

ePrep Software Release Notes					
Version	Type	Status	Features and Requirements	Bug Fixes	Date
1.40.52	Official Release	Active	<ul style="list-style-type: none">All in 1.40.50Requirement #ADO-3292. New 46ml high recovery 4P rack (Rack Class 0143)	Workflow Run: <ul style="list-style-type: none">Bugfix #ADO-3281. Workflow does not stop when the bumper is triggered while moving down to pierce a vial. (REVISITED - Safety Improvements) Database: <ul style="list-style-type: none">Bugfix #ADO-3277. System out-of-memory error when saving records to the database uSPEed: <ul style="list-style-type: none">Bugfix #ADO-3291. uSPEed cartridge is parked in the wrong position when uSPEed Load is the last task in a sequential task groupBugfix #ADO-1993. Error when starting from second position in uSPEed RackBugfix #ADO-3295. uSPEed cartridge is not picked up properly after dispensing in rack	01-Nov-24
1.40.50	Official Release	Obsolete	<ul style="list-style-type: none">All in 1.40.48Requirement #ADO-3284. Record Frequency from RFID coils in Command Log.	Workflow Run: <ul style="list-style-type: none">Bugfix #ADO-3281. Workflow does not stop when the bumper is triggered while moving down to pierce a vial. Workflow Reports: <ul style="list-style-type: none">Bugfix #ADO-3218. Viewing Report after Running Workflow takes a long time to load/show.	28-Oct-24
1.40.48	Official Release	Obsolete	<ul style="list-style-type: none">All in 1.40.45	Communications: <ul style="list-style-type: none">Bugfix #ADO-2997. Incompatible Firmware/Software version Popup showing when connecting without device connected.Bugfix #ADO-3231. Rack or Tool Scan fails reporting incorrect Adapter Plate or corrupted tool RFID. Racks: <ul style="list-style-type: none">Bugfix #ADO-3265. TWD 2 x 24 x 4mL tube rack volume formula is incorrect.Bugfix #ADO-3138. Racks covering tubes warning popup appears in every rack scan even when not running a workflow. (REVISITED)	17-Oct-24
1.40.46.1	Official Release (Hot-Fixed)	Active	<ul style="list-style-type: none">All in 1.40.45Requirement #ADO-2523. New Rack Classes for Cryovials (REVISITED)<ul style="list-style-type: none">Cryovial 34 x 1.8mL tube rack (Rack Class 007501)Waters Acquity (Cryovial) 2 x 48 x 1.3-2.2mL tube rack (RackClass 050201) with 1.3ml and 2.2ml tube types	Volume Validation: <ul style="list-style-type: none">Bugfix #ADO-3204. New Cryovial 050201 rack class fails workflow validation when vials are used as a source group Racks: <ul style="list-style-type: none">Bugfix #ADO-3193. Waters Acquity rack 2x48 1.5mL vials with caps off produce height error at beginning of workflow (REVISITED) Communications: <ul style="list-style-type: none">Bugfix #ADO-2997. Incompatible Firmware/Software version Popup showing when connecting without device connected.	23-Sep-24
1.40.45	Official Release	Obsolete	<ul style="list-style-type: none">All in 1.40.44	Vortex Mixer <ul style="list-style-type: none">Bugfix #ADO-3200. Vortex mixer rack scan not working correctly.	20-Sep-24

Version	Type	Status	Features and Requirements	Bug Fixes	Date
1.40.44	Official Release	Obsolete	<ul style="list-style-type: none"> All in 1.40.42 	<p>Volume Validation:</p> <ul style="list-style-type: none"> Bugfix #ADO-3132. Workflow validation gets stuck when validating a sequential task group using the Magnetic Stirrer Task <p>Racks:</p> <ul style="list-style-type: none"> Bugfix #ADO-3138. Racks covering tubes warning popup appears in every rack scan even when not running a workflow. Bugfix #ADO-3193. Waters Acquity rack 2x48 1.5mL vials with caps off produce height error at beginning of workflow <p>Dispense IN RACK for uSPEd:</p> <ul style="list-style-type: none"> Bugfix #ADO-3152. The bumper touches the uSPEd/SPEmx Rack when dispensing into the rack. Bugfix #ADO-3163. Unexpected error when going to the Wash Station after dispensing into the uSPEd Rack. <p>Communications:</p> <ul style="list-style-type: none"> Bugfix #ADO-2663. Comms Error: "Object reference not set to an instance of an object" when null response is not handled <p>Needle Tip Touch Off:</p> <ul style="list-style-type: none"> Bugfix #ADO-3143. Bent spSPE cartridge needle after wrong needle length measure Bugfix #ADO-2684. "Object reference not set to an instance of an object" Error during Filter Needle Tip Touch-Off 	19-Sep-24
1.40.42	Official Release	Obsolete	<ul style="list-style-type: none"> All in 1.40.41 Requirement #ADO-2826. Allow Capped and Uncapped Option for Racks Covering Tubes. 	<p>Workflow</p> <ul style="list-style-type: none"> Bugfix #ADO-3029. Workflow partially loading due to tool class issue. Bugfix #ADO-2630. Workflow is loading blank. <p>Racks</p> <ul style="list-style-type: none"> Bugfix #ADO-3123. Unexpected Error during Volume Validation for Aliquot Sample task when using Luer Tool. 	26-Aug-24
1.40.41	Official Release	Obsolete	<ul style="list-style-type: none"> All in 1.40.39 Requirement #ADO-3005. Add PTFE Free Generic syringes 	<p>Piercer Wash</p> <ul style="list-style-type: none"> Bugfix #ADO-2973. Range Error when doing Piercer Wash in the Wash Station (Revisited) <p>Import Service</p> <ul style="list-style-type: none"> Bugfix #ADO-2980. Empty User table startup error appears if previous a startup error occurred after a clean install. <p>Installer</p> <ul style="list-style-type: none"> Bugfix #ADO-3004. Access denied startup error after installing version 1.40.38 <p>PTFE Free Syringes</p> <ul style="list-style-type: none"> Bugfix #ADO-3001. Incorrect needle length for 1000 ul Syringe, PTFE free, 103mm Needle 	02-Aug-24
1.40.39	Official Release (Hot-Fixed)	Obsolete	<ul style="list-style-type: none"> All in 1.40.38 	<p>Piercer Wash</p> <ul style="list-style-type: none"> Bugfix #ADO-2973. Range Error when doing Piercer Wash in the Wash Station <p>Volume Validation</p> <ul style="list-style-type: none"> Bugfix #ADO-2974. Workflow Validation does not detect some tube index changed when dispensing the filtering holdup volume <p>Anti-Drip Procedure for Probe Dispenser</p> <ul style="list-style-type: none"> Bugfix #ADO-2996. Residual volume exception when running Make Up To Volume task with Probe Dispenser with Anti-Drip 	18-Jul-24
1.40.38	Official Release	Obsolete	<ul style="list-style-type: none"> All in 1.40.37 	<p>Database</p> <ul style="list-style-type: none"> Bugfix #ADO-2979. User table error on startup when doing a clean install 	11-Jul-24

Version	Type	Status	Features and Requirements	Bug Fixes	Date
1.40.37	Official Release	Obsolete	<ul style="list-style-type: none"> All in 1.40.36 	Export Service: <ul style="list-style-type: none"> Bugfix #ADO-2880. Software crashes when disconnecting external USB drive during data export (Workflows, Workflow Reports and Audit Trails) Installer <ul style="list-style-type: none"> Bugfix #ADO-2954. Installer fails when encountering unexpected files in INI Files. Logic for Bumper Trigger <ul style="list-style-type: none"> Bugfix #ADO-2805. Bumper counter log crashes software (wrong firmware version) (Revisited - minimum firmware version updated to 01-84) Racks <ul style="list-style-type: none"> Bugfix #ADO-2481. Shimadzu rack 105 x 1.5mL vial height is incorrect User Accounts and Access Levels <ul style="list-style-type: none"> Bugfix #ADO-2961. Default User Account lost its permissions on an update (Cannot Reproduce - Additional database integrity checks added) Database <ul style="list-style-type: none"> Bugfix #ADO-2825. Running the software as not admin shows a database error (Cannot Reproduce - improved error handling and error message) 	09-Jul-24
1.40.36	Official Release	Obsolete	<ul style="list-style-type: none"> All in 1.40.33 Pickup Routine PTFE Free Syringes Anti-Drip Procedure for Probe Dispenser Requirement #ADO-2879. Graphic changes in Export Service: select minimum required files by default. 	Accessories: <ul style="list-style-type: none"> Bugfix #ADO-2873. If an Accessory Class is updated a database error appears Tools: <ul style="list-style-type: none"> Bugfix #ADO-2911. Update not clearing redundant Tool Class properly Bugfix #ADO-2906. Wrong tool RFID serial number shows not supported Autodisconnect Tool Bugfix #ADO-2779. Corrupt RFID message Bugfix #ADO-2905. Tool Scan not scanning any tools Bugfix ADO-2911. Update not clearing redundant Tool Class properly Export Service: <ul style="list-style-type: none"> Bugfix #ADO-2821. Software crashes when disconnecting external USB drive during data export Filtering: <ul style="list-style-type: none"> Bugfix #ADO-2892. Filtering holdup waste setting allows selecting the filtering output group as waste tubes instead of collect tubes Bugfix ADO-2908. Filtering indexing error stuck in a memory Workflow Reports: <ul style="list-style-type: none"> Bugfix #ADO-2872. Viewing Workflow through Reports allows workflow to be overwritten through the Approve button Command Logs: <ul style="list-style-type: none"> Bugfix #ADO-2638. Command Log files size is too large (Revisited) Probe Dispenser: <ul style="list-style-type: none"> Bugfix ADO-2827. Probe dispenser Aspirate/Dispense default values Logic for Bumper Trigger <ul style="list-style-type: none"> Bugfix #ADO-2805. Bumper counter log crashes software (wrong firmware version) (Revisited - minimum firmware version rolled-back to 01-82) Anti-Drip Procedure for Probe Dispenser <ul style="list-style-type: none"> Bugfix #ADO-2903. Droplets on top of vials. 	27-Jun-24
1.40.33	Official Release	Obsolete	<ul style="list-style-type: none"> All in 1.40.32 	Audit Trail <ul style="list-style-type: none"> Bugfix #ADO-2869. Audit Trail - Missing Actions (Tool and Rack Scan) Accessories <ul style="list-style-type: none"> Bugfix #ADO-2873. If an Accessory Class is updated a database error appears Volume Validation <ul style="list-style-type: none"> Bugfix #ADO-2874. Wrong volume validation error when validating workflow with SPE tasks SPE Suite <ul style="list-style-type: none"> Bugfix #ADO-2875. Not all SPE cartridges are conditioned in a SPE Condition batch task even though the volume validation passes 	07-Jun-24

Version	Type	Status	Features and Requirements	Bug Fixes	Date
1.40.32	Official Release	Obsolete	<ul style="list-style-type: none"> All in 1.40.31 	Database <ul style="list-style-type: none"> Bugfix #ADO-2861. Database error during workflow run. (Revisited - Improved Concurrency Control) Workflows <ul style="list-style-type: none"> Bugfix #ADO-2866. Restore All button crashing Editor. (Button Removed) Accessories <ul style="list-style-type: none"> Bugfix #ADO-2867. Accessories configuration is lost after software update. uSPEd <ul style="list-style-type: none"> Bugfix #ADO-2870. uSPEd Crash on entering Tube Editor. Backup, Restore and Archive Service <ul style="list-style-type: none"> Bugfix #ADO-2802. Progress bar too slow for smaller record count. (Revisited) 	03-Jun-24
1.40.31	Official Release	Obsolete	<ul style="list-style-type: none"> All in 1.40.26 Requirement #ADO-2773. Update Auditor Access Level Default Permissions Configuration Requirement #ADO-2653. Auto-populate date when exporting Audi Trails Requirement #ADO-2766. Long reach syringe images Requirement #ADO-1359. Merge Init Data into Audit Trail Requirement #ADO-1360. Merge Run Data into Audit Trail Requirement #ADO-2773. Update Auditor Access Level Default Permissions Configuration Requirement #ADO-2784. Change 21 CFR Part 11 Settings Dependencies 	Syringe Conformance <ul style="list-style-type: none"> Bugfix #ADO-2747. Tool Conformance not updating Tool conformance expiry on Workflow Load. Electronic Signatures <ul style="list-style-type: none"> Bugfix #ADO-2772. Wrong keyboard label when adding additional comment in Electronic Signature User Accounts and Access Levels <ul style="list-style-type: none"> Bugfix #ADO-2774. Temporary Password it's been shown in ePrepDatabaseLog file Bugfix #ADO-2259. Unclear error message text when creating password Bugfix #ADO-2687. Expiry Date for Administrators shows as 1/Jun/5001 Bugfix #ADO-2455. Cannot remove temporary password for password reset Bugfix #ADO-2062. Change password says incorrect username Volume Validation <ul style="list-style-type: none"> Bugfix #ADO-2785. Database error during Volume Validation Bugfix #ADO-1754. Auto wash not working properly Bugfix #ADO-1942. prime volume at reagent station not considered in volume validation Bugfix #ADO-2834. Workflow validation error due to source vials not switching to the next available one. Bugfix #ADO-2865. Residual Volume is dispensed into the output vial when prime cycles for refills are zero Digital Ports in Accs Interface <ul style="list-style-type: none"> Bugfix #ADO-2755. Touch off probe not checked Graphics <ul style="list-style-type: none"> Bugfix #ADO-2769. Strange behavior when clicking around group selector Audit Trails <ul style="list-style-type: none"> Bugfix #ADO-2770. Tube references are zero based in the Audit Trail Bugfix #ADO-2771. Task deck group selection is not clear in Audit Trail Bugfix #ADO-2299. Audit trails fixes Bugfix #ADO-2793. Audit Trail Descriptions are imported every time Bugfix #ADO-2691. Report Audit Trail Event when restoring a system backup Workflows <ul style="list-style-type: none"> Bugfix #ADO-2525. Unclear message when stopping a workflow Bugfix #ADO-2630. Workflow is loading blank Bugfix #ADO-2768. Incomplete Workflow Summary (Workflow Summary has been removed) Command Logs <ul style="list-style-type: none"> Bugfix #ADO-2638. Command Log files size is too large Bugfix #ADO-2792. Old command log is not being deleted Racks <ul style="list-style-type: none"> Bugfix #ADO-2737. Bel-Art 5x6 Rack missing in Rack Choices Prompt Task <ul style="list-style-type: none"> Bugfix #ADO-1766. Issues with prompt task Deck Objects <ul style="list-style-type: none"> Bugfix #ADO-206 (8L08-787). Falcon Tubes Dead Volume Backup, Restore and Archive Service <ul style="list-style-type: none"> Bugfix #ADO-2802. Progress bar too slow for smaller record count Bugfix #ADO-2835. Silent Backup required before performing Restore. Logic for Bumper Trigger <ul style="list-style-type: none"> Bugfix #ADO-2805. Bumper counter log crashes software (wrong firmware version) Database <ul style="list-style-type: none"> Bugfix #ADO-2861. Database error during workflow run 	31-May-24

Version	Type	Status	Features and Requirements	Bug Fixes	Date
1.40.26	Official Release	Obsolete	<ul style="list-style-type: none"> All in 1.40.21 Requirement #ADO-2688. Include "System Position Check" and "Software OQ Test" workflows in installer package Requirement #ADO-2692. New vial for Supelco 36 x 6-20mL (Rack Class 0202) Requirement #ADO-2743. Update to Falcon Tube Rack Classes 0118 and 0117. 	<p>Volume Validation:</p> <ul style="list-style-type: none"> Bugfix #ADO-2715. Validation error - variable dispense - set prime to "NEVER" Bugfix #ADO-2305. Batch many to many validation error <p>User Accounts:</p> <ul style="list-style-type: none"> Bugfix #ADO-2685. New User Access Level disappear after been created Bugfix #ADO-2455. Cannot remove temporary password for password reset <p>Workflow Reports:</p> <ul style="list-style-type: none"> Bugfix #ADO-2262. Bumper fault clears workflow report content Bugfix #ADO-2263. Workflow report incorrect filename Bugfix #ADO-2137. No error information in workflow report Bugfix #ADO-2616. Wrong Tool Volume for Probe Dispenser in Workflow Report <p>Filtering:</p> <ul style="list-style-type: none"> Bugfix #ADO-2479. Misleading missing filter message <p>Workflows:</p> <ul style="list-style-type: none"> Bugfix #ADO-2219. Cannot select workflows with errors <p>Archive Service:</p> <ul style="list-style-type: none"> Bugfix #ADO-2411. Partial audit trail archive inconsistent <p>Vortex Mixer:</p> <ul style="list-style-type: none"> Bugfix #ADO-2415. System bump into a rack Bugfix #ADO-2738. Vortex mix inside Liquid Handling task allows for high RPM Bugfix #ADO-2741. RPM for Vortex Mixer reverting to default <p>Racks:</p> <ul style="list-style-type: none"> Bugfix #ADO-2737. Bel-Art 5x6 Rack missing in Rack Choices 	03-May-24
1.40.21	Official Release	Obsolete	<ul style="list-style-type: none"> All in 1.40.17 Requirement #ADO-2493. Change Archive Data default permission to YES for Administrator UAL Requirement #ADO-2523. New and Updated Rack Classes. <ul style="list-style-type: none"> - Agilent, 2 x 40 x 1.5-1.8mL tube rack (RackClass 0265) - CTC VT12 2 x 12 x 20-40mL tube rack (RackClass 0285) - 2 x Microtiter Plate, 96 well (RackClass 0090) with new 2.2ml tube type - Waters Aquity 2 x 48 x 1.3-2.2mL tube rack (RackClass 0502) with new 1.3ml and 2.2ml tube types Requirement #ADO-2521. Reduce Max and Recommended Vortex Mixer Speeds for Racks 	<p>SPE:</p> <ul style="list-style-type: none"> Bugfix #ADO-2517. Index error when running SPE workflow Bugfix #ADO-2488. Workflow stopped with SPE group across multiple racks on spSPE elute <p>User Accounts (Locking Interface):</p> <ul style="list-style-type: none"> Bugfix #ADO-2491. 21CFR11 - Prompt message allowed to continue when Interface is locked Bugfix #ADO-2489. 21CFR11 - User can ok out of error messages while interface is locked Bugfix #ADO-2071. User accounts are being deleted. Bugfix #ADO-2117. Able to interact with popup while interface locked Bugfix #ADO-2516. Software becomes un-clickable if locked out and bumper Triggered Bugfix #ADO-2406. If user is locked out while importing workflow dialog is showing, App can become un-usable Bugfix #ADO-2629. Username not reqd when unlocking Bugfix #ADO-2633. Lock Screen disappears after Prompt Task <p>Electronic Signatures:</p> <ul style="list-style-type: none"> Bugfix #ADO-2403. Can Sign workflow with errors <p>Audit Trail:</p> <ul style="list-style-type: none"> Bugfix #ADO-2299. Audit trails fixes Bugfix #ADO-2520. Audit database error, just fills database with junk <p>Deck Objects (Vials):</p> <ul style="list-style-type: none"> Bugfix #ADO-2044. 0202 Rack class update. Needle buckles in Headspace vials. Bugfix #ADO-2637. Cannot dispense to MT plate <p>Workflow Reports:</p> <ul style="list-style-type: none"> Bugfix #ADO-2524. 21CFR11: Mag Stirring task does not show in Workflow report or Audit Trail Bugfix #ADO-2134. Vortex mixer task not shown in workflow report Bugfix #ADO-2174. No pause entries in workflow reports Bugfix #ADO-2628. Tool wash and prompt not in report <p>Command Logs:</p> <ul style="list-style-type: none"> Bugfix #ADO-2265. Command Logs need additional decoding Bugfix #ADO-2474. Decoding issue for Command Log <p>Tools:</p> <ul style="list-style-type: none"> Bugfix #ADO-2261. Tool Scan images is not scaled on a tablet <p>Database Migration:</p> <ul style="list-style-type: none"> Bugfix #ADO-2548. Very long start-up time <p>Graphics:</p> <ul style="list-style-type: none"> Bugfix #ADO-2458. 21 CFR 11 beta on multiple monitors Bugfix #ADO-2431. Calibration button cut off 	08-Apr-24

Version	Type	Status	Features and Requirements	Bug Fixes	Date
1.40.17	Official Release	Obsolete	<ul style="list-style-type: none"> • All in 1.35.13.3 • 21 CFR Part 11 - User Accounts and Access Levels • 21 CFR Part 11 - Electronic Signatures • 21 CFR Part 11 - Audit Trails • 21 CFR Part 11 - Export Service (CSV and PDF) • 21 CFR Part 11 - Backup and Restore Service • 21 CFR Part 11 - Archive Service - Audit Trail and Workflow Reports • 21 CFR Part 11 - Workflow Import/Export and Approval • 21 CFR Part 11 - Workflow Reports • 21 CFR Part 11 - Targeted Updates and Import Service • 21 CFR Part 11 - Restructure Debug Commands and Raw Moves files to CSV (Command Logs) • 21 CFR Part 11 - Settings: Descriptions • 21 CFR Part 11 - System Checks and Validation - Integrity Checks • 21 CFR Part 11 - System Checks and Validation - Qualification Procedures Confirmation (IQ, OQ, PQ) • 21 CFR Part 11 - System Checks and Validation - Syringe Conformance Check • Update tracking (Settings cut-off) • Database Migration • Graphics Clean - Full Screen in 16:9 Aspect Ratio • Graphics Clean - Standarise pop-ups (C# pop-ups to look like VB pop-ups) 	<ul style="list-style-type: none"> • Bugfix #ADO-2412. Needle tip touch off is done every time a cartridge is picked up. (Revisited) • Bugfix #ADO-2393. Incomplete Workflow Reports. (Fix for 1.40) • Bugfix #ADO-2474. Decoding issue for Command Log. • Bugfix #ADO-2478. Backup service crashes the software. • Bugfix #ADO-2461. Software crash on selecting Optional Setting WFR Highest Record No. • Bugfix #ADO-2470. Software crash on selecting group that spans multiple 50mL racks. • Bugfix #ADO-2475. Pickup SPE method failing in deck group loop. 	20-Feb-24
1.40.15	Official Release	Obsolete	<ul style="list-style-type: none"> • All in 1.35.13.3 • 21 CFR Part 11 - User Accounts and Access Levels • 21 CFR Part 11 - Electronic Signatures • 21 CFR Part 11 - Audit Trails • 21 CFR Part 11 - Export Service (CSV and PDF) • 21 CFR Part 11 - Backup and Restore Service • 21 CFR Part 11 - Archive Service - Audit Trail and Workflow Reports • 21 CFR Part 11 - Workflow Import/Export and Approval • 21 CFR Part 11 - Workflow Reports • 21 CFR Part 11 - Targeted Updates and Import Service • 21 CFR Part 11 - Restructure Debug Commands and Raw Moves files to CSV (Command Logs) • 21 CFR Part 11 - Settings: Descriptions • 21 CFR Part 11 - System Checks and Validation - Integrity Checks • 21 CFR Part 11 - System Checks and Validation - Qualification Procedures Confirmation (IQ, OQ, PQ) • 21 CFR Part 11 - System Checks and Validation - Syringe Conformance Check • Update tracking (Settings cut-off) • Database Migration • Graphics Clean - Full Screen in 16:9 Aspect Ratio • Graphics Clean - Standarise pop-ups (C# pop-ups to look like VB pop-ups) 	<ul style="list-style-type: none"> • Bugfix #ADO-2012. Needle tip touch off is done every time a cartridge is picked up. • Bugfix #ADO-2229. Unused deckgroups warning not showing when "Run". • Bugfix #ADO-2024. Workflow Reports are being lost on upgrade from 1.35.6 to 1.40.6 • Bugfix #ADO-2389. Adding tube to group then selecting it breaks software - Control-Alt-Delete required. • Bugfix #ADO-2404. Supplied workflow Example #2 - uSpeed loads with errors • Bugfix #ADO-2456. Targeted Updates Inaccessible when 21 CFR Part 11 settings are not turned ON. • Bugfix #ADO-2464. Database error when importing a Workflow with empty name. • Bugfix #ADO-2465. Timestamp in CSV export of Audit trail is not human readable. • Bugfix #ADO-2469. Scanned Tools are deleted when upgrading to version 1.40. • Bugfix #ADO-2460. PTFE free syringes/tools still listed as prototypes. • Bugfix #ADO-2405. Restore backup from newer version. • Bugfix #ADO-2454. Minimum password length crash. • Bugfix #ADO-2265. Command Logs need additional decoding. • Bugfix #ADO-2300. Failed to abort rack scan. • Bugfix #ADO-2303. Targeted updates backup progress bar. • Bugfix #ADO-2332. Error not handled when Backing up. 	16-Feb-24
1.35.13.3	Official Release (Hot-Fixed)	Obsolete	<ul style="list-style-type: none"> • All in 1.35.13.1 	<ul style="list-style-type: none"> • Bugfix #ADO-2393. Incomplete Workflow Reports. 	06-Feb-24
1.35.13.1	Official Release (Hot-Fixed)	Obsolete	<ul style="list-style-type: none"> • All in 1.35.13 	<ul style="list-style-type: none"> • Bugfix #ADO-2334. Cannot vortex mix prototype rack 18 x 40ml. 	22-Jan-24
1.35.13	Official Release	Obsolete	<ul style="list-style-type: none"> • All in 1.35.12.1 	<ul style="list-style-type: none"> • Bugfix #ADO-2289. Getting tube type changed message even if the tube type was not changed. • Bugfix #ADO-2264. Bumper trigger after aspiration needle drop off. • Bugfix #ADO-2298. R Motor enabled during rack scan. 	12-Jan-24