



vControl

Revision History

Document name : 04714_01456_Revision History vControl 1.1 SP1_V1.0

Author : Laszlo Ballabas

Storage place of valid version : DHF Number 04714_01456

Corresponds to vControl Version : 3.8.12.735

1.	vControl 1.0 SP1 (Build 3.7.15.224)	3
2.	vControl 1.1 (Build 3.8.7.729)	3
2.1	Bug Fixes	3
2.2	Changes	7
3.	vControl 1.1 SP1 (Build 3.8.12.735)	8
3.1	Bug Fixes	8
3.2	Changes	8

1. vControl 1.0 SP1 (Build 3.7.15.224)

Initial Version

2. vControl 1.1 (Build 3.8.7.729)**2.1 Bug Fixes**

Bug ID	Title
123445	After last Nested FCA Tip is fetched there is an empty error message Replace Used DiTi Rack
123729	[KomodoControl] Application SW cannot be closed properly when the instrument is in simulation mode
124208	Crashdump created while vControl was closed
125050	[KomodoControl][3DSim][WorktableConversion] NAP worktable cannot be converted completely in simulation
125127	[NGS] Run is paused after AddLabware command and cannot be continued in simulation mode
125978	It is not possible to transfer a plate to the ODTC in simulation
126124	[Teaching] Device placement for devices duplicated in 3.7.x not working
126252	DiTiBox Lid doesn't add in 3DSim while "Add Labware" command is being performed for the DiTiBox with "Has Lid" option checked. Simulation Mode
126332	[vControl][OneView]: Always date most in future is visible in OneView, not the next one due for a given maintenance method
127256	[vControl][3DSim] Channel Self Test Fails in Simulation
127495	Required driver to connect to the instrument not available, when only vControl is installed
127501	Moving RGA to the home position for DeckCheck twice leads to an error
127510	Orphaned labware is not displayed, when NGS application is imported
127517	[Veya] DeckCheck producing failed test results despite worktable being set-up correctly
127519	DB update after version update shows a lot of warnings for not transferable items
127563	[Teaching][Dev] Crash when transferring robot vector coordinates for incubator
127564	Generic Error when tips are tried to pick up from a box covered with a stacking plate
127646	Controlbar Filter file is not updated during an update installation
127698	Wrong coloring for Research Use Only in about box, startup and shutdown window
127721	[vControl]: Get Tips took too long error during a pipetting script execution
127747	[DesVal]: Run Abort when trying to put plates back to source in hotel
127814	[KomodoControl]Non-informative error message is displayed in case of unreachable RGA target location.
127825	[Veya] Door remain closed after the liquid detection error and recovery not possible
127827	[Scripting] In the "On Error GoTo Group" the worktable is not updated at edit time
127842	RGA arm not powered on while teaching leads to error message
127879	Method Recovery does not work, when a loading guide script is part of the method

Bug ID	Title
127905	[vControl][NGS] Liquid Class using DispenseMode ContactEdge defaults to Tracking
127906	[vControl][NGS] Liquid Class with DispenseMode ContactEdge does not display entered and saved DispenseEdgeOffsetXmm for ContactEdge in editor
127920	It's not possible to use variables for Labware group name in process editor
127935	[vControl] FCA Move Arm command throws runtime error if well selection spans multiple columns
127938	If the tightness block with holder is on the worktable, then the initialization is not working in simulation mode
127941	[VControl] [NAP] RUP blinking blue
127945	[NAP] Doorlock does not open during pause
127949	Fatal Error During Second Incubation
127950	IVD (GP) and RUO mode not working
127969	[Compliance] Electronic signatures not triggered on reverting to a restore point
127985	Information Dialog has wrong application name - FC instead vC for Scheduler Task Handling mode
127991	[VControl] [NAP] Liquid detection error recovery lead to script end
128010	[VControl] [NAP] Doorlock does not open for user interaction
128025	[Early Access] After aborting a run, VControl froze at restarting the same script
128043	[VControl] [NAP] Liquid detection error with aspirate anyway
128081	Wrong MoveDirection in LiquidClass
128083	[Early Access] Clicking on MoveTool x,y or z did not activate any movement
128085	[Early Access] After an AspirateAnyway step, plate volume is not updated
128090	Existing LC disappear after importing of other LCs, with different names, no conflicts + Recreation of the disappeared LC with the same name not possible anymore
128093	[vControl][NGS] Run Abort due to Aspirate Error
128118	Gx mode not working due to defect in WL for washer driver
128122	[VControl] [NAP] Error recovery of a Pressure Profile error leads to script abort
128125	Cannot Continue or Abort a Paused run
128135	vControl and File Version Checker does not check the relative path of dlls
128151	NullRefException during execution when stopped and resumed
128183	[KomodoControl] - Warning Fan speed of cooling fan out of range: Expected speed 150.00, actual speed 2730.00
128198	[LiquidClass] An error appears when save changes for the duplicated LC
128238	Sample Transfer Smart Command cannot create subscript
128239	VControl froze at previewing touch tool command
128242	Duplicating a default liquid classes changes more data than expected
128289	Remove Process without references leads to App crash.
128314	[NGS] Error during recovery leads to run abort
128337	Error during aspirate command when there are two different errors and "use next when empty" active
128362	[NGS] ODTC bounding box is too small for closed ODTC door
128366	vControl crashes sporadically when user logs in at startup window

Bug ID	Title
128375	Initialization is possible with open door after instrument safety manager dialog is canceled once
128385	SW is stuck in 'Initializing Step' after running a modified method that was previously recovered
128400	Recovery Point for Loading Guide Method is not created in case another recovery point is available for another method
128403	[Robustness] FluentControl cannot properly shutdown via API and has error at next start
128407	Some carriers cannot be converted correctly when they are in the "Add Labware" command in simulation mode
128421	[Scripting] 1x32 10 mm tube runner cannot be used in Sample Transfer smart command
128429	[SampleTracking] Assigns sample ID to 2 samples on Pressure out of Range error
128444	vC crash when the run finished successfully after recovering a method
128501	Cryptic Error Message when full access of FCA channels to labware is not possible
128505	[vControl][NAP]_TouchToolsDisappearingWhileEditing
128507	Scheduled Process is aborted when (needless) Incubator commands are used in Incubation processes
128540	Context check error of potential lid-plate collision while no risk is present
128542	[Scripting] Vectors of nests changes when moving them from one segment to another
128549	[Scheduler] Pathfinder error (RGA in colliding position) because task started before other task had finished
128559	Operator User has the right to ignore a "labware could not be detected" Error
128568	Section "Decimal digits" in "Balance weight" does not function properly
128609	[vControl][NGS]Transfer Labware reports collision dependent on protocol
128620	SW crash in Edit Mode
128623	Sample Transfer Smart Command does not re-use tips even if parameter Max. Tip re-use is set to >1 (independent of liquid class selection)
128625	Error during dispense into Deep Well 0.5ml since cannot find instance (probably comma problem)
128644	Presence Check for Transfer Labware failing in 3D Sim with Channel Grippers and tip boxes
128645	Transfer From Lower Deck Segment does not work in case obstacles are removed with Remove Labware command
128661	"An Error occurred during Aspirate Command" after adding a process dynamically to a run
128684	2ml sarstedt tube position in x is often off
128737	[vControl][Service Input] Status Light issues
128747	[3DSim] Drop Tips with 5ml Tips results in error after drop
128744	[vControl][Service Input] Status light does not change color in TouchTools commands
128753	ODTC error at 'Close door' command after a method was aborted with option to abort incubation automatically in 3DSim

Bug ID	Title
128754	Wrong teaching triangle is used for reference pin on channel 8 and teaching plate on site 1 of 4 landscape 7mm nest carrier
128759	Changes to pLLD parameters in the advanced section of liquid classes are not stored.
128760	The software crashes in edit time when PMP Threshold Values are out of range in the liquid class
128776	Sometimes RGA does not move to home position before SW shutdown
128810	Fatal Error when recovery is chosen for retry with new tips and no DiTis are mounted
128815	SW got stuck while showing preparing run overlay in recovery view
128819	[vControl] Instrument connection not possible
128821	[vControl] FCA missing in move tool
128866	[vControl][Service Input] Pipetting System Check can be executed with mounted tools
128885	Dispense took place in the wrong location
128898	[vControl] Script cannot be opened when Liquid Class is missing
128968	A 61mm Nest cannot be placed on the right of a Alpaqua FLX magnet
129005	Error while editing script: simulation failed at line x
129007	[vControl] Channel Self Test failure
129046	ELISA process has to stop when Incubator drawer is not closed properly
129055	[IAM] FC crashing when only password is entered to log-in in TouchTools
129091	[IAM] Login Token expires after 24h and leads to strange behavior, e.g. Multiple lock screens appearing
129092	VControl crashed at removing one row in LC correction volume
129106	Running SVN repair in new location with 3.7 deletes database folder and doesn't show correct run output
129120	[vControl] Error with 10µL Tips on 7mm Nest
129131	[vControl][CAT Service Feedback] Error picking up FCA Channel Grippers
129144	Run finishes with warning instead of Run aborted
129147	The File Version Checker Tool shows 'Error loading version files: No Version file found'
129150	Teach worktable command must check all teaching plate triangles to decide if site is accessible or not.
129189	vControl crash after clicking Retry on Error dialog
129193	If the edit usermanagement right is assigned to a role, the user does not see all data in user management
129206	Change Severity of the LogEntry to report 'RunEnd: to IoT from Error to Info
129214	"No access to Tool" message if Operator tries to select File Menu while a Tool Editor is open
129234	Deadlock in RGA Subroutine when main script runs into an error
129235	Deadlock in FCA Subroutine when main script runs into an error
129241	Serial number contains comma in real mode
129249	Run continues with open doorlocks after LoadingID command
129258	DiTi box is not automatically refilled in 3DSim in simulation mode

Bug ID	Title
129307	[IAM] TouchTools User login does not allow input of the user name with screen keyboard
129343	[vControl] Error "TargetPosition is out of valid position range" when trying to get DiTi and get GripperFingers.
129347	Error was thrown while executing aspirate command execution because of violation of minimum distance between channels
129392	It's not possible to use IPP for an Aspirate command including x-Offset values
129398	Open door does not prevent a resume of the run
129409	[Teaching][Scan Location] Cannot select carrier by clicking and no context check for Selected Location
129426	Invalid String in notification Info on TTS
129435	[vControl] Context Check error in Pipetting Smart Commands to drop fingers even though fingers are already dropped.

2.2 Changes

ID	Title
98143	[DevBranch][UMS] User Roles and Rights need to be updated
98497	[DevBranch][UMS] No warning appears if admin tries to lock himself in the User Management
114173	[Scripting] Add Labware cannot add lid onto plate if plate is located in storage carrier
114680	[Scripting] Sample Transfer cannot select multiple runners if source labware is a variable
116043	[Robustness] FluentControl does not start causing downtime
123110	[KomodoControl][Performance] After each SW start a full instrument initialization is executed
124746	[KomodoControl][LiquidHandling] MAP automated channel placement difficult to handle with 25ml- and 600ml trough
124848	[LiquidHandling]- Support of custom LiquidProperties in vControl
124853	[KomodoControl][Scripting][DesignVal] Parser function 'SerialNumber()' is not available in vControl
124975	[vControl][Usability] User expects to find Audit Trail file (or link) within Application SW
125351	[KomodoControl][GUI] SW ignores Plate to Lid offset for placing and picking (but only in x)
125488	[ErrorRecovery] FCA doesn't fully retract before moving to the next column for DiTi pick up
125652	[Scripting] Store Well Concentration command improvement
125722	[Dev-Branch][UMS] Integrate MAP IAM (new UMS) in vC & FC
125724	[Variables][Scripting] "Replace Empty Tip Box" command to facilitate Tip Box stacking
125727	[Worklist] Load Worklist command contains "Wash Station Waste_1", which is a FluentControl item as default
125803	[LiquidHandling] Excess volume shall be dispensable back to source tubes
125883	[KomodoControl][3DSim][Custom Models] Ability to customize models in simulation
126017	[InhecoIncubator] The drawer remains opened during device initialization

ID	Title
126035	[DevBranch][UMS] MAP IAM User Migration of old UMS Users
126105	New Licence for Assurance Software for RUO
126121	Operator shall be permitted to replace inline filter and verify functionality without technical support
126223	[vControl][OneView]: Status not consistent in OneView
126312	[vControl][Teaching] Teaching sites to be created for ODTC, CPAC-PCR, (Teleshake-PCR)
126621	[OneView]: Editor for adjusting entries from config file - Avatar & lab contact
127175	[KomodoControl][Service Feedback] User shall be in the position to execute teaching command from the script
127266	[KomodoControl][ErrorText] No Error Dialog is displayed when the Target position is not reachable
127540	Push ARP Pipetting System Check results and warnings to IoT
127751	vControl Liquid Classes - Missing options
127820	[DesVal]: 3D sim config not available for 9 Grid, 8 channel FCA with RGA
127876	Lid handling (uncover plate) with absolute robot vectors does not work correctly
128159	[NAP] SW reports unnecessary SampleTracking error during EmptyTips command
128262	Non-manufacturer Carrier/labware data which is not part of the DB, shall be imported after an update installation automatically
128480	[Scripting][Usability] Variable name does not turn blue after clicking Declare button
128516	[Veya] Teleshake with 3mm orbit position adjustment
128570	[KomodoControl][LiquidHandling] Add attribute for Restricted Pipetting Area 25ml- and 600ml trough
128641	Stacked tip box transport issue
128787	[Compatibility] Update MasterMedia Contents of vControl V1.1
128796	[Veya] Change to LCs with default cLLD sensitivity set to 0.95
128915	Integrate IoT client 3.1 in vControl 1.1 and FluntControl 3.8
129195	Add WetDispense to EmptyTipsDispenseMode and Infobutton beside the DispenseMode combobox

3. **vControl 1.1 SP1 (Build 3.8.12.735)**

3.1 Bug Fixes

Bug ID	Title
129990	vControl cavity attribute "Volume" not updated correctly

3.2 Changes

ID	Title
130173	Liquid level detection sequence mode shall be set to sequential