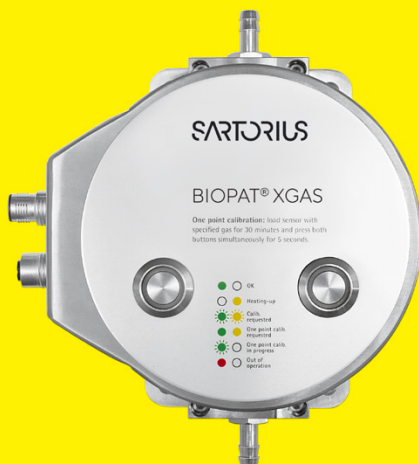


Operating Manual

BioPAT[®] Xgas Vis

Software Manual



85037-545-79



SARTORIUS

Contents

1	About This Document	5
1.1	Accompanying Documents	5
1.2	Symbols Used	6
2	Safety Information	7
2.1	General Safety Precautions	7
2.2	Intended Use and Foreseeable Misuse	7
3	Overview	8
3.1	Requirements	8
3.2	Contents of Delivery	9
4	Installation	10
4.1	Installation Steps	10
4.2	Hardware Installation	12
5	Software Installation	14
5.1	Preparation Steps	14
5.2	BioPAT® Xgas Vis Software	19
5.3	Perle Device Driver.....	24
5.4	TruePort Application	28
6	Hardware Settings	32
6.1	Checking Jumper Settings at the Perle Device	32
6.2	Resetting Perle Device.....	33

7	Configuration	34
7.1	Perle Device	34
7.2	Serial Port	41
7.2.1	Configuring Serial Port with TruePort Management Tool	41
7.2.2	Configuring Serial Port with Device Manager	47
7.3	BioPAT® Xgas Vis OPC License	49
7.4	BioPAT® Xgas Sensor	51
7.4.1	Configuring One Sensor	51
7.4.2	Configuring Multiple Sensors with one Ethernet Modbus Converter	56
7.5	Changing the Modbus Address	58
8	Installation of BioPAT® Xgas OPC Server	61
8.1	Installing MFCS/win OPC Client	61
8.2	Configuring MFCS/win Configuration Management	62
8.3	Browsing the BIOPATXGAS.OPC.1 Server with OPC Test Client	69
9	Firmware Version	72
9.1	Checking Upgrading Firmware Version of the Perle Device	72
9.2	Configuring Settings for Windows ModbusService	76
10	Troubleshooting	78

1 About This Document

All information and instructions in this operating manual have been compiled in consideration of applicable standards and guidelines, the state of technology and our many years of experience and knowledge.

These operating instructions provide you with all the information necessary to install and to configure the BioPAT® Xgas Vis Software.

In addition to the operating manual, all generally valid, legal and otherwise binding regulations for accident protection and environmental protection of the country of use must be observed.

The operating manual must always be kept at the place of use of the device.

1.1 Accompanying Documents

The operating manual describes how to install and to configure the BioPAT® Xgas Vis Software.

In addition to these operating instructions, follow the information in the operating manual "BioPAT® Xgas" and in the documentation of the other software and hardware components.

1.2 Symbols Used

NOTE

- is an indication of a function or setting on the device.
 - is an indication that caution should be exercised while working.
 - identifies useful information.
-

The following presentations will also be used:

- Texts that follow this mark are lists.
- ▶ Texts that follow this mark describe activities that must be carried out in the specified order.
- ▷ Texts that follow this mark describe the result of an action.
- “ ” Texts in quotes are references to other chapters or sections.
- [→] Texts following this symbol make reference to other chapters, sections or documents.

2 Safety Information

2.1 General Safety Precautions

The BioPAT® Xgas Vis Software may only be installed and configured after gaining familiarity with this operating manual.

Always keep the operating manual in the location where the device is in use.

In addition to the operating manual, follow all general and local regulations for accident prevention and environmental protection.

2.2 Intended Use and Foreseeable Misuse

The operational safety of the device is only ensured when it is used for its intended purpose and operated by trained personnel.

Application

The device is only intended for industrial, commercial, training, and research facilities.

The BioPAT® Xgas Vis Software is intended for use with the BioPAT® Xgas Sensor and the BioPAT® MFCS/win Software.

3 Overview

With the BioPAT® Xgas hardware, the O₂ and CO₂ concentration can be monitored and controlled for any given bioprocesses in real time.

The BioPAT® Xgas is a dual oxygen-carbon dioxide sensor featuring automatic compensation for humidity and pressure. The oxygen concentration can be measured in a range of 1–50 % and the carbon dioxide concentration in a range of 0–10 %.

The BioPAT® Xgas Sensor can be directly connected by analog signals to the DCU system. Therefore, only a cable and two free analog input ports at the DCU systems need to be available.

Alternatively, if there are no available analog input ports on the DCU system, then it is possible to connect the BioPAT® Xgas hardware using the BioPAT® Xgas Vis Software interfacing to the MFCS/win SCADA software. Herein, the direct connection of the BioPAT® Xgas Sensors to the MFCS/win software with all necessary hardware and software components is described.

3.1 Requirements

- The BioPAT® Xgas Vis Software requires MFCS/win Version 3.0
- The Firmware version of the used Perle device is minimum 4.5

Computer Configuration

General	CD ROM drive
	USB Port
Operating System	Microsoft Windows XP (32 bit)
	Microsoft Windows Server 2003 R2 (32 bit & 64 bit)
	Microsoft Windows 7 (32 bit & 64 bit)
	Microsoft Windows Server 2008 SP2 (32 bit & 64 bit)
	Microsoft Windows Server 2008 R2 Sp1 64 bit
Operating System Language	English, German

3.2 Contents of Delivery

The delivery of the BioPAT® Xgas Vis Software contains the following components:

- CD-ROM, containing all necessary software items
- Software activation Dongle
- User Manual for the BioPAT® Xgas Vis

4 Installation

4.1 Installation Steps

The following table presents an overview on the necessary steps to install and configure the BioPAT® Xgas Vis Software. Please follow the recommended order in the table step by step.

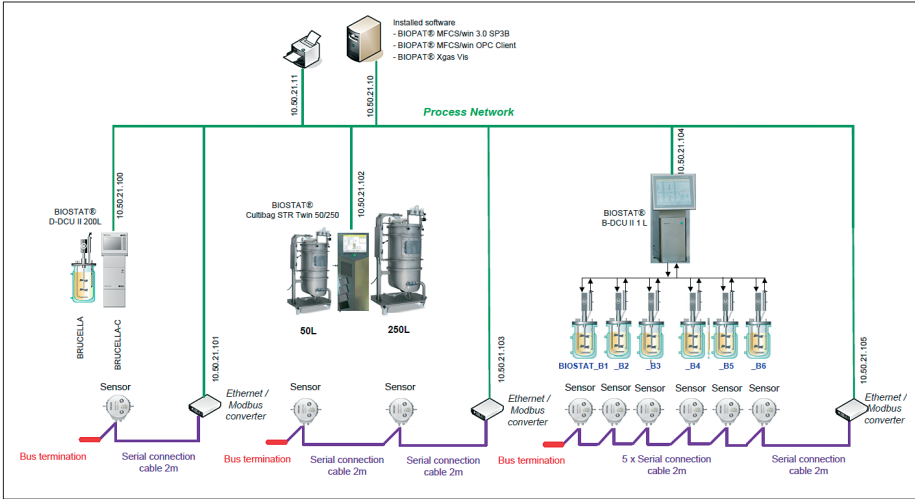
Each entry refers to a separate chapter or sub-chapter in this handbook where the procedure is described in detail.

Step	Installation Procedure	Remark
0	Hardware installation [→4.2, page 12]	
1	Preparation step for installation on Windows XP and Windows Server 2008, only [→5.1, page 14]	Only for Windows XP and Windows Server 2008
2	Software installation of BioPAT® Xgas Vis Software [→5.2, page 19]	
3	Installation and configuration of the Perle device driver [→5.3, page 24]	
4	Installation of TruePort application [→5.4, page 28]	
5	Checking hardware jumper settings at the Perle device [→6.1, page 32]	
6	Reset Perle device [→6.2, page 33]	
7	Configure Perle device [→7.1, page 34]	
8	Configure serial port with TruePort Management Tool [→7.2.1, page 41]	

Step	Installation Procedure	Remark
9	Configure serial port with Device Manager [→7.2.2, page 47]	
10	Activate the BioPAT® Xgas Vis OPC License [→7.3, page 49]	
11	Configuration of one BioPAT® Xgas sensor [→7.4.1, page 51]	
12	Configuration of multiple sensors with one Ethernet Modbus Converter [→7.4.2, page 56]	
13	How to change the Modbus Address [→7.5, page 58]	Multiple sensors only
14	Installation of BioPAT® Xgas OPC Server [→8, page 61]	
15	Installation of MFCS/win OPC Client [→8.1, page 61]	
16	Configuration work inside MFCS/win Configuration Management [→8.2, page 62]	
17	How to browse the BIOPATXGA.OPC.1 Server with OPC Test Client [→8.3, page 69]	Optional
18	How to check upgrade the firmware version of the Perle device [→9.1, page 72]	Update only necessary, if firmware < 4.5
19	Configure settings for Windows Modbus Service [→9.2, page 76]	

4.2 Hardware Installation

- ▶ Install the BioPAT® Xgas according the operating manual “BioPAT® Xgas”, delivered together with the BioPAT® Xgas hardware.
- The installation is described in chapter “4. Installation”.



Connection of BioPAT® Xgas to BioPAT® MFCs/win with BioPAT® Xgas Vis Software

Accessories

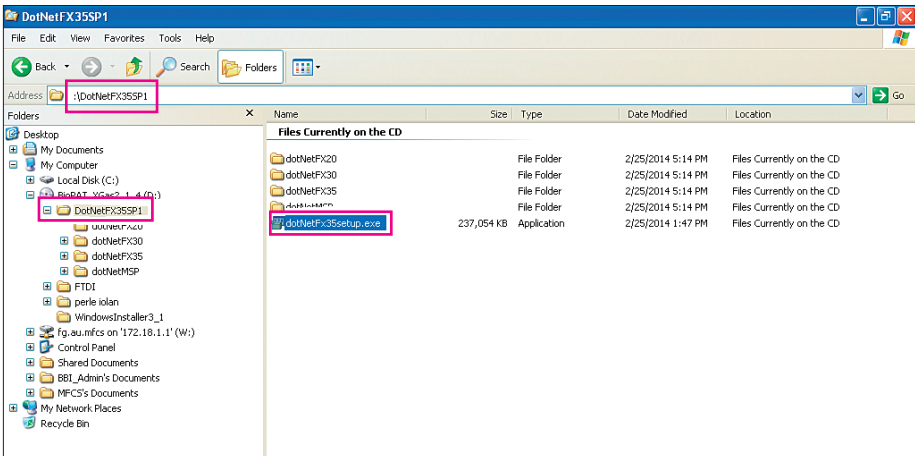
The following articles are available for the BioPAT® Xgas Vis Software:

Article Number	Description	Remark
BPX0009	BioPAT® Xgas Vis Software	Can be used with one MFCS/win server system.
BPX00(11-15)	BioPAT® Xgas Sensor	Includes power supply & cable and article number depends on vessel size and corresponding gas flow path.
BPX0018	PC connection package (includes Ethernet Modbus converter and Bus termination)	One is needed for each DCU system or independent vessel fitted with BioPAT® Xgas [see figure in chapter →6.2, page 33].
BB-34147295	Serial Modbus Connection cable 2 m	One cable (2 m or 5 m) is needed to connect each sequential BioPAT® Xgas unit.
BB-34147298	Serial Modbus Connection cable 5 m	

5 Software Installation

5.1 Preparation Steps

For Windows XP, only

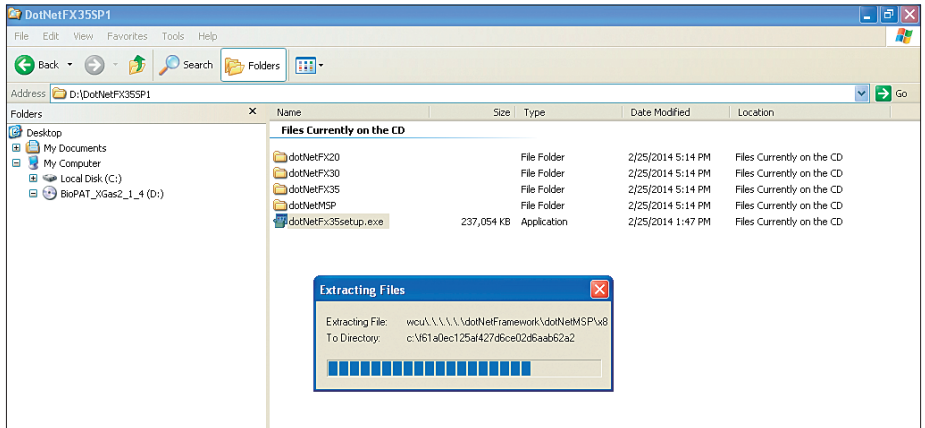


- ▶ Start “:\\DotNetFX35SP1\\dotNetFx35setup.exe” from the XGAS-CD.

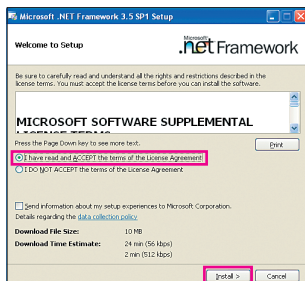
NOTE

NET Framework 3.5.1 must be installed manually from the BioPAT® Xgas VIS CD.

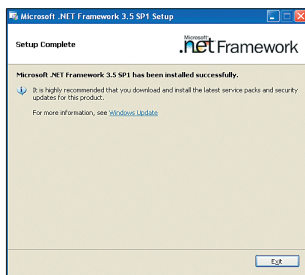
- ▶ Open Windows Explorer and browse to the CD into the directory.
- ▶ Double click the file “:\\DotNetFX35SP1\\dotNetFx35setup.exe”.



► Files will be extracted.



► Accept the terms and press “Install” button.



► Press “Exit” button.

► Continue with chapter [→5.2, page 19].

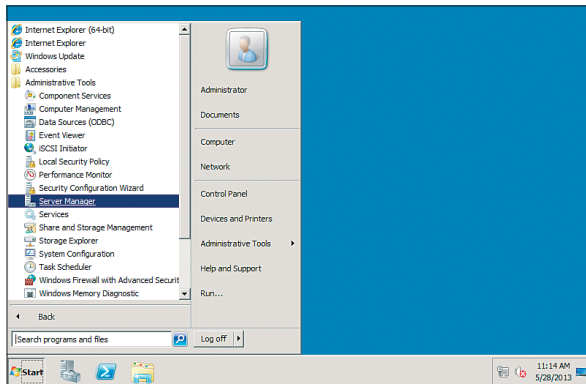
For Windows Server 2008 R2, only

NOTE

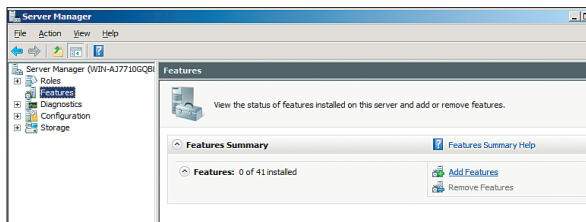
NET Framework 3.5.1 must be activated for Windows Server 2008.

- Windows Server 2008 includes the necessary .Net framework.

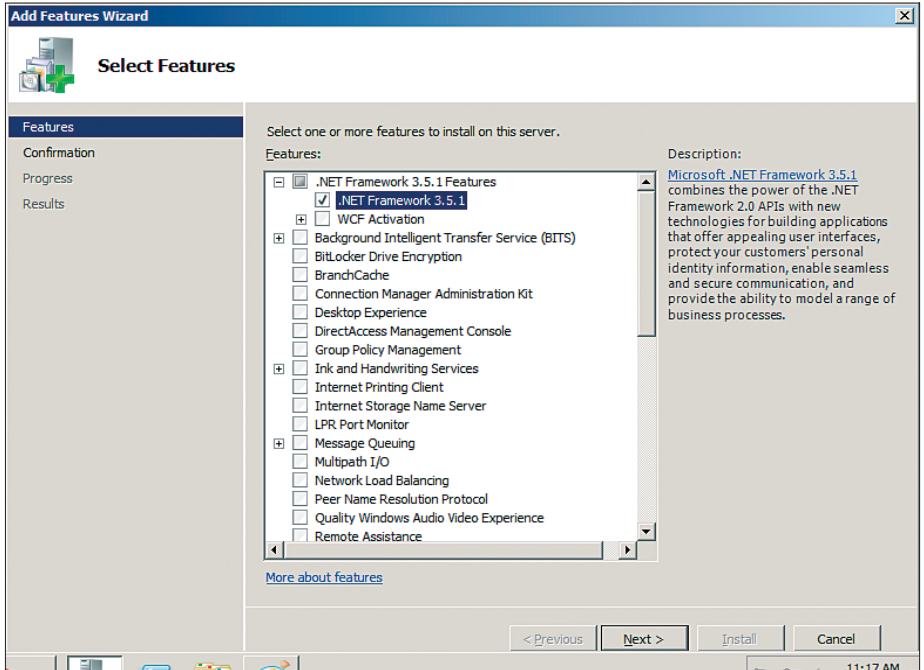
Follow the steps below in order to activate the .Net framework.



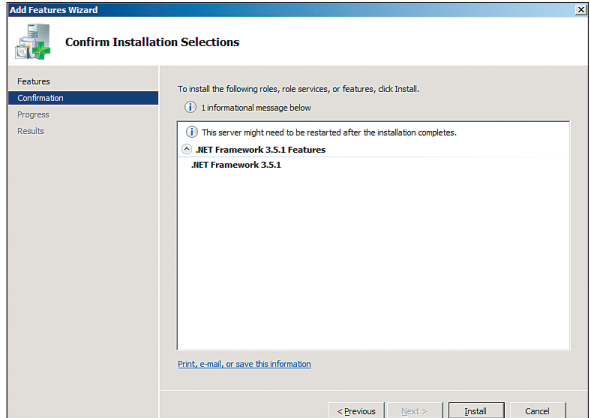
- ▶ Go to “Administrative Tools”.
- ▶ Start the “Server Manager”.



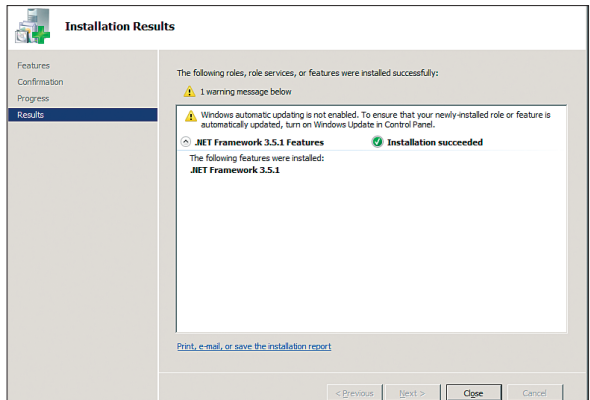
- ▶ Open “Roles / Features”.
- ▶ Click “Add Features”.



- ▶ Expand “.NET Framework 3.5.1 Features”.
- ▶ Select “.NET Framework 3.5.1”.
- ▶ Click on the “Next” button.



- ▶ “Confirm Installation Selection” opens.
- ▶ Press button “Install”.



- ▶ The installation is completed; the installation message “Installation succeeded” is displayed.
- ▶ Click “Close” button.

5.2 BioPAT® Xgas Vis Software

This chapter describes the installation of the BioPAT® Xgas Vis Software.

NOTE

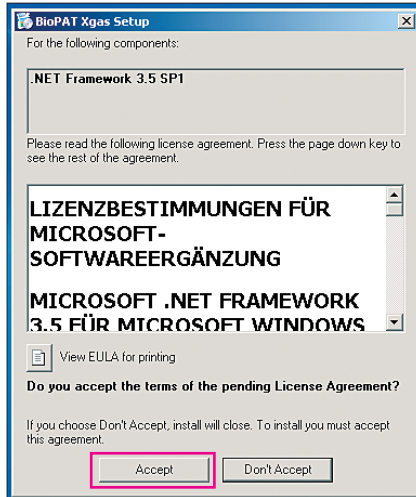
Please consider that MFCS/win must have been already installed. Please note, the installation of this software requires local administrative rights.

The CD-ROM containing the BioPAT® Xgas Vis Software is labeled as “BioPAT® Xgas Vis Software”. A separate program named SETUP.EXE performs the software installation.

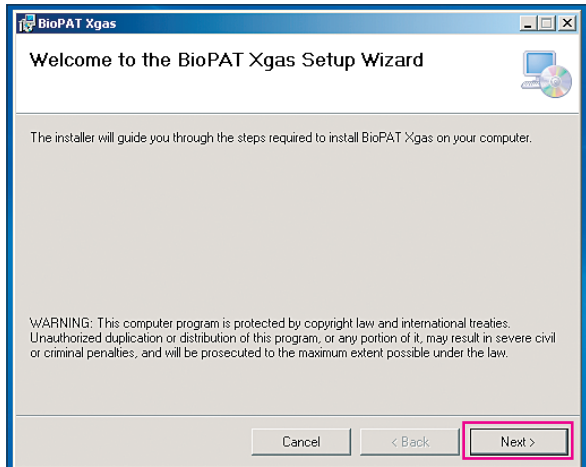
The following steps describe the installation on your computer system.

Installing the BioPAT® Xgas Vis Software

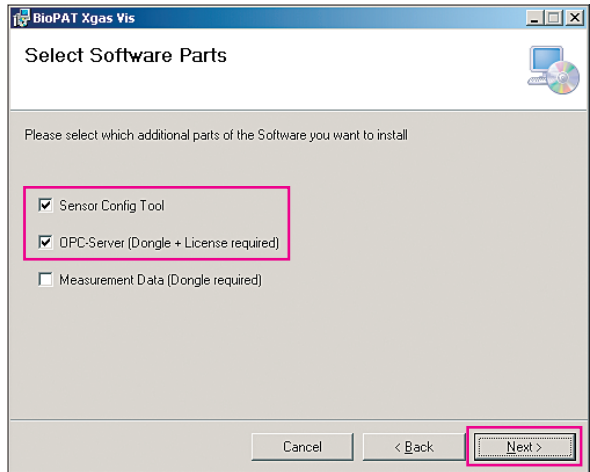
- ▶ Shutdown the MFCS/win system from the MFCS/win Shell.
- ▶ Insert the CD-ROM into the CD drive.
- ▶ Open the windows explorer and change to the CD drive.
- ▶ Double-click on file SETUP.EXE to start the installation program.
- ▶ If the “.Net Framework 3.5” is not installed, then the following dialog will be shown on the screen.



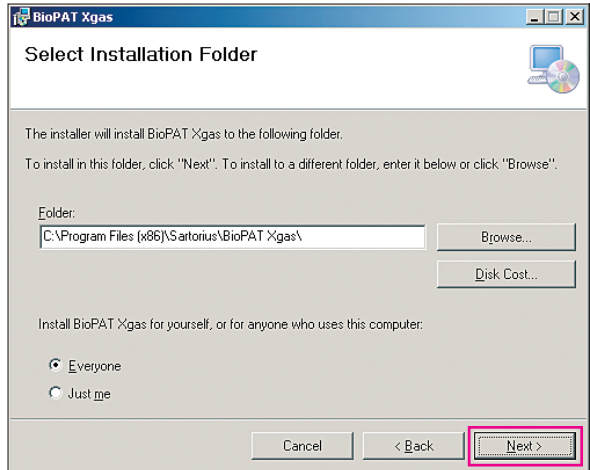
- ▶ Press the button “Accept” to continue.



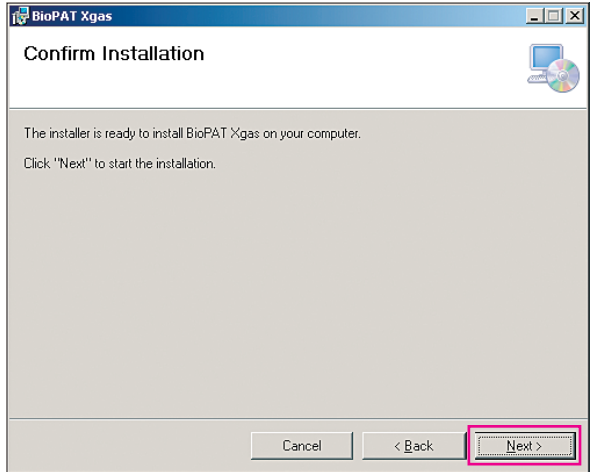
- ▶ The welcome screen is shown.
- ▶ Press button “Next”.



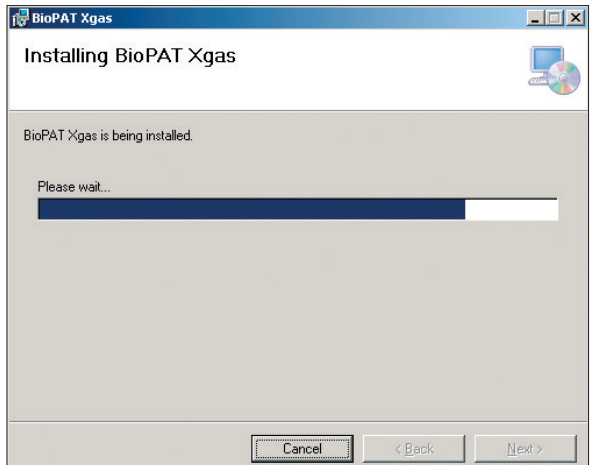
- ▶ Keep the default settings as shown in the picture.
- ▶ Press button “Next”.



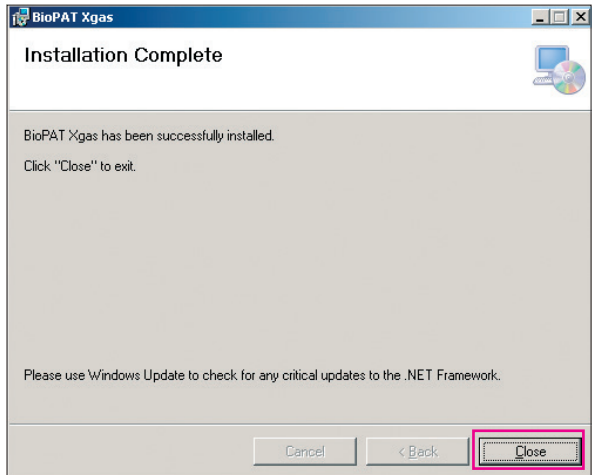
- ▶ Press button “Next” to install the software in the specified directory.



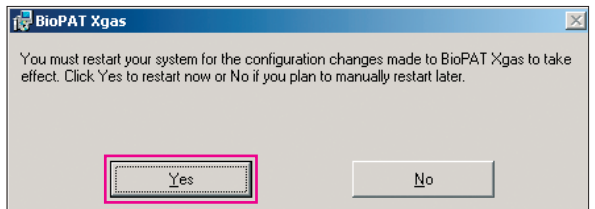
- ▶ Press button "Next" to confirm installation.



- ▶ The installation progress is shown on the screen.



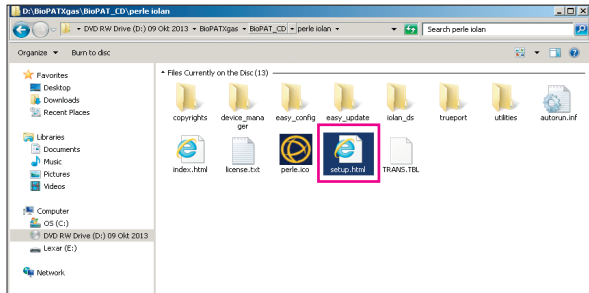
- ▷ The message "Installation Complete" appears.
- ▶ Press button "Close".



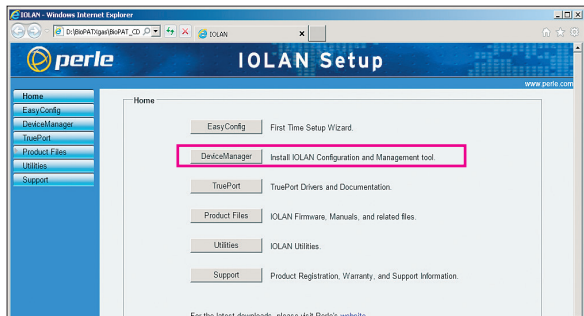
- ▷ Restart message appears.
- ▶ Press button "Yes".

5.3 Perle Device Driver

This chapter describes the installation and configuration of the Perle device driver.



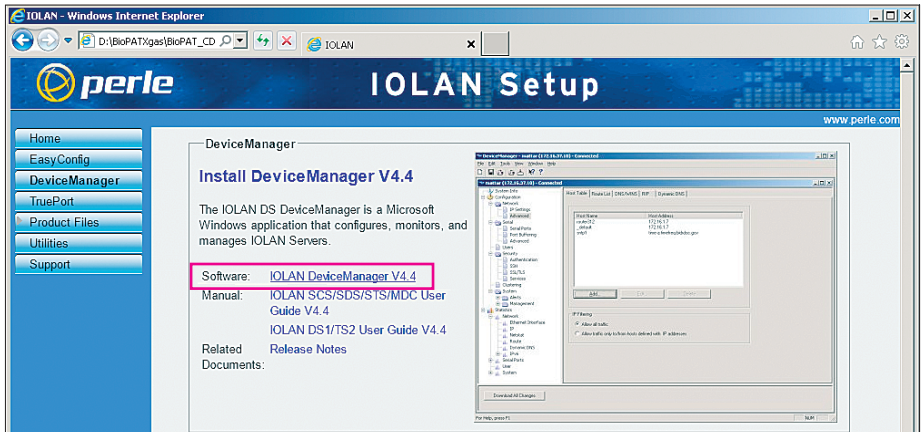
- ▶ Start the application “\perle iolan\setup.html” from the delivered CD.



- ▶ Press button “DeviceManager”.

NOTE

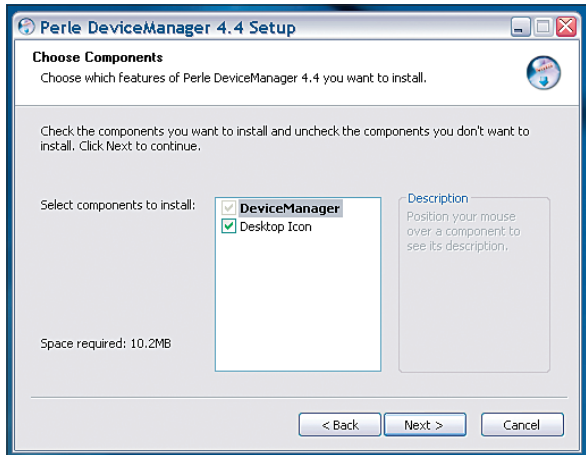
If there is a message to “allow blocked content” then enable blocked content.



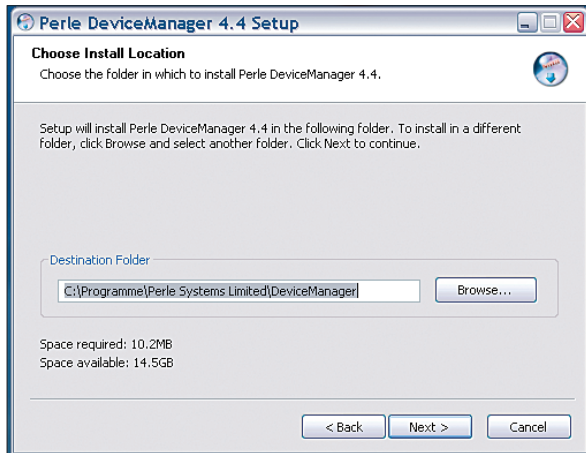
- ▶ Press link “IOLAN DeviceManager V4.4” to start Device Manager installation.



- ▶ Press “Next” button to continue with the installation.



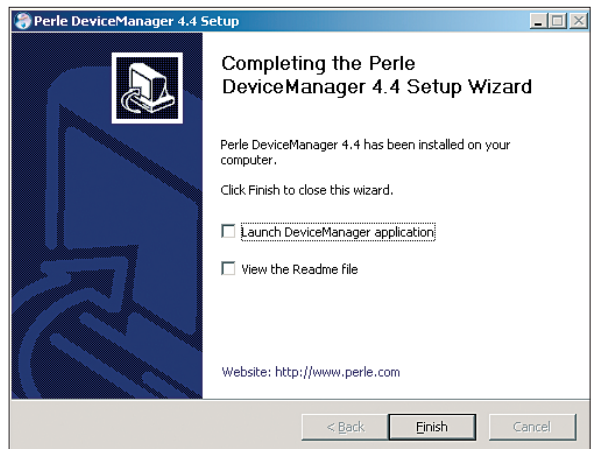
► Use default settings and press “Next” button.



► Use default settings and press “Next” button.



- Use default settings and press “Install” button.



- Uncheck both topics and press “Finish” button.

5.4 TruePort Application

The following pages describe how to install and configure the TruePort application.

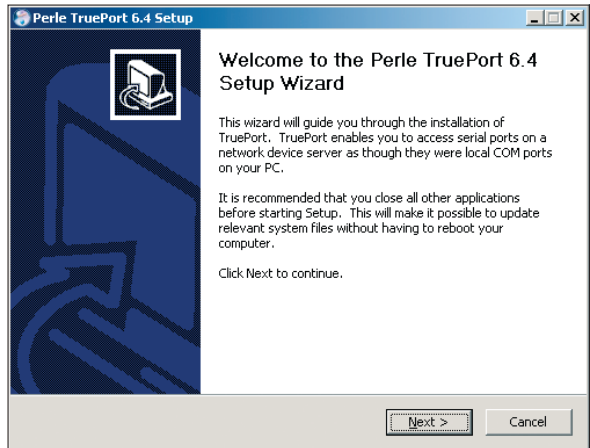
Operating System	Firmware	Documentation
Windows Server 2008 R2	6.4 (x64/AMD64)	Release Notes User Guide Sample SSL Key/Certificate (RSA)
Windows Server 2008	6.4 (x86) 6.4 (x64/AMD64)	Release Notes User Guide Sample SSL Key/Certificate (RSA)
Windows Server 2003	6.4 (x86) 6.4 (x64/AMD64) 6.4 (IA64)	Release Notes User Guide Sample SSL Key/Certificate (RSA)
Windows 7	6.4 (x86) 6.4 (x64/AMD64)	Release Notes User Guide Sample SSL Key/Certificate (RSA)
Windows Vista	6.4 (x86) 6.4 (x64/AMD64)	Release Notes User Guide Sample SSL Key/Certificate (RSA)
Windows XP	6.4 (x86) 6.4 (x64/AMD64)	Release Notes User Guide Sample SSL Key/Certificate (RSA)
Windows 2000	6.4 (x86)	Release Notes User Guide

► Click on the button “TruePort” on the left side of the IOLAB Setup.

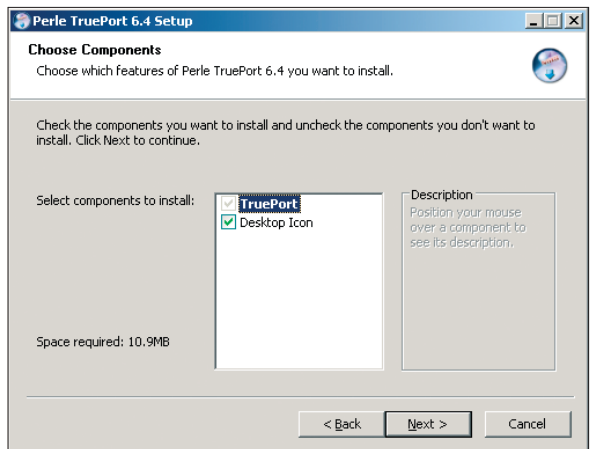
NOTE

Select the operating system which is installed at your pc.

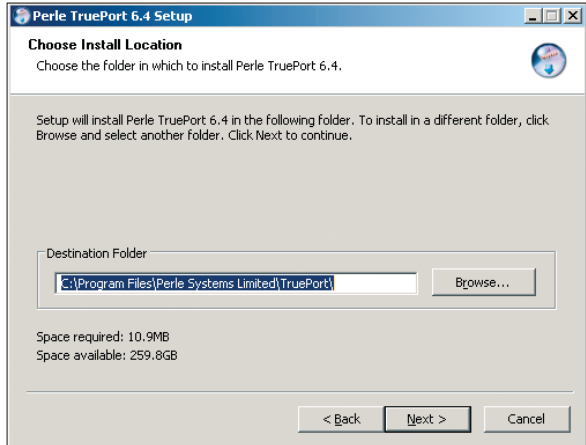
- For the 32 bit version of the operating system click to “(x86)”.
- For the 64 bit version of the operating system click to “(x64)”.



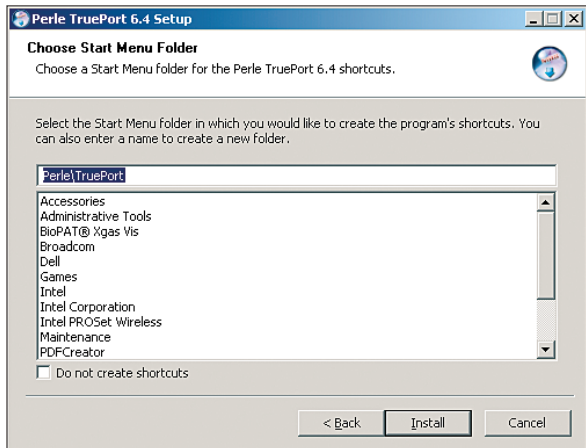
- ▶ Press "Next" button to continue.



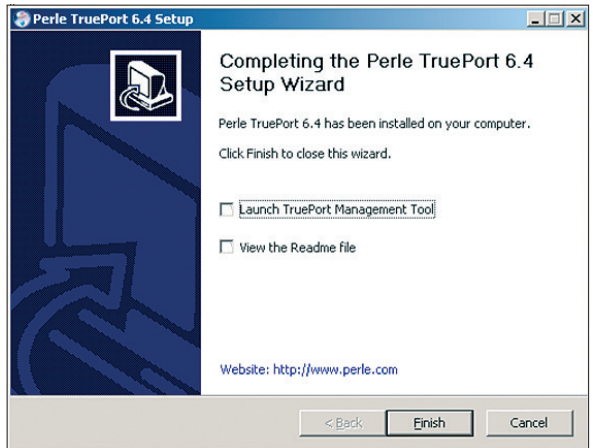
- ▶ Keep default settings and press "Next" button to continue.



- Keep default settings and press “Next” button to continue.



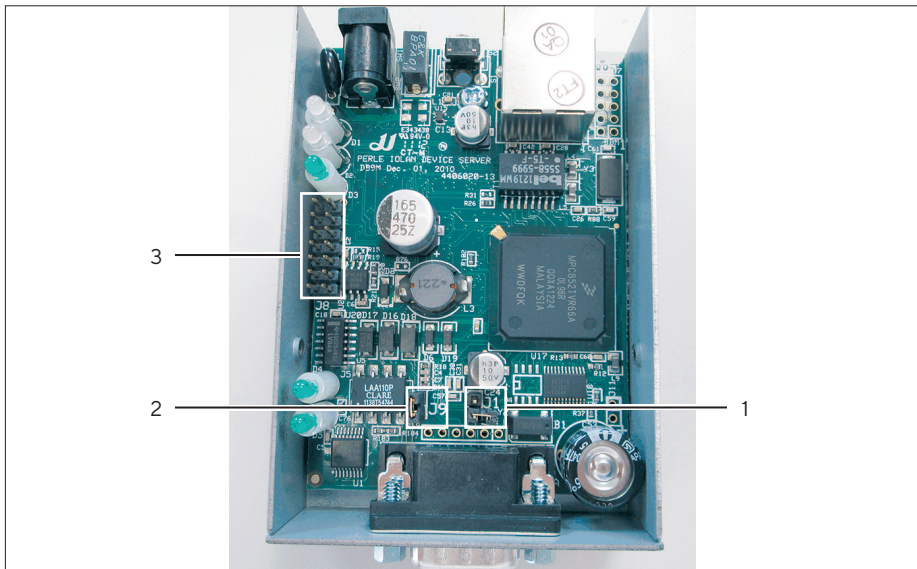
- Keep default settings and press button “Install” to continue.



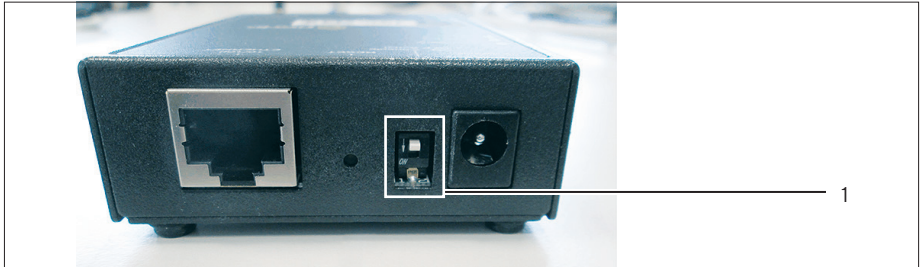
- Uncheck both topics and press “Finish” button.

6 Hardware Settings

6.1 Checking Jumper Settings at the Perle Device

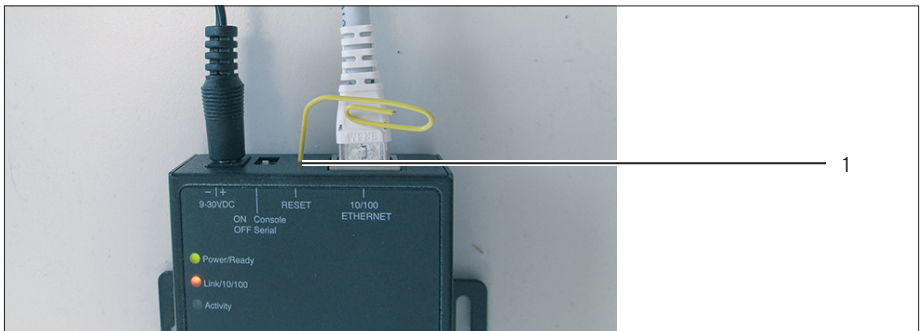


- ▶ Loosen the two screws on the side of the housing to open the Perle device.
- ▶ Set the jumpers as specified below:
 - Set jumper J1 (1) open.
 - Set jumper J8 (3) all open.
 - Set jumper J9 (2) closed.
- ▶ Close the Perle device and fasten the two screws.



- ▶ Set the switch (1) to “OFF” (upper position).
- ▷ The perle device is set to serial mode.

6.2 Resetting Perle Device



- ▶ Press the “RESET” button with a paper clip to reset the Perle device.
- ▷ After 5 s the lamp “Power/Ready” changes color from green to orange.

NOTE

After “Reset” please wait at least 2 min to let the Perle device finish the reset.

7 Configuration

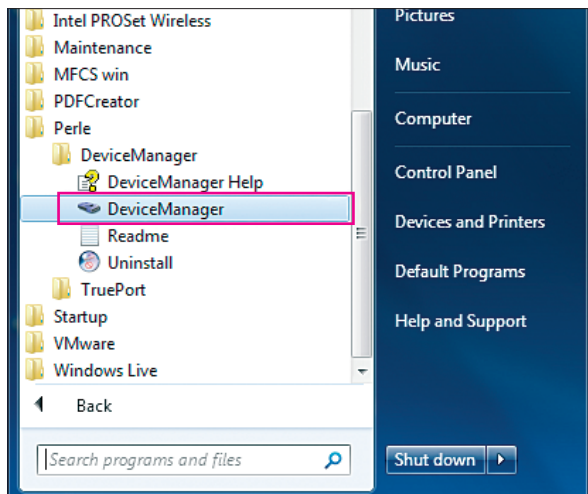
7.1 Perle Device

NOTE

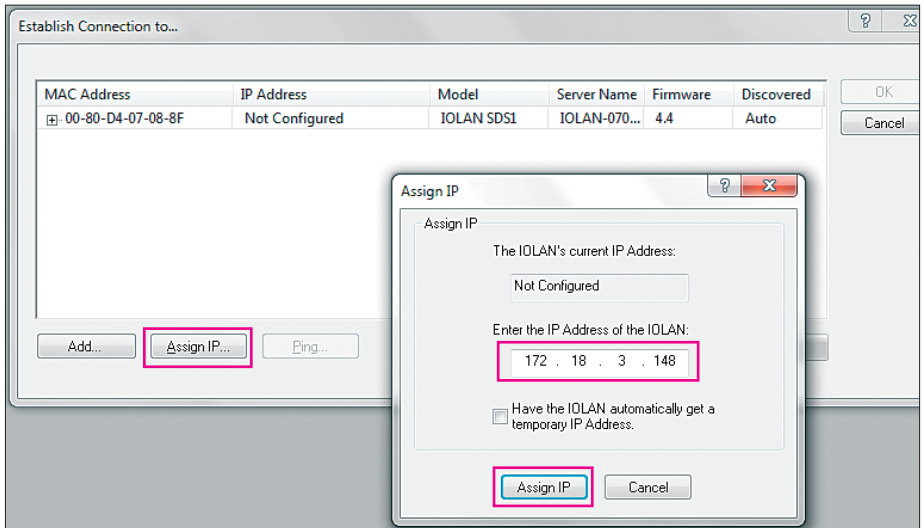
Double check, whether the Windows firewall is disabled or the Perle device manager is enabled in the firewall. In Windows firewall open all ports especially for the “devicemanager.exe” (UDP) and for the “tftp_svr.exe” (TCP) or temporarily disable the firewall.

NOTE

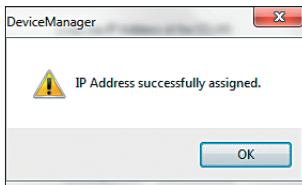
For server OS deactivate or stop “DHCPClient” service before starting the configuration. Otherwise the device will receive an IP address automatically.



► Start Perle “Device Manager”.

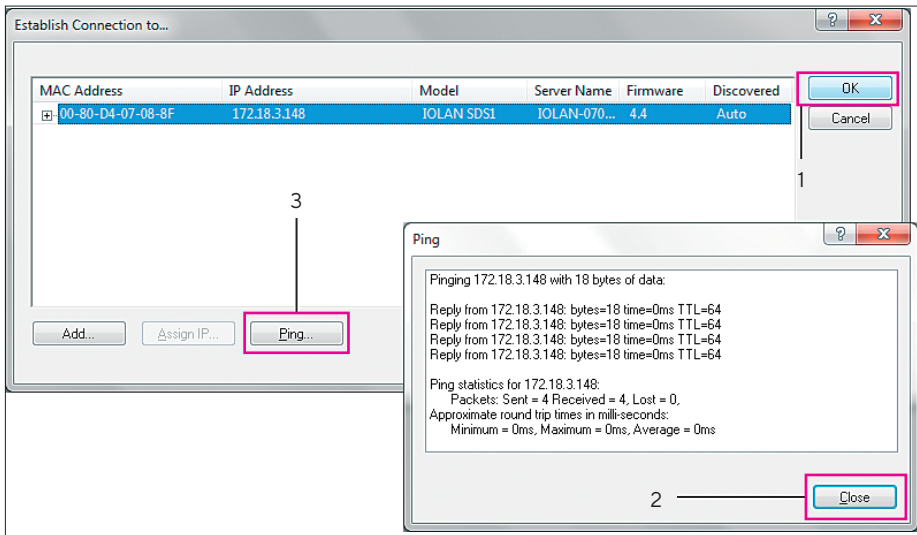


- ▶ Press button "Assign IP".
- ▶ Enter your IP address for the Perle device.
- ▶ Press button "Assign IP".
- ▶ The following message is shown.

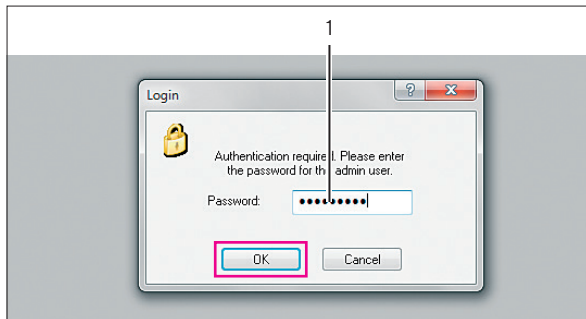


- ▶ Press "OK" button.

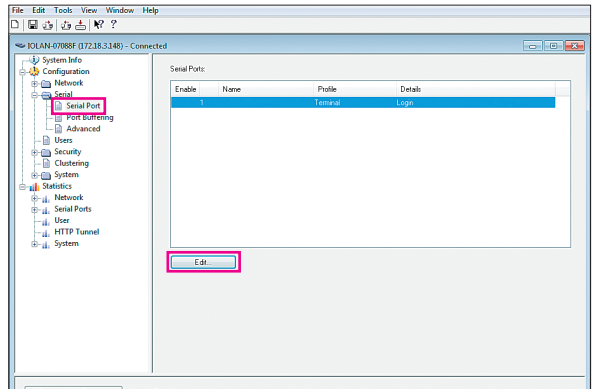
Test connection



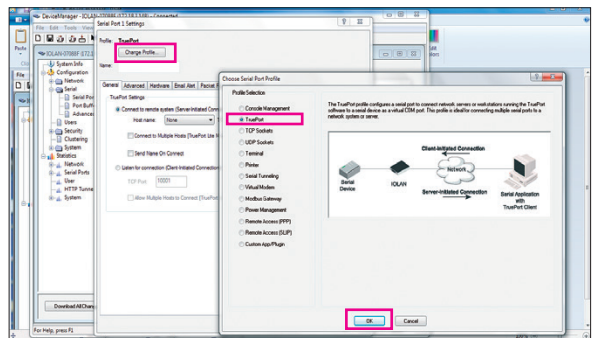
- ▶ Press “Ping” (3) button to test connection.
- ▶ Close Ping window (2).
- ▶ Press “OK” button (1) to connect to Perle device.



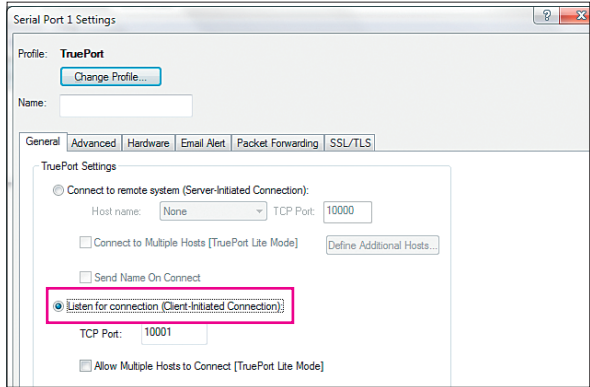
- ▶ Enter the default password “superuser” (1).
- ▶ Press the “OK” button.



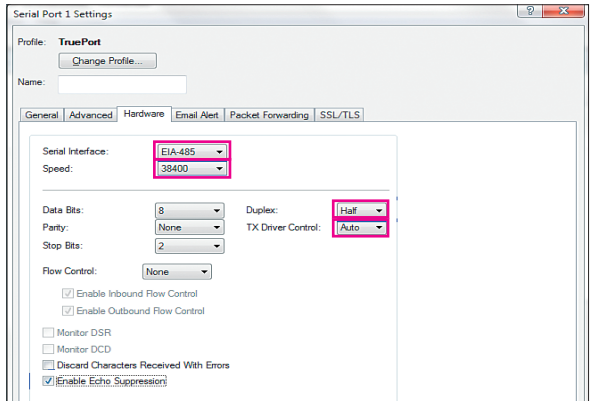
- ▶ Open menu “Configuration / Serial / Serial Port”.
- ▶ Press “Edit” button.



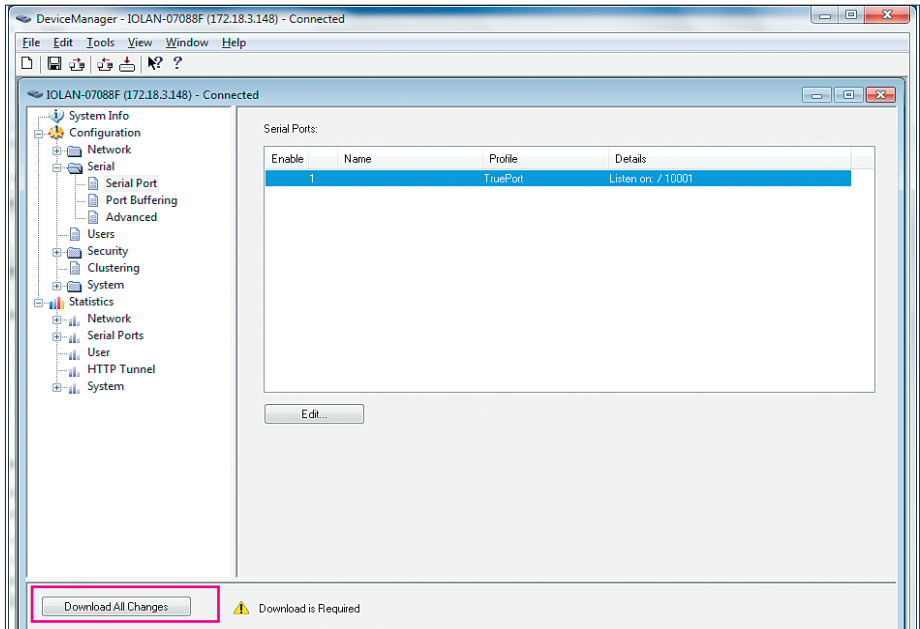
- ▶ Press button “Change Profile”.
- ▶ Select option “TruePort” and press “OK” button.



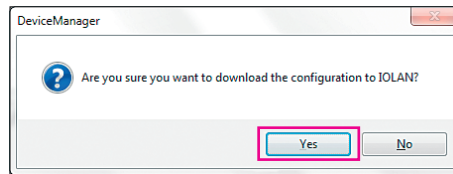
- ▶ Select option “Listen for connection” under tab “General”.
- ▶ Press “OK” button.



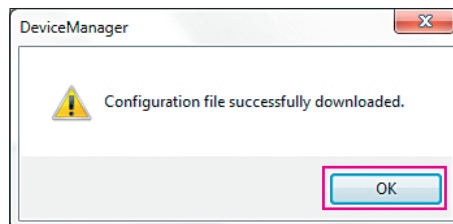
- ▶ Change to tab “Hardware”.
- ▶ Make the selections as shown in the picture.
- ▶ Press “OK” button.



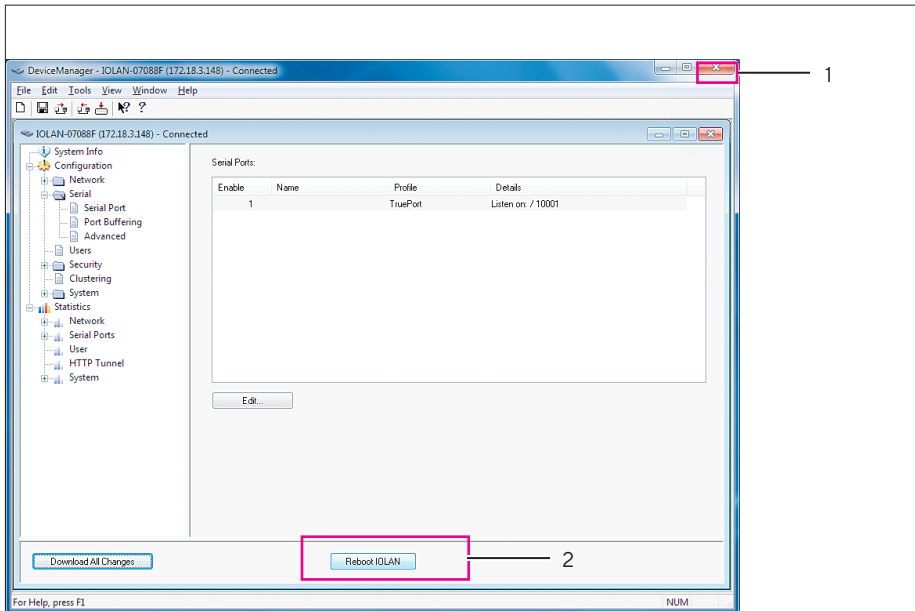
- ▶ Press button "Download All Changes".



- ▶ Press button "Yes" to confirm.



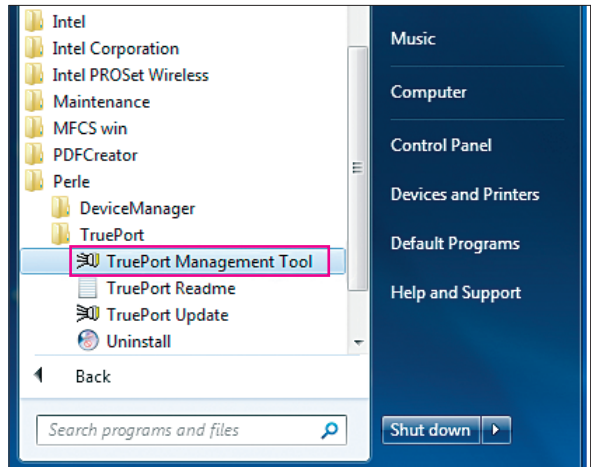
- ▶ Press "OK" button.



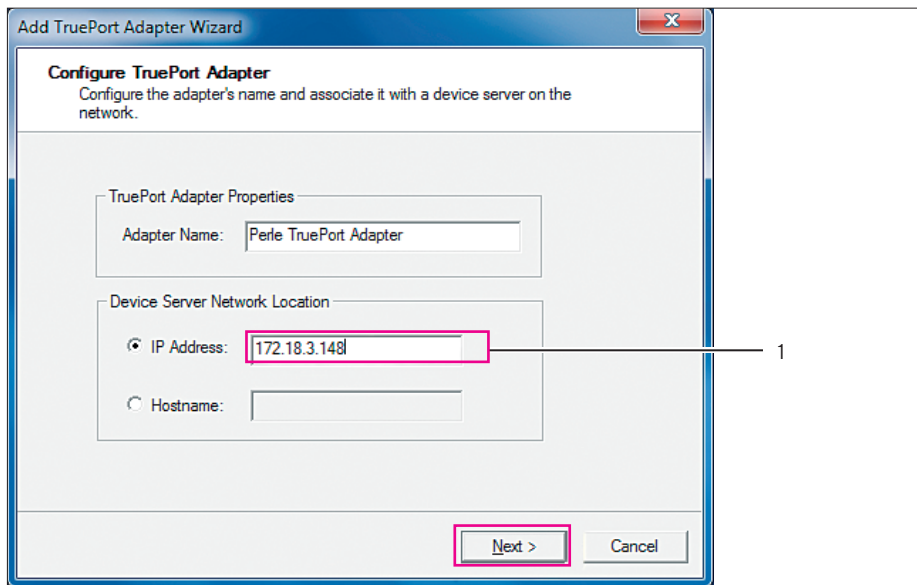
- ▶ Press button “Reboot IOLAN” (2).
- ▶ Close the DeviceManager (1).

7.2 Serial Port

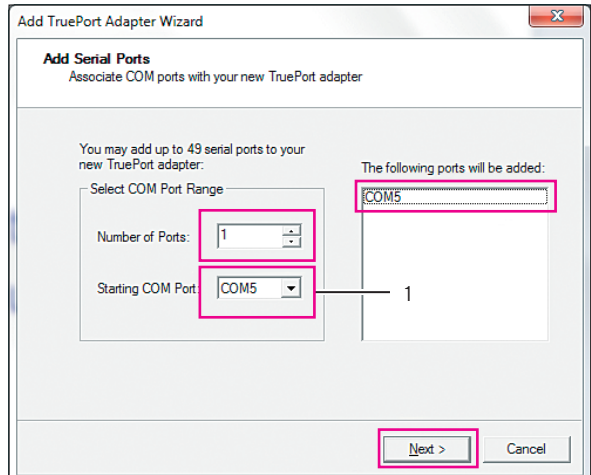
7.2.1 Configuring Serial Port with TruePort Management Tool



- ▶ Start the “TruePort ManagementTool” to configure the serial port.



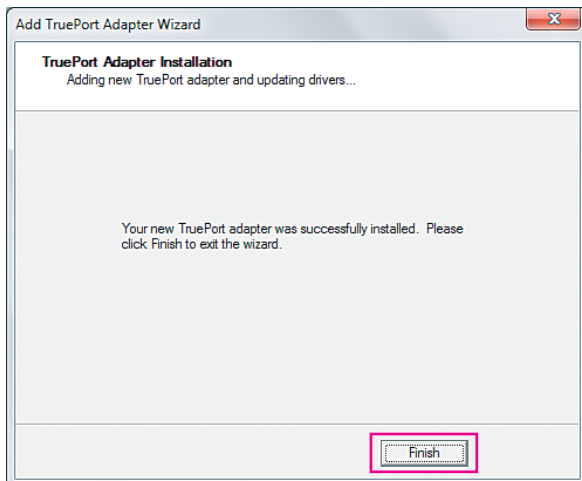
- ▶ Enter the IP address of the Perle device to the field "IP Address" (1)
- ▶ Press "Next" button.



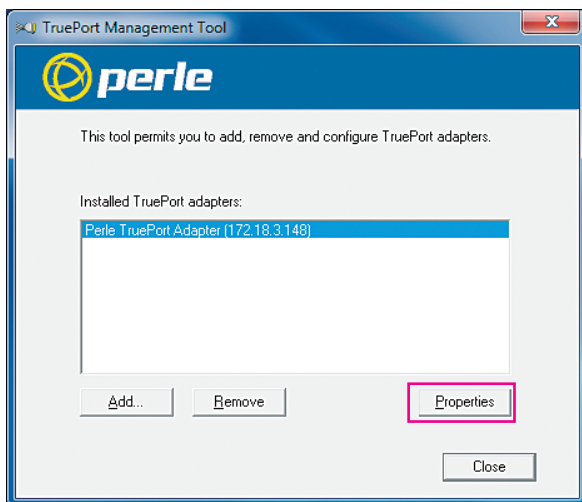
- ▶ Select an available port from the drop down list “Starting COM Port” (1).
- ▶ Press “Next” button.

NOTE

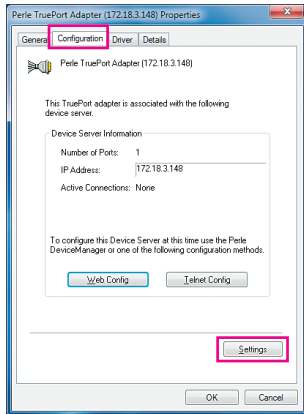
It is recommended to start with COM5, because COM5 will not be used (and therefore available) on most of the systems. Continue with COM6, COM7 if more Perle devices will be connected.



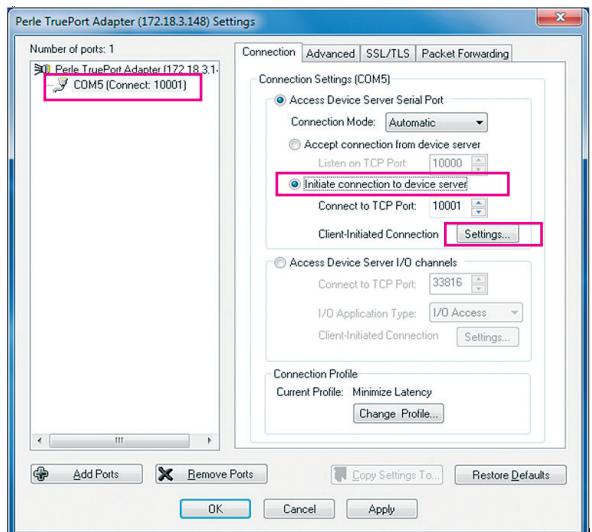
► Press button “Finish”.



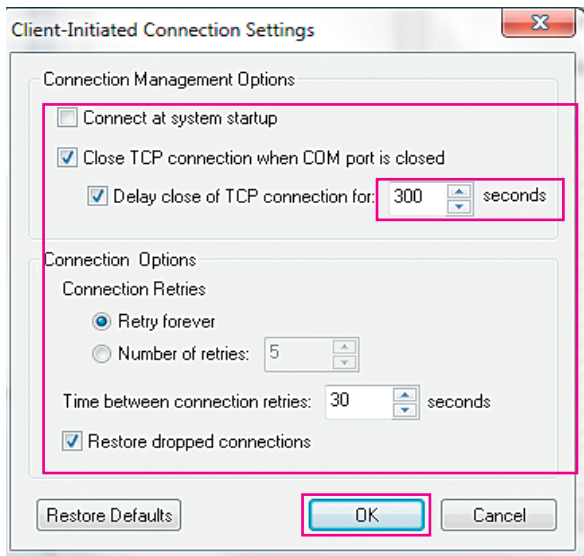
► Select the recently configured device and press button “Properties”.



- ▶ Change to register card “Configuration”.
- ▶ Press button “Settings”.

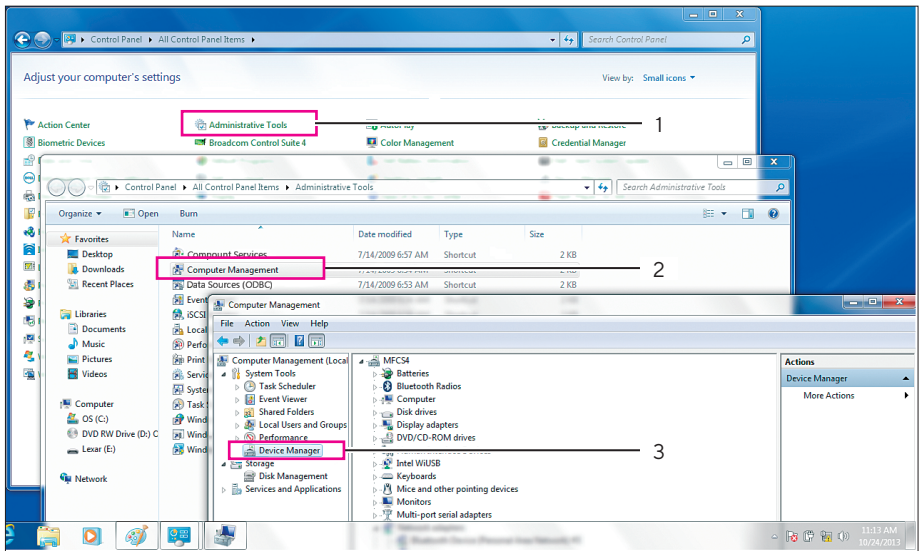


- ▶ Select the COM port.
- ▶ Activate “Initiate connection to device server”.
- ▶ Press button “Settings”.

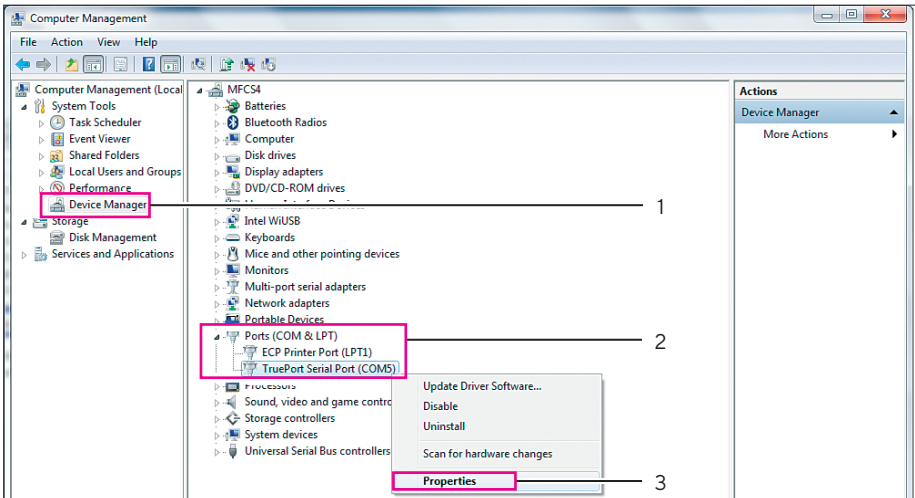


- ▶ Change setting “Delay close of TCP connection for” to 300 seconds.
- ▶ Press button “OK” and close all the remaining windows.

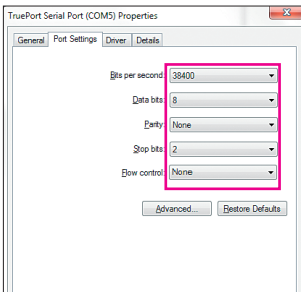
7.2.2 Configuring Serial Port with Device Manager



- ▶ Open the Windows "Control panel".
- ▶ Select "Administrative Tools" (1).
- ▶ Select "Computer Management" (2).
- ▶ Select Device Manager (3).



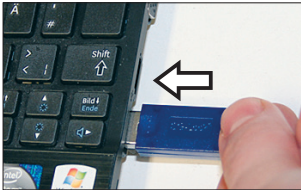
- ▶ Select Ports (COM & LPT) (2).
- ▶ Right mouse click to "Trueport Serial Port (COM5)" (2).
- ▶ Select "Properties" (3).



- ▶ Change to register card "Port Settings".
- ▶ Perform the settings shown in the picture above.
- ▶ Press "OK" button and close the Windows device manager.

7.3 BioPAT® Xgas Vis OPC License

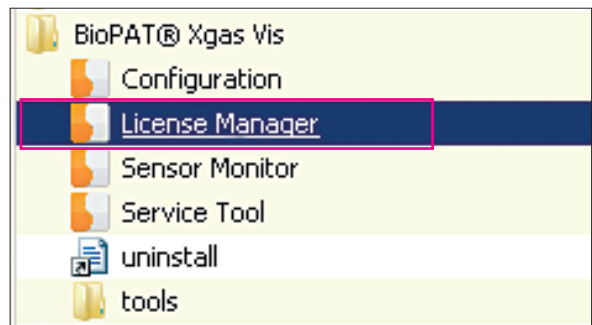
This chapter describes how to activate the BioPAT® Xgas Vis OPC License.



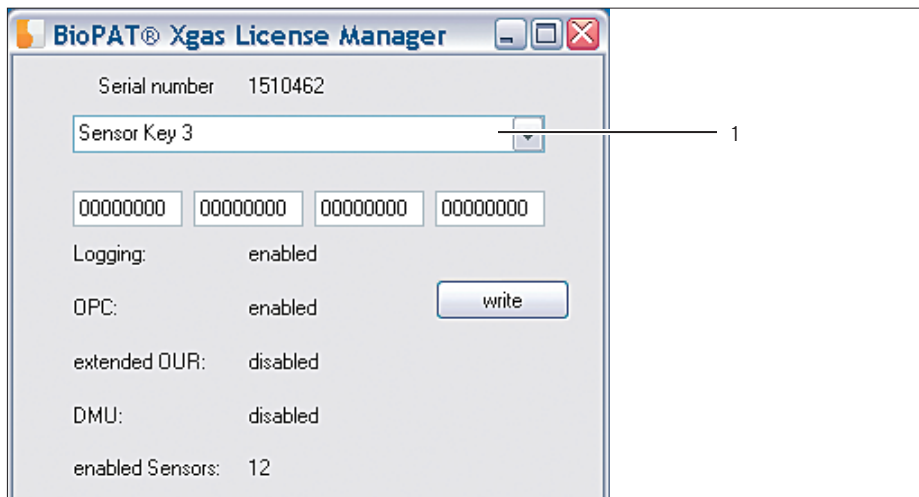
- ▶ Plug the blue license dongle (SG-Lock) into a free USB port of the MFCS/win computer.

NOTE

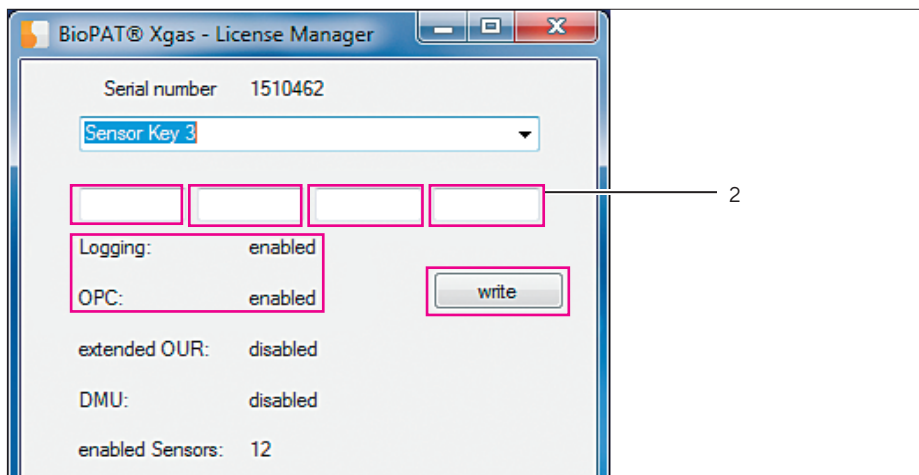
The dongle must stay in the USB port all the time if you want to use all features. If the application without the dongle is used, then only the measurement of one analyzer in the BioPAT® Xgas Vis software (without logging) can be used in the Service Tool for the calibration and the changing the MODBUS ID.



- ▶ Start the “License Manager” to activate the BioPAT® Xgas VIS OPC license.



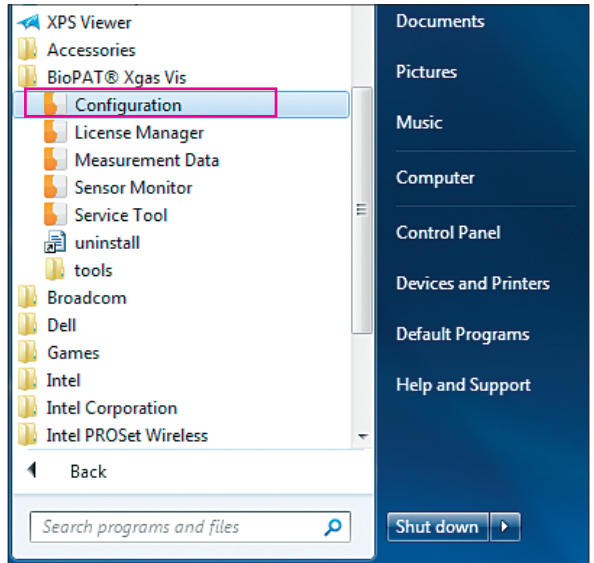
- ▶ Select “Sensor Key 3” (1) from drop down menu.



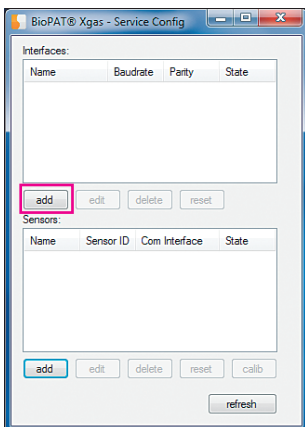
- ▶ Enter the corresponding key in the four input fields (2).
- ▶ Press the button “write”.
- ▷ The unlocked features and the number of enabled sensors is shown.
- ▶ Close the “License Manager”.

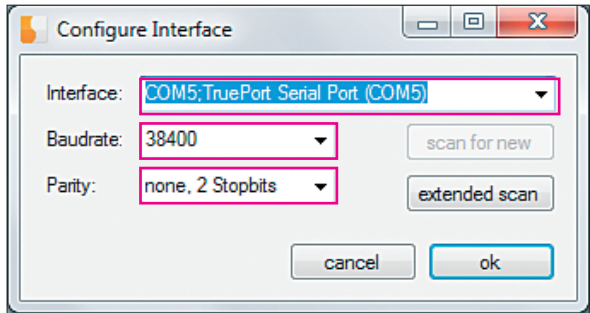
7.4 BioPAT® Xgas Sensor

7.4.1 Configuring One Sensor

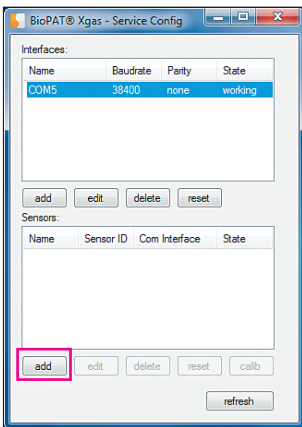


- ▶ Start the “Configuration” tool for BioPAT® Xgas to configure one XGAS Sensor.
- ▶ Press the “add” button for interfaces to add a new “Interface”.

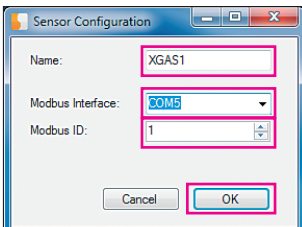




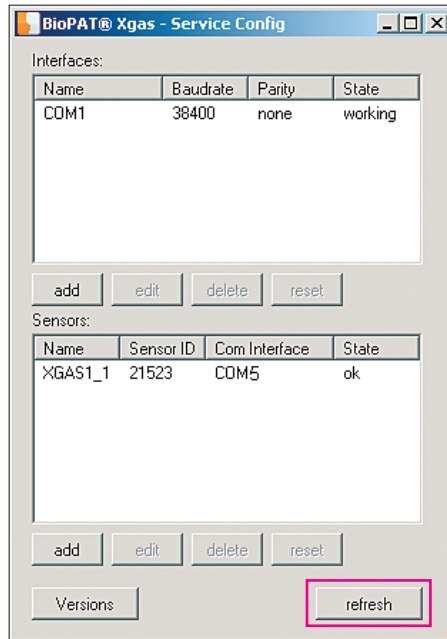
- ▶ Select
 - the correct Interface.
 - Baudrate "38400".
 - Parity "none, 2 Stopbits".
- ▶ Press "OK" button.



- ▶ Press the "add" button for sensors to add a new sensor.



- ▶ Enter a name, the Modbus interface and the Modbus ID 1.
- ▶ Press the "OK" button.



- ▷ After the sensor has been added, it is shown in the Sensor list.
- ▶ If state "ok" is not shown in the Sensor list, press the "refresh" button to refresh the list.

NOTE

For sensors naming conventions refer to examples on the following pages.

7.4.1.1 Naming conventions for XGAS Sensors

Example 1:

Only one BioPAT® Xgas Sensor with a Biostat® fermentor

	Name	Modbus Interface	Modbus ID	Description
1. Sensor	XGAS1_1	COM5	1	Single

Example 2:

Two BioPAT® Xgas Sensors with a twin fermentor system

	Name	Modbus Interface	Modbus ID	Description
1. Sensor	XGAS1_1	COM5	1	Twin A
2. Sensor	XGAS1_2	COM5	2	Twin B

Example 3:

Four BioPAT® Xgas Sensors with a B-DCU II 6-fold fermentor system

	Name	Modbus Interface	Modbus ID	Description
1. Sensor	XGAS1_1	COM5	1	B6 A
2. Sensor	XGAS1_2	COM5	2	B6 B
3. Sensor	XGAS1_3	COM5	3	B6 C
4. Sensor	XGAS1_4	COM5	4	B6 D
5. Sensor	XGAS1_5	COM5	5	B6 E
6. Sensor	XGAS1_6	COM5	6	B6 F

Example 4:

Nine BioPAT® Xgas Sensors with:

- One Biostat® B single fermentor system
- One Biostat® B twin fermentor system
- One Biostat® B-DCU II 6-fold fermentor system

	Name	Modbus Interface	Modbus ID	Description
1. Sensor	XGAS1_1	COM5	1	Single
2. Sensor	XGAS2_1	COM6	1	Twin A
3. Sensor	XGAS2_2	COM6	2	Twin B
4. Sensor	XGAS3_1	COM7	1	B6 A
5. Sensor	XGAS3_2	COM7	2	B6 B
6. Sensor	XGAS3_3	COM7	3	B6 C
7. Sensor	XGAS3_4	COM7	4	B6 D
8. Sensor	XGAS3_5	COM7	5	B6 D
9. Sensor	XGAS3_6	COM7	6	B6 D

7.4.2 Configuring Multiple Sensors with one Ethernet Modbus Converter

If more than one BioPAT® Xgas Sensor is to be connected to one Perle device (Ethernet Modbus converter), then additional configuration steps are needed, because the default delivery Modbus address of each BioPAT® Xgas Sensor is "1".

► Please perform the following configuration steps one after each other:

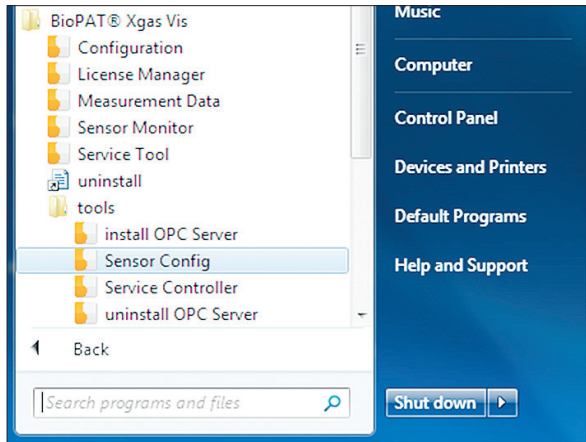
Configuration Step	Description
1	Connect Sensor #1. Verify that it already has Modbus address #1.
2	Label Sensor #1 with the following information: Modbus ID : 1.
3	Disconnect Sensor #1.
4	Connect Sensor #2. Change the Modbus address for #1 to #2. Please refer to chapter [→7.5, page 58].
5	Label Sensor #2 with the following information: Modbus ID : 2.
6	Disconnect Sensor #2.
7	Connect Sensor #3. Change the Modbus address for #1 to #3. [→7.1, page 34].
8	Label Sensor #3 with the following information: Modbus ID : 3.
9	Disconnect Sensor #3. Do the same with all the others, which will be connected to the same Perle device.

Configuration Step	Description
10	Connect Sensor #n. Change the Modbus address for #1 to#n [→7.5, page 58].
11	Label Sensor #n with the following information: Modbus ID : n.
12	Disconnect Sensor #n.

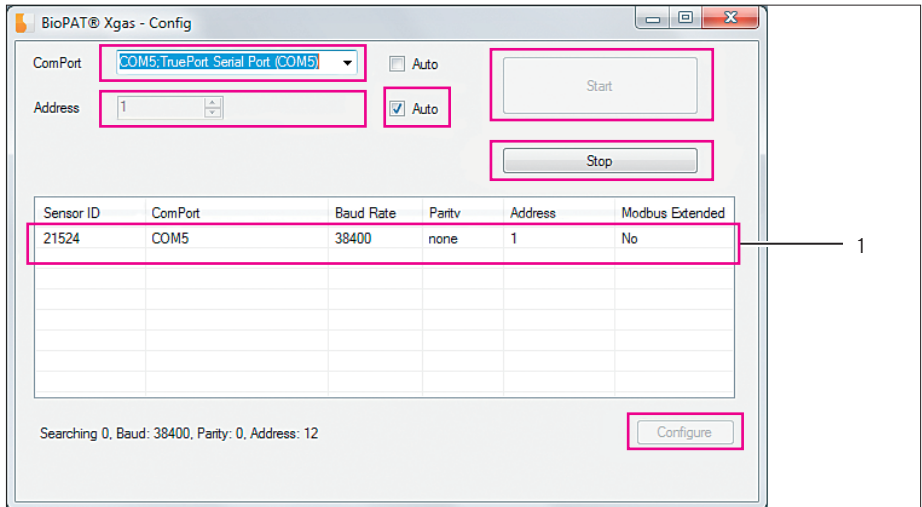
After the Modbus addresses of all BioPAT® Xgas Sensors has been set correctly, perform the following steps:

- ▶ Connect all BioPAT® Xgas Sensors with the delivered Modbus cables.
- ▶ Connect the Modbus termination to the last sensor as shown in figure [→see chapter 4.2, page 12].

7.5 Changing the Modbus Address



- ▶ Start the tool “BioPAT® Xgas Vis / Tools / Sensor Config” to change the Modbus address.

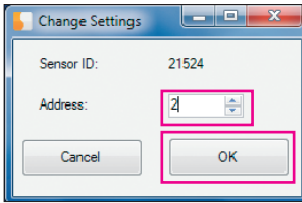


- ▶ Select the ComPort, where the sensor is connected.
- ▶ Check the field (Address Auto).
- ▶ Press “Start” button.
- ▶ Wait until the sensor appears (1).
- ▶ Press “Stop” button.
- ▶ Select the sensor from the list.
- ▶ Press “Configure” button.

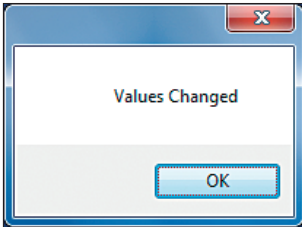
NOTE

A few seconds after pressing the “Start” button, the message “Searching” must appear in the upper left corner of the “Sensor Config” window.

- If this message is not shown, press “Start” button again.



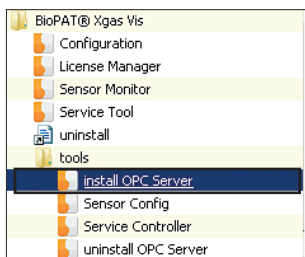
- ▶ Enter the new address for the sensor.
- ▶ Press "OK" button.



- ▶ The success message is shown.
- ▶ Press "OK" button.

8 Installation of BioPAT® Xgas OPC Server

Start the program “install OPC server” to install the BioPAT® Xgas OPC Server.



- ▶ Open “BioPAT® Xgas Vis / tools menu / install OPC server”.



- ▶ Press button “OK” to finish the installation.

8.1 Installing MFCS/win OPC Client

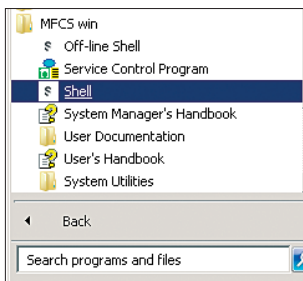
If the MFCS/win OPC DA Client is already installed, then continue with next chapter.

- ▶ If the MFCS/win OPC DA Client is not installed, install the MFCS/win OPC DA Client according the delivered documentation “MFCS/win OPC DA Client Handbook (MFCSOPCC)”.

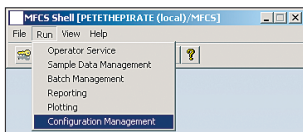
8.2 Configuring MFCS/win Configuration Management

Inside the MFCS/win SCADA software first a control unit of type “OPC” must be configured.

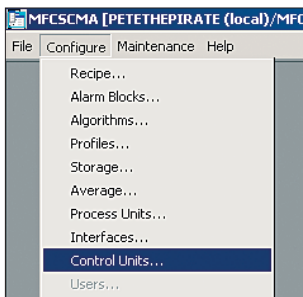
Then the variables for O₂ and CO₂ of the different fermentors must be configured.



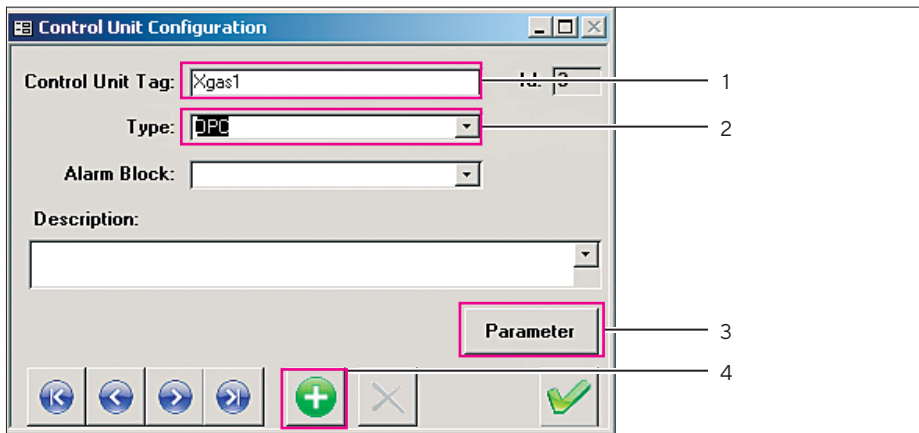
▶ Start MFCS/win Shell.



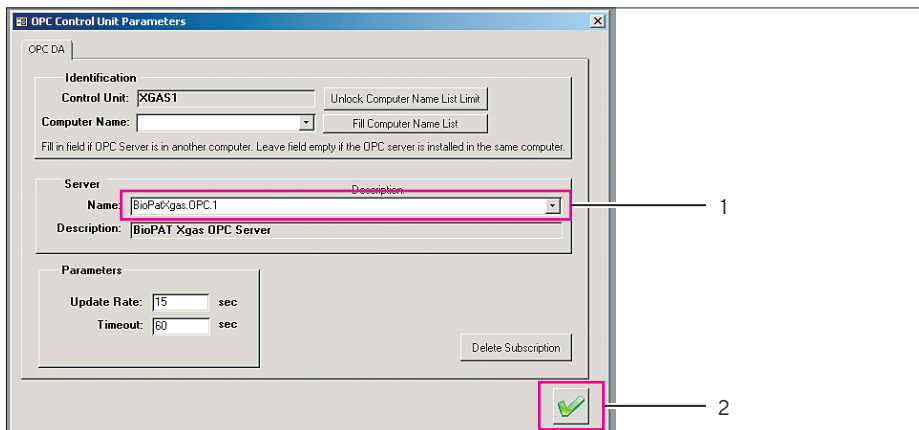
▶ Start “Configuration Management”.



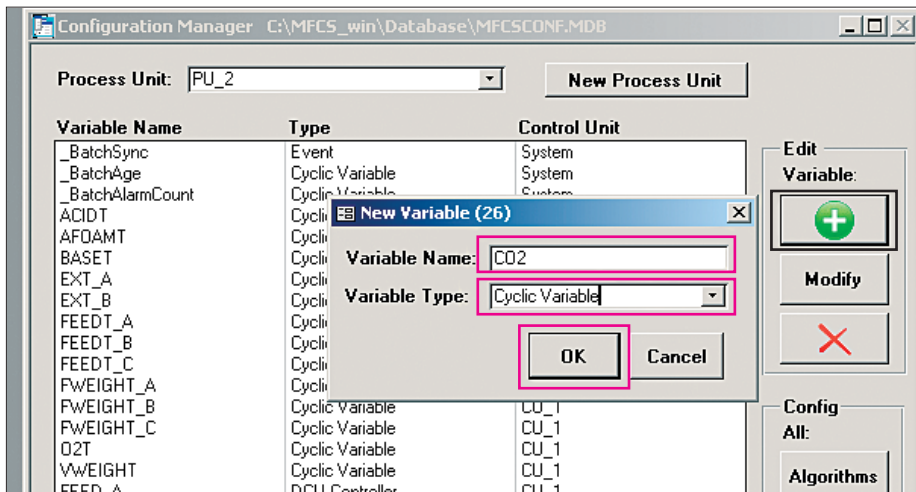
▶ Select “Configure / Control Units”.



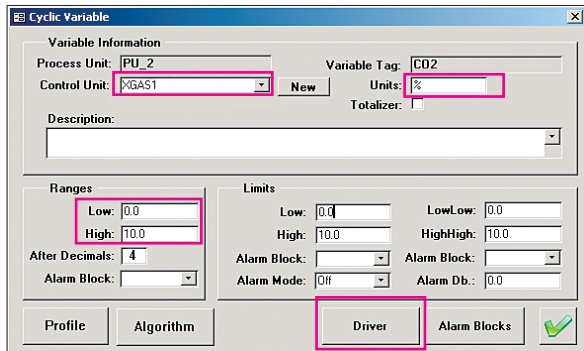
- ▶ Create a new Control Unit of type “OPC” and enter a name (e.g. Xgas1):
 - Press “New” button (4).
 - Select Control Unit name (1).
 - Select type „OPC” (2).
 - Press button “Parameter” (3).



- ▶ Select the “BioPatXgas.OPC.1” server from the drop down list (1) and confirm (2).



- ▶ Create a new cyclic variable "CO₂".
- ▶ Create a new cyclic variable "O₂".



- ▶ Select control unit "XGAS1".
- ▶ For CO₂ configure the range 0-10 %
- ▶ For O₂ configure the range 1-50 %.
- ▶ Press "Driver" button.
- ▶ The "OPC Item Configuration" opens (see following picture).

OPC Item Configuration

Process Unit: PU_2

Variable: CO2

Control Unit: XGAS1

Process Value Tag

Value: XGAS1_1.Values.CO2

Range: 0 (Low) 10 (High)

Controller Tags

Setpoint: Range: (Low) (High)

Output: Range: (Low) (High)

- ▶ If the sensor name is “XGAS1_1”, enter the OPC Tag “XGAS1_1.Values.CO2”.
- ▶ Confirm with “OK” button.

OPC Item Configuration

Process Unit: PU_2

Variable: O2

Control Unit: XGAS1

Process Value Tag

Value: XGAS1_1.Values.O2

Range: 1 (Low) 50 (High)

Controller Tags

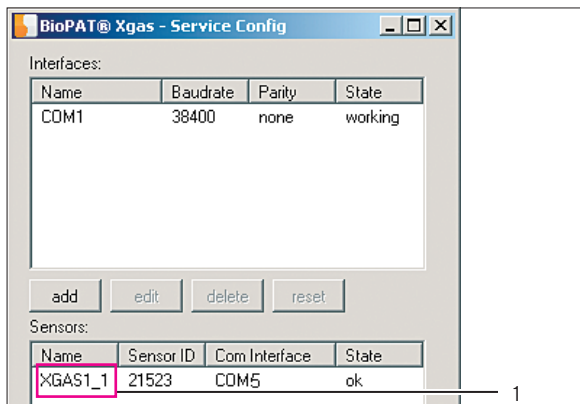
Setpoint: Range: (Low) (High)

Output: Range: (Low) (High)

Mode: Range: (Low) (High)

- ▶ If the sensor name is “XGAS1_1”, enter the OPC Tag “XGAS1_1.Values.O2”.
- ▶ Confirm with “OK” button.

Sensor name



NOTE

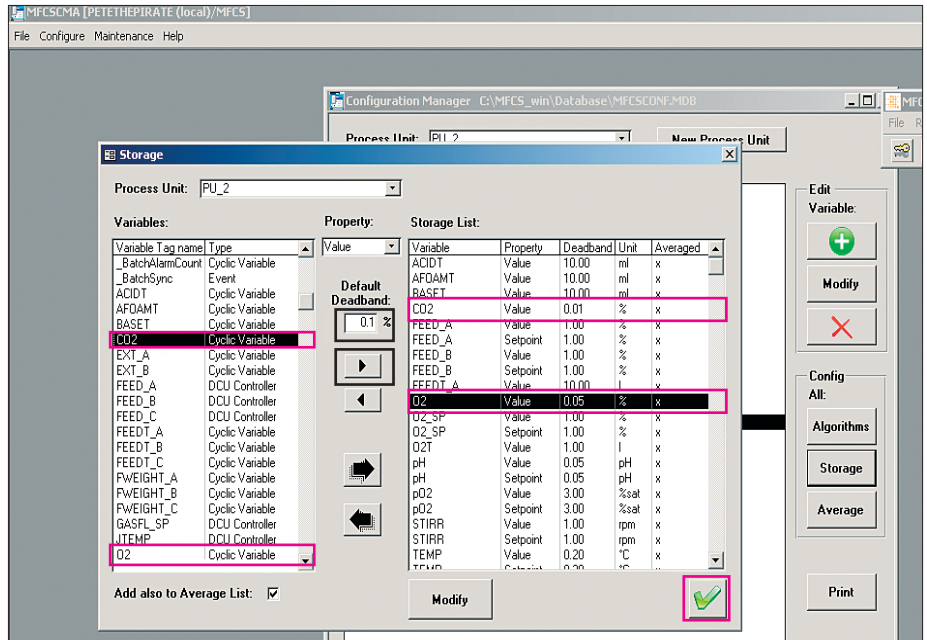
The sensor name (1) can be looked up in the Sensor configuration as described in chapter [→7.4, page 51].

NOTE

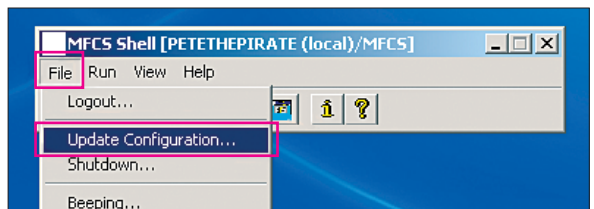
Instead of directly filling in the OPC Tags, an OPC Test Client can be used to browse and then the OPC tags can be filled in with copy & paste.
 – Refer to chapter [→8.3, page 69].

Adding new variables to Storage list

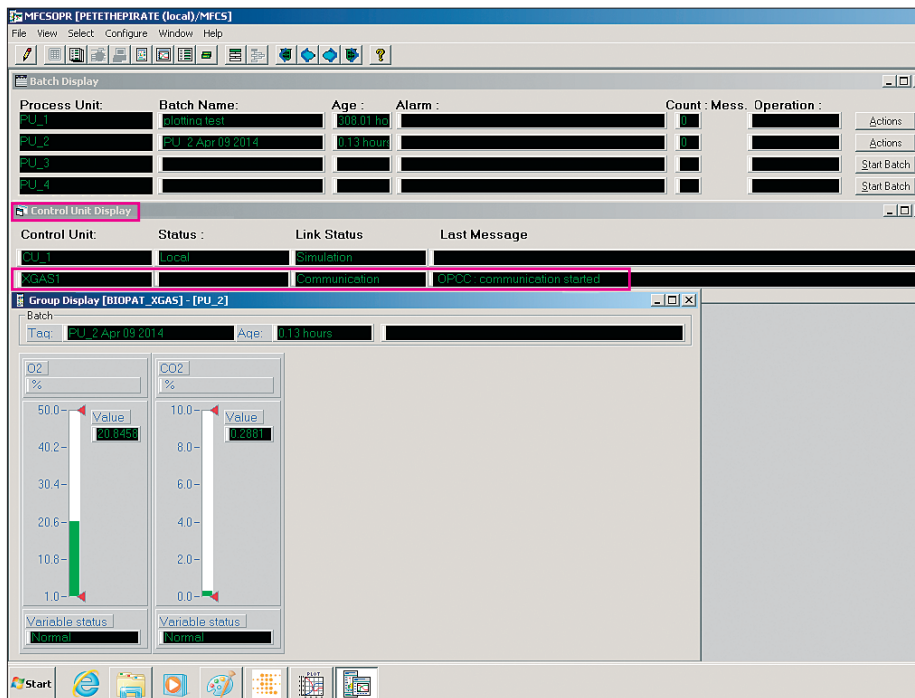
After the interface configuration is finished, the new variables must be added to the Storage list.



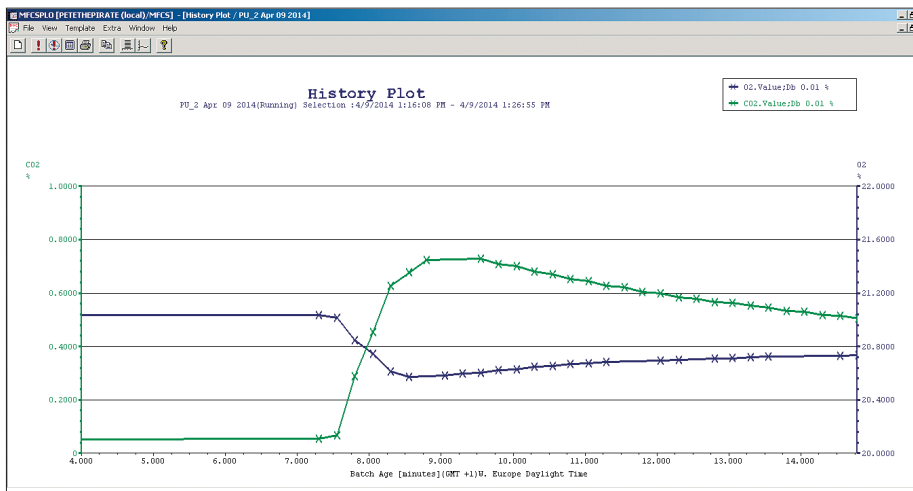
- ▶ Add the new variables O₂ and CO₂ to the Storage list as shown on the picture above.



- ▶ Perform "Update Configuration" from the "MFC5/win Shell to finalize the MFC5/win configuration".



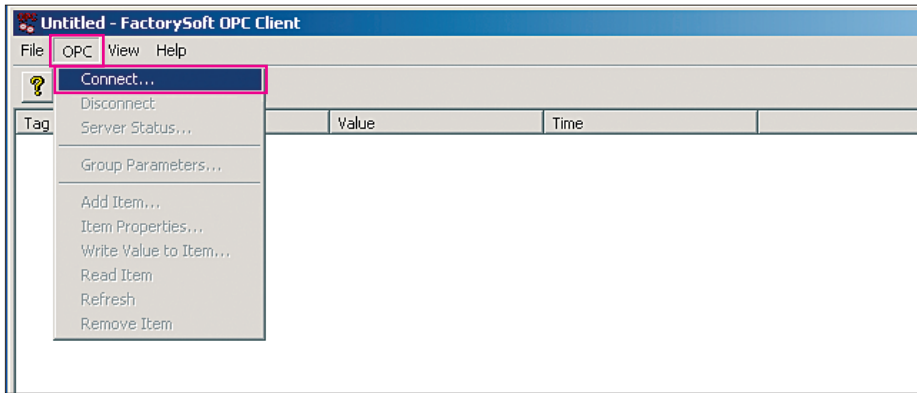
Verification of the communication between MFCS/win and BioPAT® Xgas from the MFCS/win Operator Service Program



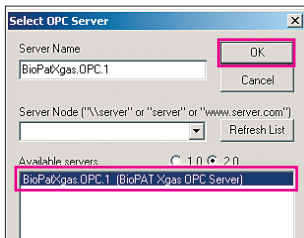
Verification of the communication between MFCS/win and BioPAT® Xgas with the MFCS/win Plotting

8.3 Browsing the BIOPATXGAS.OPC.1 Server with OPC Test Client

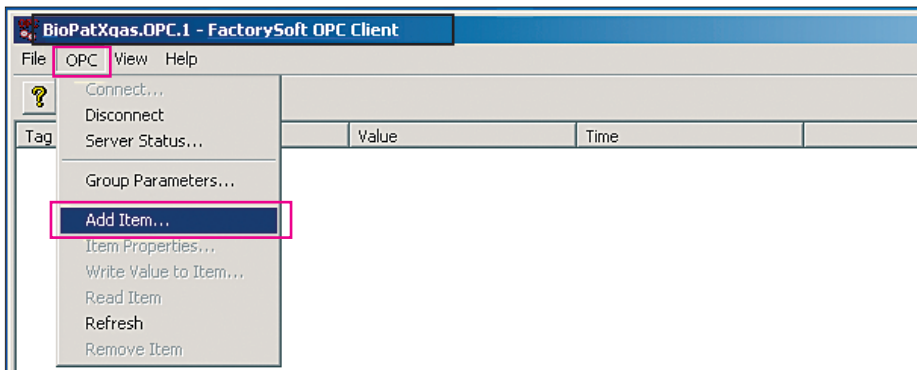
- ▶ Copy the tool "OPC_Client.exe" to the MFCS/win server computer into directory ":\MFCS_win\Tasks\OPC\".
- This tool is available at the BioPAT MFCS/win OPC DA Client CD in directory "<CD-ROM>:\SupportTools\".
- ▶ Start the file "OPC_Client.EXE" at the MFCS/win Server computer.
- ▷ Now the connection to the OPC server can be done, by doing the following steps.



► Connect to OPC Server with “OPC / Connect”.

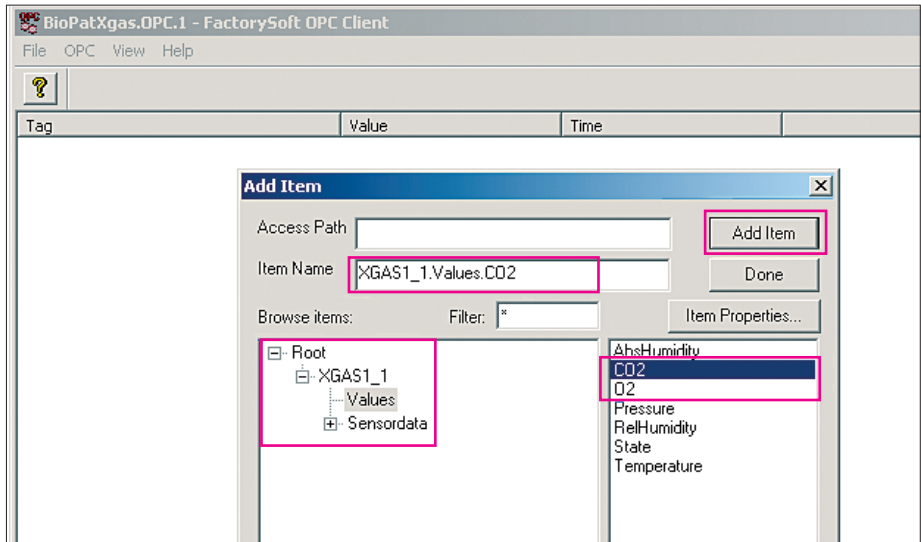


► Select the “BioPatXgas.OPC.1” server from the list.



After the Test Client is connected to the BioPATXgas.OPC.1 server, add items.

► Click “OPC / Add Item...”.



- ▶ Browse for O₂ and CO₂.
- ▶ Click "Add Item".

NOTE

If the CO₂ variable is selected in the list, then the OPC Tag will be shown in the field "Item Name".

This OPC Tag can be transferred with cut & paste to the MFCS/win configuration as shown in section [→page 66].

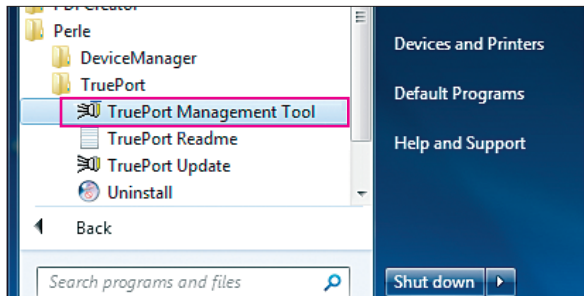
O₂ and CO₂ values in OPC Test Client:

Tag	Value	Time
XGAS1_1.Values.CO2	0.0730031	04/09/14 14:49:05
XGAS1_1.Values.O2	21.0131	04/09/14 14:49:05

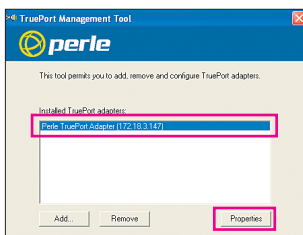
9 Firmware Version

9.1 Checking | Upgrading Firmware Version of the Perle Device

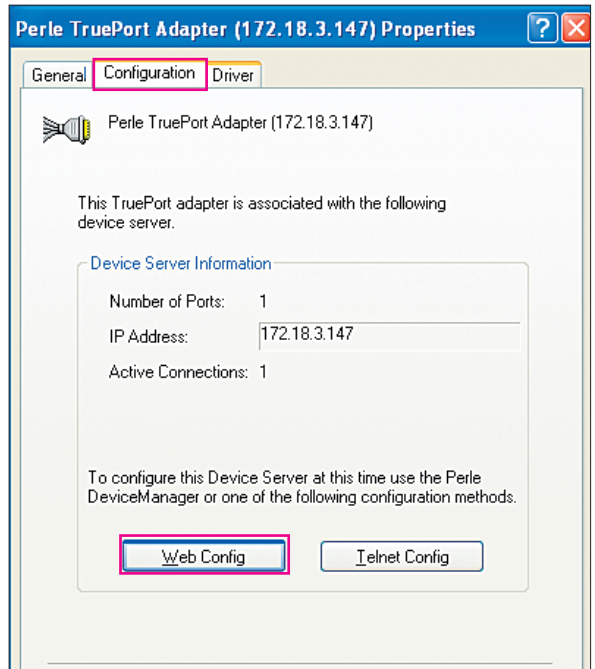
For trouble free operation it is necessary, that the firmware version 4.5 is available in the Perle device. How to check the installed firmware version of the Perle device is described in this chapter.



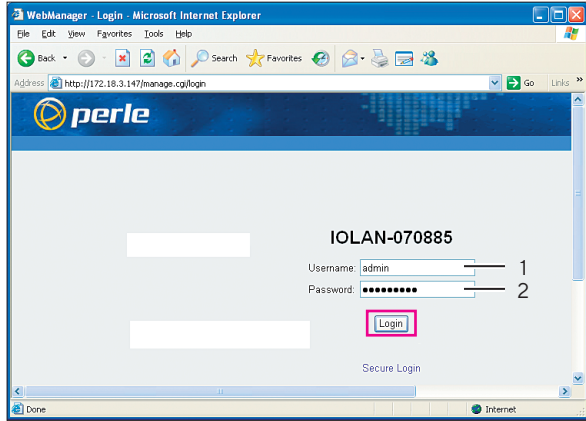
- ▶ Start the “TruePort ManagementTool” for verification.



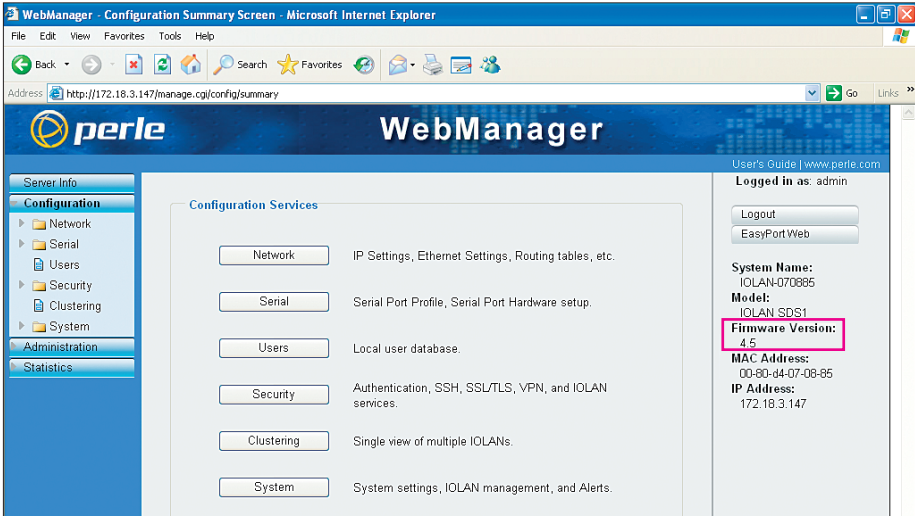
- ▶ Select the Perle TruePort Adapter in question.
- ▶ Press the “Properties” button.



- ▶ Change to register card "Configuration".
- ▶ Press button "Web Config".

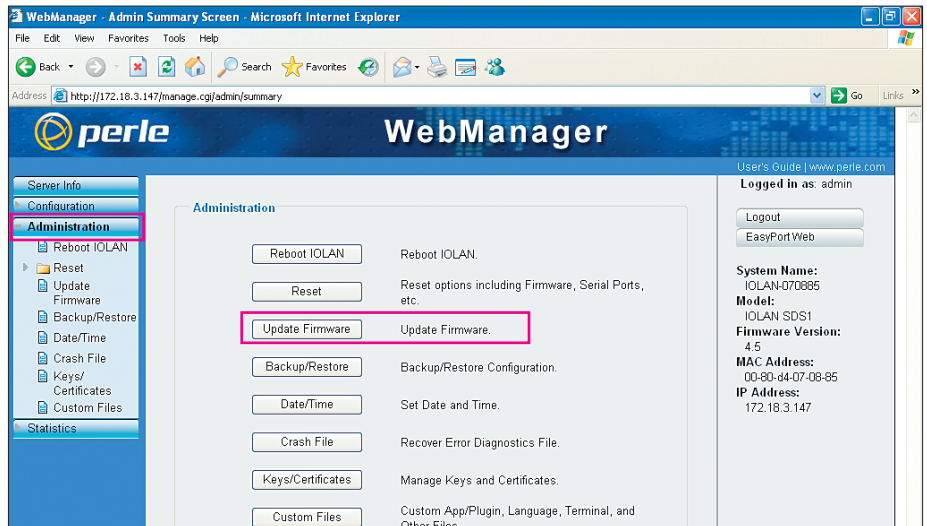


- ▶ Enter username “admin” (1) and password “superuser” (2).
- ▶ Press button “Login”.



- ▷ Informations about the used Perle device are shown on the right side including the installed firmware version.

Updating firmware

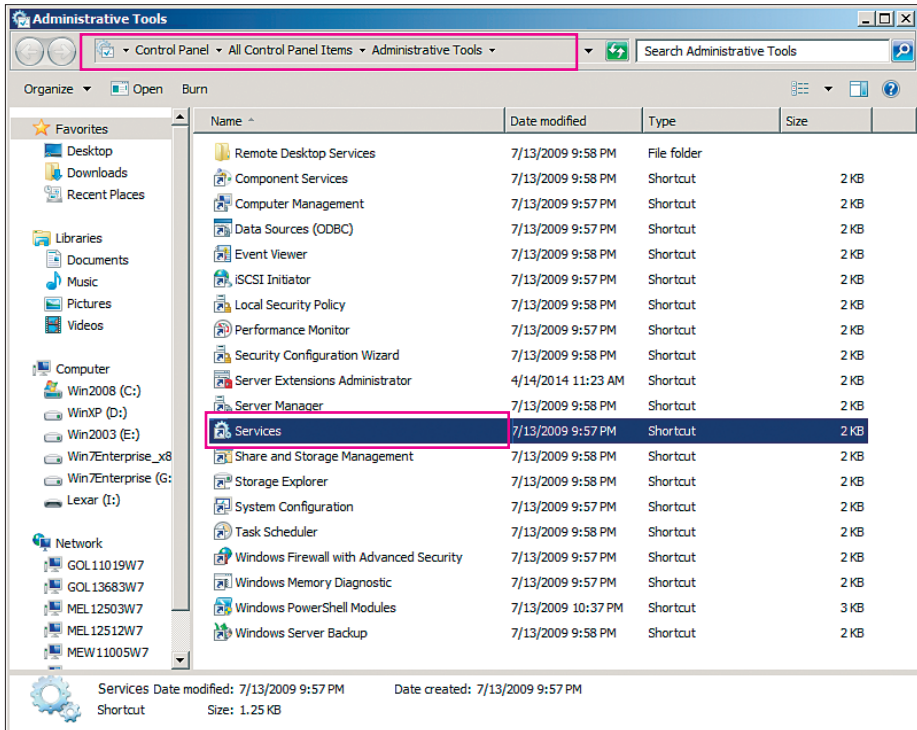


- ▶ If the installed firmware is below version 4.5, upgrade the firmware version of the Perle device.
- The newest firmware version can always be downloaded from the internet www.perle.com/downloads.

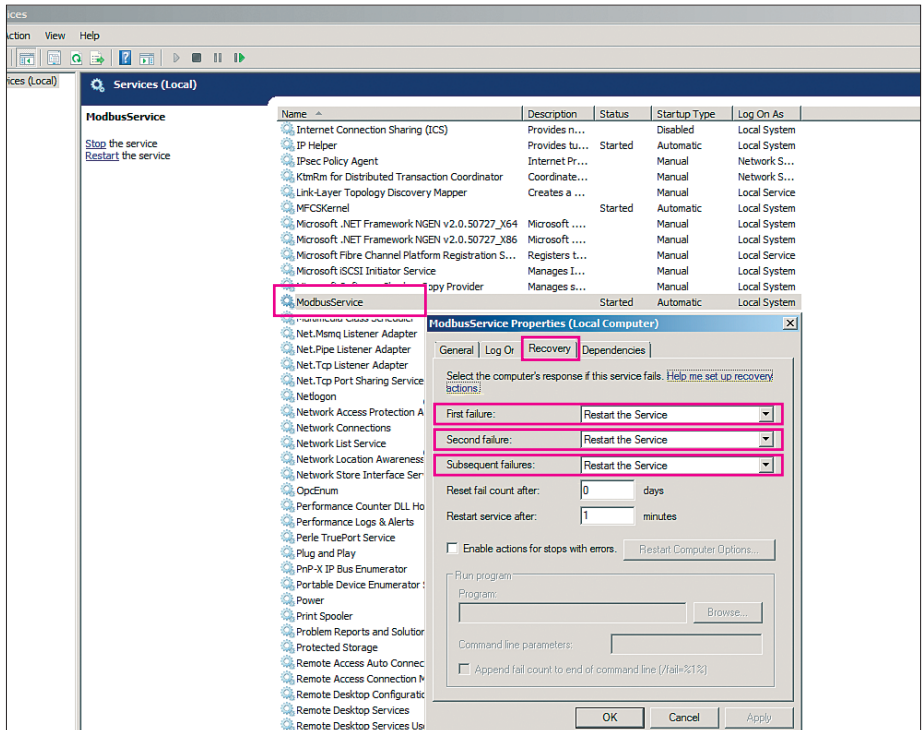
NOTE

How the firmware will be installed at the Perle device is described in chapter “Downloading IOLAN Firmware” on page 348 in the manual “IOLAN_SCS-SDS-ST5_UG_v4.4.pdf”, available on the BioPAT® Xgas VIS CD.

9.2 Configuring Settings for Windows ModbusService



- ▶ Open “Windows Control Panel / All Control Panel Items / Administrative Tools”.
- ▶ Double click “Services”.



- ▶ Scroll down the list until “ModbusService” appears in the list.
- ▶ Double click “ModbusService” and select register card “Recovery”.
- ▶ Select the setting “Restart the Service” from the drop down list for the three items “First failure”, “Second failure” and “Subsequent failure”.

10 Troubleshooting

If you cannot correct a fault independently, contact:

Sartorius Stedim Systems GmbH
Robert-Bosch-Straße 5-7
34302 Guxhagen, Germany

Tel.: +49 5665 407 0
www.sartorius.com

e-mail: mfcs.service@sartorius-stedim.com

Sartorius Stedim Systems GmbH
Robert-Bosch-Strasse 5-7
34302 Guxhagen, Germany

Phone: +49 5665 407 0
www.sartorius.com

The information and figures contained in these instructions correspond to the version date specified below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

Masculine or feminine forms are used to facilitate legibility in these instructions and always simultaneously denote the other gender as well.

Copyright notice:

This instruction manual, including all of its components, is protected by copyright.

Any use beyond the limits of the copyright law is not permitted without our approval.

This applies in particular to reprinting, translation and editing irrespective of the type of media used.

© Sartorius Germany

Last updated:
06 | 2021

Printed in the EU on paper bleached
without chlorine. | AM
Publication No.: SBI6018-e210602