

vControl

Revision History

Document name : 04714_01578_Revision History vControl 1.2_V1.0

Author : Laszlo Ballabas

Storage place of valid version : DHF Number 04714_01578

Corresponds to vControl Version : 3.9.4.1178

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
4. vControl 1.2 (Build 3.9.4.1178)	8
4.1 Bug Fixes.....	8
4.2 Changes	11

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
127905	[vControl][NGS] Liquid Class using DispenseMode ContactEdge defaults to Tracking

Bug ID	Title
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
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

Bug ID	Title
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
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

Bug ID	Title
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
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

Bug ID	Title
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
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

ID	Title
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

4. vControl 1.2 (Build 3.9.4.1178)

4.1 Bug Fixes

Bug ID	Title
124535	vControl switches into GP color scheme, although being in demo mode
125356	[3DSim] Collision with tilted vial runners during aspiration
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

Bug ID	Title
127280	[API] FluentControl becomes not responding when try to open Settings in Menu
127937	FluentControl crashes when starting a method just after it finished executing one
128387	Recovered Run cannot be finished. 'Ending run is currently not allowed'
128606	Disposable tip status is not correctly displayed or reset when restarting a method
129050	[Robustness][Scheduler] Runtime errors caused by memory leak
129105	Doorlocks are not opened when LoadingID runner needs to be reloaded after a system pause
129178	[Compliance] Audit trail does not properly display Chinese method names
129236	Deadlock on Startup
129300	[vControl] Status light not active during run
129390	Aspirate failed error during execution of Nap simulation script
129412	[vControl] Error after Teach Worktable Command Channel cannot be positioned inside Ranges
129478	Well Offset in Sample Transfer does not work for runners if more than one runner is needed
129494	[Teaching] Placement issues after upgrade from 3.5 to 3.7 SP1
129507	Deadlock in FCA/RGA Subroutine when pausing and resuming the run during pathfinding
129514	Run is not aborted, when the incubator door does not close when the plate is brought to the incubator
129546	Move To in Labware Editor not functional in RC3
129553	Aborting a method triggers an unclear error "The script was automatically aborted by the system"
129557	Pressing F1 in user administration shows File not found
129559	Block UI until 3D Sim is connected
129588	[Compliance] Failed to open the Data Audit Tool
129620	Vial access on "2x12 3ml to 4ml Tilted Vial Runner" denied
129627	Instrument connection is not established on first attempt quite often
129631	9 Position Base Plate cannot be saved after editing
129717	VControl crashed at take reference scene
129732	Move to does not work when the channels are moved up too high
129826	Error message when you try to open the 3rd Party license file from about box
129832	After an update installation, the worktable conversion in simulation is not working properly
129863	Not possible to aspirate from a 6 Well Flat Plate
129865	Modifications in Dimensions section in my carrier are not persisted after successful saving, closing and opening again - occurs sporadically
129896	Complete Filter Change command has no context check error for missing tightness block
129906	Step by Step execution and breakpoints are not working for immediate finished commands
129911	Alpaqua Magnum FLX cannot be converted in simulation mode
129915	Add Labware does not work in combination with eppendorf tubes and 5 Pos 1.5ml tube adapter in sim mode
129923	It is not possible to transport a plate to an alpaqua magnum flx carrier in simulation mode

Bug ID	Title
129930	Error "Target Position is out of valid position range" and led to run abort
129934	Unable to re-initialize and restart of instrument was needed when error "Grip distance violated between channels" occurred
129969	vControl crashes after Copy/Paste of SystemTraceView
129973	Tip selection has no effect on well selection after reloading a script
129977	TouchTools config file missing after installation in some cases
130024	Stacking Plate cannot be transferred on a 1000µl DiTi box
130059	Sample Transfer well selection changes itself after re-open script (beta version at customer site under NRSW agreement)
130067	Liquid Classes not in dependencies of script export
130070	200 µl Filtered SBS box not detected at 7 mm nest after transfer from another 200 µl Filtered SBS box
130096	Many subfolders are not shown at Save as
130137	vControl got stuck in the loading screen
130142	9Grid_XCA_4Channel_XGA Instrument configurations has different module version than other instrument configurations
130154	First click on a move button in move tool is not registered, when Switch Zero G Button is used
130155	Timeout when the move tool button to move 1mm in x is used with modulated speed
130164	Warning "System.Security.Cryptography.CryptographicException: The system cannot find the file specified." is overloading the log
130166	Connection to instrument cannot be established, when the startup is blocked by a temporary not running audit trail service
130168	A warning message pops up when recovering a method with Tips mounted
130222	[vControl]: Reagent distribution command shows needle instead of DiTi
130244	vControl crashes after upgrade
130250	The "Reset all" button in the disposable tips dialog does not reset the position of the stacking plate
130266	Unnecessary Move after LLD error handling
130351	[3DSim] Labware missing MapTemplateName
130370	Misalignment of the paddle station
130377	Move Tool - Small 'Up' arrow button moves the arm down (and vice versa)
130395	Aspirate with well offset does not work when it has to switch to the next plate in a labware group
130415	It is not possible to change the color for FCA and RGA commands
130430	After successful initialization the run is not started showing an error as not initialized
130472	[Dev] Memory leak in ContextCheckController (Dev branch)
130508	While executing dispense command in script, error message shown only in system trace view, vcontrol status paused and it's frozen
130512	[IAM] Multiple lock screens are appearing sporadically
130540	Start Loading command text not readable in light mode
130551	One or more environmental sensors are not functional
130552	vControl freezing in the middle of a run

Bug ID	Title
130553	New value instead of old value is transferred into subroutine with Register Event Handler command
130558	Some items cannot be converted
130561	Selected tips are not reflected in well selection properly
130572	[Performance] Importing zeia files containing external files causes internal import error
130659	vControl IVD and non IVD graphics are shown both at the same time
130690	After pressing pause at the beginning of a command, it is not possible to stop the run
130700	[Dev] Erroneous "Data set out of sync" error reported at startup
130754	Imported Script is automatically approved after import
130764	Dialogue to validate checksum is not showing up
130792	While dropping 5 mL tips, vControl always handles Channel 8 separately
130802	Starting vControl after an update installation triggers the installer again
130803	There is no automatic restart after the DUF import
130804	It is not possible to change the well selection when several tips use the same well
130871	[vControl] Unhandled exception and application crash after instrument communication loss
130956	vControl cavity attribute "Volume" not updated correctly
130994	[vControl]: ArgumentException on Empty Tips Command Due to Invalid heightInCavity Parameter
130174	UI issue with tube runner in Sample transfer command
130696	The system tries to initialize with gripped plate when method is recovered after a collision
130754	Imported Script is automatically approved after import
131046	Data out of sync error

4.2 Changes

ID	Title
115771	[TipHandling] On Empty trigger scenario with Hotel is being occupied.
117680	[Cybersecurity] Weak hash is used for checksum: Protection Commands
123006	[Performance] Replace nuget package Microsoft.Office.Interop.Excel with better performing Package (DocumentFormat.OpenXML)
124831	Harmonize UI of vControl
125068	[FCA] Remove workaround for unreachable source position due to minimum channel distance violation
125544	[Performance] SetVariableStatement performance issue
125652	[Scripting] Store Well Concentration command improvement
127163	[Driver][Compatibility] Support driver development for drivers with x64
127244	[Compatibility] Update MasterMedia Contents of FluentControl 3.8
127480	[Visualizer][Scripting] INSTRUMENTDoorOpenedError in Simulation
127959	[Compatibility][SampleTracking] FluentControl and vControl shall install latest Sample Tracking
128299	[ReleaseBranch][Performance] Improve usability Import workflow
128422	[vControl] cLLD curves shall be loggable

ID	Title
128479	[SnapshotTool] Troubleshooting data related to MAP IAM shall be automatically collected
128647	New FCA command "Tip Adapter Check" to Implement execution of ventilation/cone tightness test
128679	[vControl][NGS] Improvements for mixing: definition of 2 volumes for mixing
128682	[vControl][NGS] Improvements for mixing: splitting of final dispense volume and adding delays
128710	[Performance] Improve Edit Time
128712	[Performance] Improve FluentControl start up time
128787	[Compatibility] Update MasterMedia Contents of vControl V1.1
128867	[vControl][Service Input] Pipetting System Check shall return success message
128876	[Field Feedback]: Allow going to >1 plates in parallel
128878	[Early Feedback]: Ability to query temperature, humidity and ambient pressure with a function in order to produce a report
128904	Bottom detection liquid class
128980	[Veya] Channel self check shall be replaced with inline filter test
128981	New FCA command "Complete filter change"
129033	[Robustness] FluentControl cannot start because of problem with file ChangeLog.xml in attic folder
129200	[Cybersecurity] FluentControl can be blocked by IT security software because it is not signed as a verified software
129264	Alpaqua FLX Magnet is updated to the latest version.
129269	Allow Tip- touching in Liquid Class for dispense
129270	Add Method progress list to OneView
129272	Add humidity chart for OneView
129273	[Usability] Calculate runtime based on command parametrization
129321	[Troubleshooting] vControl shall write installation log by default
129324	[Scripting] Create new Register Event Handler command
129339	[Troubleshooting][Performance] Switch default for Log Level to Debug
129341	[Database] Update 1.5ml eppendorf tube definition to include the correct cap
129402	Update vControl Icon
129599	Default screen for TouchTools for Veya shall be No.1
129605	[Neptune][vControl][XLS] Integrate XLS Functionality in the Application
129653	[Database] New stacking plate adapter definition
129653	[Database] New stacking plate adapter definition
129666	[Compatibility] Update MasterMedia Contents of vControl V1.2
129700	[KomodoControl][3DSim][WorktableConversion] Integrate missing 3D Sim components for application packages - Part 1
129827	Add "MountOffset" custom attribute for XCA
129891	MapTemplateName is directly linked to the 3D Sim worktable conversion, these should be independent from another.
129929	vControl IVD intended use shall be implemented the same way as the Fluent IVD intended use
130017	[Usability] Zero-G Button state is not updated correctly when Zero-G is activated or deactivated
130098	[xRuntime] Upgrade to .Net 8

ID	Title
130230	Liquid level detection sequence mode shall be set to sequential
130361	[ETON] Allow definition of Tip types defined in MAP
130385	[Performance] Improve Edit Time
130394	[Performance] Improve performance of Context Check
130421	[Database] New 6 Grid Base worktable containing sites for hotel in the back
130470	[Dev][Performance] Improve usability Import workflow
130870	Veya] Sequential liquid level detection to removed