

98-70009

ePrep Software Version 1.45.32 Release Notes

ePrep Software Version 1.45.32 Release Notes

Introduction

This document contains lists of new features, minor updates and bug fixes included in the ePrep Software version 1.45.32. For the changes included in previous versions, please refer to the Release Notes of that release.

ePrep Software Version Details

- **Version Number:** 1.45.32
- **Release Type:** Version Release
- **Release Date:** 21-Aug-2025

Installation

Please follow the instructions in the **98-65009 MAINTENANCE PROCEDURE, SOFTWARE UPDATE** to install the ePrep Software.

Feature Updates

The following is the list of new features included in the ePrep Software version 1.45.32.

Feature	Description
Workflow Editor Rack Selection	This function enhances Workflow Adapter Plate and Rack selection within the Workflow Editor by enabling users to search by rack names, types, and sizes, while also allowing them to select and view favourites for streamlined workflow creation.
Syringe Calibration Function	<p>The syringe calibration feature enables users to validate and/or calibrate syringes for use on the ePrep ONE. The calibration function uses a single-point calibration, where the user selects a target volume and takes 10 gravimetric aliquots to calculate the average measured volume, which is then used to determine a specific syringe calibration.</p> <p>The software automatically calculates a correction factor that can be applied to all plunger movements for dispensed volume. Expiry dates can also be set for recalibration.</p>
Advanced Workflow Runtime Estimation	This function provides a comprehensive runtime estimation for ePrep ONE workflows through Workflow Validation and Run Screens by executing a complete background simulation of the workflow, delivering time estimates that are typically accurate within $\pm 10\%$ of actual runtime.
Probe Dispenser Anti-Drip Function	Probe Dispenser operations can occasionally cause dripping after vial removal. The new Anti-Drip Function eliminates this issue by withdrawing a small volume from the Probe Dispenser Needle tip before removal of the needle from the vial. The withdrawal volume is automatically compensated on subsequent dispenses.

ePrep Software Version 1.45.32 Release Notes

Feature	Description
Enable Workflow Registered Tools	<p>Verified application-specific workflows are available for ePrep ONE, such as EPA 3511 and 1633. ePrep have created these workflows to use specific "Registered" syringes allocated to tasks for speed and cross-contamination purposes.</p> <p>This syringe reallocation function enables users to import external verified workflows and manage the controlled reassignment of "registered" syringes to syringes available in the user's inventory. The system prevents the workflow from executing until all required syringes have been properly reallocated.</p>
Allow Multiple 4P and 2P Vortex Mixers	<p>The ePrep ONE software now enables placement of multiple 2P and 4P Vortex mixers on the deck, with each unit independently controlled through software Workflow Tasks.</p> <p>This feature extends ePrep's capabilities for applications such as EPA methods, where large sample batches are needed.</p>

Minor Updates and Bug Fixes

ePrep Software version 1.45.32 includes the following **minor updates** from version **1.44.28**.

ID	ePrep Software Area	Description
3551	Workflow Volume Validation	Improve volume validation calculations.
4295	System Error Handling	Add additional initialisation error information.
4250	PFAS Free Syringes	Update PFAS free syringe images.
4340	Set Sample Volume Task	Enable Set Sample Volume for 20-40mL Enviro Vials in 56 x 4-Pin Rack.

ePrep Software version 1.45.32 includes the following **bug fixes** from version **1.44.28**.

ID	ePrep Software Area	Description
4220	SPEmx Task	Remove "Rinse Needle" wash from SPEmx Elute.
4257	Workflow Editor	Improve Workflow name titles when importing XML files.
4237	Workflow Editor	Improve Group Names after opening a workflow.
4182	Vortex Mixer Task	Improve multiple deck accessory communications.
4140	Vortex Mixer Task	Correct Vortex Mix Task times are reset to 0s after a software upgrade.

ePrep Software Version 1.45.32 Release Notes

ID	ePrep Software Area	Description
4303	System Error Handling	Correct system recovery logic during foot movements.
4232	System Error Handling	Improve User error messages.
4312	Probe Dispenser Tasks	Robustness of probe dispenser wash port selection.
4343	Probe Dispenser Tasks	Correct Probe Dispenser multi-dispense Anti-Drip function.
4318	Probe Dispenser Tasks	Correct Probe Dispenser needle tip rinse.
4304	Piercer Wash	Remove Piercer Wash from uSPEed Syringe operation.
4328	Needle Position	Change vial extraction of syringe needles to avoid move errors.
4339	Extra Tool Wash Tasks	Improve Post Extra Tool Wash for reliability.
4354	μSPEed Tasks	μSPEed Task logic to avoid misuse of Aspiration Needle.

Other minor bug fixes and system reliability improvements have been included in this version.

ePrep Software Version 1.45.32 Release Notes

Revision history

Rev No	Date	Changes
01	21/Aug/2025	Initial Release.