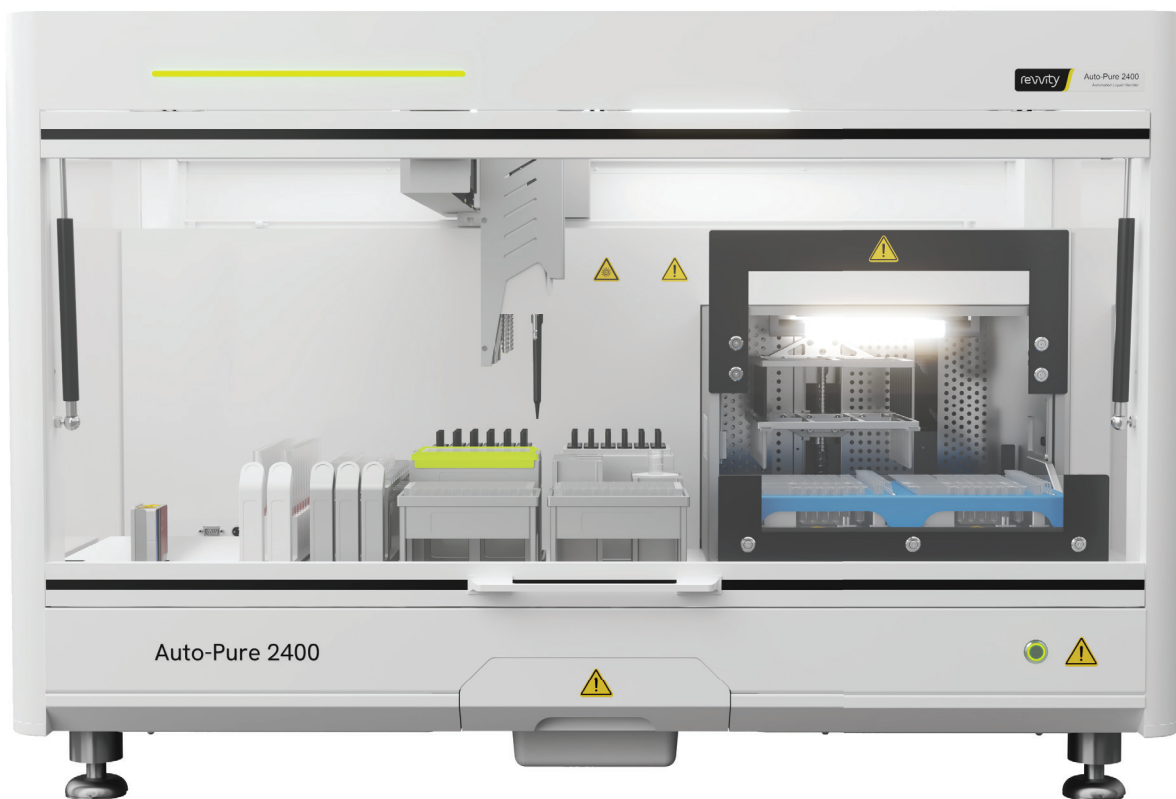




revvity

# Auto-Pure 2400

Automated liquid handler with integrated magnetic isolation



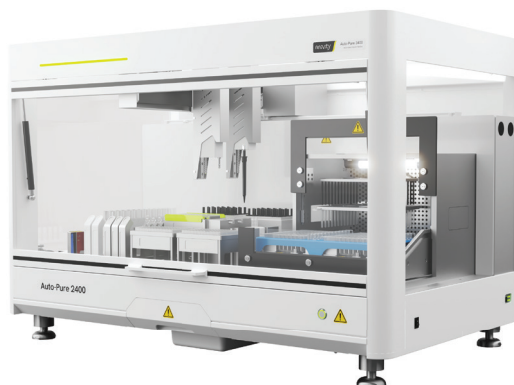
PRODUCT SPECIFICATION SHEET

# Auto-Pure 2400

For use with T-SPOT™.TB

Designed for testing up to 24 samples per run, the Auto-Pure 2400 is an automated liquid handling platform with integrated magnetic cell isolation technology.

The Auto-Pure 2400 can perform day 1 T-SPOT.TB workflows in under 3.5 hours with a single mid-run user interaction.



**Less than 10 mins hands-on time** once the Auto-Pure 2400 workflow has commenced \*

Model	Auto-Pure 2400	
Product order details	Auto-Pure 2400 (incl. PC and Monitor)	Order code: AUTOPURE 2400
Format	Automated liquid handler with integrated magnetic isolation technology. 4 000 µL pipettor and 1 000 µL pipettors.	
Sample run size	Up to 24 (1-24 T-SPOT.TB tests)	
Run time	Approx. 3 hrs per 24 sample T-SPOT.TB run (excluding count time)	
Count preparation support	Cellaca MX (2 × 12 format) and single tube (@1:4 dilution)	
Worklist file support	Cellaca MX and ELISPOT readers	
Barcode support	CODE 128, Code 39, Codabar, ITF 2 of 5, UPC-A, EAN-8, EAN-13	
Operating System	Windows 10/11, 64 bit	
Dimensions	1 090 mm × 713 mm × 734 mm (L × W × H)	
Power	Standby 30 W, maximum 600 W	
Weight	130 kg	
Consumables	Product code	Pack size
AP2400_12 Tip Comb	AP2400-12TC	200
AP2400_Tube Strip	AP2400-TS	500
AP2400_1000uL Tip	AP2400-TIP1000	1 536
AP2400_4000uL Tip	AP2400-TIP4000	576
AP2400_2mL Microcentrifuge Tube	AP2400-2MLT-E	4 000
AP2400_5mL Sample Tube	AP2400-5MLT	120
AP2400_100mL Reagent Tank	AP2400-100MLRT	200

\* Mid run user intervention with the Auto-Pure 2400 workflow and the Cellaca MX.

Products may not be licensed in accordance with the laws in all countries. Please check with your local representative for availability.



revvity