

revvity

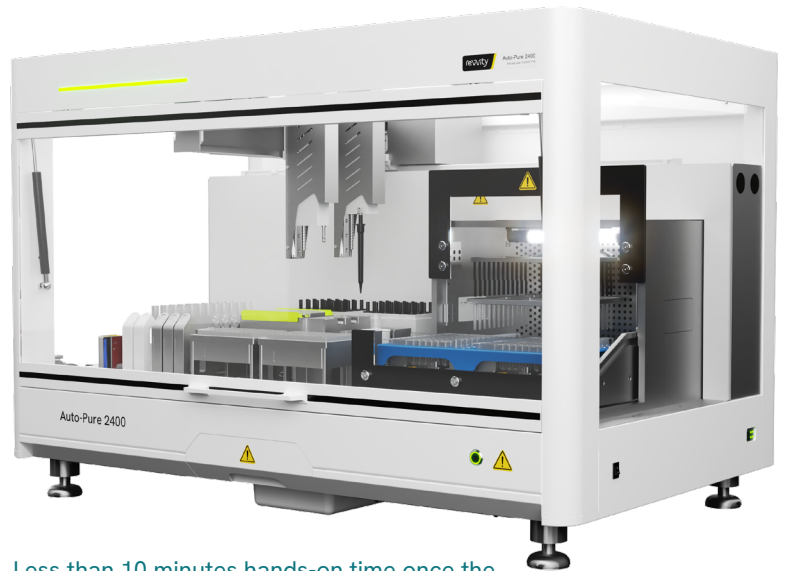
# Efficient workflow, more time for you.

## Auto-Pure 2400

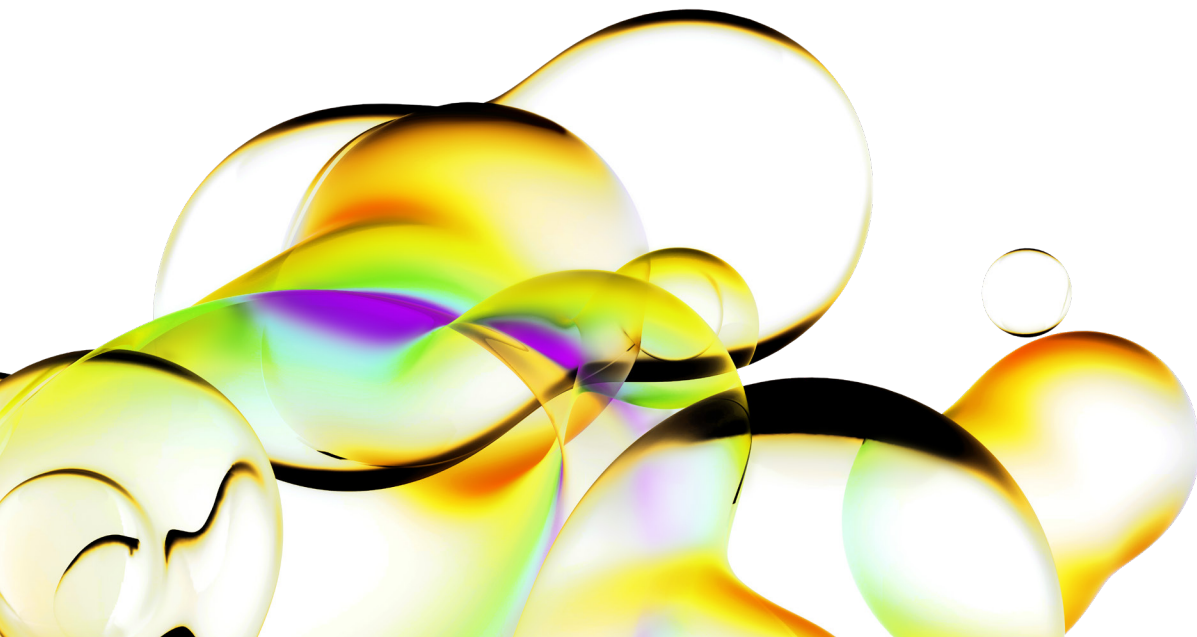
For use with T-SPOT™.TB

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 minutes hands-on time once the Auto-Pure 2400 workflow has commenced\*



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

# Easy to use, easy to get right.

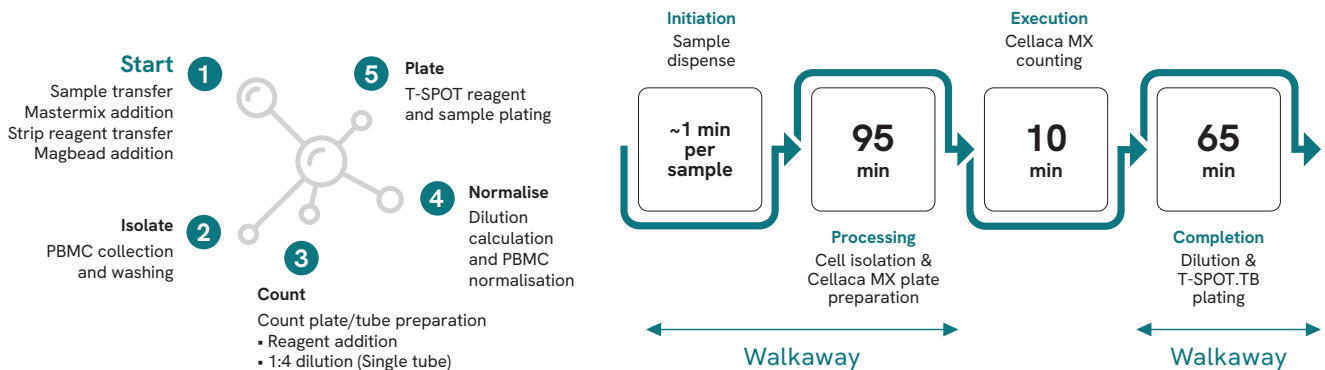
## Simplified workflow allowing for increased productivity.



### Auto-Pure 2400 key features:

- Auto-Pure magnetic cell isolation technology for efficient peripheral blood mononuclear cell (PBMC) extraction
- Compact: 1090 mm x 713 mm x 734 mm (L x W x H)
- Automated worklist generation for cell counting and spot visualisation hardware
- Cellaca™ MX cell counter file support
- User access controls
- Post run reporting for traceability
- Efficient use of consumables
- Liquid level sensing tip technology

### Automated system processes and timing for T-SPOT.TB



For more information on T-SPOT.TB testing with the Auto-Pure 2400, please contact your local Revvity representative or visit [www.revvity.com/contact-us/sales](http://www.revvity.com/contact-us/sales).

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

revvity