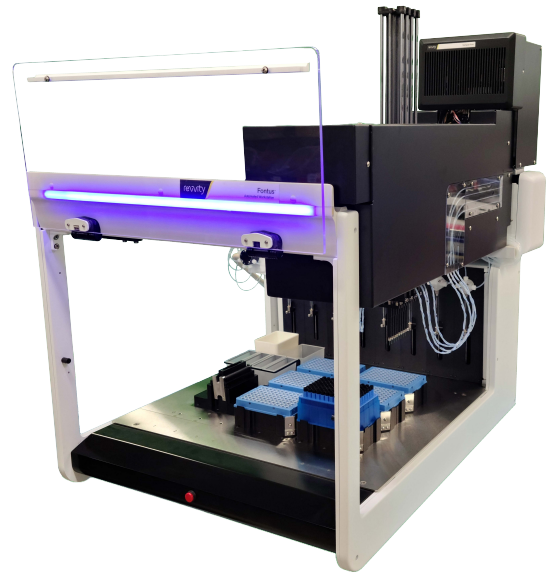




Small footprint. Big impact.

Fontus™ 8i Mini workstation - affordable & compact
Expanding the Fontus liquid handling family

Geared to support volumes ranging from 1 µL to 5000 µL with an 8 variable spanning pipetting module with independent Z-axis (up and down) movement control and 12 deck locations, the Fontus 8i Mini workstation is a fully capable and user-friendly automated liquid handling instrument.



Cornerstone of next generation automation, the Fontus 8i Mini liquid handler provides:

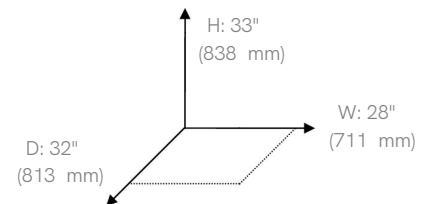
- a small footprint with no enclosure,
- integration or standalone capability,
- easy robotic access with completely open side and front access,
- outstanding flexibility such as disposable and fixed pipetting tips and liquid level detection,
- ease-of-use and efficiency with Revvity's WinPREP™ software and Application Assistant.

This system can easily go inside a system enclosure, bio-safety cabinet hood or a HEPA filtration system.

More than ever before, researchers need their instruments to run unattended, enhancing laboratory productivity.

Pipetting specifications

Varispan	Volume	Performance
20 µL disposable tip	1 µL	CV < 4% Distilled water, 500 µL syringe
20 µL disposable tip	5 µL	CV < 2% Distilled water, 500 µL syringe
Fixed tip	50 µL	CV < 1% Distilled water, 500 µL syringe



Certifications CE, CSA

Simplicity with the power of reliability

Flexibility without complexity - easily adapts to any labware or tips

- Compatible with tubes, vials, standard and high-density microplates, the Fontus 8i Mini workstation is adaptable to any application labware.
- Its Varispan™ span-8 pipetting provides independent Z, liquid level sensing and variable probe spacing for multi-tip throughput regardless of labware utilized.
- Its VersaTip™ adapter enables fixed washable, or disposable tips, or both, usable in the same protocol.
- The Varispan™ Gripper option enables labware movement automation, space optimization, walkaway time, and protocol development.

Software features - intuitive user interface

- The Fontus 8i Mini workstation with WinPREP™ software and Application Assistant simplifies setup, method creation, and protocol execution. It enables real-time research monitoring, sample tracking, and device integration for automated workflows.
- Designed to adapt to labs' evolving research needs, easy-to-use WinPREP simplifies modifying existing protocols or creating new ones without needing a programmer, reducing downtime and costs.
- The performance file library offers optimized pipetting parameters and the ability to modify pump speeds, air gaps, waste and blowout volumes, and liquid delays.

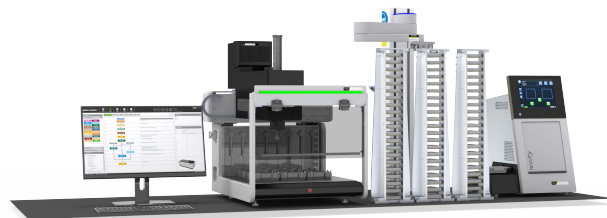
Add extended functionality - for complete automated applications solutions

From cellular screening and imaging applications to high-throughput screening and genomics-based applications, custom explorer™ G3 integrated workstations provide innovative application-focused lab automation solutions to simplify microplate handling, liquid handling, and detection.



Varispan Arm options

- Fixed tips: regular and low volume and VersaTip sampling tip
- Disposable tips: 20, 25, 50, 175, 200, 900, 1000, 5000 µL tips available.
- Conductive (black) and clear, filter, non-filter and pre-sterilized formats available
- Syringe volumes: 250, 500, 1000, 2500, 5000 µL



explorer™ G3 integrated workstations

Increase your lab's throughput and efficiency even further by integrating other Revvity instrumentation.



Release the power of Fontus automation in your workflows.

[Learn more](#)

revvity