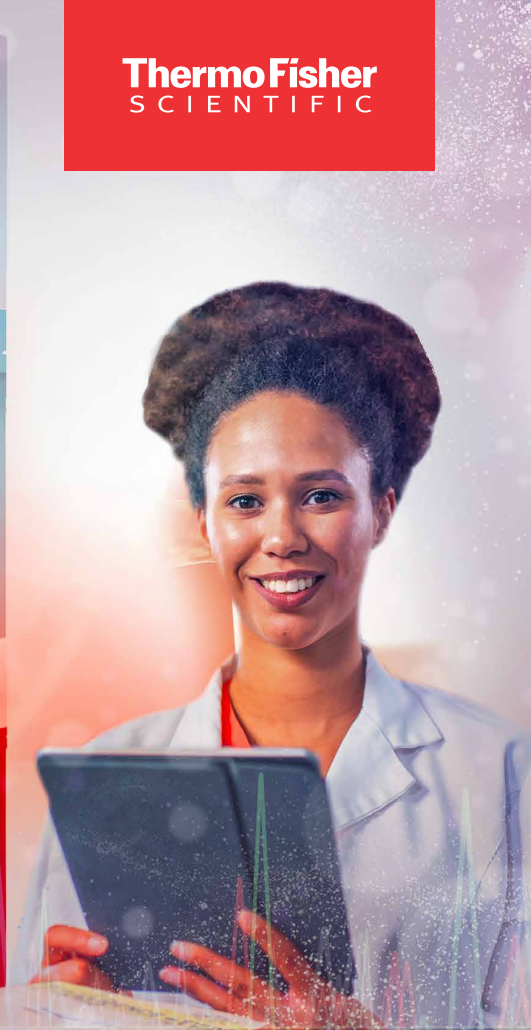




Automated cell-free DNA purification



## Automate your cfDNA isolation workflows with a high-throughput integrated solution

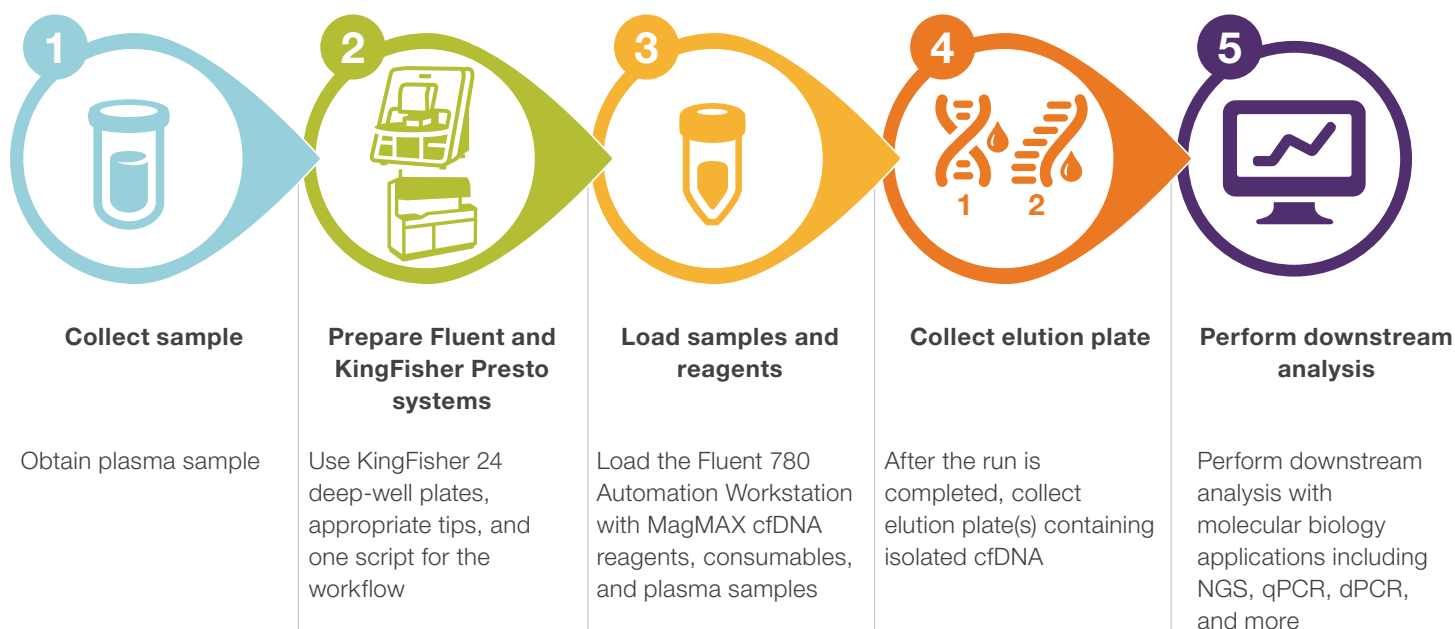
Isolate high-quality cfDNA to help save time by streamlining your workflow and reducing multiple manual steps

Integrate the Thermo Scientific™ KingFisher™ Presto Purification System into a Tecan® Fluent® 780 Automation Workstation to achieve automated workflows without the need to prepare plates and reagents for sample preparation (Figure 1). Reduce operator hands-on time with workflows optimized to be consistent and reproducible without sacrificing sample quality.

The Applied Biosystems™ MagMAX™ Cell-Free DNA Isolation Kit can be leveraged on a high-throughput integrated solution, i.e., the Fluent 780 Automation Workstation with the KingFisher Presto instrument, for isolation of cell-free DNA (cfDNA). Optimized and qualified scripts are available to assist you in setting up your high-throughput workflow.



Fluent 780 Automation Workstation and  
MagMAX Cell-Free DNA Isolation Kit



**Figure 1. Basic workflow of high-throughput cfDNA isolation.** The KingFisher Presto system is integrated into the Fluent 780 Automation Workstation and used with the MagMAX Cell-Free DNA Isolation Kit (Cat. No. A29319).

## Chemistry specifications

Description	
Qualified sample prep kit	MagMAX Cell-Free DNA Isolation Kit
Throughput	24–96
Sample type	Plasma (4 mL)
Sample input format	KingFisher 24 deep-well plate (8)
Sample output format	KingFisher 24 deep-well plate (4)

## Instrument specifications

Description	
Automation workstation	Fluent 780 system
System integration	KingFisher Presto system (2)
Required arm(s)	Flexible Channel Arm™ (FCA), 8-channel Robotic Gripper Arm™ (RGA)
Dimensions (D x W x H)	78.5 x 165 x 123.6 cm (30.9 x 65.0 x 48.6 in.)
Dimensions (D x W x H) with Kingfisher Presto system integration	107.2 x 165 x 123.6 cm (42.2 x 65.0 x 48.6 in.)
Weight (not including the KingFisher Presto system)	~308.6 lb (140 kg)
Warranty	1 year included, extended warranties available

## Operation

Description	
Power input	100–240 V AC, 50–60 Hz, 5 A
Operating temperature	15–32°C (59–90°F)
Relative humidity	30–80% noncondensing

## Computer\* and software specifications

Description	
Operating system	Microsoft™ Windows™ 10 Enterprise LTSC 2019 (version 1809)
Processor	Intel™ Core™ i7 processor or equivalent
Memory	32 GB RAM
Hard drive	Solid-state hard drive with 512 GB free space
Graphics card	NVIDIA™ GPU with minimum 5 GB RAM and CUDA™ node support
Network card	1–2 GigE network cards with Intel™ chipsets
Communication	1 USB port
Monitor	Vertical resolution of 1,024 pixels

\* Minimum specification.

## Performance

Flexible Channel Arm—air displacement pipetting										
Gravimetric measurement using deionized water in single pipetting mode under optimized conditions										
Tecan® Disposable Tip (DiTi) size	DiTi 10		DiTi 50		DiTi 200		DiTi 350		DiTi 1,000	
Volume (μL)	CV	ACC	CV	ACC	CV	ACC	CV	ACC	CV	ACC
0.5	≤6%	±9.5%	–	–	–	–	–	–	–	–
1	–	–	≤4%	±8%	–	–	–	–	–	–
10	≤1%	±2%	≤0.5%	±1%	≤2%	±2%	≤2%	±2%	≤1.2%	±2%
100	–	–	–	–	–	–	–	–	≤0.3%	±0.5%
200	–	–	–	–	≤0.2%	0.5%	0.2%	0.5%	–	–

The combination of the Fluent 780 Automation Workstation and KingFisher Presto system, developed through a collaboration between Tecan and Thermo Fisher Scientific, is a comprehensive and unified solution. The combination workstation optimizes automated nucleic acid extractions from biological samples like blood, plasma, saliva, and plasmids in bacteria. Utilizing magnetic

bead-based processing in both 24- and 96-well formats, the combined workstation of the Fluent 780 and KingFisher Presto systems can easily be adapted to evolving lab requirements while simultaneously delivering high-quality results, all within a compact footprint—excellent for preserving valuable laboratory space.

## Ordering information

Description	Quantity	Cat. No.
KingFisher Presto Purification System, 96 deep-well head	1 each	5400830
KingFisher Presto Purification System, 24 deep-well head	1 each	5400840
MagMAX Cell-Free DNA Isolation Kit	1 kit	A29319
KingFisher 96 deep-well plate	Case of 50	95040450
KingFisher 96 tip comb for deep-well magnets	Case of 100	97002534
KingFisher 24 deep-well plate	Case of 50	95040470
KingFisher 24 deep-well tip comb and plate	Case of 50	97002610

 For more information about Tecan's Fluent Automation Workstation, visit [lifesciences.tecan.com/fluent-laboratory-automation-workstation](https://lifesciences.tecan.com/fluent-laboratory-automation-workstation)

 Learn more at [thermofisher.com/magmaxcfdna](https://thermofisher.com/magmaxcfdna)