

Software release note



File Name : SRN PSS0052-04 Archimedes Software.docx

Page: 1 of 5

1. Version Control

Part	Description	New Issue	Previous Issue
PSS0052	Archimedes Software	1.21	1.20

2. Release Control

ECN Number:		
Compatibility:	Windows 7 32-bit, Windows 7-64 bit	
Media:	DVD and Extranet Download	
Operating Systems Tested On:	Partially (e.g. installed, measurement taken, printed)	Full alpha test
	Windows 7-64 bit and 32 bit	

Software release note



File Name : SRN PSS0052-04 Archimedes Software.docx

Page: 2 of 5

3. Installation Instructions

3.1. Installing the Software/Firmware

It is assumed that you have authority to install or update software within your company's SOPs. If you do not have this authority please consult with your I.T. support department before proceeding.

It is assumed that you have Administrator rights for the computer. This is required by the installation process.

It is recommended that any previous Archimedes Software version be uninstalled before version 1.20 is installed. This should be done by following the uninstall instructions in section 3.3.

Do not restart the computer until the entire install process is complete – select "Restart later" when prompted.

3.2. Connecting the Archimedes to the PC

With the software installed, the instrument should be connected via the USB port, and the system switched on. Once the connection is detected, the Windows system will automatically start to install the drivers for the Archimedes hardware.

Once the driver installation is completed, Windows will indicate that the new device driver software is installed and ready to use.

3.3. Uninstall procedure

It is important to uninstall all previous versions of the software prior to attempting to install the new Archimedes software on any computer.

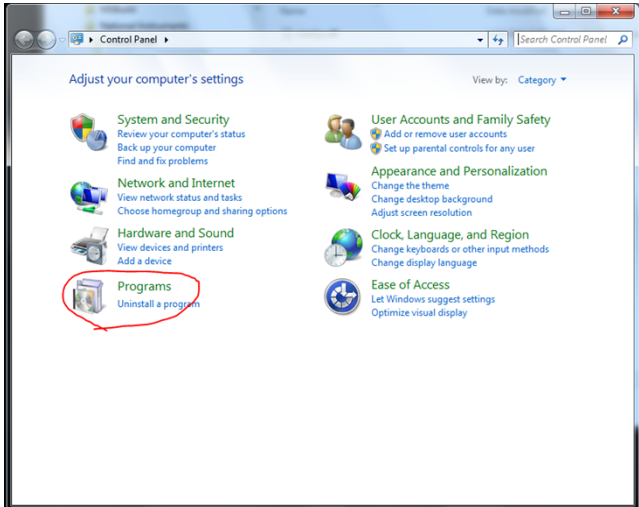
Previous versions of the "Archimedes" software were called "Particle Lab" software and it can be removed by using the standard Add/Remove feature in the Windows 'Control panel'.

Some of the software is based on algorithms completed with National Instruments software and it is important this is also removed prior to attempting any new installation. Details of how to remove this software can be found in the section, "Uninstalling National Instruments Software".

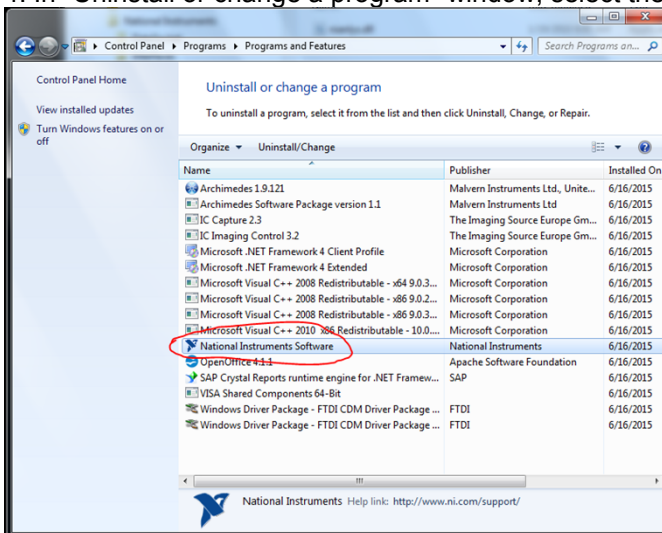
Uninstalling National Instruments software

1. Restart the computer to make sure that no application or relevant driver is actively running while doing installation.
2. Go to the Windows "Control Panel" (Start -> Control panel);
3. In the "Control Panel" select the "Uninstall a program" to open the "Uninstall" window;

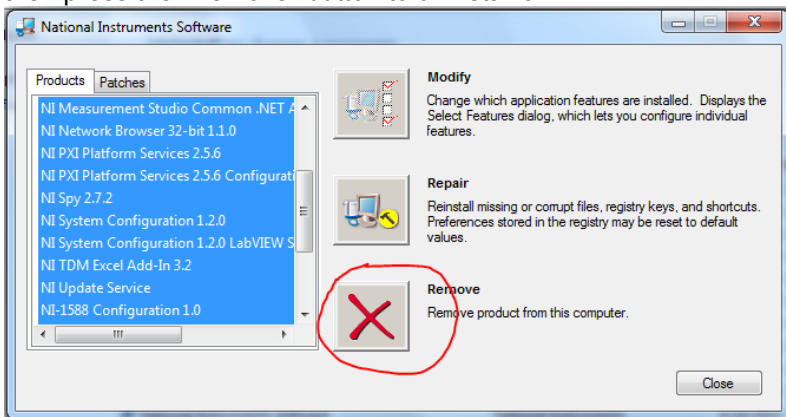
Software release note



4. In "Uninstall or change a program" window, select the "National Instrument Software" to uninstall;



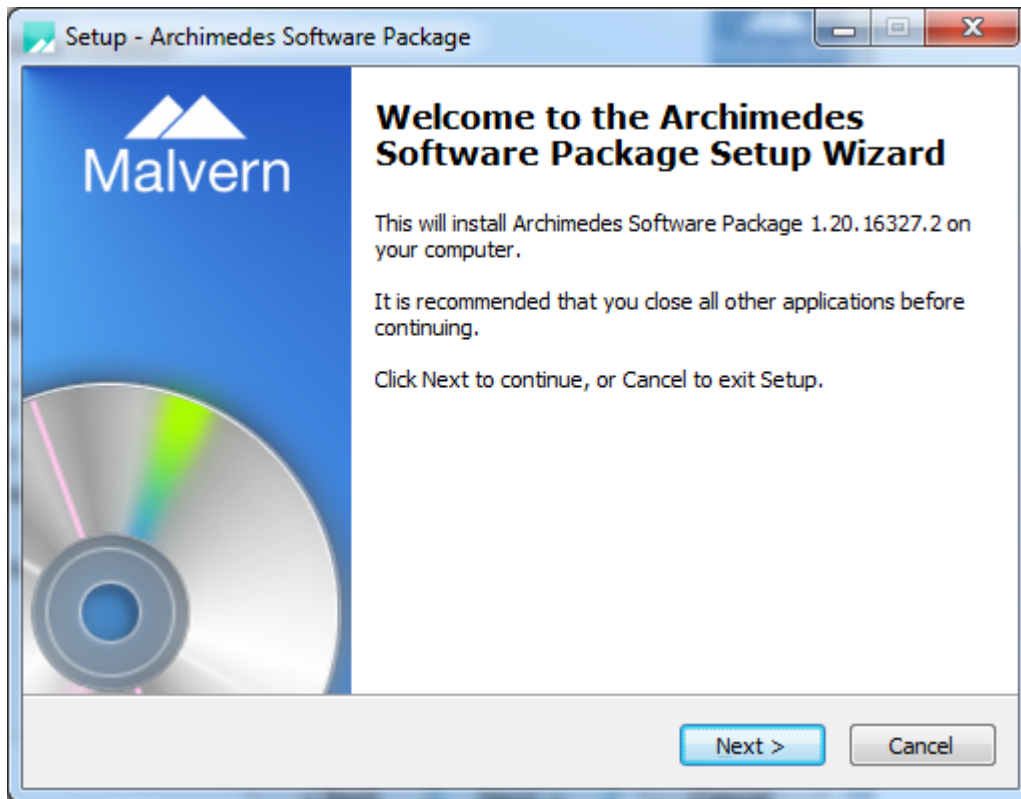
5. In the following dialog box for the "National Instrument Software", select all of the items in the "Products" tab, then press the "Remove" button to uninstall all.



Software release note

- When the uninstallation of the "National Instrument Software" is completed, follow the prompt message to restart the computer again.
- After the machine is restarted, install the Archimedes software.

Do not restart the computer until the entire install process is complete – select "Restart later" when prompted.



- When the installation is completed, follow the prompt message to restart the computer again.

4. Known Issues

Reference(s)	Issue	Workaround
N/A	Issue where when user pressed "stop" button during the measurement, all buttons turned gray.	Restart the software
N/A	When trying to program the board, were unable to disable the tracking whereas it could be done on the old version of Particle Lab on a different PC	Close the software, right click on the icon and choose 'Run as administrator'
N/A	Trouble installing the NI components during installation of Archimedes Software – seen on PCs with prior versions of Particle Lab or Archimedes software installed.	Restart the computer, follow the steps necessary to uninstall 'National Instrument Software' through the Windows 'Control panel,' restart the computer and install Archimedes Software.

Software release note



File Name : SRN PSS0052-04 Archimedes Software.docx

Page: 5 of 5

Bug 47652	After changing sensor, the tune plot and Q value for the previous sensor persists	Optimizing the amplitude and tuning the sensor are required when using a new sensor.
NA	Software is grey after starting despite all hardware being attached, log file states: InvalidOperationException: Cannot load Counter Name data because an invalid index '' was read from the registry.	Click Start, type cmd right click cmd.exe, and select Run as administrator. At the prompt, type lodctr /r and press ENTER.

5. Change Control

5.1. Changes made via PBI or Bug Report (TFS)

ID	Work Item Type	Title
73013	Bug	Build date incorrect in 'About Archimedes' window
73019	Bug	Copyright out of date in 'About Archimedes' window
73050	Bug	Calibration experiment 'select' button cut off on the right-hand side
73107	Bug	'Edit Calibration source' window can be maximised
74812	Bug	If a report shows positive buoyancy data, any additional analysis data will not show "Limit of Detection"
74880	Bug	Software cannot connect to DMx cameras
75053	Bug	Exporting positive buoyancy particle data has blank LOD
76241	Product Backlog Item	Changes to support Pico sensors