



Auto-Pure 10B

Automated cell isolation system

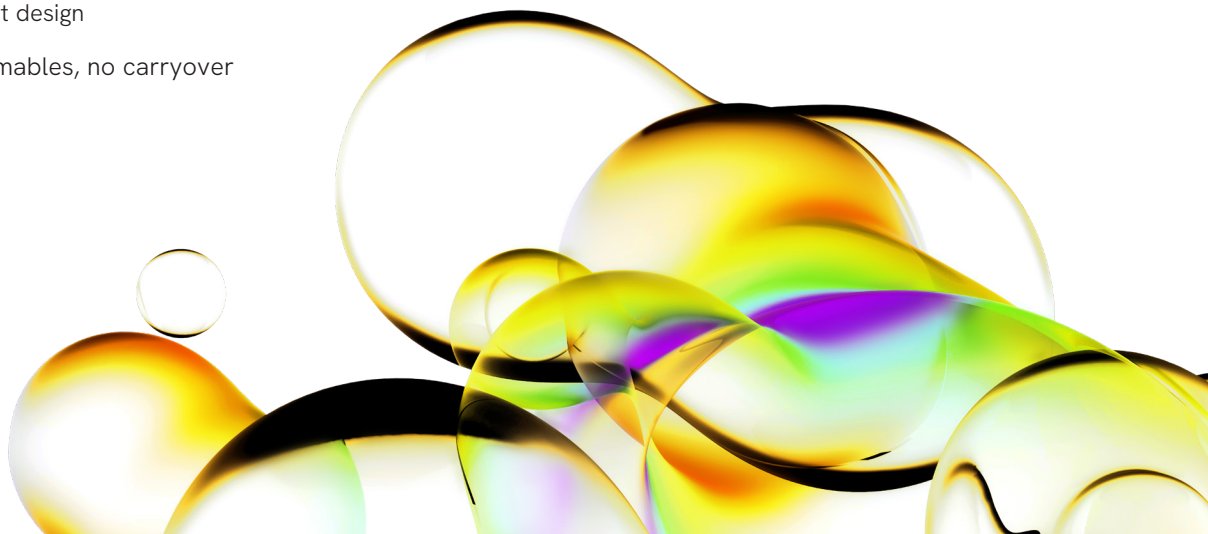
The Auto-Pure 10B serves as a versatile automated cell isolation platform for the T-SPOT™.TB test. It's an ideal solution for laboratories with lower throughput requirements, while also providing valuable backup support in high-volume testing environments.



A fully 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 10B uses magnetic rods to transport PBMCs through the various purification phases; mixing, binding, washing and elution.

Built in flexibility

The Auto-Pure 10B cell isolation system utilises a strip based workflow, bringing enhanced operational efficiency to PBMC isolation. This allows users to deal with runs of variable sizes, optimising consumable usage.

- Affordable, easy to use with minimal maintenance
- Simple touch screen interface
- Efficient hands-free processing; handles up to 10 samples in under 45 minutes
- Pre-loaded validated protocol
- Internal lighting and UV sterilization function
- Bench-top, compact design
- Disposable consumables, no carryover or contamination



Model	Auto-Pure 10B
Features	AUTOPURE 10B
Processing volume	3.5 mL*
Processing time	Less than 45 minutes
Number of samples	1-10
Protocol storage	Pre-loaded cell isolation protocol
External interfaces	Standard USB, 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)	340 × 350 × 410 mm
Weight	18 kg
Regulatory status	CE IVD
Consumables	Order code
AP20B/10B_Tube Strips	TUBESTRIPS10/20B 
AP20B/10B_Magnetic Rod Tips	AUTOPURETIPCOMB 

*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.

