



Auto-Pure 24

Automated cell isolation system

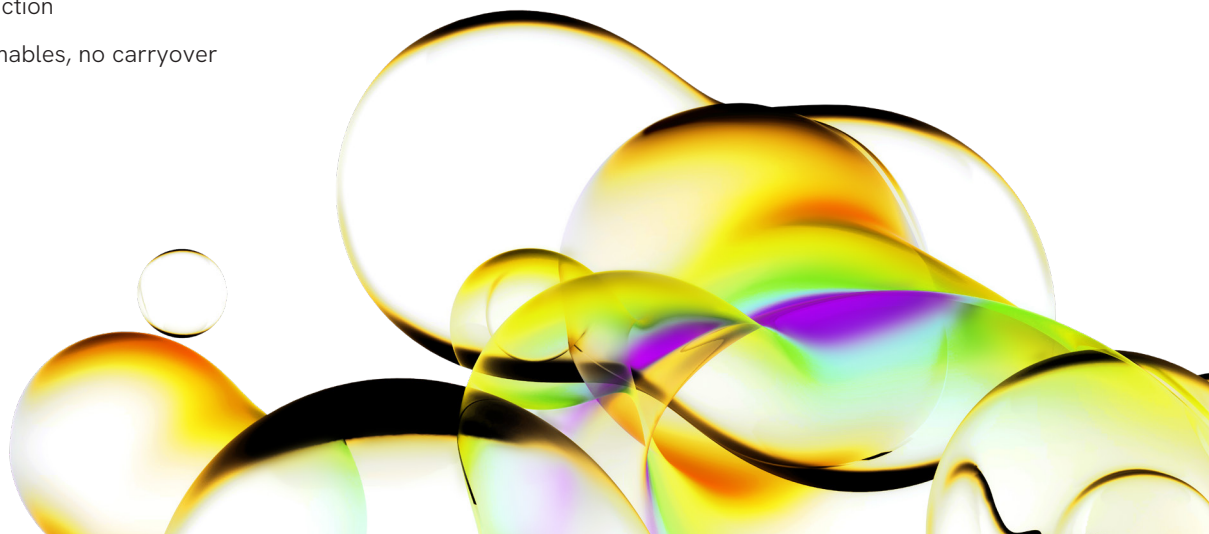
The Allsheng Auto-Pure 24 is a versatile cell isolation platform designed for mid to high-throughput laboratories. It offers dual functionality, seamlessly integrating with fully automated liquid handling systems while also capable of standalone operation for independent use. This adaptability makes the Auto-Pure 24 an ideal solution for various laboratory setups and workflow requirements, providing flexibility in cell isolation processes.

Designed for automation

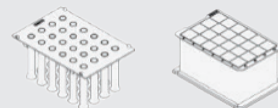
The plate-based design allows users to conveniently process up to 24 samples per run with the aid of a pre-loaded protocol. These plates are compatible with liquid handling platforms, enabling users to take advantage of the increased walk-away time automation brings.

A validated protocol developed by Revvity enables users to quickly and efficiently automate the important peripheral blood mononuclear cell (PBMC) isolation stage of the **T-SPOT™.TB** test. The Auto-Pure 24 utilises a tip-comb and 24 deep-well plates to process samples. Using positive selection, the magnetic rods collect the cell-antibody-bead complex, which is then washed and eluted.

- Deep-well plate format allows integration into fully automated liquid handler workflows
- Pre-loaded validated protocol with optimised mixing and agitation for efficient cell extraction
- Efficient hands-free processing: handles up to 24 samples in under 45 minutes
- Easy to use with minimal maintenance
- Simple touch screen interface
- Rotary table design with large access door for easy loading/unloading
- UV sterilization function
- Disposable consumables, no carryover or contamination



Model	Auto-Pure 24
Features	AUTOPURE 24
Processing volume	3.5 mL*
Processing time	Up to 45 minutes
Number of samples	1-24
Protocol storage	Pre-loaded cell isolation protocol
External interfaces	4 standard USB port, ethernet port, Wi-Fi, built-in SD card
Operating conditions	Ambient temperature: 10 - 35 °C Relative humidity: ≤ 70 % Input voltage: AC 100 - 240 V, 50 Hz/60 Hz
Service information	2 levels of service contract available (Gold/Silver)
Power supply	300 W
Dimensions (WxDxH)	560 × 620 × 500 mm
Weight	54 kg
Regulatory status	CE IVD and FDA approved
Consumables	Order code
AP24_14.6mL DWPs and Magnetic Rod Tip (1:4) (10 pcs)	AP24-DWP4-TC1



*Could vary depending on age and immune status

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

