Description
30042430	SET FINGER REGULAR MP ECCENTRIC	Standard eccentric RGA fingers
30042431	SET FINGER REGULAR MP CENTRIC	Standard centric RGA fingers
30042432	SET FINGER REGULAR TUBE STD	Standard RGA tube fingers
30042433	SET FINGER REGULAR MP ECC LONG	Long eccentric RGA fingers
30042425	BARCODE READER MODULE	Optional Barcode reader module for mounting on regular gripper head.



Finger Exchange System™ (FES)

The Finger Exchange System is an optional head for the RGA that allows to exchange fingers 'on the fly'.

- Finger Exchange System - option for RGA
- Extends range of labware and devices to be handled in a process
- Extends walkaway time and scope of process
- Each finger type uniquely identified
- Custom fingers possible via Tecan Labwerx™

Part number	Part name	Description
30042415	FINGER EXCHANGE SYSTEM ECCENTRIC	ECCENTRIC fingers
30042416	FINGER EXCHANGE SYSTEM CENTRIC	CENTRIC fingers
30042417	FINGER EXCHANGE SYSTEM TUBE STD. 43MM	Tube fingers
30042418	FINGER EXCHANGE SYSTEM ECC. LONG	ECCENTRIC long fingers. Required for 4 slot MIO cell plate incubator
30042437	DOCKING STATION - WIDE	Wide docking station is for e.g. Eccentric 30042415 or Eccentric long 30042418 FES fingers.
30042438	DOCKING STATION - NARROW	Narrow docking station is for e.g. Centric 30042416 or Tube 30042417 FES fingers.



30042415



30042416



30042417



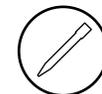
30042418



30042437

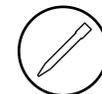


30042438



Deck segments

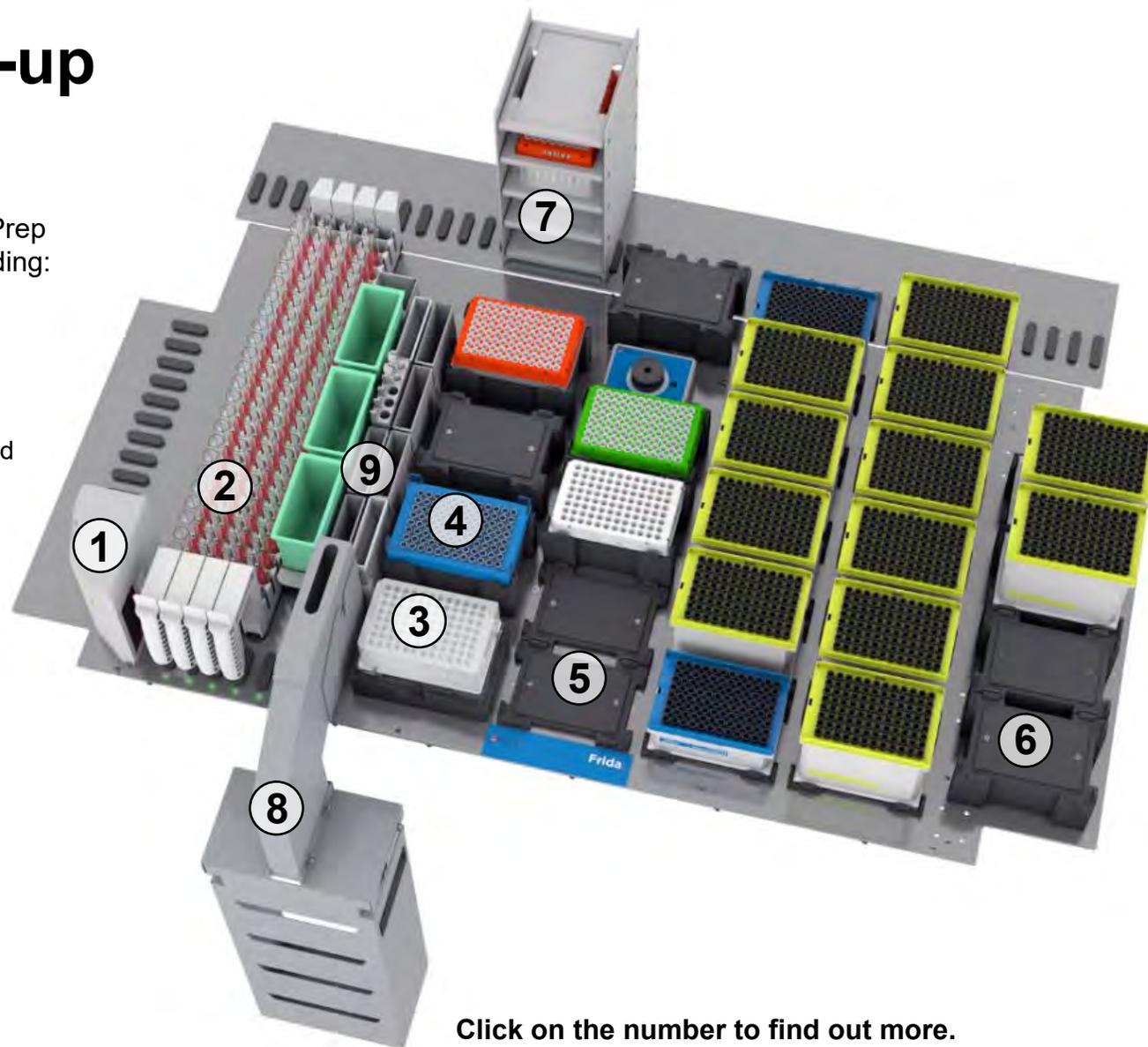
Wash and waste segments	17
Trough segments	18
Microplate segments	19-20
Plates and tubes adapters	21
Fluent tube runners	22
Fluent ID tube runners	23
Trough runners	24
Deck segments with grid pins	25



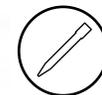
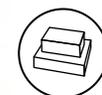
Example of a worktable set-up

The following image showcases the worktable for the Fluent DreamPrep NAP workstation including some deck segments and modules, including:

- 1 Fluent ID™ for automated barcode scanning on samples, handheld barcode scanner for reagent barcode tracking
- 2 Tube ID tube runners
- 3 Q Instruments Bioshake D-30 T-ELM shaker
- 4 Alpaqua Magnum FLX® Enhanced Universal Magnet Plate
- 5 Frida Reader™ on-deck module
- 6 Carriers for disposable tips and 96-well plates
- 7 Hotel
- 8 Waste chute
- 9 Trough segments



Click on the number to find out more.



Wash and waste segments

Fluent deck segments support wash stations for fixed tips and waste liquid, and waste chutes for disposable tips and labware. For cabinet-mounted Fluents the liquid and waste containers can be placed in the cabinet underneath the system.

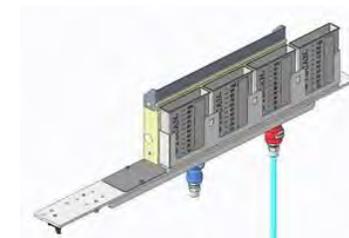
Part number	Part name	Description
30042795	WASTE CHUTE FRONT 2 GRID	DiTi Waste chute with cover for bench top installation.
30042749	SEGMENT ACTIVE WASH WASTE TROUGH 100ML	Deck Segment for Fluent liquid FCA. Active wash and waste station for low volume fixed tips.
30042777	HOLDER BAG CHUTE THRU 2 GRID	For 2 grid waste.
30042745	SEGMENT WASH WASTE THRU TROUGH 100ML	DiTi Waste chute with cover. (*)
30042523	SEGMENT WASTE THRU TROUGH 3*320ML	Includes disposable tip waste chute with cover and 3 reagent trough positions 320 mL. (*)
30042524	SEGMENT WASTE THRU TROUGH 8*100ML	Includes a disposable tip waste chute with cover and 8 reagent trough positions 100 mL. (*)
30042737	SEGMENT NEST 61MM LAND.4 WASTE THRU	Waste with four 61 mm nests for microplates. (*, **)
30042738	SEGMENT NEST 7MM LAND.4 WASTE THRU	Waste with four 7 mm nests for microplates. (*, **)
30042741	SEGMENT NEST FCA DITI TRAY 4 WASTE THRU	Waste with four DiTis in trays (not SBS DiTis). (**)

(*) Can be used only with a cabinet.

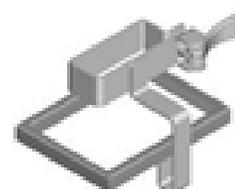
(**) Nest positions can be removed; an extra one can be added.



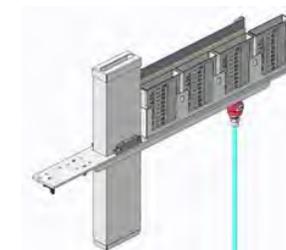
30042795



30042749



30042777



30042745



30042524



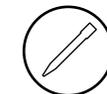
30042737



30042738

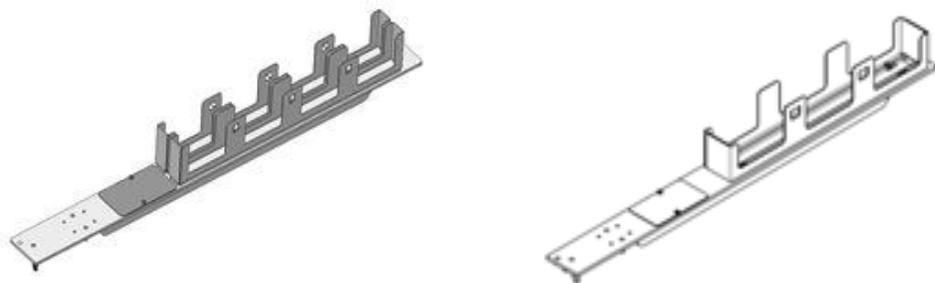


30042741

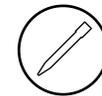


Trough segments

Fluent deck segments support 100 mL and 320 mL reagent troughs. The segments can be easily positioned on the deck and locked in place during use. The reagent positions can be loaded manually prior to use.

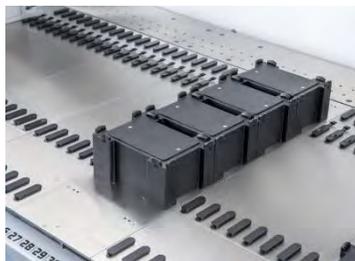


Part number	Part name	Description
30042521	SEGMENT TROUGH8*100ML	8 reagent troughs for 100 mL.
30042522	SEGMENT TROUGH 3*320ML	3 reagent troughs for 320 mL.

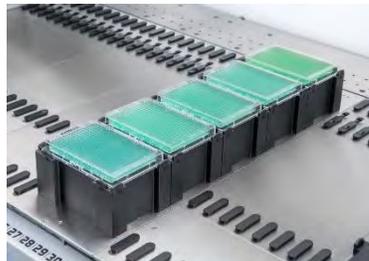


Microplate nest segments

Fluent microplate nest segments provide support for up to 6 ANSI/SBS/SLAS format pieces of labware, such as microplates, deep-well plates or tip boxes. Segments can be easily exchanged between methods to optimize deck usage for each protocol.



30042720



30042722



30042724



30042721

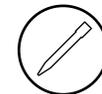


30042723



30042725

Part number	Part name	Description
30042720		4 positions, landscape orientation
30042722	SEGMENT NEST 61MM	5 positions, landscape orientation
30042724		6 positions, landscape orientation
30042721		4 positions, landscape orientation
30042723	SEGMENT NEST 7MM	5 positions, landscape orientation
30042725		6 positions, landscape orientation



Microplate and disposable tips nests

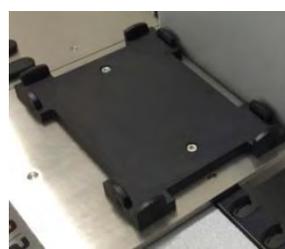
Each ANSI/SBS/SLAS nest position on a deck can be manually exchanged to optimally configure the Fluent deck layout for disposable tip stacks or trays, microplates or similar labware.



30042780



30042781



30042782



30042784



30042787



30042785



30042783

Part number	Part name	Description
30042780	NEST MP/DITI 7MM	7 mm nests for micro plates or standalone DiTi boxes. Includes adapter (30042784).
30042781	NEST MP/DITI 61MM	61 mm nests for micro plates or standalone DiTi boxes. Includes adapter (30042784).
30042782	NEST MP/DITI 7MM BASE	7 mm nests for micro plates or standalone DiTi boxes. Does not include adapter (30042784).
30042784	ADAPTER NEST MP STANDAR D	Can be used with Nests 30042782 and 30042783
30042787	NEST MP/DITI 6MM SPACER	Can be used with Nests 30042782 and 30042783
30042785	NEST FCA DITI TRAY	Can be used with Nests 30042782 and 30042783
30042783	NEST MP/DITI 61MM BASE	61 mm nests for micro plates or standalone DiTi boxes. Does not include adapter nest.

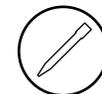


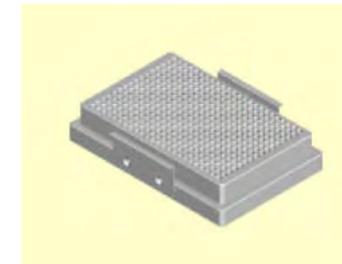
Plate and tube adapters

Each microplate nest can be adapted to hold other items such as tubes, by use of ANSI/SBS/SLAS-format adapters. These include:

Part number	Part name	Description
10762022	ADAPTER MP PCR 96 WELL	Plate Adapter, PCR, 96-well, for Standard Microplate Nest (10613038)
30012982	ADAPTER PCR 384WELL	Plate Adapter (has flat bottom), PCR, 384-well, Temperature Conductive, fits on Microplate Nests (exclusive for Tecan US: 30011871, 30011674).
10762021	ADAPTER MP CRYOVIAL 0.5ML	Tube Adapter, 0.5 mL Microfuge with Hinged Caps, 24-Position (4 x 6), Microplate Format. Has holes for cap storage.
10762017	ADAPTER FOR MP 1.5ML EPPENDORF 24POS	Tube Adapter, 1.5 mL or 2ml with hinged caps, 24-Position (4 x 6), Microplate Format, Aluminum. Have holes for cap storage.
30012613	ADAPTER BLOCK EPPENDORF 1.5/0.5ML	Tube Adapter, 1.5 mL Microfuge with Hinged Cap, 12-Position (4 x 3), and 0.5 mL Microfuge with Hinged cap, 12-Position (4 x 3), Microplate Format, Aluminium.
30014924	RACK ADAPTER 96 POS. MICRONIC TUBE	Tube Adapter, Minitube, 96-Position, Microplate Format, Aluminium. Tested with Micronic and Matrix Minitubes.
10762020	ADAPTER MP CORNINGTUBES 13MM 48POS	Tube Adapter, 13 mm Tubes, 48-Position, Microplate Format, Aluminium.
30012856	ADAPTER 60 POS. WHEATON VIAL 2ML CPL.	Vial Adapter, 13 mm (2 mL Wheaton Vial), 60-Position (6 x 10), microplate format, Aluminium.



10762022



30012982



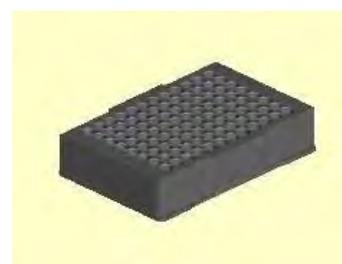
10762021



10762017



30012613



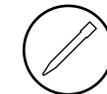
30014924



10762020



30012856



Fluent tube runners

Fluent tube runners allow tube collections to be slid onto the worktable on a Fluent grid segment.

Part number	Part name	Description
30042762	RUNNER TUBE 10MM 1*24	1 x 24 10 mm tubes
30042763	RUNNER TUBE 13MM 1*24	1 x 24 13 mm tubes
30042764	RUNNER TUBE 16MM 1*16	1 x 16 15 mm tubes
30042766	RUNNER TUBE 13MM 4*24	4 x 24 13 mm tubes
30042767	RUNNER TUBE 16MM 6*16	6 x 16 16 mm tubes
30042769	RUNNER TUBE EPPENDORF 1.5-2ML 1*24	1 x 24 1.5 - 2.0 mL Eppendorf tubes
30042770	RUNNER TUBE FALCON 15ML 1*16	1 x 16 15 mL Falcon tubes
30042771	RUNNER TUBE FALCON 50ML 1*10	1 x 10 50 mL Falcon tubes



30042762



30042763



30042764



30042766



30042767



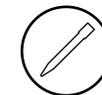
30042769



30042770



30042771



Fluent ID™ tube runners

Fluent ID runners include barcoded positions for barcoded tubes and are designed to be used with the Fluent ID module(s). The runners can also be placed onto standard grid segments for systems (or positions) without Fluent ID module(s).

Part number	Part name	Description
30042506	RUNNER TUBE 10MM 1*32 Fluent ID	1 x 32 10 mm tubes
30042507	RUNNER TUBE 13MM 1*32 Fluent ID	1 x 32 13 mm tubes
30042508	RUNNER TUBE 16MM 1*26 Fluent ID	1 x 26 16 mm tubes
30042499	RUNNER TUBE 17MM 1*24 Fluent ID	1 x 24 17 mm tubes
30042509	RUNNER TUBE EPPENDORF 2ML 1*32 Fluent ID	1 x 32 2 mL tubes



30042506



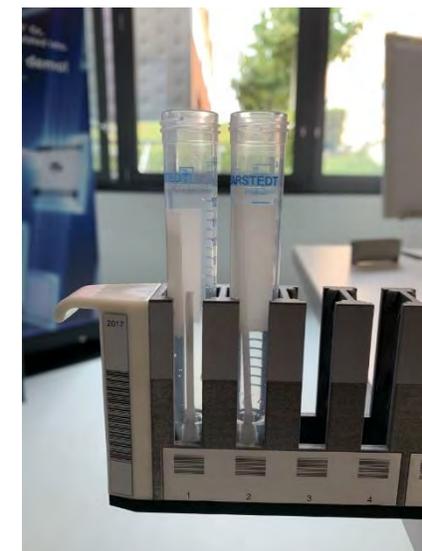
30042507



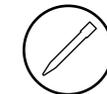
30042508



30042509



30042499

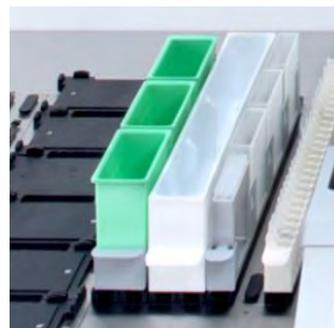


Trough runners

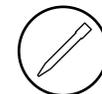
Trough runners allow reagents to be loaded into the runner outside the instrument and loaded into place on a Fluent Grid Segment



30042773

30042774 (green troughs)
30042775 (white trough)

Part number	Part name	Description
30042773	RUNNER TROUGH 100ML 1*4	Occupies 1 grid
30042774	RUNNER TROUGH 320ML 1*3	Occupies 2 grids
30042775	RUNNER TROUGH 1000ML 1*1	Occupies 2 grids



Deck segments with grid pins (required for runners)

Fluent grid segments support the positioning of tube runners, troughs, plate hotels and devices. The segments are mounted on the Fluent deck and can easily be exchanged for other segment types allowing the deck layout to be optimized for each protocol.

Grid segments are available to support Freedom EVO® carriers and others can also provide access points for the Robotic Gripper Arm to reach below the deck (for device integration etc.).

Part number	Description
30042700	8 grids
30042701	6 grids
30042702	3 grids
30042703	2 grids
30042705	9 grids with access whole that is 7 grids wide
30042707	8 grids, required for Freedom EVO carriers on Fluent deck
30042708	6 grids, required for lower deck segments



30042700



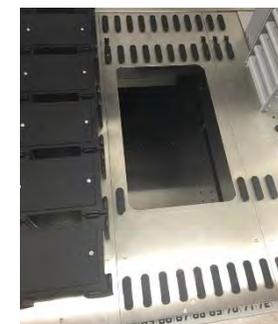
30042701



30042702



30042703



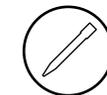
30042705



30042707



30042708



Modules

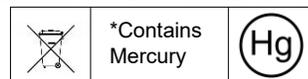
Containment	27-28
Barcode scanning	29-30
Logistics and storage	31-33
Heating, cooling, shaking	34-36
Separation and purification	37-39
Frida Reader	40
Mix and Pierce Workstation (tube rotator)	41
Incubators	42
Microbiology modules	43-46
Microplate sealer and seal remover	47



HEPA Hood

Containment solution manufactured by Bigneat to fit all Fluent sizes. Delivers clean air, according to clean room class ISO 5, to maintain sterile environments. Application examples include genomic workflows, cell-culturing and cell-based assays.

- Clean air and UVC protection
- Laminar air flow
- BS EN ISO 146644-1/1999, ISO Class 5
- Maintains compact instrument footprint
- Integrated via touch screen
- UVC light*; UVC resistant panel sets
- Standard and long-Z Robotic Gripper Arm
- Available for each Fluent instrument size
- HEPA on cabinet & Long-Z gripper = 2,900 mm



Custom BSLII & III Enclosures

Containment solutions for laboratories with higher biosafety levels (or non-biological hazards such as chemicals) can be offered through our customized solutions service, Tecan Labwerx™. Labwerx can adapt systems to fit into containment solutions from other manufacturers and upgrade access doors with door locks to ensure operator safety while the instrument is running.



Speak with your local sales representative to receive a quote for a BSLII or III or chemical safety enclosure.

Part number	Description		
30042948			Fluent 480
30042949	HEPA hood with touch screen and standard Z-Standoff	110 volt	Fluent 780
30042950			Fluent 1080
30042945			Fluent 480
30042946	HEPA hood with touch screen and standard Z-Standoff	240 volt	Fluent 780
30042947			Fluent 1080



Dust covers

The dust cover eliminates dust from falling onto the Fluent worktable and is suitable for applications that do not require a laminar flow of clean air, or operator protection. The dust cover features two ventilation ducts with built-in dust filters to allow heat to escape from the worktable. An optional UV light cover is available to provide UV sterilization for the system.

Part number	Description	
30042592		Fluent 480
30042593	Metal Top Cover	Fluent 780
30042594		Fluent 1080
30042530		Fluent 480
30042530	UV Light Option	Fluent 780
30042531		Fluent 1080



Fluent ID™

Fluent ID barcode scanning module for tubes

- Scanning 1D barcodes while loading (less than 30s for 96 samples)
- Immediate feedback on every tube runner on touchscreen
- Immediate error handling
- Loading is simplified by customizable lights and runner stickers
- 1 Fluent ID can have up to 6 runners (up to 32 tubes per runner),
- Max. 4 Fluent ID for Fluent 780 and 1080 (768 tubes / batch) and 3 Fluent ID for Fluent 480 (576 tubes / batch)
- Unlimited batches per run (i.e. reloading during a run is possible)
- Innovative module complements an array of other barcode reading options – including arm-mounted and 2D ANSI/SBS/SLAS rack scanners – to maximize process security



Check out
Fluent ID tube runners



30042504



30042505

Part number	Description
30042504	SEGMENT Fluent ID 1D LEFT
30042505	SEGMENT Fluent ID 1D MIDDLE

Each Fluent ID segment requires the placement of a reflector on the grid immediately to the right of the Fluent ID segment. Two reflector options are supplied with every segment Fluent ID:

- a reflector tab to be mounted on a segment trough (30042522, 30042523, 30042521 or 30042524) or a segment with grid pins (30042700, 30042701, 30042702, 30042703, 30042705)
- a reflector sticker to be affixed on the left wall of a front waste chute (30042522, 30042748, 30042749 or 30042521 with 2-grid waste chute 30042795 or 30042720, 30042721, 30042739, or 30042786 with 6-grid waste chute 30042794)



Ziath DataPaq™ Express Rack Reader

The Ziath DataPaq Express Rack Reader is a 2D data matrix tube camera reader with an image capture time of only one second. DataPaq Express is only marginally larger than the ANSI/SBS/SLAS nest itself, making it ideal for robotic integration.

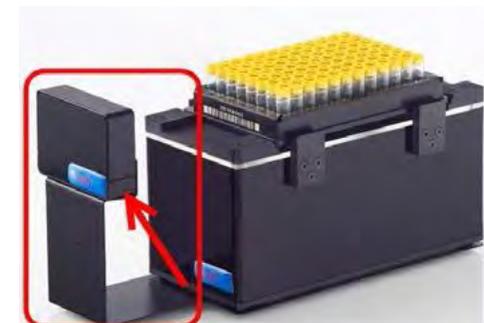
Compatible with all 2D barcoded tube racks in ANSI/SBS/SLAS format, the DataPaq Express is the smallest rack scanner currently available. Its camera-based scanning technology delivers rapid imaging and decoding.

Key Features:

- Scan all makes of 2D barcoded tubes
- Export data to Excel, text file, XML, JSON
- Essential tool for sample tracking



30099754



30099756
(option to read the 1D code on the side / edge of racks)

Part number	Description
30099754	BCR ZIATH 2D 96POS. DATAPAQ EXPRESS
30099756	BCR ZIATH 1D SCANNER FOR DATAPAQ EXPR.



Fluent Stacker™

Serial access storage for plates & tips on right and / or left of Fluent.

- Stores up to 50 microplates or 40 tip racks per stack for fast transfer to the worktable
- Up to 4 stacks per stacker (maximum of 200 microplates or 160 tip racks)
- Allows pipetting directly into plates on its transfer station for added speed
- Used tip boxes or plates discarded directly to a waste bin
- Benchtop or cabinet mounted



Part number	Description
30070071	Fluent Stacker

Fluent Carousel™

Random access storage for plates and or disposable tip boxes on the right and / or left of the Fluent.

- Capacity: 210 standard microplates
- Average plate delivery time: 15 seconds (incl. BC reading)
- Linear transfer stations at 103, 253 & 383 mm for presenting microplates/tips into Fluent work area
- Plate presence sensor
- Includes 10 freely selectable stackers
- 7 shelf pitches from 19 mm (1536-well MP) – 131 mm (5000 µL tips)
- Benchtop or cabinet mounted



Part number	Description
30070070	Fluent Carousel



Hotels

Hotels are a simple solution to extend your storage space on the work deck. Accessing plates in a hotel requires a robotic gripper arm (either standard RGA or RGA with FES) with eccentric gripper fingers. Hotels can be fitted with a barcode scanner to register barcodes during plate movements.



Example pictures of hotels



Barcode reader for hotel

Part number	Description
30042868	Static hotel for 9 micro plates
30042869	Static hotel for 5 deep well plates
30042863	Hotel for micro plates
30042864	Hotel for cell plates
30042865	Hotel for deep well plates
30042867	Barcode reader for hotel

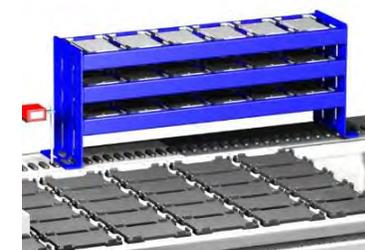
Cooled Hotel

Hotel tower with 2 shelves, each for 6 MP/PCR plates. Each tower can be equipped with up to 4 shelves (maximum capacity of 24 MP/PCR plates). Accessible with gripper using eccentric fingers (standard RGA or RGA with FES).

Requires an external recirculating cooling/heating system. Depending on the cooling performance, a 400 W cooling unit may be required for per shelf.

Key features

- Stores liquids in micro plates from 6-50 degrees Celsius
- Dimensions: L=701 mm, W=180 mm, H=260 mm
- Positioned in the back of the Fluent, allowing for 5 x ANSI/SBS/SLAS positions in front of the hot-plate
- Flexible setup of the hotel
- No additional ambient heat generated



Part number	Description
30174517	Hotel Cooling/Heating 12 MTP/PCR Plates



Cabinets

- All Fluent systems can be bench-top or free standing
- The cabinet allows use of the space under the Fluent for integration of devices such as microplate readers, washers and shakers etc.
- The cabinet allows access to larger floor mounted devices through the side of the cabinet, such as the Cytomat™ C10*
- Liquid containers and disposables waste can be fitted into pull-out carts for easy installation and removal
- The cabinet has adjustable feet that allow the Fluent to be rolled into position and locked in place, allowing for easy movement around the laboratory
- Door locks can be equipped on each cabinet door for operator safety



Standard cabinet for Fluent extension

Part number	Description
30042015	Standard Cabinet for Fluent 480
30042025	Standard Cabinet for Fluent 780
30042035	Standard Cabinet for Fluent 1080
30042050	Standard Cabinet for Fluent Extension

Azenta® IntelliXcap™

Azenta's IntelliXcap™ is a next-generation automated sample tube capper/decapper, engineered for increased throughput, ease of use and high system reliability. This device is ideal for any laboratory that manages compound libraries or biological sample stores.

The IntelliXcap™ is fast and is compatible with multiple 96-well, 48-well and 24-well plates from manufacturers such as FluidX, Matrix, Thermo-Nunc, Micronic, LVL and Greiner

Automated high-speed capping and decapping reduces sample processing time, increasing the throughput of a workflow. The IntelliXCap™ is an Ideal solution for medium to high throughput laboratories managing compound libraries, biobanks or other biological samples



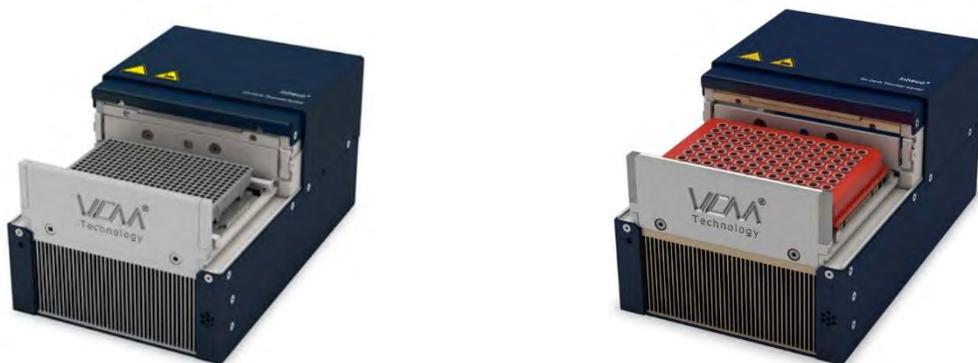
Speak with your local sales representative to receive a quote.



INHECO® On Deck Thermal Cycler®

The On Deck Thermal Cycler (ODTC) is the most compact end-point thermal cycler for use with automated liquid handling workstations. It delivers unrivalled thermal performance with the new VCM technology.

- Temperature range:
 - thermal block: +4 °C to +99 °C
 - heated lid: +30 °C to +115 °C
- Freely programmable temperature profiles and lid temperature
- Horizontally opening heated lid
- Automated adjustment of lid pressure for different disposables and sealing options



Part number	Description	
30112557	384 well plates	Ventilation on the back
30112558		Ventilation on the left side
30101851	96 well plates	Ventilation on the back
30101852		Ventilation on the left side

Visit [Inheco ODTC](#) website for more information

Alpaqua® POGO®

The POGO is a 24-position tube block with integrated Spring Cushion Technology. Constructed from solid aluminium to facilitate temperature control when placed on a temperature-controlled microplate nest position.

Each of the 24 positions is independently equipped with a spring-loaded tube support, capable of moving a full 3 mm down to compensate for variations in Z-positioning of pipette channels.

Each position of the POGO has a diameter of 10.90 mm and a depth of 35.25 mm when the supporting spring is fully compressed.



Part number	Description
30166199	ADAPTER ALPAQUA POGO 24 POS

Visit POGO [Alpaqua](#) website for more information



Qinstruments® BioShake® Series

The BioShake microplate shaker series provides high-speed shaking, heating and cooling functionality in a footprint that is only marginally larger than a ANSI/SBS/SLAS microplate position. The 2-dimensional shaking axis thoroughly mixes samples in a short period of time and the constant orbit reduces the need to centrifuge plates after shaking. The patented Edge Locking Mechanism (ELM) secures the plate during shaking actions and precisely positions the plate afterwards, to allow liquid handling operations in microplates directly upon the Bioshake.

- Options with the ability to heat, cool and/or shake (model-dependent)
- Adjustable shaking speeds (model-dependent)
- Temperature control between -20° and 99.9°C (model-dependent)
- Easily exchangeable plate adapters for optimal heat transfer
- Robust build quality with a housing made of aluminium



BioShake Q1



BioShake 3000 ELM, BioShake 3000 ELM DWP



BioShake 3000-T ELM, BioShake D30-T ELM



ColdPlate

Part number	Description	
30187662	BioShake Q1	Cooling & heating & mixing tubes, microplates and deep well plates
30127732	BioShake 3000-T ELM	Heating & mixing microplates & deep well plates
30127733	BioShake 3000 ELM DWP	Mixing deep well plates
30127731	BioShake 3000 ELM	Mixing microplates
30125516	BioShake D30-T ELM	Heating & mixing microplates & deep well plates
30165954	ColdPlate	Heating & cooling



Visit [Qinstruments](https://www.qinstruments.com) website for more information

Inheco® CPLC Kit

The Inheco CPLC kit is a ready-to-install deck segment with three temperature-controlled ANSI/SBS/SLAS microplate positions. Each CPLC unit can be equipped with a large variety of labware adapters to ensure optimal heat transfer. The CPLC units are liquid cooled with an external heat exchanger unit, reducing heat emissions on the worktable.

- Dimensions W x D x H: 88 mm x 128 mm x 80 mm
- Temperature range from +4°C to 70°C.
- Provided with an electronic module to be mounted in an Inheco TEC controller unit. Tecan standard cable length is 1.60 m (63 inches)
- Different adapters available depending on the choice of labware



Part number	Description
30187625	SET Fluent 3 CPLC INHECO COOL/HEA

Inheco® CPAC

The Inheco CPAC Ultraflat HT 2-TEC unit is a compact all-in-one heating/cooling device. The CPAC can be equipped with a large variety of adapters to suit labware such as microplates, tubes or reservoirs/troughs

- Dimensions W x D x H: 88 mm x 128 mm x 80 mm
- Temperature range from +4°C to 110°C.
- Provided with an electronic module to be mounted in an Inheco TEC controller unit. Tecan standard cable length is 1.60 m (63 inches)
- Accuracy at 37°C target temperature = $\pm 0.3^{\circ}\text{C}$
- Uniformity at 37°C target temperature = $\pm 0.5^{\circ}\text{C}$
- Different adapters available depending on the choice of labware



Part number	Description
30053221	POSITION CPAC ULTRAFLAT COOLING/HEATING



Te-VacS™

The Te-VacS module is an option for automating high-throughput cartridge-based filtration and purification with two independently controlled vacuum positions.

Several vacuum blocks, spacers and adapters are available to make sure the cartridge plate and elution plate can be correctly positioned.

The Te-VacS module can be placed on a grid segment together with two adapters, or alternatively on the Freedom EVO 8-grid deck segment.

The Te-VacS module includes the vacuum pump control tower, a 4 L waste bottle with liquid level sensor, tubing and a user manual. Plate bases, vacuum blocks, adapters, and spacers are sold separately.



Part number	Description
10760680	Option Te-Vacs

Te-Chrom™

Te-Chrom is a module to automate Repligen's OPUS® RoboColumns® for small-scale column chromatography. The Te-Chrom can process 1-96 RoboColumns in a batch and has an integrated plate shuttle that allows for automated fraction collection. Te-Chrom requires the use of a L-FCA, equipped with specialized stainless-steel tips.

- Fully automated column chromatography
- Capable of collecting fractions
- FluentControl® smart command for Te-Chrom available with FC 3.3 and higher (no license required)
- Downholder designed to hold and align short and long columns

Order information

- Te-Chrom is fitted to a 7-grid segment. Three 61 mm Nests can be placed directly behind the Te-Chrom module
- The Te-Chrom segment is equipped with Freedom EVO pin holes, therefore a Te-VacS can be installed behind the Te-Chrom without the need for "Fluent-to-EVO" grid pin adapters
- The plate shuttle spans 20 grids and covers the first nest position



Part number	Description
30201015	Segment Te-Chrom incl. Shuttle III and downholder
30125527	Tip std. stainless steel conical end
30200229	Strip column downholder Te-Chrom
30045303	Spacer 29.9 mm Te-Chrom
30131588	Option pump waste Te-Chrom
30136973	Driver for waste pump (only for FluentControl 3.2 and lower)



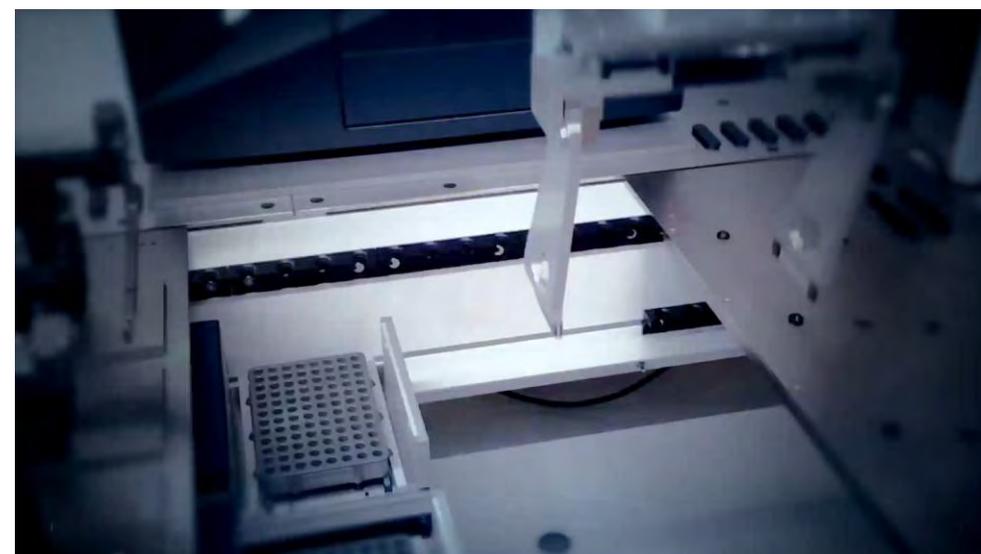
Column base plate and OPUS RoboColumns order directly from [Repligen](#) and their partners.



Hettich® Centrifuge Rotana® 460 Robotic

The Rotanta 460 Robotic is a powerful high-capacity centrifuge that can reach reciprocal centrifugal forces (g-force) of up to 6446 RCF. The Rotanta 460 can hold up to four microplates or up to 80 blood tubes per run. The centrifuge features an integrated refrigerator for cooling of samples during centrifugation. When working with tubes, Tecan Labwerx™ offers automation-friendly centrifugation buckets that can be moved using the Robotic Gripper Arm™ (RGA) of the Fluent liquid handling platform.

- MAX. RCF: 6,446
- MAX. RPM: 6,200 min⁻¹
- WEIGHT: approx. 159 kg
- REFRIGERATION: Yes
- DIMENSIONS (W x D x H): 554 mm x 697 mm x 684 mm



Part number	Description
30126978	CENTRIFUGE HETTICH PLATES 2017 230V EU
30126979	CENTRIFUGE HETTICH PLATES 2017 115V US
30162918	HANGERS CENTRIFUGE 4488-R/4



Frida Reader™

The Frida Reader is an add-on module for the Fluent Automation Workstation that provides UV-based concentration and purity measurements for nucleic acids, from a 'hanging drop' of liquid. This unique, patented method completely avoids sample loss, as the hanging drop is aspirated back into the tip and can be used for further processing (e.g., normalization) or returned to the source location. No additional preparation steps, labware or reagents are required for the measurement, which offers precision and accuracy comparable to a reference reader over a range of 2 ng/μL - 1,000 ng/μL (1 μg / μL).

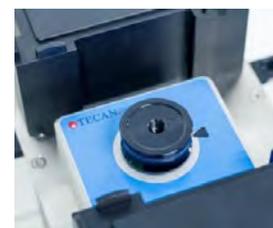
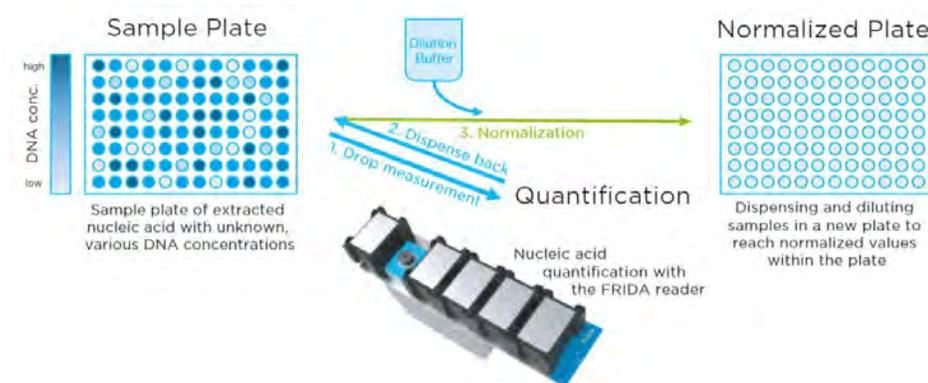
Key features

- No loss of sample
- Only one single ANSI/SLAS/SBS position required on deck
- Eliminate consumables (no extra DiTi or UV measurement plate required)
- No RGA, Fluent cabinet or deck extension for reader integration required
- Can provide quality control after extraction and become the linking element between the Fluent DreamPrep NAP and Fluent DreamPrep NGS
- FluentControl template scripts are available that contain all relevant parameters for a user-friendly quantitation and normalization workflow

Part number	Description
30168273	MODULE FRIDA READER
30042230	SEGMENT NEST FRIDA READER
30042231	DECK TRAY FRIDA READER

Integrated quantification and normalization

The combination of the Frida Reader and Fluent Automation Workstation allows walkaway quantitation and normalization of 96 samples (including blank and blank buffer measurement) in around 40 minutes, quantitation alone for 96 samples in 30 minutes, and quantitation of a single sample in only a few seconds. The illustration below describes the workflow:



MODULE
FRIDA
READER



SEGMENT NEST WITH
MODULE FRIDA
READER



DECK TRAY
FRIDA
READER



Mix and Pierce Workstation (Tube Rotator™ & Tecan Piercing Tips™)

The Mix and Pierce modules allow capped tubes to be handled without removing the cap. Samples can be aliquoted from and/or reagents can be added to tubes and the tubes can then be inverted, rocked, and aliquoted after mixing.

Key advantages:

- Easy operation, even for less experienced users, with TouchTools™
- ~1 hour 15 minutes walkaway time during a 120-sample run
- Automated 1D barcode scanning for full sample traceability
- Optimal extraction condition by oscillating mixing using the Tube Rotator™
- Integrated liquid level detection for high pipetting precision
- Standardized extraction conditions provide reliable results

Unique use case: The Mix and Pierce Workstation enables users to develop workflows for forensics. Crime scene samples are unique, irreplaceable and may be used in legal proceedings. The Mix and Pierce Workstation reduces the risk of sample loss by adding extraction solvents directly to the sample collection tubes and mixing in a controlled manner.

Part number	Description
30042215	Safety Shield Liquid FCA
30042213	Protection tip piercing 100
30042214	Tool tip piercing removal 1 pce
30042217	Kit FCA Channel Piercing Tip
30042210	Tip piercing with lock nut
30042552	Rotator tube 12 grid



Monitored Incubation Option (MIO™)

Precise control of microplate incubation

Our range of incubator options is designed to address the varying environmental conditions and capacity requirements of a wide variety of life-science and diagnostic applications. The MIO allow incubation of either four or six microplates (model dependent) on the work deck, with temperature control from room temperature +5 °C, up to 60 °C. The MIO is available with or without shaking (common to all positions). It is ideally suited to ELISA and cell-based assay applications, and several incubator modules can be used in parallel to provide different temperatures and environmental conditions.

Temperature-controlled carriers are also available for reagent and media reservoirs, tubes and microplates, ensuring consistent assay conditions.

Key features of the MIO

- 4-position Incubator (for 28 mm cell-culture plates)
- 6-position Incubator (for standard microplates)
- Shaking option (model-dependent)
- Temperature range:
 - Ambient +5 °C – 53 °C (Calibrated at 37 °C or 46 °C)
 - Ambient +5 °C – 63 °C (Calibrated at 37 °C or 60 °C)
- Optional fan (for room temperature incubation with shaking)



Part number	Description	
30027295	Calibrated to 37 °C	
30027296	Calibrated to 46 °C	Position for 6 microplates Shaking
30027297	Calibrated to 60 °C	
30027292	Calibrated to 37 °C	
30027293	Calibrated to 46 °C	Position for 6 microplates No shaking
30027294	Calibrated to 60 °C	
30052119	Calibrated to 37 °C	Shaking
30052130	Calibrated to 60 °C	No Shaking
10750752	Optional Fan for RT incubation with MIO2 6 positions	



SciRobotics Pickolo™ Colony Picker

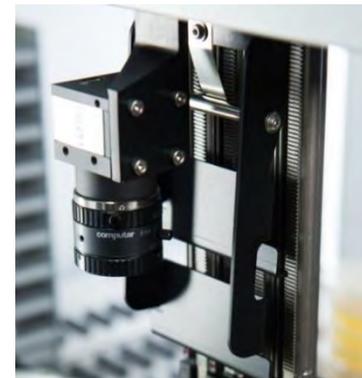
SciRobotics' Pickolo Colony Picker is an add-on for Tecan liquid handling platforms. The Pickolo comes with a 10 megapixel camera that is attached to the Flexible Channel Arm, as well as a backlight illumination table, and powerful software. Add Pickolo to a Fluent and transform it into a powerful and fully-automated colony-picking robot. It can pick:

- Bacteria
- Fungi
- Algae
- Phage
- Mammalian cell colonies
- Hybridoma

... and almost any other small object that can be detected by eye and fits through a standard pipette tip, like zebra fish egg or plant seeds.

- Colony Plate format: 90 mm Petri Dish & ANSI/SBS/SLAS Plates
- Throughput: Fluent system with A-FCA = up to 800 colonies per hour
- A-FCA with 20 μL / 350 μL tips recommended
- Backlight table requires 6 grids
- Three USB slots are needed (camera, light table and software dongle)

Part number	Description
30149954	SciRobotics Pickolo upgrade kit for Fluent
30147067	Adapter Petri dish finger (optional)
30149956	PetReader – barcode reader for Petri dish
30084808	Petri dish stacker 4x15 (capacity = 45 Petri dish, one stack is empty)
30084804	Petri dish stacker 2x15 (capacity = 15 Petri dish, one stack is empty, can be placed in the back or side)



30149954



SciRobotics PetriPlater™

SciRobotics' PetriPlater is a spiral plating and streaking add-on device for Tecan's liquid handling platforms. Microbial plating and streaking is used in many laboratories for either colony counting or colony-picking applications. When preparing serial sample dilutions for plating, the functions of liquid handling and microbial plating integrate seamlessly.

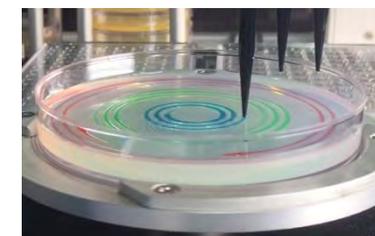
Combine plating with liquid handling for a flexible, fully automated high-throughput solution.

- Process up to 60 plates per hour, unattended
- Use pre-defined flexible plating patterns
- Provide multi-segment plating – saving space, time and money
- Continuous plate loading with the integration of a SciRobotics PetriSel™
- PetriPlater requires 6 grids
- Two USB slots are required (PetriPlater and software dongle)

Part number	Description
30126611	SciRobotics Petri Plater upgrade kit for Fluent and Freedom EVO
30147067	Petri dish gripper adapters for Fluent (optional)
30149956	PetReader – barcode reader for Petri dish
30084808	Petri dish stacker 4x15 (capacity = 45 Petri dish, one stack is empty)
30084804	Petri dish stacker 2x15 (capacity = 15 Petri dish, one stack is empty, can be placed in the back or side)



30126611



SciRobotics PetriSel™

PetriSel is a Petri dish storage add-on for Tecan liquid handling automation platforms, providing an automated and easily-integrated storage solution for up to 180 Petri dishes. The PetriSel™ includes 12 stackers containing up to 1 Petri dishes each.

The PetriSel includes an ultrasonic sensor that automatically detects the quantity Petri dishes in each tower. The robotic arm (with eccentric grippers) can access the top-most stacker position for either returning or removing a Petri dish to / from the stack.

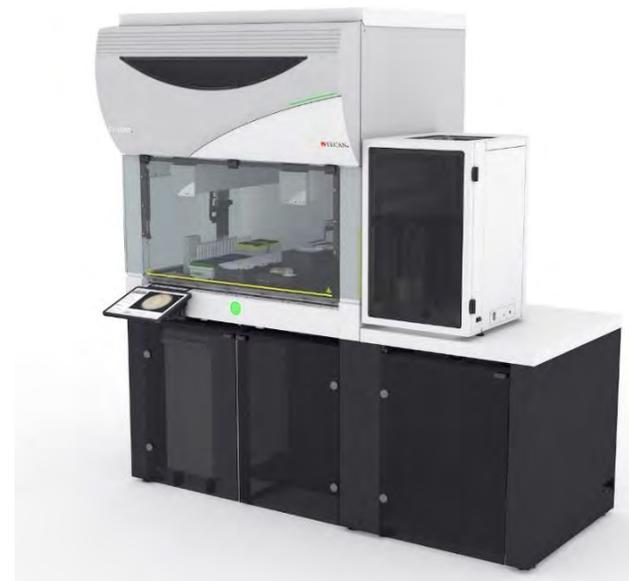
Features

- Up to 180 Petri dishes stacked in 12 towers
- Automatic plate height detection via ultrasonic sensor
- RGA finger adapters designed to handle Petri dishes
- Easy integration with FluentControl
- No random access (only the top-most position can be accessed)

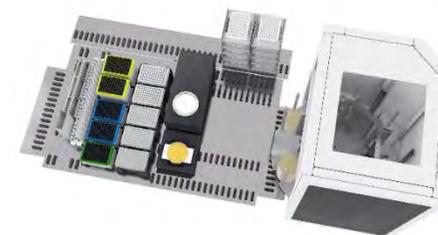
Integration information

- PetriSel cannot be placed directly on an extension deck of a standard cabinet
- The feet of PetriSel require elevation by 103.5 mm above the Fluent feet
- If the Fluent is placed on a standard cabinet with an extension cabinet (Cabinet 340), a plinth of 46.55 mm is required
- PLEASE contact Tecan Labwerx™ with a URD for customization

Visit [SciRobotics PetriSel](#) website for more information



30149955



Part number	Description
30149955	KIT UPGRADE CAROUSEL PETRI-DISH PETRISEL
30147067	Petri dish gripper adapters for Fluent (optional)
30042050	Standard cabinet for Fluent Extension 340
30042667	Deck extension 340
30042711	Segment deck side wall right 87mm



SciRobotics accessories

Multiple accessories are available for picking and plating with SciRobotics's Pickolo™ colony picker and PetriPlater™

Gripper finger adapter

- The transport of Petri dish requires special adapters for the RGA eccentric fingers
- Adapters can be fitted on FES or fixed eccentric fingers

Petri dish storage

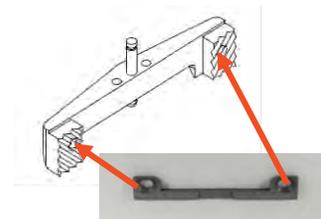
Petri dish passive stacker without random access. One stack of 15 Petri dishes must be empty for returning plates after use. E.g., max capacity of 60 Petri dish stacker = 45 Petri dishes

- 30 Petri dish stacker (2x15) can be placed on the side or in the back of the Fluent deck.
- 60 Petri dish stacker (4x 15) can be placed on the Fluent deck

PetReader™ barcode reader

- 2D barcode reader can be fixed in a laying or sitting position to read barcodes on the bottom or side of a Petri dish
- Can be fixed as stand alone on a Fluent deck, or attached to Pickolo or PetriPlater module

Visit [SciRobotics](https://www.scirobotics.com) website for more information



30147067



30084804



30084808



30149956



Part number	Description
30147067	Adapter Petri dish finger
30084804	30 Petri dish stacker 2x15
30084808	60 Petri dish stacker 4x15
30149956	PetReader – 2D barcode reader for Petri dish



Agilent® PlateLoc thermal microplate sealer

For fast, easy, reliable microplate sealing, the Agilent PlateLoc Thermal Microplate Sealer has distinguished itself as the premier fully automatic thermal sealer. The PlateLoc offers a well-established platform, robust and compact design, flexibility, and versatility with an easy-to-use interface. The PlateLoc Sealer automatically accommodates deep-well, assay, PCR, and compound-storage microplates, and is available with a variety of seals for different applications.

- Fast – Fast cycle times with no cool-down period enable fast plate sealing
- Flexible – Automatically adjusts plate stage to accommodate a wide range of microplates. A total of four types of plate stage inserts are included to provide support for even the most challenging microplate types
- Energy efficient – Sleep mode shuts off the screen and hotplate after a designated time of inactivity to save energy
- Integration – Integrated on the side of a Fluent, this device will not occupy space on the work deck

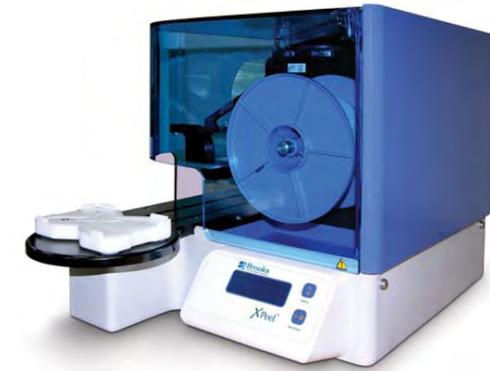


Speak with your local sales representative to receive a quote.

Visit [Agilent PlateLoc](#) website for more information

Azenta® XPeel automated plate seal remover

- The Azenta XPeel automated plate seal remover is compatible with an extensive range of microplate and seal types
- The patented XTape removal medium eliminates the need for mechanical removal mechanisms, which are often prone to failure



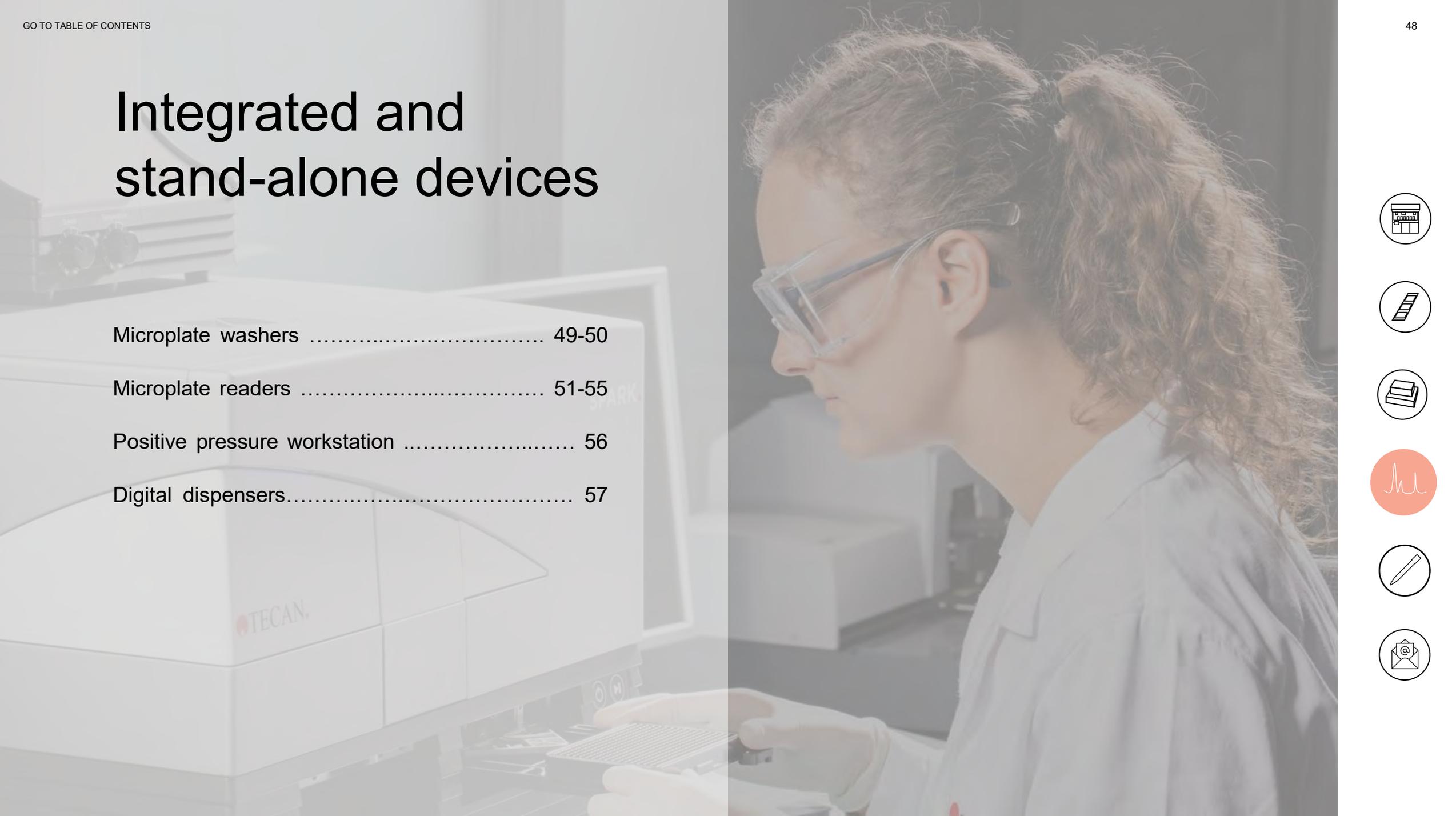
Speak with your local sales representative to receive a quote.

Visit [Brooks XPeel](#) website for more information



Integrated and stand-alone devices

Microplate washers	49-50
Microplate readers	51-55
Positive pressure workstation	56
Digital dispensers.....	57



HydroFlex™

Cells. ELISAs in 96-well plate format

The HydroFlex microplate washer is a truly flexible and upgradeable platform for automated microplate-strip washing, magnetic-bead washing and vacuum filtration in 96-well plate formats. It is an ideal solution for a wide range of cell-, bead- and ELISA-based applications, reflecting over 30 years of Tecan expertise in advanced liquid handling.



Every HydroFlex washer comes with microplate shaking and automated buffer switching (when two or more input channels are selected) included, ready for ELISA washing and gentle washing of cells. In addition, optional user-interchangeable plate carriers are available for easy switching between washing, bead separation, and filtration applications.

Features and benefits

- Ready for automated washing of ELISAs, cells and protein arrays
- Magnetic-bead washing using powerful neodymium magnets
- Automated system for low-level vacuum filtration using vacuum sensor
- Control of wash head positions and speed settings for gentle cell washing
- Up to two aspiration points per well for minimal residual volumes
- Automated buffer switching (when at least two input channels are configured) and plate shaking capabilities
- Controls up to four wash buffers
- Ready for robotic integration
- Designed to meet the IVDR (EU) 2017/746 (HydroFlex Plus)
- Supports FDA 21 CFR Part 11 if used in combination with HydroControl™ software

Part number	Part name
30025471	HYDROFLEX
30190374	HYDROFLEX PLUS

Speak to your sales rep about configuration of this device.

For more information visit our [website](#).



HydroSpeed™

For cells, beads and ELISAs in 96 & 384-well plates

The HydroSpeed is Tecan's advanced microplate washer, designed for quick, parallel washing of entire 96- and 384-well plates. Washing all wells in parallel eliminates variation between wells for uniform results. The HydroSpeed provides full control of wash power (aspiration rate and dispense speed), as well as customized head positioning, for optimized washing of cells, beads and ELISAs.



The HydroSpeed plate washer is ready for ELISA washing, as well as gentle washing of adherent and weakly-adherent cells. It supports soak steps with microplate shaking, as well as buffer switching. With full on-board operation via the easy-to-use touchscreen interface, the HydroSpeed offers incredible flexibility and control for cell, bead and ELISA washing applications.

Features and benefits

- Adjustable wash head positions, speed settings and vacuum level for extra gentle cell washing
- Anti-clogging - active prevention of needle clogging
- Easy X-change wash heads for intense cleaning in ultrasonic bath
- Patented design with two rare-earth magnets per well enables fast bead settling and high bead recovery rates
- Automated buffer switching of four liquids and plate shaking capability
- Up to four aspiration points per well for minimal residual volumes
- Automated system for vacuum filtration using vacuum sensor
- Individual vacuum settings for each filtration step within a protocol
- pLLD for maintenance-free liquid level detection in the waste bottle
- Robotic integration requires "OPTION ROBOTIC HYDROSPEED" (30060646), ordered separately
- Plate assistant function for easy teaching of new plates
- Supports FDA 21 CFR part 11 if used in combination with HydroControl™ software

Part number	Part name
30063691	HYDROSPEED

Speak to your sales rep about configuration of this device.

For more information visit our [website](#).



Spark®

The Spark® multimode microplate reader - your research partner.

The Spark multimode microplate reader offers solutions for virtually any life science research or drug discovery applications. Spark is a modular platform that is built to specification, but is also able to be upgraded at a later time. The Spark is an instrument that can meet the customer's needs both today and in the future.



Part number	Part name
30086261	SPARK

Speak to your sales rep about configuration of this device.

For more information visit our [website](#).

Spark microplate readers are for research use only. Not for use in diagnostic procedures.

Capabilities

- High-speed monochromators for free wavelength selection and fast wavelength scanning (from 200 nm to 1,000 nm)
- Monochromator- or filter-based fluorescence measurements, or
- Unique Fusion optics combines the flexibility of monochromators with the sensitivity of filters – for all fluorescence measurements
- Bright-field cell-imaging option for counting of a broad range of cells in suspension, using Tecan's Cell Chip™
- Luminescence scanning and multicolor luminescence measurements
- Fully-integrated Gas Control Module (GCM™) with CO₂ and O₂ regulation for cell kinetic studies
- Humidity cassette for evaporation protection, and integrated lid lifter system
- Patented Te-Cool™ module for active temperature regulation from 18°C to 42°C
- Te-Cool empowers optimal performance of enzymatic assays, Alpha Technologies® assays, or kinetic assays using live organisms
- Automated conditional workflows minimize hands-on time and help to increase reproducibility for long term experiments
- Supports 6- to 1,536-well plates with enhanced detection modules
- Bright-field imaging option supports 6- to 384-well plates



Spark® Cyto

Plate reader with live cell imaging and real-time cytometry.

Spark Cyto is a multi-mode microplate reader with fluorescence imaging and cytometry capabilities, unlocking new possibilities for cell-based research. By combining live-cell imaging with industry-leading detection technologies, the Spark provides the ability to unite qualitative and quantitative information into unique multi-parameter datasets.

Spark Cyto works in real-time, using parallel data acquisition and analysis to deliver meaningful insights faster.



Part number	Part name
30167885	SPARK CYTO

Speak to your sales rep about configuration of this device.

Capabilities

Imaging applications

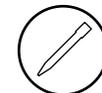
- Label-free cell counting
- Nuclei counting
- Transfection efficiency
- Cell viability
- Apoptosis
- Confluence assessment

Detection modes

- Fluorescence imaging (blue, green, red, far red)
- Bright field imaging
- Digital phase contrast-imaging
- Absorbance – incl. UV/VIS
- Fluorescence top and bottom

For more information visit our [website](#).

Spark Cyto is for Research Use Only. Not for use in diagnostic procedures.



Infinite® 200 PRO Series

The Infinite 200 PRO is an easy-to-use multimode plate reader family that offers affordable, high performance detection solutions, referenced in more than 1800 peer reviewed publications.

Based on this customer experience and feedback, the Infinite 200 PRO microplate plate reader has been continuously engineered for excellence. This proven platform has now evolved into a budget-friendly reader family consisting of six application-focused configurations, delivering performance and value to every lab, every day, for every assay.



Part number	Part name
30108056	INFINITE FILTER 200
30111058	INFINITE MONO 200

Speak to your sales rep about configuration of this device.

For more information visit our [website](#).

Capabilities

- Quad4 Monochromators for free wavelength selection
- High wavenlength accuracy and reproducibility
- Spectral scanning in both absorbance and fluorescence reading modes
- Optimized detectors for absorbance, fluorescence and luminescence
- Glow, flash and dual-color luminescence with NanoBRET™ filters
- z-focussing as integrated background correction
- Field-upgradeable injector module
- 3-in-1 reader for microplates (6-384 well), NanoQuant™ plate and cuvettes

Applications:

- Nucleic acid quantification and purity checks
- Protein quantification
- ELISA
- Cell based assays such as MTT-assay and LDH-release
- Growth curves and proliferation studies with microorganisms
- HTRF®, DELFIA®, LanthaScreen
- DLR®
- BRET (including NanoBRET™)



Infinite® F50

The Infinite F50 and Infinite F50 Robotic absorbance plate readers, together with Magellan™ data analysis software, are ideal solutions for photometric measurements of ELISAs and kinetic assays in the 400 nm to 750 nm wavelength range. The systems' innovative design combines eight measurement channels with long-life LED lights for fast parallel reading and minimal maintenance costs.



Features and benefits

- Ideal for absorbance-based ELISAs, kinetic and multi-label reading
- Ready for qualitative and quantitative data analysis using Magellan data analysis software
- Magellan's example files and tutorials enable you to get started easily
- 405 nm, 450 nm, 492 nm and 620 nm filters included (additional absorbance filters can be purchased separately)
- Eight-channel optics for fast reading + reference channel for optimized light intensity
- Long-life LED lights offer a ten-fold increase in maintenance-free operation, compared to conventional halogen lamps
- Two year warranty and tested reliability of >100,000 plates
- Linear plate shaking
- Designed to meet IVDR (EU) 2017/746 (Infinite F50 Plus)
- Provides FDA 21 CFR Part 11 functionality for regulated environments via Magellan Tracker software
- MultiCheck™ QC plate for IQ/OQ



Speak to your sales rep about configuration of this device.

For more information visit our [website](#).

Part number	Part name
30051599	INFINITE F50
30183570	INSTRUMENT INFINITE F50 PLUS

Sunrise™

Advanced. Reliable. Validated.

The Sunrise absorbance reader, complemented by Tecan's Magellan™ universal reader control and data analysis software, is an ideal solution for a range of applications, such as ELISAs, enzyme kinetics, cell viability studies, veterinary assays, and many more. It supports a broad wavelength range – from 340 nm to 750 nm – and can measure an entire 96-well plate in less than six seconds, ensuring it is ready to meet the needs of routine diagnostic and research labs.



For more information visit our [website](#).

Part number	Part name
30001687	SUNRISE

Features and benefits

- Ideal for colorimetric ELISAs, kinetic assays and agglutination measurements
- Temperature control option for enzyme kinetics
- Tunable wavelength option for greater flexibility and wavelength scanning
- Powered by Magellan data analysis software
- Advanced curve fitting and QC functions, plus data export to your LIS
- Advanced 12-channel optics for fast results + additional reference channel for optimizing light intensity
- Prolonged lamp lifetime via auto switch-off function when not in use
- Microplate shaking capability
- Ready for robotic integration
- Three-year warranty, demonstrating Tecan's commitment to quality
- IQ/OQ tools and procedures available
- Provides FDA 21 CFR Part 11 in combination with Magellan Tracker software



Speak to your sales rep about configuration of this device.

Resolvex® A200

The Resolvex A200 is a compact, industry-leading, that automates positive pressure and solid phase extraction (SPE) with exceptional efficiency. It demands minimal manual intervention while granting complete workflow automation. The spectrum of applications it facilitates is vast, encompassing protein desalting, nucleic acid extraction, and beyond. The adoption of positive pressure workflows yields a cascade of benefits: cleaner samples, heightened accuracy, precision, and method robustness, and an extended lifespan of your analytical instrument.

The system effortlessly takes on two distinct roles: a stand-alone semi-automated processing solution, or when harmonized with Tecan's liquid handlers, a fully automated powerhouse. The choice of a 96-channel or a 24-channel configuration amplifies its capabilities, rendering the process of 1 ml SPE columns and 96-well plates as effortless as handling larger volumes within 3 ml, 6 ml, and 24-well SPE labware.

Venturing further, the Resolvex Kit Upgrade Omics integrate a HEPA filter air system, effectively curbing particulate contamination – a boon for applications sensitive to airborne pollutants such as keratin.

Part number	Part name
253-1160	Resolvex A200 (96-channell format)
30199839	Resolvex A200 Omics (96-channel with HEPA filter)
30219820	Resolvex A200-24 (24-channel format)
30189741	Resolvex Kit Upgrade Omics (Hepa filter air system)

Applications

- Bioanalysis
- Analyte clean-up and enrichment for therapeutic drug monitoring, drug screening and other clinical chemistry workflows
- Forensic and toxicology testing
- Nucleic acid screening and purification
- Life sciences research
- Food safety testing
- Environmental monitoring



Key features of the A200

- Automated liquid dispensing of up to 11 connected solvents
- Positive pressure manifold using pressurized N2 gas or compressed air
- Intuitive touchscreen user interface
- Light curtain ensuring operator safety
- High-capacity waste reservoir with direct drain to waste
- Sensor controlled movements to ensure precision solvent dispensing
- Several spacers available to accommodate many SPE labware formats
- Standalone system or integrated into liquid handler

Visit our [website](#) for more information. Full catalogue with SPE labware and instrumentation available upon request.

Tecan digital dispenser D300e

The Tecan digital dispenser D300e is a bench-top device to simplify the generation of dose-response curves, enzyme profiles, synergy experiments, and PCR reactions. Using HP's direct digital dispensing technology, it provides picoliter to microliter non-contact dispensing of liquids directly into the assay plate, saving time, minimizing consumption of valuable samples and accelerating research.

Key features

- Precise dispensing: Dispense liquids in picoliter to microliter volumes directly into assay plates, eliminating the need for serial dilutions.
- Time-saving: Generate dose-response curves directly from stock solutions, reducing the time spent on manual processes.
- Resource optimization: Lower consumption of precious samples and accelerate research processes.

Part number	Part name
30100152	DIGITAL DISPENSER D300E
30097370	CASSETTE DISPENSEHEAD HP T8 PLUS 20PC
30097371	CASSETTE DISPENSEHEAD HP D4 PLUS 20PC

Tecan digital dispenser D300e:



Dispensehead cassettes:



30097370

30097371

Visit our [e-shop](#) to order and our [website](#) for more information.

For research use only. Not for use in diagnostic procedures. Tecan D300e Digital Dispenser is a product of HP Inc. CA, USA



Uno Single Cell Dispenser™

Tecan's Uno Single Cell Dispenser is an automated benchtop instrument designed for ease of use, efficiency, and precision in single-cell workflows. Uno empowers a wide range of single-cell applications across the “-omics” landscape, including MS-based single-cell analysis, iPSC libraries, 3D cell research, cell-line development, and beyond.

With Uno, you can achieve accurate, viable, and precise single-cell dispensing, managing reagent volumes from picoliters to microliters within minutes.



C1a – Single Cell
Dispensehead
Cassette



D1 sf / D1 Reagent
Dispensehead
Cassette



T1 sf / T1 Reagent
Dispensehead
Cassette

Part number	Part name
30230840	INSTRUMENT DISPENSER SINGLE CELL UNO
30230841	CASSETTE DISPENSEHEAD C1A 20PC
30230842	CASSETTE DISPENSEHEAD T1 30PC
30230843	CASSETTE DISPENSEHEAD D1 30PC
30230844	CASSETTE DISPENSEHEAD T1 SF 20PC
30230845	CASSETTE DISPENSEHEAD D1 SF 20PC

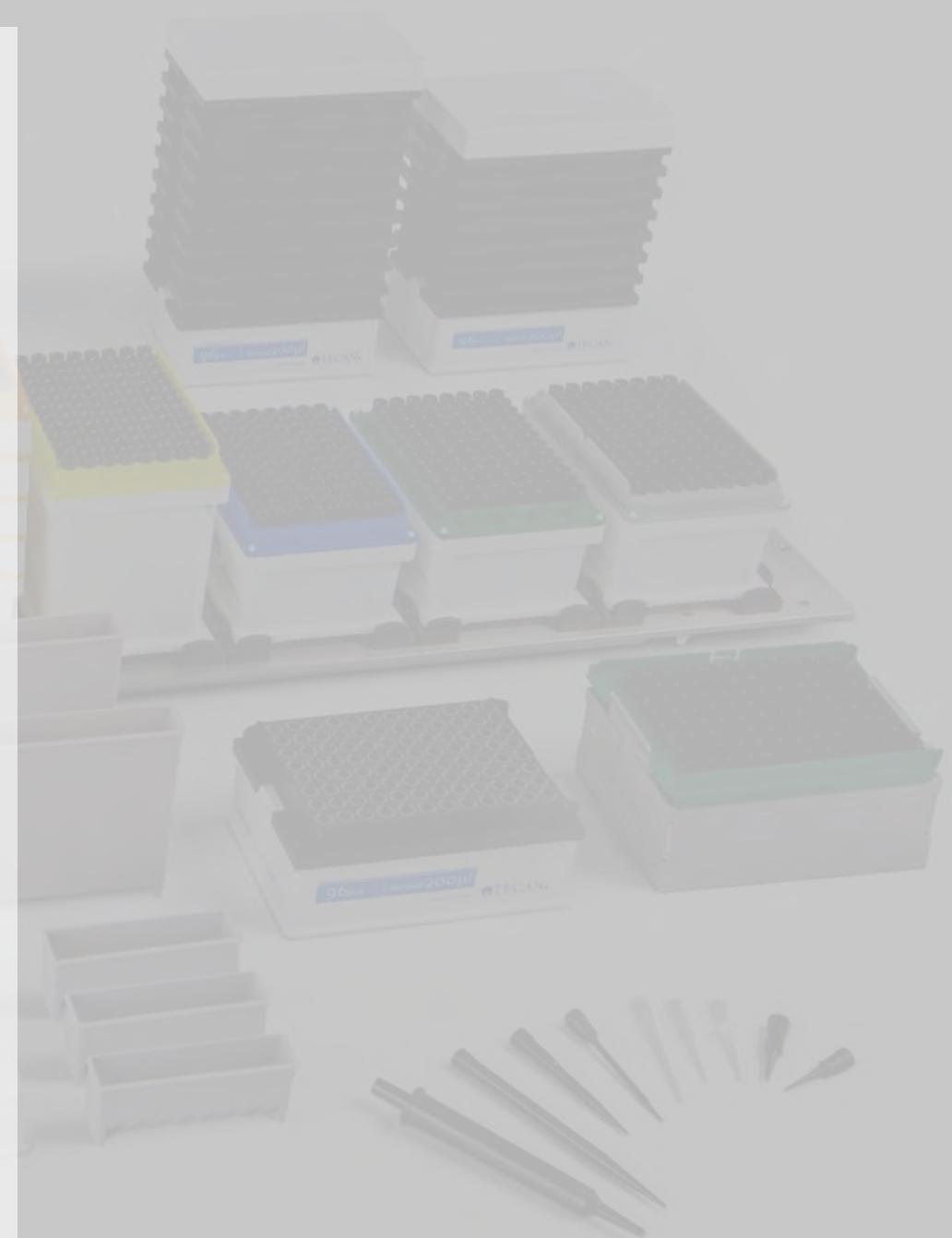
Visit our [e-shop](#) or our [website](#) for more information.

For research use only. Not for use in diagnostic procedures. Uno single cell dispenser is a product of HP Inc. CA, USA



Plastic consumables

Disposable tips	60
LiHa tips	60-62
MCA384 tips	63
MCA96 tips	63
Labware	64
Troughs	64



Liquid handling (LiHa) tips

Carefully designed, produced and validated to ensure reliable performance using Liquid Handling (LiHa) Arms and Flexible Channel (FCA) Arms on Tecan platforms.

Available in a variety of tip volumes, with or without filters, and made of conductive polypropylene to allow capacitive liquid level detection. Available in three purity levels (Tecan Standard, Tecan Pure and Tecan Sterile), tailored towards specific applications (and price points).



Hanging / Standard tray

Part number	Part name	Purity
30104803	DITI LIHA 10µL CONDU. 2304 PCE.	Pure
30104804	DITI LIHA 10µL CONDU.FIL. 2304 PCE.	Pure
30000627	DITI LIHA 200µL CONDU.2304 PCE.	Pure
30000629	DITI LIHA 200µL CONDU.FIL. 2304 PCE.	Pure
30000630	DITI LIHA 1000µL CONDU.2304 PCE.	Pure
30000631	DITI LIHA 1000µL CONDU.FIL.2304 PCE.	Pure
30200712	DITI LIHA 50µL CONDU.FIL. 2304 PCE.	Pure
30032114	DITI LIHA 50µL CONDU.FIL. 2304 PCE.	Pure
30032115	DITI LIHA 50µL CONDU.2304 PCE.	Pure
30115239	DITI LIHA 1000UL COND.FIL. 960 PCE WIDE	Pure



Combi Tray tips

Part number	Part name	Purity
30201339	DITI COMBI 50µL CONDU.FIL. 2304 PCE.	Pure
30201341	DITI COMBI 200µL CONDU.FIL. 2304 PCE.	Pure
30199162	DITI COMBI 1000µL CONDU.FIL. 2304 PCE.	Pure
30201344	DITI COMBI 1000µL COND.FIL.WB. 2304 PCE.	Pure

Combi Tray™ tips are compatible with the following arms:
LiHa, FCA; Multiple Channel 96.

Not compatible with: Multi Channel Arm 96, Multiple Channel Arm 384,
Multiple Channel Arm 384 w/ EVA 96 adapter

Also available in Tecan standard purity – however for application security its preferred to recommend Pure or Sterile products



Liquid handling (LiHa) tips

ANSI/SBS/SLAS refill – box is refilled with tray of tips



ANSI/SBS/SLAS tray in blister

ANSI/SBS/SLAS box

Part number	Part name	Purity
30057811	DITI LIHA 50µL CONDU. 2304 PCE. SBS	Pure
30057813	DITI LIHA 50µL CONDU.FIL. 2304 PCE. SBS	Pure
30057814	DITI LIHA 200µL CONDU. 2304 PCE. SBS	Pure
30057815	DITI LIHA 200µL CONDU.FIL. 2304 PCE. SBS	Pure
30057816	DITI LIHA 1000µL CONDU. 2304 PCE. SBS	Pure
30057817	DITI LIHA 1000µL CONDU.FIL. 2304PCE. SBS	Pure
30058506	LIHA DITI SBS BOX REFILL SMALL 10PCE.	Pure
30058507	LIHA DITI SBS BOX REFILL LARGE 10PCE.	Pure
30104973	DITI LIHA 10µL COND. 2304 PCE. SBS	Pure
30104974	DITI LIHA 10µL COND.FIL. 2304 PCE. SBS	Pure

ANSI/SBS/SLAS Sterile

Comes with tray and tip box combined due to higher purity requirements



Part number	Part name	Purity
30057818	DITI LIHA 50µL CONDU. STE. 3840 PCE.	Sterile
30057819	DITI LIHA 50µL CONDU.FIL.STE. 3840 PCE.	Sterile
30057820	DITI LIHA 200µL CONDU.STE. 3840 PCE.	Sterile
30057821	DITI LIHA 200µL CONDU.FIL.STE. 3840 PCE.	Sterile
30057822	DITI LIHA 1000µL CONDU.STE. 3840 PCE.	Sterile
30057823	DITI LIHA 1000µL CONDU.FIL.STE. 3840 PCE	Sterile
30104976	DITI LIHA 10µL COND.FIL.STE. 3840 PCE.	Sterile
30104975	DITI LIHA 10µL COND. STE. 3840 PCE.	Sterile

ANSI/SBS/SLAS Nested



Part number	Part name	Purity
30083400	DITI LIHA 350µL COND.NESTED 7680 PCE	Pure
30083401	DITI LIHA 350µL COND.STE.NESTED 7680 PCE	Sterile
30104977	DITI LIHA 10µL COND.NESTED 7680 PCE	Pure
30104978	DITI LIHA 10µL COND.NESTED FIL.7680 PCE	Pure
30104979	DITI LIHA 10µL COND.NESTED.FIL.STE.7680	Sterile

5 mL Large volume tip



Part number	Part name	Purity
30059897	DITI LIHA 5000µL CONDU. 240 PCE.	Pure
30059898	DITI LIHA 5000µL CONDU.FIL.STE. 240 PCE.	Sterile
30065423	DITI LIHA 5000µL CONDU.FIL. 240 PCE.	Pure



Liquid handling (LiHa) tips

Disposable Tip (DiTi) Boxes, empty, for use with refill insert tips, ANSI/SBS/SLAS format. Tecan Pure purity level – tested and certified to be free of human DNA, Rnase, Dnase and PCR inhibitors. For use with the Freedom EVO Liquid Handling Arm (LiHa), Fluent Flexible Channel Arm (FCA) or Fluent Multiple Channel Arm 96.

Large ANSI/SBS/SLAS box



Part number	Part name	Purity
30058507	LIHA DITI SBS BOX REFILL LARGE 10PCE.	Pure

For ANSI/SBS/SLAS refill tip sizes:

1000µL

1000µL WB

Small ANSI/SBS/SLAS box



Part number	Part name	Purity
30058506	LIHA DITI SBS BOX REFILL SMALL 10PCE.	Pure

For ANSI/SBS/SLAS refill tip sizes:

10µL

50µL

200µL



MCA 384 tips

Proprietary sealing system

The advanced design of MCA 384 disposable tips incorporates multiple stabilization guides and flexible sealing rings to provide excellent seal integrity, without the use of O-rings. This proprietary sealing system ensures accurate and reproducible low-volume pipetting across a broad volume range, offering a reliable solution to maximize productivity, while reducing maintenance requirements.



Part number	Part name	Purity
30138080	DITI 15UL 40X384P MCA384	Pure
30138085	DITI 15UL STE. 40X384P MCA384	Sterile
30138086	DITI 15UL STE. FIL. 40X384P MCA384	Sterile
30138087	DITI 50UL 40X384P MCA384	Pure
30138088	DITI 50UL STE. 40X384P MCA384	Sterile
30138089	DITI 50UL STE. FIL. 40X384P MCA384	Sterile
30138090	DITI 125UL 40X384P MCA384	Pure
30138091	DITI 125UL STE. 40X384P MCA384	Sterile
30138092	DITI 125UL STE. FIL. 40X384P MCA384	Sterile

MCA 96 tips

Designed, validated and manufactured to ensure optimal performance for the MultiChannel Arm 96. Available in four volumes – 50 µL, 100 µL, 200 µL, and 500 µL. Supplied in ANSI/SBS/SLAS-format racks, allowing up to 768 tips per nest (stack of 8 trays) on the worktable, or 3,840 tips in a Te-Stack™ column.



Part number	Part name	Purity
30038606	DITI 50µL 3840 PCE. MCA96 SBS	Pure
30038607	DITI 50µL STE. 3840 PCE. MCA96 SBS	Sterile
30038608	DITI 50µL STE.FIL. 3840 PCE. MCA96 SBS	Sterile
30038609	DITI 50µL 3840 PCE. MCA96 NESTED	Pure
30038610	DITI 50µL 7680 PCE. MCA96 TE-STACK SBS	Pure
30038611	DITI 100µL 3840 PCE. MCA96 SBS	Pure
30038612	DITI 100µL STE. 3840 PCE. MCA96 SBS	Sterile
30038613	DITI 100µL STE.FIL. 3840 PCE. MCA96 SBS	Sterile
30038614	DITI 100µL 3840 PCE. MCA96 NESTED	Pure
30038616	DITI 200µL 3840 PCE. MCA96 SBS	Pure
30038617	DITI 200µL STE. 3840 PCE. MCA96 SBS	Sterile
30038618	DITI 150µL STE.FIL. 3840 PCE. MCA96 SBS	Sterile
30038619	DITI 200µL 3840 PCE.MCA96 NESTED	Pure
30038620	DITI 200µL 7680 PCE. MCA96 TE-STACK SBS	Pure
30046341	DITI 500µL 4800 PCE. MCA96 SBS	Pure
30046342	DITI 500µL STE.FIL. 4800 PCE. MCA96 SBS	Sterile
30048822	DITI 50µL STE. 3840 PCE. MCA96 NESTED	Sterile
30048823	DITI 100µL STE. 3840 PCE. MCA96 NESTED	Sterile
30048824	DITI 200µL STE. 3840 PCE. MCA96 NESTED	Sterile
30050348	DITI 200µL 3840 PCE. MCA96 WIDE BORE SBS	Pure
30050349	DITI 150µL 3840 PCE.FIL. MCA96 WIDE BORE	Sterile



Troughs

Tecan offers a collection of reagent troughs of varying volumes and shapes that are designed for the liquid handling arms on Fluent and Freedom EVO Liquid Handling Platforms.

Compatibility with strict application needs is guaranteed by our certified purity levels.



Part number	Part name	Purity
10613048	TROUGH DISPOSABLE 100ML PP TRA. 108 PCE.	Standard
10613049	TROUGH DISPOSABLE 100ML PP GREY 108 PCE.	Sterile
30055743	TROUGH DISPOSABLE 25ML PP 120PCE.	Pure



Ordering information

There are three ways to get in contact with us:

[Connect with Tecan: Your Next Step in Lab Excellence](#)

If you do not have a local Tecan sales representative yet, fill out the form and we will contact you shortly.

Visit Tecan e-shop

ORDER
ONLINE



You can buy a selection of devices and consumables directly on the Tecan e-shop :
<https://shop.tecan.com/>

Already have a contact? Contact your sales representative directly.



This Tecan product catalog contains a collection of Fluent modules, deck segments and devices. If something additional is needed, please contact us.



Tecan – Who we are

At Tecan we are driven to improve people's lives and health. We do this by empowering our customers to scale healthcare innovation globally from life science to the clinic. Tecan is a pioneer and global leader in laboratory automation. As an original equipment manufacturer (OEM), Tecan is also a leader in developing and manufacturing OEM instruments, components and medical devices that are then distributed by partner companies. Founded in Switzerland in 1980, the company has more than 3,300 employees, with manufacturing, research and development sites in Europe, North America and Asia, and maintains a sales and service network in over 70 countries. Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TECN; ISIN CH0012100191).

Tecan Group Ltd. makes every effort to include accurate and up-to-date information, 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. Changes 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 please contact your Tecan representative. This may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

© Tecan Trading AG, Switzerland,
all rights reserved.

www.tecan.com

Australia

+61 39 647 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

402731 V1.0, 2023-09

