

JANUS G3 Automated Workstations

Versatile JANUS G3 Automated Liquid Handling Platforms



Description

JANUS® G3 Automation Workstations deliver real-time and future adaptability in throughput, capacity, and dynamic volume range from 0.5 µl to 5000 µl for consistent and reproducible sample preparation with a choice of pipetting heads, gripper options, and application accessories for complete, walk-away automation. JANUS G3 WinPREP user friendly software with Application Assistant interface provides “drag-and-drop” applications programming without complexity.

The JANUS instrument family includes four workstations, the Mini, Standard, Integrator, and Expanded platforms, providing complete flexibility to best suit your particular application and deck capacity requirements; additionally, NGS and BioTx application specific workstations designed with customer input for optimal automated NGS library preparation and small scale purification, respectively.

Product Features

- The unique JANUS G3 platform design allows you to add up to two (2) pipetting arm(s) and a plate movement arm to enhance your workstation's overall functionality. Standard pipetting volumes range from 0.5 µl to 5000 µl, and with the nanohead pipetting arm option, the low-end volume range is 20 nL. (Available for all models except JANUS G3 NGS Express and BioTx Workstations).
- JANUS G3 platform MDT arm provides interchangeable and automated head functionality for 96- and 384-channel dispense heads in less than 25 seconds using a powerful and fail-safe interlocking mechanism. With 'on the fly' head swapping you can easily dispense into 96-, 384- and 1536-well shallow and deep well microplates, without user intervention. (Available for all models except JANUS G3 NGS Express and BioTx Workstations).
- Optional JANUS G3 enclosure helps to minimize contamination (available for all models except JANUS G3 BioTx Workstation).
- Ability to automate plate movement and access off-deck plates without the need for an additional robotic arm.
- VersaTip® provides a unique sampling tip design to deliver the benefits of both a fixed washable tip and disposable tip adapter.
- Varispan 4- or 8-tip arm provides independent motion and liquid level sensing in Z direction to access test tubes, microplates, deep well plates, and vials. Sampling probes offer variable spacing from 9 mm to 20 mm for 4-tip arm or 9 mm to 40 mm for 8-tip arm.
- Accusense™ our patented independent liquid level sensing technology efficiently monitors your sample volumes, with the ability to detect samples as low as 50 ul in microplates and 100 ul in test tubes. Low or non-ionic polar solutions and organic solvents such as DMSO or methanol are also detectable.
- Active Identification provides automatic recognition of the MDT head type.
- Carry-over elimination via a peristaltic pump for high volume/high throughput sample tip washing for carry-over to less than 1:1,000,000 with appropriate washing or use of disposable tips for improved elimination of carry-over.

