

Thermo Scientific

KingFisher Presto

Integration Guide

Cat. No. N17647 Rev 1.0 2016

Thermo
SCIENTIFIC

Thermo Fisher Scientific Inc. provides this document to its customers with a product purchase to use in the product operation. This document is copyright protected and any reproduction of the whole or any part of this document is strictly prohibited, except with the written authorization of Thermo Fisher Scientific Inc.

The contents of this document are subject to change without notice. All technical information in this document is for reference purposes only. System configurations and specifications in this document supersede all previous information received by the purchaser.

Thermo Fisher Scientific Inc. makes no representations that this document is complete, accurate or error-free and assumes no responsibility and will not be liable for any errors, omissions, damage or loss that might result from any use of this document, even if the information in the document is followed properly.

This document is not part of any sales contract between Thermo Fisher Scientific Inc. and a purchaser. This document shall in no way govern or modify any Terms and Conditions of Sale, which Terms and Conditions of Sale shall govern all conflicting information between the two documents.

Release history:

For Research Use Only. Not for use in diagnostic procedures.

Introduction

This guide includes all documentation related to integrating the KingFisher Presto instrument to an automation environment.

The following documents are included in this guide:

- *KingFisher Presto Developer's Guide*
- *KingFisher Presto Interface Specification*
- *KFModule.dll Interface Specification*
- *ThermoUSB.dll Interface Specification*
- *ThermoLAN.dll Interface Specification*
- *ThermoCOM.dll Interface Specification*

Related Documentation

In addition to this guide, Thermo Fisher Scientific provides the following documents for KingFisher Presto:

- *Thermo Scientific™ KingFisher™ Presto User Manual* (Cat.no. N17413)
- *Thermo Scientific™ BindIt™ Software User Manual* (Cat. no. N07974)

Safety and Special Notices

Make sure you follow the precautionary statements presented in this guide. The safety and other special notices appear in boxes.

Safety and special notices include the following:



CAUTION Highlights hazards to humans, property, or the environment. Each CAUTION notice is accompanied by an appropriate CAUTION symbol.

IMPORTANT Highlights information necessary to prevent damage to software, loss of data, or invalid test results; or might contain information that is critical for optimal performance of the system.

Note Highlights information of general interest.

Tip Highlights helpful information that can make a task easier.

Contacting Us

For the latest information on products and services, visit our website at:

www.thermofisher.com/kingfisher

Thermo Scientific

KingFisher Presto

Developer's Guide

Contents

1	KingFisher Presto Developer's Guide	1
1.1	Contents	1
1.2	Confidential	1
1.3	Introduction	2
2	System Description	3
3	Interfacing to KFModule.dll	7
3.1	Building the Integration Sample	7
3.2	Running Tests with KingFisher Presto Simulator	8
4	Creating and Uploading Protocols	9
4.1	BindIt Protocol Editor	10
4.2	Creating "My Test Protocol"	10
4.3	Uploading "My Test Protocol"	11
4.4	Uploading and Downloading Protocols	11
5	Protocol Execution Methods	13
5.1	Step-by-Step Execution	13
5.2	Event-Based Execution	15
5.3	Executing Multiple Overlapping Protocols	17
5.4	Precautions for Heater Control	22
6	Error Handling	25
6.1	Communication Interfaces	25
6.2	Communication Protocol	25
6.3	KingFisher Presto Protocol Integrity	26

Chapter 1

KingFisher Presto Developer's Guide

1.1 Contents

- [Introduction](#)
- [System Description](#)
- [Interfacing to KFModule.dll](#)
- [Creating and Uploading Protocols](#)
- [Protocol Execution Methods](#)
- [Error Handling](#)

1.2 Confidential

This document has been prepared by Thermo Fisher Scientific Oy to be used solely for the purposes defined by Thermo Fisher Scientific Oy. Use for other purposes is not authorized.

Please note that any and all information contained in this document is the property of Thermo Fisher Scientific Oy. This confidential information ("Confidential Information") shall not be reproduced in whole part or disclosed to any third party without the prior written approval of Thermo Fisher Scientific Oy. The receiving party shall ensure that its employees, officers, representatives and agents shall not disclose to third parties any Confidential Information.

Upon written request from Thermo Fisher Scientific Oy, the receiving party shall promptly return all Confidential Information or destroy all Confidential Information.

1.3 Introduction

This document is intended for software designers writing computer programs for controlling the instrument. It contains information necessary to know in order to be able to write such a program. It is assumed that the reader of this document is familiar with the function of the instrument.

Chapter 2

System Description

Figure 3-1 shows the basic architecture of a system using the KingFisher instrument. The system is constructed into three layers: **Application layer**, **Communication layer** and **Hardware layer**. Components using the layers are **Automation System**, **BindIt Software** and **KingFisher Presto Instrument**.

System Components

Automation System refers to a process management system (PMS) software module provided by an automation integrator and running on a computer with a MS Windows operating system.

BindIt Software is used to create KingFisher Presto protocols. It can also upload and download protocols to and from the KingFisher Presto instrument using KFModule.dll, but it is not part of the automation system and it is not used to run the protocols in the automation system.

KingFisher Presto instrument provides communication interfaces for transferring and executing KingFisher Presto protocols created in BindIt software.

Application Services Layer

KFModule.dll dynamic link library provides easy access to the KingFisher Presto instrument interface through communication libraries. The physical connection between a computer and the instrument can be USB, RS232 or LAN.

Communication Services Layer

This is a middle layer between the KingFisher Presto instrument and the KFModule.dll and can be ignored by the Automation System. Depending on a physical connection, one of the communication libraries will be used: **ThermoUSB.dll**, **ThermoCOM.dll** or **ThermoLAN.dll**.

Hardware Interfaces

The KingFisher Presto instrument has three alternate communication ports for connecting to a computer: RS232, USB and LAN. All connections can be accessed either through higher level service layers or directly through physical ports.

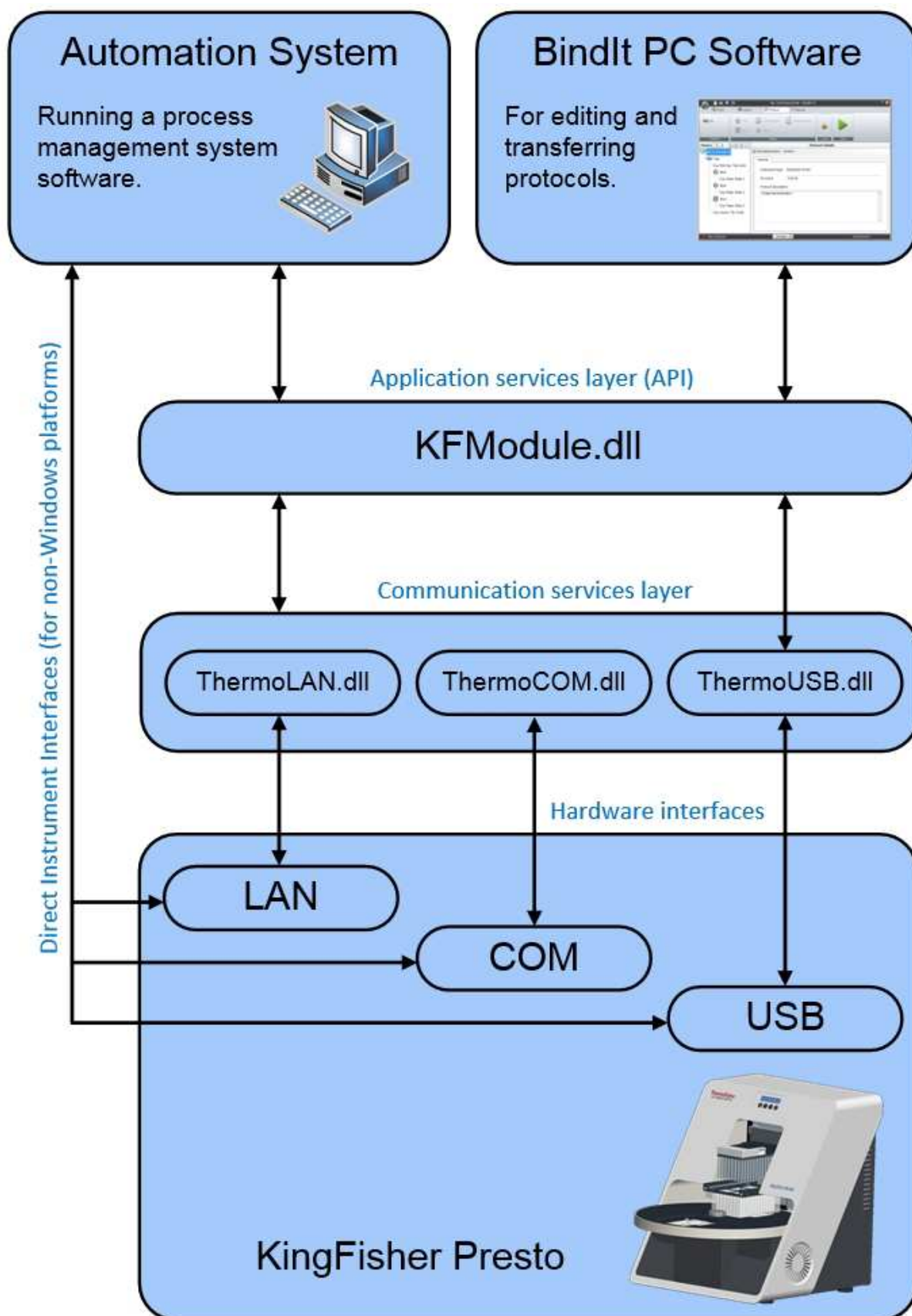


Figure 2.1: System description

Related Documents

BindIt User Manual: Thermo Scientific BindIt Software for KingFisher Instruments User Manual (Cat. No. N07974)

[KingFisher Presto Developer's Guide](#)

[KingFisher Presto Interface Specification](#)

[KFModule.dll Interface Specification](#)

[ThermoUSB.dll Interface Specification](#)

[ThermoLAN.dll Interface Specification](#)

[ThermoCOM.dll Interface Specification](#)

Chapter 3

Interfacing to KFModule.dll

This chapter introduces an integration sample (C#) which demonstrates how to use the functions of the KFModule.dll.



Note. The sample solution is created with Microsoft Visual Studio 2012 and provided as is, it has not been fully tested for stability and may malfunction when used in unexpected way.

For detailed information of the KFModule.dll exported functions and instrument commands, see: [KFModule.dll Interface Specification](#) and [KingFisher Presto Interface Specification](#).

3.1 Building the Integration Sample

Look for a folder named **IntegrationSample** provided together with this document. From that folder, open the **Sample.sln** with Microsoft Visual Studio 2012 or later.

Select Debug configuration and build the solution. You should get an output listing similar to one below.

```
1>----- Build started: Project: KFModuleWrapperLibrary, Configuration: Debug Any CPU -----
1>  KFModuleWrapperLibrary
   -> C:\temp\IntegrationSample\KFModuleWrapperLibrary\bin\Debug\KFModuleWrapperLibrary.dll
2>----- Build started: Project: Sample, Configuration: Debug Any CPU -----
2>  Sample -> C:\temp\IntegrationSample\Sample\bin\Debug\Sample.dll
2>          1 file(s) copied.
2>          1 file(s) copied.
2>          1 file(s) copied.
===== Build: 2 succeeded, 0 failed, 0 up-to-date, 0 skipped =====
```

Note the project named **KFModuleWrapperLibrary**. It provides an interface to the KFModule.dll and can be used as a starting point when designing your own interface library.

The project **Sample** contains test methods demonstrating how to communicate with a KingFisher Presto Simulator. The difference between connecting to the simulator or to a real instrument is very small. You just choose to call a function `KFModule_OpenSimulator()` instead of `KFModule_Open()`. The simulator is included in the **IntegrationSample**.

3.2 Running Tests with KingFisher Presto Simulator

Open Test Explorer from the Visual Studio and you should see these four unit tests:

1. TEST1_SimulatorStarted
2. TEST2_UploadProtocol
3. TEST3_RunProtocol
4. TEST4_RemoveProtocol

The KingFisher Presto simulator **KFPresto.exe** needs to be started before running the tests. See also **StartSimulator.bat** in the IntegrationSample package. It can be used to start the simulator with logging enabled. Then all communication messages between IntegrationSample and simulator are written to a file named **kfm.log**.

After simulator is started, run the tests in the given order and you should succeed.

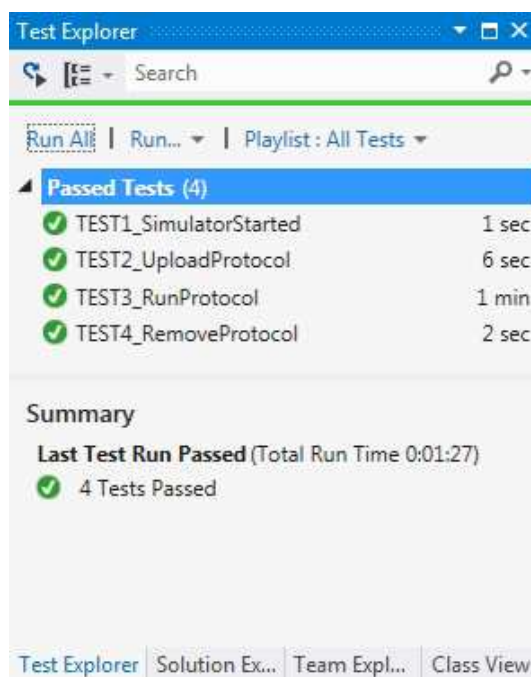


Figure 3.1: Passed tests

Chapter 4

Creating and Uploading Protocols

This chapter contains instructions how to create KingFisher Presto protocols and upload them into the instrument using BindIt PC software or KFModule.dll API. An example protocol named "My Test Protocol" is introduced. The example protocol is addressed later on in this document, when it is explained how to execute uploaded protocols in the KingFisher Presto instrument.

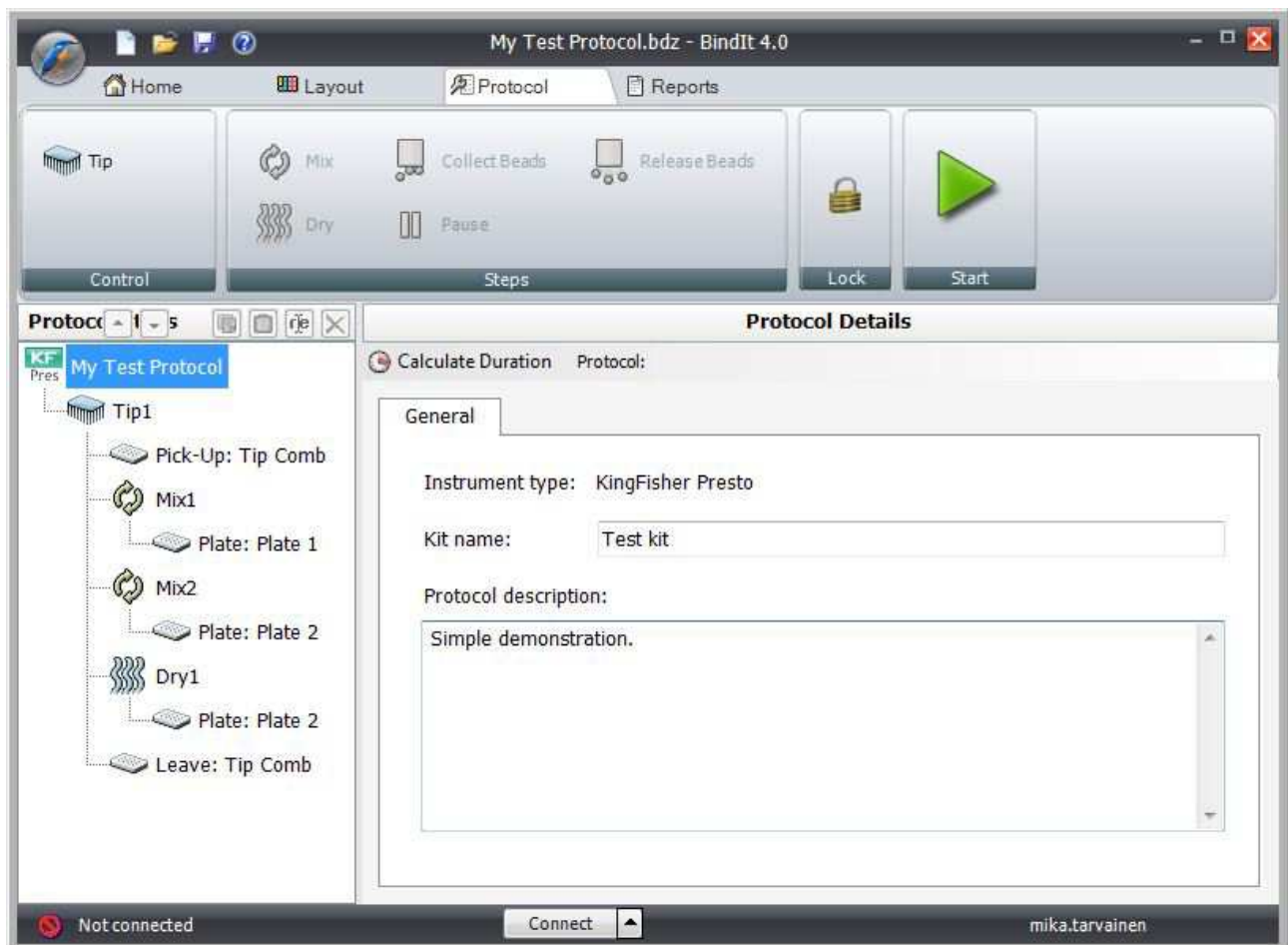


Figure 4.1: BindIt protocol editor

4.1 BindIt Protocol Editor

KingFisher Presto protocols are created with **BindIt** PC software. It can also upload and download the protocols to and from the KingFisher Presto instrument but it is not part of the Automation System. It can be used to run protocols "manually", meaning that a human person needs to handle all plate load/remove events.

4.2 Creating "My Test Protocol"

Here are listed the steps needed to create the "My Test Protocol" example protocol. The outcome is a protocol structure with few plates and steps as shown in figure 5.2. Detailed instructions of how to create protocols can be found in the Thermo Scientific BindIt Software for KingFisher Instruments User Manual.

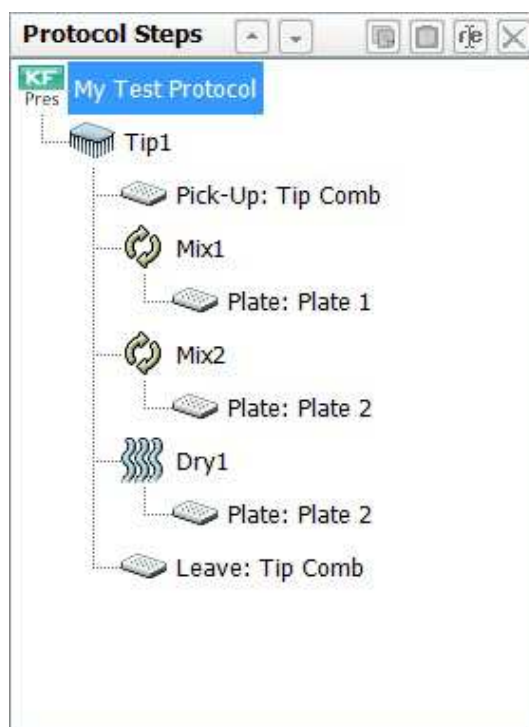


Figure 4.2: My Test Protocol

1. Open BindIt and select "Home" tab.
2. Press button "New" from the "Protocol" button group.
3. Select KingFisher Presto from the opened list.
4. Press button "New" from the "Plate" button group in "Layout" tab.
5. Select "96 DW plate"
6. Type "Tip Comb" to the "Plate name:" text box.
7. Add a dummy reagent by typing "None" to the reagent name row with guide text "Type a name to add new...". Default 50 µl volume is ok.
8. Create two more plates in a similar way, but name them "Plate 1" and "Plate 2".
9. Select "Protocol" tab.

10. Select "Tip1" from the "Protocol Steps" list and then "96 DW tip comb" from the "Tip:" dropdown list on the right on "General" tab.
11. Select "Pick-Up" from the "Protocol Steps" list and then "Tip Comb" from the "Plate:" dropdown list on the right on "Details" tab.
12. Select "Leave" from the "Protocol Steps" list and then "Tip Comb" from the "Plate:" dropdown list on the right on "Details" tab.
13. Select "Tip1" from the "Protocol Steps" list.
14. Press button "Mix" from the "Steps" button group in "Protocol" tab to create a mix step. Default parameters are ok.
15. Select "Plate" under the "Mix1" in the "Protocol Steps" list.
16. Select "Plate 1" from the from the "Plate:" dropdown list on the right on "Details" tab.
17. Create another Mix step in a similar way, but Select "Plate 2" for it.
18. Create a Dry step in a similar way and select "Plate 2" for it.
19. Save the protocol as "My Test Protocol" by pressing the save symbol on top and left of the window.
20. Verify that you created identical step list as in figure 5.2

4.3 Uploading "My Test Protocol"

The easiest way to upload and test a protocol with a KingFisher Presto instrument is by connecting the instrument to BindIt and pressing a green Start symbol. However, the protocol is not saved permanently to the memory of the instrument and is available only during the execution of the protocol in BindIt software.

The green Start symbol can be found from the "Start" group, which is visible in all tabs and enabled after a protocol is created and saved. See quick guide steps below.

1. Open BindIt.
2. Power on a KingFisher Presto instrument and attach it to the computer running the BindIt with USB cable.
3. Connect BindIt to the instrument, see Thermo Scientific BindIt Software for KingFisher Instruments User Manual.
4. Open "My TestProtocol".
5. Press Start symbol from the "Start" group.

4.4 Uploading and Downloading Protocols

Connect to a KingFisher Presto instrument and open protocol transfer view in BindIt Software by pressing the "Transfer" symbol in the "Instrument" group of the "Home" tab. In this view it is possible to upload, download and remove protocols to/from the instrument. Note that the "Instrument" group is enabled only after the KingFisher Presto instrument is connected to the BindIt Software. Notice the button "Connect" or "Disconnect" in the status bar in the bottom of the window. See Thermo Scientific BindIt Software for KingFisher Instruments User Manual for more details about connecting to the instrument.

Protocols can be also saved to the file system folders of the computer running the BindIt Software. The file format of these files is "BindIt export data file" and the file extension is ".bdz". BindIt protocols can be uploaded and downloaded to/from the KingFisher Presto instrument through KFFModule.dll API functions `KFFModule_UploadProtocol()` and `KFFModule_DownloadProtocol()`. See [KFFModule.dll Interface Specification](#) for more details.

Chapter 5

Protocol Execution Methods

The automation interface offers two distinct ways to execute protocols on the KingFisher Presto instrument. The integrator can decide which option suits his/her needs the best. The first **Step-by-step** approach leaves the total control of the process to the automation system. The instrument control can be seen as a master-slave relationship between automation system and instrument. Alternatively, KingFisher Presto protocols can be executed using an **event-based** approach where the automation system has to react to events sent by the instrument. KingFisher Presto supports also overlapped multi-protocol executions as an application of Step-by-step method.



Caution. The validities of the KingFisher Presto protocols provided by ThermoFisher Scientific are verified using [Event-Based Execution](#) method. It is up to the integrator to ensure that the protocols are executed correctly when using [Step-by-Step Execution](#) method or when [Executing Multiple Overlapping Protocols](#).

5.1 Step-by-Step Execution

The automation can be build to run sequential step execution scripts where the turntable rotation commands are placed between the steps. The system does not need to react to the events coming from the instrument, it can poll the status of the execution instead.

KingFisher Presto protocols are created with BindIt protocol editor and uploaded to the instrument through the communication interface of the instrument. Protocol steps are executed by addressing them by name. As said before, turntable rotation is handled by the automation or the process management system. Understanding of the content of the protocol is not required for an integrator. Only the order of the steps needs to be understood. The execution of the "My Test Protocol" example protocol is descibed below. It assumes that the protocol is already uploaded, see [Creating and Uploading Protocols](#).



Note. The example demonstrates how the KingFisher Presto instrument replies to the commands and how the automation system needs to poll the instrument status before sending new commands. Not all of these repeatedly needed commands and responses are shown in order to keep the example readable. Furthermore, all event messages are stripped off. See [KingFisher Presto Interface Specification](#) for more details.

1. Rotate turntable to a startup position, meaning that nest 1 will be in the processing position. If the command can be executed, then the KingFisher Presto instrument replies immediately with "ok" response and starts rotating the turntable.

```
<Cmd name="Rotate" nest="1" position="1" />
```

```
<Res name="Rotate" ok="true" />
```

2. Read the status of the instrument until the reply from the instrument is "Idle" instead of "Busy".


```
<Cmd name="GetStatus" />
<Res name="GetStatus" ok="true"><Status>Busy</Status></Res>
<Cmd name="GetStatus" />
<Res name="GetStatus" ok="true"><Status>Busy</Status></Res>
...
...
...
<Cmd name="GetStatus" />
<Res name="GetStatus" ok="true"><Status>Idle</Status></Res>
```
3. Load Tip Comb plate to the nest 2, which is in the load position, and rotate it to the processing position. Poll again for the instrument to return to the "Idle" state.


```
<Cmd name="Rotate" nest="2" position="1" />
<Res name="Rotate" ok="true" />
<Cmd name="GetStatus" />
<Res name="GetStatus" ok="true"><Status>Idle</Status></Res>
```
4. Start the "Pick-Up" step. Again, wait for the "Idle" status after the instrument has first replied with "ok" to the start command.


```
<Cmd name="StartProtocol" protocol="My Test Protocol" tip="Tip1" step="Pick-Up" />
<Res name="StartProtocol" ok="true" />
<Cmd name="GetStatus" />
<Res name="GetStatus" ok="true"><Status>Busy</Status></Res>
<Cmd name="GetStatus" />
<Res name="GetStatus" ok="true"><Status>Busy</Status></Res>
...
...
...
<Cmd name="GetStatus" />
<Res name="GetStatus" ok="true"><Status>Idle</Status></Res>
```
5. Place "Plate 1" to the nest 1, which is now in the load position and rotate it to the processing position.

Note. The plate could have been loaded also during the time when the instrument was running the "Pick-Up" step.

```
<Cmd name="Rotate" nest="1" position="1" />
```
6. Unload "Tip Comb" plate from the nest 2 in the load/unload position of the instrument and start the first mix step "Mix1" for the "Plate 1", which is now in the processing position.


```
<Cmd name="StartProtocol" protocol="My Test Protocol" tip="Tip1" step="Mix1" />
```
7. Place "Plate 2" to the nest 2, which is now in the load position and rotate it to the processing position.


```
<Cmd name="Rotate" nest="2" position="1" />
```
8. Unload "Plate 1" from the nest 1 and start step "Mix2" for the "Plate 2" in the processing position.


```
<Cmd name="StartProtocol" protocol="My Test Protocol" tip="Tip1" step="Mix2" />
```
9. Start step "Dry1" for the "Plate 2" in the processing position.


```
<Cmd name="StartProtocol" protocol="My Test Protocol" tip="Tip1" step="Dry1" />
```
10. Load Tip Comb plate to the nest 1, which is in the load position, and rotate it to the processing position. "Plate 2" will be rotated to the load/unload position.

```
<Cmd name="Rotate" nest="1" position="1" />
```

11. Unload Plate 2 from the load/unload position of the instrument and start step "Leave" to drop off the tips to the "Tip Comb" plate.

```
<Cmd name="StartProtocol" protocol="My Test Protocol" tip="Tip1" step="Leave" />
```

12. Rotate the turntable one more time to get the "Tip Comb" plate to the load/unload position. **Note.** This specific command example addresses the position 2, which is the load/unload position. Command "rotate nest 2 to the position 1" would have the same effect.

```
<Cmd name="Rotate" nest="1" position="2" />
```

13. Unload Tip Comb plate from the load/unload position after the instrument has finished the final rotation.

```
<Cmd name="GetStatus" />
```

```
<Res name="GetStatus" ok="true"><Status>Idle</Status></Res>
```



Caution. Minimize delay times between steps.

Tips/magnets of the KingFisher Presto are always lifted up from a microwell plate and thus from the liquids within the plate when a step execution is finished. It is very important to minimize the delay before the next step so that the tips do not dry between the steps. Otherwise some applications may suffer. In practice, all the plates must be prepared beforehand. It is also recommended that the plate changes are designed in the way that a next plate is loaded to the load position during the time when a previous plate is being processed.

5.2 Event-Based Execution

KingFisher Presto protocols can be executed in a mode where the instrument handles turntable rotation by itself and sends events to the automation system whenever it requires attention. See example below demonstrating messaging between the instrument and an automation system running the "My Test Protocol" execution in event mode. The example is explained in detail after the listing. Note that not all events are listed in order to keep the listing readable. These events include **StepStarted** **ProtocolTimeLeft** and **Temperature**. See [KingFisher Presto Interface Specification](#) for more details about events, commands and responses.

```
<Cmd name="StartProtocol" protocol="My Test Protocol" /> (1)
<Res name="StartProtocol" ok="true" />
```

```
<Evt name="LoadPlate" plate="Tip Comb" optional="false" /> (2)
<Cmd name="Acknowledge" />
<Res name="Acknowledge" ok="true" />
```

```
<Evt name="LoadPlate" plate="Plate 1" optional="true" /> (3)
<Evt name="LoadPlate" plate="Plate 1" optional="false" />
<Cmd name="Acknowledge" />
<Res name="Acknowledge" ok="true" />
```

```
<Evt name="ChangePlate" optional="true" > (4)
  <Evt name="RemovePlate" plate="Tip Comb" optional="true" />
  <Evt name="LoadPlate" plate="Plate 2" optional="true" />
</Evt>
<Cmd name="Acknowledge" />
<Res name="Acknowledge" ok="true" />
```

```

<Evt name="ChangePlate" optional="true" >                                     (5)
  <Evt name="RemovePlate" plate="Plate 1" optional="true" />
  <Evt name="LoadPlate" plate="Tip Comb" optional="true" />
</Evt>
<Cmd name="Acknowledge" />
<Res name="Acknowledge" ok="true" />

```

```

<Evt name="RemovePlate" plate="Plate 2" optional="true" />                                     (6)
<Cmd name="Acknowledge" />
<Res name="Acknowledge" ok="true" />

```

```

<Evt name="RemovePlate" plate="Tip Comb" optional="false" />                                     (7)
<Cmd name="Acknowledge" />
<Res name="Acknowledge" ok="true" />

```

```

<Evt name="Ready" />                                                         (8)

```

1. Protocol is started by sending command "StartProtocol" to which the instrument immediately replies.
2. The first "LoadPlate" event is sent by the instrument when it is requesting "Tip Comb" plate to be inserted to the load position of the turntable. Value "false" of the attribute "optional" means that the instrument does not continue its operation until the automation system acknowledges the event. After a command "Acknowledge" is received and replied, the instrument rotates the turntable and starts processing the plate, which in this example means picking up the tip comb.
3. The second "LoadPlate" event is sent immediately after the processing of the tip comb is started. Attribute "optional" is now "true" indicating that the "Plate 1" can be placed to the load position but it is not obligatory. This event is ignored by the automation system in this example and so the instrument continues processing. After finished with the "Tip Comb" plate, the instrument sends another event requiring the "Plate 1" to be loaded. Attribute "optional" is now "false" and the plate needs to be placed to the load position. The instrument continues after it receives command "Acknowledge" and sends a reply to it. Again, the turntable is rotated and processing of the "Plate 1" starts.
4. There are plates in both nests of the turntable now. Processing of the "Plate 1" is under progress and the next plate needed is "Plate 2". Instrument sends a "ChangePlate" event to the automation system, telling that now it is a good time to remove "Tip Comb" from the load position and replace it with "Plate 2". Automation system does that and acknowledges. Instrument replies and sends no further events of this particular need, but rotates the turntable automatically after processing of the "Plate 1" is finished and starts processing "Plate 2" without any delay.
5. See previous. Event "ChangePlate" is sent in order to remove "Plate 1" and load "Tip Comb".
6. At this phase of the protocol execution the instrument has rotated the turntable again and is starting to drop off the tip comb to the "Tip Comb" plate. Before that, it sends a "RemovePlate" event suggesting that the "Plate 2" could be removed. Note that attribute "optional" is "true" meaning that it is just a suggestion, automation system could also remove the plate later. In this example the plate is removed by this time and command "Acknowledge" is sent and replied.
7. The instrument is finished with the "Tip Comb" plate and is now requesting it to be removed. Event "RemovePlate" is sent with attribute "optional" value "false". Automation system removes the last plate and sends a final "Acknowledge" command. Instrument replies.
8. Event "Ready" is sent by the instrument when the protocol execution is completed. This event needs not to be acknowledged.

5.3 Executing Multiple Overlapping Protocols

KingFisher Presto supports overlapped multi-protocol executions.



Caution. If a protocol contains **Mix** steps with pre-heating option enabled, then there is some differences in heater control between straightforward protocol execution and overlapped multi-protocol executions. See [Precautions for Heater Control](#) for more details.

See figure 6.1 below for an example protocol created in the BindIt protocol editor. The name of the protocol is "Protocol A" and it uses two microplates: one for a tip comb and another for mixing. Actual applications would never use just one plate for processing, but this way the following discussion is not too exhaustive.

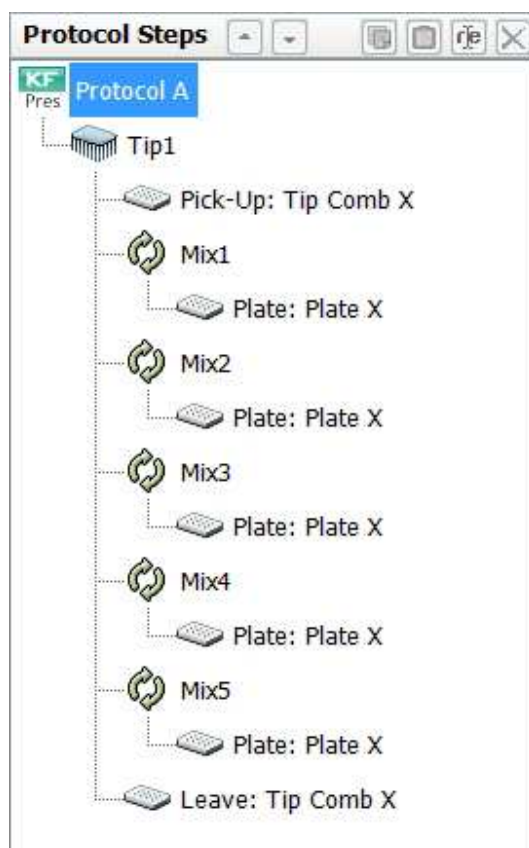


Figure 5.1: Protocol A

Probably the most likely scenario for multiprotocol execution is to run the protocols sequentially. See Figure 6.2 for example. However, some automation system integrators may find it useful that [Step-by-Step Execution](#) enables multi-protocol execution in an overlapped way, meaning that an automating system can execute several instances of a same protocol or different protocols by simply switching between the step lists. See Figure 6.3 for example.

It should be noted that typically every separate execution uses different tips and plates than the other executions, meaning that the context switch between executions requires "**Pick-Up**" and "**Leave**" steps around the actual step to be executed. Conceptually, this means that the context switch is actually a plastics switch where plastics are microwell plates and tip combs for the magnets. Need for a plastics switch is especially clear in Figure 6.3 in comparison with straightforward repeated execution as in the Figure 6.2 below.

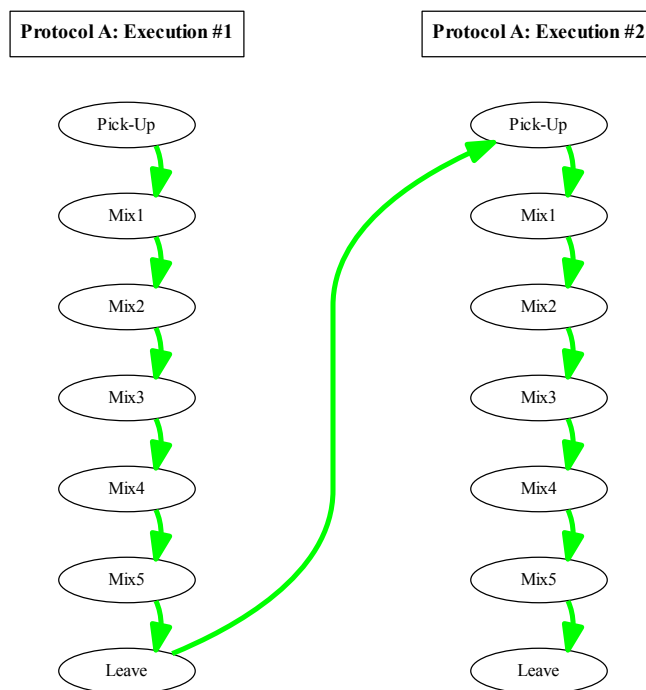


Figure 6.2: Consecutive executions of Protocol A

Figure 6.2 above and the command listing on the next page illustrate how a **Protocol A** is executed twice using Step-by-step execution method. The example protocol executions contain tips in a "Tip Comb" plates 1 and 2 and the actual plates to be processed are named as "Plate 1" and "Plate 2".

Note that the execution status polling and command reply messages are not listed in order to keep the example more readable. See [Step-by-Step Execution](#) for more details on those.

1. `<Cmd name="Rotate" nest="1" position="2" />`
2. Place "Tip Comb 1" plate to the turntable.
3. `<Cmd name="Rotate" nest="2" position="2" />`
4. Place "Plate 1" to the turntable.
5. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Pick-Up" />`
6. `<Cmd name="Rotate" nest="2" position="1" />`
7. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix1" />`
8. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix2" />`
9. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix3" />`

10. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix4" />`
11. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix5" />`
12. `<Cmd name="Rotate" nest="2" position="2" />`
13. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Leave" />`
14. Remove "Plate 1" from the turntable.
15. `<Cmd name="Rotate" nest="1" position="2" />`
16. Remove "Tip Comb 1" plate from the turntable.
17. Place "Tip Comb 2" plate to the turntable.
18. `<Cmd name="Rotate" nest="2" position="2" />`
19. Place "Plate 2" to the turntable.
20. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Pick-Up" />`
21. `<Cmd name="Rotate" nest="2" position="1" />`
22. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix1" />`
23. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix2" />`
24. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix3" />`
25. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix4" />`
26. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix5" />`
27. `<Cmd name="Rotate" nest="2" position="2" />`
28. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Leave" />`
29. Remove "Plate 2" from the turntable.
30. `<Cmd name="Rotate" nest="1" position="2" />`
31. Remove "Tip Comb 2" plate from the turntable.

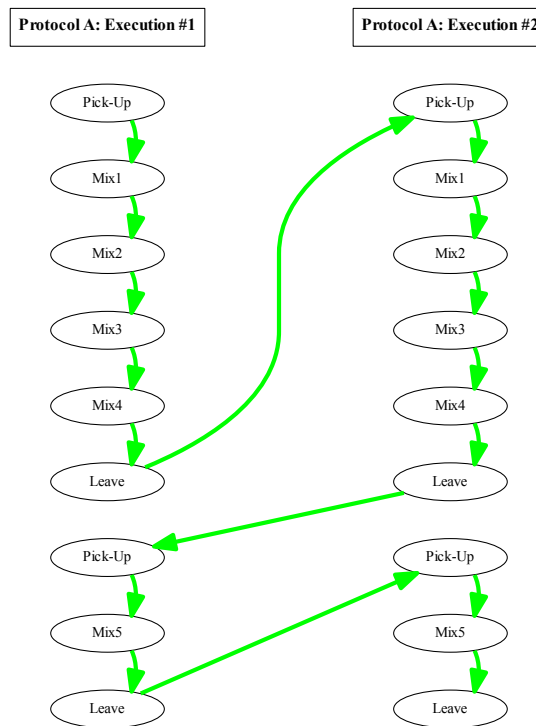


Figure 6.3: Context switch between two executions of Protocol A

Figure 6.3 above and the command listing below demonstrates switching of plastics when two instances or executions of a **Protocol A** are executed in an overlapped fashion.

The example protocol executions contain tips in a "Tip Comb" plates 1 and 2 and the actual plates to be processed are named as "Plate 1" and "Plate 2". Note that the execution status polling and command reply messages are not listed in order to keep the example more readable. See [Step-by-Step Execution](#) for more details on those.

1. `<Cmd name="Rotate" nest="1" position="2" />`
2. Place "Tip Comb 1" plate to the turntable.
3. `<Cmd name="Rotate" nest="2" position="2" />`
4. Place "Plate 1" to the turntable.
5. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Pick-Up" />`
6. `<Cmd name="Rotate" nest="2" position="1" />`
7. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix1" />`
8. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix2" />`
9. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix3" />`
10. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix4" />`
11. `<Cmd name="Rotate" nest="1" position="1" />`
12. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Leave" />`

13. Remove "Plate 1" from the turntable.
14. Place "Tip Comb 2" plate to the turntable.
15. `<Cmd name="Rotate" nest="2" position="1" />`
16. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Pick-Up" />`
17. Remove "Tip Comb 1" from to the turntable.
18. Place "Plate 2" to the turntable.
19. `<Cmd name="Rotate" nest="1" position="1" />`
20. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix1" />`
21. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix2" />`
22. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix3" />`
23. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix4" />`
24. `<Cmd name="Rotate" nest="2" position="1" />`
25. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Leave" />`
26. Remove "Plate 2" from the turntable.
27. Place "Tip Comb 1" plate to the turntable.
28. `<Cmd name="Rotate" nest="1" position="1" />`
29. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Pick-Up" />`
30. Remove "Tip Comb 2" from to the turntable.
31. Place "Plate 1" to the turntable.
32. `<Cmd name="Rotate" nest="2" position="1" />`
33. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix5" />`
34. `<Cmd name="Rotate" nest="1" position="1" />`
35. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Leave" />`
36. Remove "Plate 1" from the turntable.
37. Place "Tip Comb 2" plate to the turntable.
38. `<Cmd name="Rotate" nest="2" position="1" />`
39. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Pick-Up" />`
40. Remove "Tip Comb 1" from to the turntable.
41. Place "Plate 2" to the turntable.
42. `<Cmd name="Rotate" nest="1" position="1" />`
43. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Mix5" />`
44. `<Cmd name="Rotate" nest="2" position="1" />`
45. `<Cmd name="StartProtocol" protocol="Protocol A" tip="Tip1" step="Leave" />`
46. Remove "Plate 2" from the turntable.
47. `<Cmd name="Rotate" nest="2" position="2" />`
48. Remove "Tip Comb 2" from to the turntable.

5.4 Precautions for Heater Control

Mix steps in KingFisher Presto protocols may have heating option enabled, meaning that a heating block is heated and lifted up against a plate during mixing. Furthermore, a preheat option can be enabled to ensure that the heating block will reach the target temperature just before the actual step where the heating is required. The heating block is positioned some distance below the plate during preheating.

There are some differences in heater control between straightforward protocol execution and overlapped multi-protocol executions when the preheat option is enabled. See Figure 6.4 for an example where pre-heat is enabled for the step **Mix5**. When the protocol is executed using [Event-Based Execution](#) method, then the KingFisher Presto instrument calculates the moment when the preheating needs to be started and then turns the heater on automatically. The moment arises in our example during step **Mix4**.

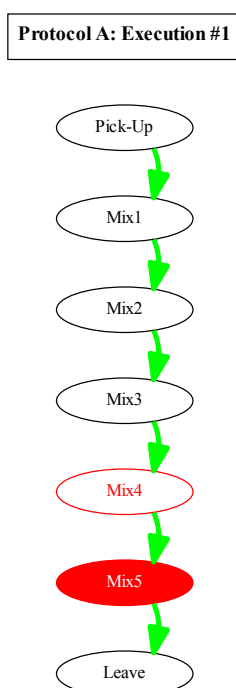


Figure 6.4: Protocol A with preheat enabled in Mix5

If [Step-by-Step Execution](#) method is used or when [Executing Multiple Overlapping Protocols](#), then the KingFisherPresto follows the same automatic preheat control logic as when using [Event-Based Execution](#) method but with some restrictions:

1. If preheat is on and the step to be executed is not from the same protocol as previous executed step, then the preheat is turned off unless:
 - Step to be executed is "Pick-Up" or "Leave".
2. If preheat is on and the step to be executed is from the same protocol as previous executed step then the preheat is turned off unless:
 - Step to be executed is "Pick-Up" or "Leave".
 - Step to be executed is same as previous executed step.
 - Step to be executed is next in the protocol after the previous executed step.

3. Preheat is turned off automatically if a new step execution is not started within 10 minutes from the end of the previous step execution.
4. Preheat is never turned on during "Pick-Up" or "Leave" steps when [Step-by-Step Execution](#) method is used or when [Executing Multiple Overlapping Protocols](#). This may cause a small delay to the preheating when compared to [Event-Based Execution](#) method.
5. If the step to be executed is the first step in a protocol, then the heater will be turned off at the beginning of the step execution.
6. If the step to be executed is the last step in a protocol and the step to be executed is from the same protocol as previous executed step, then the heater will be turned off at the end of the step execution.



Note. If preheat is turned off, because of the above mentioned restrictions, then a warning event is sent to the automation system. Same warning is also included to the reply message of instrument status query command as long as the step is being executed.



Caution. If a **Mix** step with heating and preheating contains long **End of step** actions e.q. **Postmix** or/and **Collect beads** and is switched between multiple execution instances, then the preheating is compromised. That is because the heater is always turned off after **Mixing** (e.q. before **Postmix**) and the heating block will be cooled down during the **Postmix**.

Chapter 6

Error Handling

There are several different error handling mechanisms in the communication interfaces of the KingFisher Presto instrument. Some of them are implemented in DLL level while some others are features of the instrument itself. See related interface specifications for detailed information:

- [KFModule.dll Interface Specification](#)
- [KingFisher Presto Interface Specification](#)

6.1 Communication Interfaces

Low level error handling is based on the mechanisms of the underlying communication protocol, for example on guaranteed data integrity of the USB protocol. Some connection errors can be captured in DLL level, whereas some others require polling of the instrument status.

KFModule.dll is able to detect when USB connection is lost. Function **KFModule_AttachEvent()** can be used to subscribe **KFMODULE_ERROR** events. If the connection is lost, then event **KFM_ERROR_DISCONNECTED** will be sent.

There is no detection for failed connection when using RS232 interface. Then the user application e.g. an automation system needs to poll the instrument status with **GetStatus** command in order to be sure that the instrument is still online.

It is recommended to use the above mentioned **GetStatus** method also when using LAN interface, because the KFModule.dll can only detect a failed connection after some timeout when attempting to send commands to the instrument. Meaning that any spontaneous events or some replies are lost if the physical connection is disconnected.

6.2 Communication Protocol

KFModule.dll function calls all return error codes if they do not succeed.

Responses to instrument commands use "ok" attribute for replying if a command is accepted or rejected. Responses may also contain error and/or warning codes.

Reply to **GetStatus** instrument status query command may indicate "In error" status and contain error and/or warning codes. The controlling automation system needs to send an error acknowledge command to the instrument in order to get the instrument out of the error state.

Failed protocol and step executions and other spontaneous errors are reported with error events which may require that the controlling system sends an error acknowledge command before the instrument continues its operation.

Command set includes **Stop** and **Abort** commands for stopping and disconnecting the instrument. Abort command can be generated also from the physical maintenance interface by pressing the red stop symbol.

6.3 KingFisher Presto Protocol Integrity

Protocol upload command of the KingFisher Presto includes a 32-bit cyclic redundancy check (CRC) error-detecting code from the protocol data to be transferred. The instrument verifies that the CRC matches to the CRC of the received data. The CRC value is saved to the internal memory of the instrument together with the protocol data. The KFModule.dll API function KFModule_UploadProtocol() calculates CRC codes automatically.

Furthermore, the instrument calculates CRC from a saved protocol every time a protocol or a step from a protocol is about to be started. An error is generated if the calculated CRC does not match to the CRC value that was saved when the protocol in question was uploaded to the instrument.

Thermo Scientific

KingFisher Presto

Interface Specification

Contents

1	KingFisher Presto Interface Specification	1
1.1	Contents	1
1.2	Confidential	1
1.3	Introduction	2
2	Hardware Requirements	3
2.1	RS232	3
2.1.1	Abort	3
2.2	USB	3
2.2.1	Sending commands to the instrument	4
2.2.2	Receiving responses from the instrument	4
2.2.3	Abort	4
2.2.4	Flow Control	4
2.3	LAN	4
2.3.1	Abort	5
3	Communication Protocol	7
3.1	General	7
3.2	Character Encoding	7
3.3	Maximum Line Width	7
3.4	XML Format	7
3.5	Attributes and Tag Data	7
3.6	Root Tag Types	8
3.6.1	<Cmd> - Commands	8
3.6.2	<Res> - Responses	8
3.6.3	<Evt> - Events	8
3.7	Error Handling	9
3.8	Error Codes	10
3.9	Warning Codes	10
4	Commands and Responses	11
4.1	Abort	12
4.2	Acknowledge	14
4.3	Connect	15
4.4	Disconnect	17
4.5	DownloadProtocol	18
4.6	ErrorAcknowledge	19
4.7	GetProtocolDuration	20
4.8	GetProtocolTimeLeft	22
4.9	GetStatus	23
4.10	ListProtocols	24
4.11	RemoveProtocol	25
4.12	Rotate	26
4.13	SetTemperatureReporting	27
4.14	StartProtocol	28
4.15	Stop	29

4.16 UploadProtocol	30
5 Events	31
5.1 Aborted	32
5.2 ChangeMagnets	33
5.3 ChangePlate	34
5.4 Error	35
5.5 LoadPlate	36
5.6 Pause	37
5.7 ProtocolTimeLeft	38
5.8 Ready	39
5.9 RemovePlate	40
5.10 StepStarted	41
5.11 Temperature	42
6 Appendix - XML path syntax	43

Chapter 1

KingFisher Presto Interface Specification

1.1 Contents

- [Introduction](#)
- [Hardware Requirements](#)
- [Communication Protocol](#)
- [Commands and Responses](#)
- [Events](#)

1.2 Confidential

This document has been prepared by Thermo Fisher Scientific Oy to be used solely for the purposes defined by Thermo Fisher Scientific Oy. Use for other purposes is not authorized.

Please note that any and all information contained in this document is the property of Thermo Fisher Scientific Oy. This confidential information ("Confidential Information") shall not be reproduced in whole part or disclosed to any third party without the prior written approval of Thermo Fisher Scientific Oy. The receiving party shall ensure that its employees, officers, representatives and agents shall not disclose to third parties any Confidential Information.

Upon written request from Thermo Fisher Scientific Oy, the receiving party shall promptly return all Confidential Information or destroy all Confidential Information.

1.3 Introduction

This document is intended for software designers writing computer programs for controlling the instrument. It contains information necessary to know in order to be able to write such a program. This document can also be used for black box testing of the instrument. It is assumed that the reader of this document is familiar with the function of the instrument.

Syntaxes of commands, responses and events are documented using a style similar to the `Xpath` syntax. For examples and more detailed description of the usage, see [Appendix - XML path syntax](#).

Chapter 2

Hardware Requirements

The KingFisher Presto instrument has three alternate communication ports for connecting to a computer: RS232, USB and LAN.

2.1 RS232

Serial port parameters are fixed to the following values: baud rate: 115200, 1 start bit, 8 data bits, 1 stop bit and no parity.

The serial connector on the instrument is a 9 pin male D connector. Reception is through pin 3, transmission through pin 2 and signal ground is at pin 7.

XON/XOFF flow control is used. No hardware handshaking is used. The XON character is 0x11 and the XOFF character 0x13. No frames or checksums are used in the RS232 protocol. The instrument ignores any NUL bytes (0x00) it receives.

Maximum length of a command line is limited by the receive buffer size, which is 512 bytes. Note however that XOFF will be sent when the buffer is half full (256 characters).

2.1.1 Abort

To send the Abort command to the instrument, the computer must first flush it's transmit buffer and force transmit flow control to enabled state. Only then is the Abort command sent.

Abort command is a single character, the escape control code, which is decimal 27 or hex 0x1B.

2.2 USB

The instrument has a standard USB series "B" receptacle for connecting to a PC. The USB interface complies with the USB 2.0 full speed device specification. The device class of the instrument is HID (Human Interface Device), and it has only one configuration which is numbered 1. In this configuration there is one interface with an interrupt IN endpoint (0x81) and an interrupt OUT endpoint (0x01).

The HID Report descriptor of the instrument defines three reports: A 64 byte Input report, a 64 byte Output report and a 2-byte Feature report. As there is only one type of each report, no report ID:s are used. The usage of all the reports is vendor specific. Read the USB Device Class Definition for Human Interface Devices for more information of the HID device interface.

The following identification information is returned during enumeration:

Vendor id: 0x0AB6

Product id: 0x02C9

Manufacturer: Thermo Fisher Scientific Oy

Serial number: The serial number string of the instrument.

2.2.1 Sending commands to the instrument

Commands are sent using the 64 byte HID class Output report. The first byte of the report must be the number of actual data (command) bytes in the report. The command data starts from the second byte of the report. If the command data does not fill up the whole report, the instrument discards the remaining bytes of the report.

2.2.2 Receiving responses from the instrument

The PC software receives the instrument responses in a 64 byte HID class Input report. The first byte of the report is the number of actual response bytes in the report. The response data starts from the second byte of the report. If the response data does not fill up the whole report, the PC software must ignore the remaining bytes of the report.

2.2.3 Abort

Because there is no way the PC application can force an Abort to the interrupt OUT endpoint past previous commands, aborting must be performed in two phases.

First the PC must send the two-byte Feature report through the control endpoint. The first byte of the report must be nonzero and the second zero. This causes the instrument to discard all commands it may already have received and also discard all commands it receives until the Abort command is received.

Then the PC must send the Abort character through the interrupt OUT endpoint just like any other command.

2.2.4 Flow Control

When the receive buffer of the instrument becomes full, it simply stops reading the Output reports. This means that any command writes the PC software makes do not complete until the instrument resumes reading the reports.

If the PC application cannot process the Input reports at the same rate as the instrument sends them, it should send a Feature report to the instrument to request it stop sending the Input reports. The first byte of the Feature report must be zero and the second nonzero. When the PC software can again receive more Input reports, it should send a Feature report with both bytes zero.

2.3 LAN

The instrument can optionally be connected through a Local Area Network. The interface is IEEE 802.3 compliant and has a RJ-45 connector for connection to a 10BASE-T network.

Before the LAN interface can be used, the Ethernet interface MAC address and the TCP port number must be programmed using the PAR 38 and PAR 39 commands, respectively.

The instrument gets an IP address dynamically from a DHCP server. Unless the server is configured to give a fixed IP address to the instrument, a PC software wishing to connect to the instrument through LAN must use the WS-Discovery protocol to find the transport address of the instrument.

There are three strings the instrument uses for matching the Types of a WS-Discovery Probe message: ThermoDevice, KingFisherPresto and SN_*, where * is the serial number string of the instrument. An example Probe looking for KingFisherPresto instrument serial number 12345:

```
<?xml version="1.0" encoding="UTF-8"?>
<s:Envelope
  xmlns:s="http://www.w3.org/2003/05/soap-envelope">
```

```

xmlns:a="http://www.w3.org/2005/08/addressing">
xmlns:d="http://docs.oasis-open.org/ws-dd/ns/discovery/2009/01">
  <s:Header>
    <a:Action>http://docs.oasis-open.org/ws-dd/ns/discovery/2009/01/Probe</a:Action>
    <a:MessageID>urn:uuid:dc280767-49b0-4c5e-97f3-95f961df0053</a:MessageID>
    <a:ReplyTo>
      <a:Address>http://www.w3.org/2005/08/addressing/anonymous</a:Address>
    </a:ReplyTo>
    <a:To>urn:docs-oasis-open-org:ws-dd:ns:discovery:2009:01</a:To>
  </s:Header>
  <s:Body>
    <Probe xmlns="http://docs.oasis-open.org/ws-dd/ns/discovery/2009/01">
      <d:Types>ThermoDevice KingFisherPresto SN_12345</Types>
    </Probe>
  </s:Body>
</s:Envelope>

```

If the instrument is connected to LAN, it responds with a ProbeMatches message. The <d:Xaddrs> field of the response is the transport address to use to communicate with the instrument. The transport address consists of the instrument IP address and the TCP port number the instrument is listening.

Although the instrument always responds to WS-Discovery messages, it only lets one client at a time to connect to the TCP port. The client should close the port when done communicating with the instrument to allow another client to connect.

2.3.1 Abort

Sending the Abort character to the instrument is a two phase process.

First the PC must send an UDP message containing text "Abort" (without quotes or newline) to the same UDP port number as is used for the TCP and wait a moment for an identical UDP response from the instrument. The UDP response is sent to the same UDP port as the PC used for sending the UDP message. If no UDP response, retry at least two times using short timeout.

Then, when an UDP response is received or no response after retries, send Abort character (0x1B) to the TCP port.

Chapter 3

Communication Protocol

3.1 General

This document describes the commands and the responses as they are seen by the controlling software and the instrument without the framing added by the data transport layer.

3.2 Character Encoding

All commands and responses are ASCII encoded except values, which are encoded using UTF-8 character encoding.

3.3 Maximum Line Width

Communication is based on CR or/and LF terminated lines.



Note. The maximum length of a input line including carriage return and line feed characters is **200**.

3.4 XML Format

All commands, responses and events are XML elements. Elements may or may not contain line feed and carriage return characters. Both start-tag-end-tag and empty-element-tag formats are supported. Tag and parameter names are not case sensitive.

Syntaxes of commands, responses and events are documented using a style similar to the [XPath](#) syntax. For examples and more detailed description of the usage, see [Appendix - XML path syntax](#).

3.5 Attributes and Tag Data

Attribute values can be booleans, integers, floating point numbers or strings. Each command and response contains different amount of attributes and tags and the order of those is not fixed.

3.6 Root Tag Types

See table below for the three different kind of root tags used in the communication protocol between the KingFisher Presto instrument and a controlling system.

Tag	Description
Cmd	Command from the controlling system to the KingFisher Presto instrument
Res	Response to a command
Evt	Event message from the KingFisher Presto instrument to the controlling system

3.6.1 <Cmd> - Commands

Commands are sent from a controlling system to the KingFisher Presto instrument. They are used for example to start a protocol execution.

Examples:

```
<Cmd name="MyCommand" parameter1="Something">
    <Something>More complex value</Something>
</Cmd>

<Cmd name="MyOtherCommand" parameter1="Nonething" />
```

See [Commands and Responses](#) for the complete list of available commands.

3.6.2 <Res> - Responses

Responses to the commands contain attribute named "ok", which can be used to quickly check if a command was successful.

Examples:

```
<Res name="MyCommand" ok="false">
    <Thing>Lots of things</Thing>
</Res>

<Res name="MyOtherCommand" ok="true" />
```

See [Commands and Responses](#) for detailed information.

3.6.3 <Evt> - Events

Events are used to indicate data periodically or in the case of events occurring. Some events require acknowledgement from the controlling system, see command [Acknowledge](#) for more information.

Examples:

```
<Evt name="Temperature" value="23"/>

<Evt name="LoadPlate" plate="Plate 1"/>

<Evt name="LoadPlate" plate="Plate 2"/>
```

```
<Evt name="ChangePlate">
  <Evt name="RemovePlate" plate="Plate 2"/>
  <Evt name="LoadPlate" plate="Plate 3"/>
</Evt>

<Evt name="RemovePlate" plate="Plate 3"/>
```

See [Events](#) for the complete list of events sent by the KingFisher Presto instrument.

3.7 Error Handling

Low level error handling is based on the mechanisms of the underlying communication protocol, for example on guaranteed data integrity of the USB protocol.

[Commands and Responses](#) use "ok" attribute for replying if a command is accepted or rejected, see [<Res> - Responses](#). Responses to commands may also contain [Error Codes](#) and/or [Warning Codes](#).

Failed protocol and step executions and other spontaneous errors are reported with [Error](#) events which may require that the controlling system sends an [ErrorAcknowledge](#) command before the instrument continues its operation.

Examples

```
<Cmd name="StartProtocol" protocol="KF Blood 12 DW"/>

<Res name="StartProtocol" ok="true" />

<Cmd name="StartProtocol" protocol="KF Blood 12 DW"/>

<Res name="StartProtocol" ok="false">
  <Error code="124">Protocol already running.</Error>
</Res>

<Evt name="Error" ack="true">
  <Error code="321">Execution failed</Error>
</Evt>
```

3.8 Error Codes

Code	Description
2	Received an unknown command.
3	Already connected to another port.
4	Head position error.
5	Magnets position error.
6	Turntable position error.
7	Heater unit position error.
8	Lock position error.
11	Invalid command argument.
13	Protocol memory error.
14	Protocol memory is full.
15	No protocols found from the protocols memory.
16	Protocol was not found from the protocols memory.
17	Given tip name was not found from the protocol.
18	Given step name was not found from the given tip of the protocol.
19	A name of a step to start was not given.
20	A name of a tip where to start the step was not given.
23	Protocol name is invalid. Maximum length of the name is 100 bytes e.g. 100 ASCII characters.
24	Invalid protocol file.
25	Protocol is not executable.
27	Protocol is too large and can't be loaded.
28	Instrument is executing, please wait.
32	No protocol is currently running.
33	Data transmit to USB port failed (timed out).
34	Cannot run magnets down without tips.
35	Magnetic head is missing.
38	Plate not detected in processing position.
39	Plate detected in processing position.
40	Plate not detected in load position.
41	Plate detected in load position.
43	Protocol is not for this instrument.

3.9 Warning Codes

Code	Description
101	Instrument is already connected.
102	Previous command was incomplete.
103	Date/time string is invalid, instrument time was not set.
104	Protocol execution was aborted by the user.
105	Step execution was aborted by the user.
106	Existing protocol was overwritten.
107	Heater preheat was turned off between single step executions.
108	Heater was turned off after 10 minutes since last single step execution.

Chapter 4

Commands and Responses



Note. Remember to terminate commands with a new line character (ASCII 10).

- [Abort](#)
- [Acknowledge](#)
- [Connect](#)
- [Disconnect](#)
- [DownloadProtocol](#)
- [ErrorAcknowledge](#)
- [GetProtocolDuration](#)
- [GetProtocolTimeLeft](#)
- [GetStatus](#)
- [ListProtocols](#)
- [RemoveProtocol](#)
- [Rotate](#)
- [SetTemperatureReporting](#)
- [StartProtocol](#)
- [Stop](#)
- [UploadProtocol](#)

4.1 Abort

Aborts any ongoing operation in the instrument.

Syntax

The Abort command is an exception to the normal command syntax. For example for RS232 interface, it is a single character, the escape control code, which is decimal 27 or hex 0x1B. When the Abort character is received, the execution of a possible ongoing protocol, step or some other process is stopped. Heating is turned off and all the motors are stopped. Heating block is driven down if possible. Response to the Abort command is replied immediately and event [Ready](#) is sent after motors are stopped/driven. In case of errors during driving, event [Error](#) is sent.

Abort command differs from the command [Stop](#) in the way, that the message buffers of the communication port are flushed and an existing communication connection is disconnect. Event [Aborted](#) is sent before disconnection to a connected communication port if the Abort command was received from some other communication port.

Immediate response to the Abort command follows the same syntax as the other [Commands and Responses](#) use.

Syntax

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "Abort"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if the command was accepted

Details

For the instrument to react to the Abort command immediately it must reach the instrument promptly after it is sent. This means that the Abort command must override all previous commands queued to the instrument. This is not trivial and is done differently for different interfaces.

Serial port Before sending the Abort character, the flow control must be forced to XON state so that the Abort character will actually be sent.

USB port Because there is no way the PC application can force an Abort to the interrupt OUT endpoint past previous commands, aborting must be performed in two phases.

First the PC must send the two-byte Feature report through the control endpoint. The first byte of the report must be nonzero and the second zero. This causes the instrument to discard all commands it may already have received and also discard all commands it receives until the Abort command is received.

LAN port First the PC must send an UDP message containing text "Abort" (without quotes or newline) to the same UDP port number as is used for the TCP and wait a moment for an identical UDP response from the instrument. The UDP response is sent to the same UDP port as the PC used for sending the UDP message.

If no UDP response, retry at least two times using short timeout.

When an UDP response is received or no response after retries, send Abort character (0x1B) to the TCP port.

Using KFModule.dll By far the easiest way to send Abort to the instrument is to use the `KFModule_Abort()` function of the `KFModule.dll` library. It will automatically select the right Abort procedure depending on which interface is used.

Example

```
<Res name="Abort" ok="true"/>
```

4.2 Acknowledge

General notification message for acknowledging the KingFisher Presto instrument.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "Acknowledge"	Name of the command

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "Acknowledge"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was succesfull

Details

This command is used to acknowledge the instrument in various different kind of situations, for example after [LoadPlate](#) event.

Example

```
<Cmd name="Acknowledge"/>
```

```
<Res name="Acknowledge" ok="true"/>
```

4.3 Connect

Establish a connection between a controlling system and a KingFisher Presto instrument.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String:"Connect"	Name of the command
Cmd@setTime	Date/time string in format: "YYYY-MM-DD hh:mm:ss"	Optional attribute for setting the instrument date and time

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "Connect"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was succesfull
Res/Instrument/text()	String: Max 100 chars	Name/type of the instrument
Res/Version/text()	String: Max 100 chars	Firmware version string
Res/Serial/text()	String: Max 100 chars	Instrument serial number

Details

The instrument does not accept any other commands before this command is succesfully received. Instrument serial number, type and firware version number is returned in the reply message.

Optional attribute "setTime" can be used to set the date and time of the instrument. There is no battery backup for the calendar of the instrument, so the date and time needs to be re-set every time the instrument powers on. Note that for example KingFisherDII Windows driver does this automatically.

Response message may contain warning and error codes depending on the state of the instrument. See command [Get-Status](#). See also examples below for most typical use cases.

Example 1 - Succesfully connected

```
<Cmd name="Connect"/>

<Res name="Connect" ok="true">
  <Instrument>KingFisher Presto</Instrument>
  <Version>0.0.0</Version>
  <Serial>123-456</Serial>
</Res>
```

Example 2 - Connect with date/time setting

```
<Cmd name="Connect" setTime="2015-09-04 09:29:59"/>

<Res name="Connect" ok="true">
  <Instrument>KingFisher Presto</Instrument>
  <Version>0.0.0</Version>
  <Serial>123-456</Serial>
</Res>
```

Example 3 - Already connected

```
<Cmd name="Connect" />

<Res name="Connect" ok="true">
  <Warning code="101">Instrument is already connected.</Warning>
  <Instrument>KingFisher Presto</Instrument>
  <Version>0.0.0</Version>
  <Serial>123-456</Serial>
</Res>
```

Example 4 - Connection is reserved for another communication port

```
<Cmd name="Connect" />

<Res name="Connect" ok="false">
  <Error code="3">Already connected to another port.</Error>
</Res>
```

Example 5 - Connection is established but the instrument is in error state

```
<Cmd name="Connect" />

<Res name="Connect" ok="true">
  <Error code="4">Head position error.</Error>
  <Instrument>KingFisherPresto</Instrument>
  <Version>0.0.1</Version>
  <Serial>123-456</Serial>
</Res>
```

4.4 Disconnect

Disconnect an established connection between a controlling system and a KingFisher Presto instrument.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "Disconnect"	Name of the command

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "Disconnect"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was succesfull

Details

Note that the instrument does not reply to this command if the connection is already closed. See command [Connect](#) for more details.

Example

```
<Cmd name="Disconnect"/>
```

```
<Res name="Disconnect" ok="true"/>
```

4.5 DownloadProtocol

Download a protocol from the internal memory of the KingFisher Presto instrument.

Note! It's much easier to use `KFModule_DownloadProtocol()` API function from `KFModule.dll` than this command directly. `KFModule_DownloadProtocol` handles base64 decoding and outputs BindIt .bdz files.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "DownloadProtocol"	Name of the command
Cmd@protocol	String: Max 100 bytes	Name of the protocol to be transferred, case sensitive.

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "DownloadProtocol"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was successful
Res/CDATA	Base64 string/strings	XML character data section containing base64 encoded protocol file

Details

Protocol is received as base64 encoded binary data in a CDATA section of the response message.

Example

```
<Cmd name="DownloadProtocol" protocol="KingFisher Presto Blood 24 DW"/>

<Res name="DownloadProtocol" ok="true">
  <![CDATA[
    PHN0ZXA+DQpoZXJlIGlzIGFu
    IGV4YW1wbGUNCmZvciB0aGUg
    YmFzZTY0IGVuY29kZXINCjwv
    c3RlcD4=
  ]]>
</Res>
```

4.6 ErrorAcknowledge

This command must be used to clear instrument errors.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "ErrorAcknowledge"	Name of the command

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "ErrorAcknowledge"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was succesfull

Details

When an error is detected during execution, it is reported using [Error](#) event. This command must be send in order to clear the error state of the instrument. See also command [GetStatus](#).

Example

```
<Cmd name="ErrorAcknowledge"/>
```

```
<Res name="ErrorAcknowledge" ok="true"/>
```

4.7 GetProtocolDuration

Get KingFisher Presto protocol duration.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "GetProtocolDuration"	Name of the command
Cmd@protocol	String: Max 100 chars	Name of a protocol, case sensitive.

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "GetProtocolDuration"	Name of the command.
Res@ok	Boolean string: "true" or "false"	Error status quick peek.
Res/Init@protocol	String	Name of the protocol.
Res/Init/TimeStamp[1,2]@at	Time string: hh:mm:ss	Start or stop time of the protocol initialization.
Res/Init/TimeStamp[1,2]@type	Type string of the timestamp: "Start" or "Stop"	Init step uses only "Start" and "Stop" types.
Res/Init/TimeStamp[1,2]@step	String "Init"	Fixed step name.
Res/Init/TimeStamp[2]@duration	XML Duration data type	Duration of the protocol initialization phase.
Res/Tip/TimeStamp[n]@at	Time string: hh:mm:ss	Timestamp referenced to the start time of the protocol initialization.
Res/Tip/TimeStamp[n]@type	Type string of the timestamp: "Start", "Event" or "Stop"	Time of step start/stop or an pause event. See command GetProtocolTimeLeft .
Res/Tip/TimeStamp[n]@step	String	Name of a step, fixed "pseudo" step or user defined, see event StepStarted .
Res/Tip/TimeStamp[n]@duration	XML Duration data type	Duration of a step.
Res/Tip/TimeStamp[n]@msg	String	Message string for "Event" type.
Res/Tip/TimeStamp[n]@plate	String	Name of a plate for "Event" type.
Res/Tip/TimeStamp[n]@remove	String	Name of a plate to remove for "ChangePlate" event type.
Res/Tip/TimeStamp[n]@load	String	Name of a plate to load for "ChangePlate" event type.
Res/Init@protocol	String	Name of the protocol.
Res/Finish/TimeStamp[1,2]@at	Time string: hh:mm:ss	Start or stop time of the protocol finalization.
Res/Finish/TimeStamp[1,2]@type	Type string of the timestamp: "Start" or "Stop"	Finish step uses only "Start" and "Stop" types.
Res/Finish/TimeStamp[1,2]@step	String "Init"	Fixed step name.
Res/Finish/TimeStamp[2]@duration	XML Duration data type	Duration of the protocol initialization phase.
Res/Total@duration	XML Duration data type	Total duration of the protocol.

Details

This command returns the full time structure of a KingFisher Presto protocol as a step by step list of timestamps in addition to the total duration.

Examples

```

<Cmd name="GetProtocolDuration" protocol="My Test Protocol" />

<Res name="GetProtocolDuration" ok="true">
  <Init protocol="My Test Protocol">
    <TimeStamp at="00:00:00" type="Start" step="Init"/>
    <TimeStamp at="00:00:00" type="Stop" step="Init" duration="PT0S"/>
  </Init>
  <Tip name="Tip1">
    <TimeStamp at="00:00:00" type="Start" step="Pick-Up"/>
    <TimeStamp at="00:00:00" type="Event" msg="Load plate" plate="Tip Comb"/>
    <TimeStamp at="00:00:08" type="Stop" step="Pick-Up" duration="PT8S"/>
    <TimeStamp at="00:00:08" type="Start" step="Mix1"/>
    <TimeStamp at="00:00:08" type="Event" msg="Load plate" plate="Plate 1"/>
    <TimeStamp at="00:01:41" type="Stop" step="Mix1" duration="PT1M33S"/>
    <TimeStamp at="00:01:41" type="Start" step="Mix2"/>
    <TimeStamp at="00:01:41" type="Event" msg="Change plate" remove="Tip Comb" load="Plate 2"/>
    <TimeStamp at="00:03:15" type="Stop" step="Mix2" duration="PT1M34S"/>
    <TimeStamp at="00:03:15" type="Start" step="Dry1"/>
    <TimeStamp at="00:04:17" type="Stop" step="Dry1" duration="PT1M2S"/>
    <TimeStamp at="00:04:17" type="Start" step="Leave"/>
    <TimeStamp at="00:04:17" type="Event" msg="Change plate" remove="Plate 1" load="Tip Comb"/>
    <TimeStamp at="00:04:24" type="Stop" step="Leave" duration="PT7S"/>
    <TimeStamp at="00:04:24" type="Start" step="Unload"/>
    <TimeStamp at="00:04:24" type="Event" msg="Remove plate" plate="Plate 2"/>
    <TimeStamp at="00:04:24" type="Event" msg="Remove plate" plate="Tip Comb"/>
    <TimeStamp at="00:04:26" type="Stop" step="Unload" duration="PT2S"/>
  </Tip>
  <Finish protocol="My Test Protocol">
    <TimeStamp at="00:04:26" type="Start" step="Finish"/>
    <TimeStamp at="00:04:26" type="Stop" step="Finish" duration="PT0S"/>
  </Finish>
  <Total duration="PT4M26S"/>
</Res>

```

4.8 GetProtocolTimeLeft

Get time left of a KingFisher Presto protocol or single step execution.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "GetProtocolTimeLeft"	Name of the command
Cmd@protocol	String: Max 100 chars	Name of a protocol, case sensitive.

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "GetProtocolTimeLeft"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek.
Res/TimeToPause@value	XML Duration data type	Time to next plate load/change/remove or pause event
Res/TimeLeft@value	XML Duration data type	Total time left of the protocol/step execution

Details

This command returns time left to a next pause event and total time left of a protocol execution.

Note! Element **TimeToPause** is not included when running single step execution. Note also that then the value of **TimeLeft** refers to the total execution time of a step being executed.

A pause event is an event which needs to be acknowledged with [Acknowledge](#) command. Non-optional [LoadPlate](#), [ChangePlate](#), [RemovePlate](#) and [ChangeMagnets](#) events are also pause events in addition to obvious [Pause](#) event.

Note that [Error](#) is not a pause event and it needs to be acknowledged with a special [ErrorAcknowledge](#) command.

See also event [ProtocolTimeLeft](#).

Examples

```
<Cmd name="StartProtocol" protocol="My Test Protocol"/>
  <Res name="StartProtocol" ok="true"/>
  <Cmd name="GetProtocolTimeLeft"/>

  <Res name="GetProtocolTimeLeft" ok="true">
    <TimeToPause value="PT0S"/>
    <TimeLeft value="PT30M53S"/>
  </Res>

  <Cmd name="StartProtocol" protocol="My Test Protocol" tip="Tip1" step="Mix1"/>
  <Res name="StartProtocol" ok="true"/>
  <Cmd name="GetProtocolTimeLeft"/>

  <Res name="GetProtocolTimeLeft" ok="true">
    <TimeLeft value="PT2M42S"/>
  </Res>
```

4.9 GetStatus

Get instrument status: **Idle**, **Busy** or **In error**.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "GetStatus"	Name of the command

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "GetStatus"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was succesfull and the instrument is not in an error state.
Res/Status/text()	String: "Idle", "Busy" or "In error"	Status of the instrument. Idle: waiting for a new commmand, Busy: instrument is executing a command, In error: instrument is in an error state.
Res/Error@code	uint16_t: see Error Codes	Error code number
Res/Error/text()	String: Max 100 characters	Description of an error

Details

This command is used to query the status or state of the instrument. When the instrument is executing [StartProtocol](#) or [Rotate](#) command, it will reply with **Busy** status. When an execution is finished or after succesfull power on initialization, the instrument will reply with **Idle** status. If an error occurs during an execution or initialization, then the instrument will transition to an error state and reply with **In error** status to the [GetStatus](#) command. An [ErrorAcknowledge](#) command must be send in order to clear the error. No other commands are accepted until the error is cleared. Expect commands [Disconnect](#) and [Connect](#). Note that [Error](#) event is also sent if and an error occurs during execution.

Examples

```
<Cmd name="GetStatus"/>

<Res name="GetStatus" ok="true">
  <Status>Idle</Status>
</Res>

<Res name="GetStatus" ok="true">
  <Status>Busy</Status>
</Res>

<Res name="GetStatus" ok="false">
  <Error code="5">Turntable position error.</Error>
  <Status>In error</Status>
</Res>
```

4.10 ListProtocols

List protocols in the internal memory of the KingFisher Presto instrument.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "ListProtocols"	Name of the command

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "ListProtocols"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was successful
Res/Protocols/Protocol[n]	String: Max 100 chars	Unique name of a protocol "n"
Res/MemoryUsed@value	uint16_t: 0 - 100	Protocol memory usage percentage

Details

This command returns a list of all available protocols in the instrument and a protocol memory usage percentage.

Example

```
<Cmd name="ListProtocols"/>

<Res name="ListProtocols" ok="true">
  <Protocols>
    <Protocol>KingFisher Presto Blood 24 DW</Protocol>
    <Protocol>KingFisher Presto Cells 96</Protocol>
    <Protocol>KingFisher Presto Demo 24 DW</Protocol>
  </Protocols>
  <MemoryUsed value="35"/>
</Res>
```

4.11 RemoveProtocol

Remove a protocol from the internal memory of the KingFisher Presto instrument.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "RemoveProtocol"	Name of the command
Cmd@protocol	String: Max 100 bytes	Name of the protocol to be removed, case sensitive.

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "RemoveProtocol"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was succesfull

Details

Protocol is removed permanently from the internal memory of the instrument.

Example

```
<Cmd name="RemoveProtocol" protocol="KingFisher Presto Blood 24 DW"/>
```

```
<Res name="RemoveProtocol" ok="true"/>
```

4.12 Rotate

Rotate the turntable.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "Rotate"	Name of the command
Cmd@nest	uint16_t: 1, 2	Nest of the turntable to be positioned
Cmd@position	uint16_t: 1, 2	The position where the nest is to be rotated. Position 1 is the processeing position and position 2 is the load/unload position.

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "Rotate"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was succesfull

Details

Response to the command is replied immediately and event [Ready](#) is sent by the instrument after the rotation is completed. If rotation fails, then an [Error](#) event will be sent instead. See also command [GetStatus](#).

Example

```
<Cmd name="Rotate" nest="1" position="2" />
```

```
<Res name="Rotate" ok="true"/>
```

4.13 SetTemperatureReporting

Enable/disable temperature events with desired interval.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "SetTemperatureReporting"	Name of the command
Cmd@interval	XML Duration data type	Interval of the temperature event. Min 1 second.

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "SetTemperatureReporting"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was succesfull and the instrument is not in an error state.

Details

This command can be used to trigger [Temperature](#) events with a minimum interval of one second. If the interval is set to "PT0S" or zero, then the events are disabled.

Note The interval is fixed to five seconds in KingFisher Presto Windows simulator and it can vary depending on the the operating system load and version.

Examples

```
<Cmd name="SetTemperatureReporting" interval="PT3S"/>

<Res name="SetTemperatureReporting" ok="true" />

<Evt name="Temperature">
  <Ambient value="22.1"/>
  <Heater value="37.7"/>
</Evt>

<Evt name="Temperature">
  <Ambient value="22.2"/>
  <Heater value="37.6"/>
</Evt>
```

4.14 StartProtocol

Start executing a protocol or a step from a protocol.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "StartProtocol"	Name of the command
Cmd@protocol	String: Max 100 chars	Name of the protocol to be started, case sensitive.
Cmd@tip	String: Max 100 chars	Optional, name of a tip containing the step to be started.
Cmd@step	String: Max 100 chars	Optional, name of the step to be started.

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "StartProtocol"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was successful

Details

This command is accepted if the instrument is not already running a step or a protocol. Event [Ready](#) is sent by the instrument after the execution. If attributes "tip" and "step" are given, then a single step execution is started instead of a full protocol execution. If execution fails, then an [Error](#) event will be sent instead. See also command [GetStatus](#).

Note. The protocol needs to be uploaded to the internal memory of the instrument beforehand. See commands [List-Protocols](#) and [UploadProtocol](#).

Note. If this command is used to start single step execution, then the controlling system needs to handle turntable rotations by sending separate [Rotate](#) commands.

Event [StepStarted](#) is sent for every step in the protocol being executed.

Examples

```
<Cmd name="StartProtocol" protocol="KingFisher Presto Blood 96 DW"/>
<Res name="StartProtocol" ok="true"/>

<Cmd name="StartProtocol" protocol="KingFisher Presto Blood 96 DW" tip="Tip1" step="Mix1"/>
<Res name="StartProtocol" ok="true"/>
```

4.15 Stop

Stop execution.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "Stop"	Name of the command

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "Stop"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was successful

Details

The Stop command is used to stop an ongoing execution of protocol, step or some other process. Response to the command is replied immediately and event [Ready](#) is sent after all the motors of the instrument are stopped. In case of errors during stopping, event [Error](#) is sent.

Note that in addition to the Stop command there is also the `isUsbAbort` procedure for stopping the instrument. The `isUsbAbort` procedure differs from the normal command-reply message exchange, see `isPageUsbInterface` for more detailed information.

Example

```
<Cmd name="Stop" />
```

```
<Res name="Stop" ok="true" />
```

4.16 UploadProtocol

Transfer a protocol to the internal memory of the KingFisher Presto instrument.

Note! It's much easier to use `KFModule_UploadProtocol()` API function from `KFModule.dll` than this command directly. `KFModule_UploadProtocol` takes `BindIt .bdz` files as input parameter and handles both CRC calculation and base64 encoding.

Syntax

Tag / Attribute	Data type and range/limits	Description
Cmd@name	String: "UploadProtocol"	Name of the command
Cmd@protocol	String: Max 100 bytes	Name of the protocol to be transferred, case sensitive.
Cmd@crc	uint32_t	CRC value of the BindIt protocol file data.
Cmd[CDATA]	Base64 string/strings	XML character data section containing base64 encoded protocol file

Reply

Tag / Attribute	Data type and range/limits	Description
Res@name	String: "UploadProtocol"	Name of the command
Res@ok	Boolean string: "true" or "false"	Error status quick peek, true if command was successful

Details

Existing protocol is overwritten. Possible error codes are listed in the response message.

Note! Lines in the CDATA section must be n x 4 characters of **base64** encoded data. Except the last line.

Example

```
<Cmd name="UploadProtocol" protocol="KingFisher Presto Blood 24 DW" crc="123456" >
  <![CDATA[
    PHN0ZXA+DQpoZXJlIGlzIGFu
    IGV4YWlwbGUNCmZvcjB0aGUg
    YmFzZTY0IGVuY29kZXINCjwv
    c3RlcD4=
  ]]>
</Cmd>

<Res name="UploadProtocol" ok="true"/>
```

Chapter 5

Events

- [Aborted](#)
- [ChangeMagnets](#)
- [ChangePlate](#)
- [Error](#)
- [LoadPlate](#)
- [Pause](#)
- [ProtocolTimeLeft](#)
- [Ready](#)
- [RemovePlate](#)
- [StepStarted](#)
- [Temperature](#)

5.1 Aborted

Instrument aborted event.

Syntax

Tag / Attribute	Data type and range/limits	Description
Evt@name	String: "Aborted"	Name of the event

Details

This event is sent to a connected communication port if the KingFisher Presto is aborted for some other communication port. See command [Abort](#). More specifically, see [Abort](#) for USB, [Abort](#) for LAN and [Abort](#) for RS232.

Examples

```
<Evt name="Aborted"/>
```

5.2 ChangeMagnets

KingFisher Presto is requesting a changing of the magnet heads.

Syntax

Tag / Attribute	Data type and range/limits	Description
Evt@name	String : "ChangeMagnets"	Name of the event
Evt@tips	String: Max 100 chars	Name of the tips for the magnet heads

Details

This event is sent by the KingFisher Presto instrument when a protocol with a multiple types of magnet heads is requiring a change of the magnet heads.

Note that the changing of the magnet heads is a manual operation.

Instrument continues the execution of the protocol after the command [Acknowledge](#) is sent by the controlling system.

Example

```
<Evt name="ChangeMagnets" tips="Tip 24 DW"/>
```

5.3 ChangePlate

KingFisher Presto is requesting a plate change.

Syntax

Tag / Attribute	Data type and range/limits	Description
Evt@name	String: "ChangePlate"	Name of the event
Evt@optional	Boolean string: "true" or "false"	If true, then the plate change is possible but not required. If false, then the instrument will not continue operation until the plate is changed.
Evt/Evt[1]@name	String: "RemovePlate"	Name of the wrapped event
Evt/Evt[1]@plate	String: Max 100 chars	Plate name
Evt/Evt[1]@optional	Boolean string: "true" or "false"	See Evt/@optional
Evt/Evt[2]@name	String: "LoadPlate"	Name of the wrapped event
Evt/Evt[2]@plate	String: Max 100 chars	Plate name
Evt/Evt[2]@optional	Boolean string: "true" or "false"	See Evt/@optional

Details

The instrument sends this event to a controlling system, when a plate can be changed in the load position. Event parameter "optional" may be set to "true" to indicate that the plate change is not yet obligatory. The controlling system can perform the change operation at this point and notify the instrument using [Acknowledge](#) command. The protocol execution time can be minimized by changing the plate immediately after this first event, but the controlling system can also choose to wait when the plate change is actually needed to be done.

If the event with the "optional" parameter is neglected and not acknowledged by the controlling system, then the instrument will send another event with the parameter "optional" set to "false". Now the controlling system must change the plate and notify the instrument using [Acknowledge](#) command so that the protocol under execution can continue.

The change plate event is structured in the way, that it contains [RemovePlate](#) and [LoadPlate](#) events. Note that only one [Acknowledge](#) command is required though.

Example

```
<Evt name="ChangePlate" optional="true">
  <Evt name="RemovePlate" plate="Plate 1" optional="true"/>
  <Evt name="LoadPlate" plate="Plate 2" optional="true"/>
</Evt>

<Evt name="ChangePlate" optional="false">
  <Evt name="RemovePlate" plate="Plate 1" optional="false"/>
  <Evt name="LoadPlate" plate="Plate 2" optional="false"/>
</Evt>
```

5.4 Error

Spontaneous error message from the KingFisher Presto instrument.

Syntax

Tag / Attribute	Data type and range/limits	Description
Evt@name	String: "Error"	Name of the event
Evt/Error@code	uint16_t: see Error Codes	Error code number
Evt/Error/text()	String: Max 100 characters	Description of the error

Details

This event is sent for example after failed turntable rotation, see command [Rotate](#). The instrument will transition to an error state and an [ErrorAcknowledge](#) command must be send in order to clear the error.

Examples

```
<Evt name="Error">
  <Error code="5">Turntable position error.</Error>
</Evt>
```

5.5 LoadPlate

KingFisher Presto is requesting a plate to be inserted to the load position.

Syntax

Tag / Attribute	Data type and range/limits	Description
Evt@name	String : "LoadPlate"	Name of the event
Evt@plate	String: Max 100 chars	Plate name
Evt@optional	Boolean string: "true" or "false"	If true, then the loading of the plate is possible but not required. If false, then the instrument will not continue operation until the plate is loaded.

Details

When the instrument is ready for a new plate to be loaded during protocol execution, then it sends this event to the controlling system. Event parameter "optional" may be set to "true" to indicate that the plate can be loaded but it is not yet needed. The controlling system can perform the load process at this point and notify the instrument using [Acknowledge](#) command. The controlling system can minimize the protocol execution time by loading the plate immediately after this first event, but it can also choose to wait when the plate is actually required.

If the event with the "optional" parameter is neglected and not acknowledged by the controlling system, then the instrument will send another event with the parameter "optional" set to "false". Now the controlling system must load the plate and notify the instrument using [Acknowledge](#) command so that the protocol under execution can continue.

Example

```
<Evt name="LoadPlate" plate="Plate 1" optional="true"/>
<Evt name="LoadPlate" plate="Plate 1" optional="false"/>
```

5.6 Pause

Event from a Pause step. User is requested to perform actions to a plate.

Syntax

Tag / Attribute	Data type and range/limits	Description
Evt@name	String: "Pause"	Name of the event
Evt@plate	String: Max 100 chars	Plate name
Evt/Message[text()]	String: Max ??? chars	Message for the user

Details

Instrument must be notified using [Acknowledge](#) command so that the pause step can continue.

Example

```
<Evt name="Pause" plate="Plate 1" >  
  <Message>Dispense 10 µl of ethanol to each well.</Message>  
</Evt>
```

5.7 ProtocolTimeLeft

Spontaneous time left event from a KingFisher Presto protocol or single step execution.

Syntax

Tag / Attribute	Data type and range/limits	Description
Evt@name	String: "ProtocolTimeLeft"	Name of the event
Evt/TimeToPause@value	XML Duration data type	Time to next plate load/change/remove or pause event
Evt/TimeLeft@value	XML Duration data type	Total time left of the protocol/step execution

Details

This event is sent automatically after every [StepStarted](#) event. See command [GetProtocolTimeLeft](#) for more details.

Note! Element **TimeToPause** is not included when running single step execution. Note also that then the value of **TimeLeft** refers to the total execution time of a step being executed.

Examples

```
<Evt name="ProtocolTimeLeft">
  <TimeToPause value="PT0S"/>
  <TimeLeft value="PT30M53S"/>
</Evt>
```

```
<Evt name="ProtocolTimeLeft">
  <TimeLeft value="PT2M42S"/>
</Evt>
```

5.8 Ready

Step or protocol execution completed.

Syntax

Tag / Attribute	Data type and range/limits	Description
Evt@name	String: "Ready"	Name of the event

Details

This event is sent by the instrument after execution of a step or a protocol started with command [StartProtocol](#). See also [Rotate](#).

Examples

```
<Evt name="Ready" />
```

5.9 RemovePlate

KingFisher Presto is requesting a plate to be removed from the load position.

Syntax

Tag / Attribute	Data type and range/limits	Description
Evt@name	String: "RemovePlate"	Name of the event
Evt@plate	String: Max 100 chars	Plate name
Evt@optional	Boolean string: "true" or "false"	If true, then the removing of the plate is possible but not required. If false, then the instrument will not continue operation until the plate is removed.

Details

The instrument sends this event to a controlling system, when a plate can be removed from the load position. Event parameter "optional" may be set to "true" to indicate that the plate can be removed but it is not obligatory. The controlling system can perform the removal operation at this point and notify the instrument using [Acknowledge](#) command. The protocol execution time can be minimized by removing the plate immediately after this first event, but the controlling system can also choose to wait when the plate is actually needed to be removed.

If the event with the "optional" parameter is neglected and not acknowledged by the controlling system, then the instrument will send another event with the parameter "optional" set to "false". Now the controlling system must remove the plate and notify the instrument using [Acknowledge](#) command so that the protocol under execution can continue.

Example

```
<Evt name="RemovePlate" plate="Plate 1" optional="true"/>
```

```
<Evt name="RemovePlate" plate="Plate 1" optional="false"/>
```

5.10 StepStarted

Step start notification from the KingFisher Presto instrument.

Syntax

Tag / Attribute	Data type and range/limits	Description
Evt@name	String: "StepStarted"	Name of the event
Evt/Step@name	String	Name of the step
Evt/Step@pseudo	Boolean string: "true" or "false"	If this attribute is "false", then the step is a normal BindIt Protocol step e.q. Mix, Dry, Pause, Collect or Release. If attribute is "true" then the step is a fixed pseudo step like Init, Pick-Up, Leave and Unload.
Evt/Step@tip	String	Name of the tip group under the step is in the BindIt protocol
Evt/Step@plate	String	Name of a plate used in the step

Details

This event is sent when a protocol step is started. See also command [StartProtocol](#).

Examples

```
<Evt name="StepStarted">
  <Step name="Init" pseudo="true"/>
</Evt>

<Evt name="StepStarted">
  <Step name="Pick-Up" pseudo="true" tip="Tip1" plate="Tip Comb" />
</Evt>

<Evt name="StepStarted">
  <Step name="Mix1" pseudo="false" tip="Tip1" plate="Plate 1"/>
</Evt>

<Evt name="StepStarted">
  <Step name="Unload" pseudo="true" tip="Tip1" />
</Evt>
```

5.11 Temperature

Temperature measurements event.

Syntax

Tag / Attribute	Data type and range/limits	Description
Evt@name	String: "Temperature"	Name of the event
Evt/Ambient@value	float	Ambient temperature in Celciuss degrees
Evt/Heater@value	float	Heater block temperature in Celciuss degrees

Details

This event reports ambient and heater block temperatures. See command [SetTemperatureReporting](#).

Examples

```
<Evt name="Temperature">
  <Ambient value="22.1"/>
  <Heater value="37.7"/>
</Evt>
```

Chapter 6

Appendix - XML path syntax

Syntaxes of [Events](#), [Commands and Responses](#) are documented using a style similar to the [XPath](#) syntax. Here is a few examples of the usage. Note that also the actual syntaxes are explained using detailed examples.

Example XML block:

```
<T1>
  <T2>
    <T3 a1="123">
      This is a sentence.
    </T3>
    <T4>
      <T5 a2="321"/>
      <T5 a2="432">
      <T5 a2="543">
      <T5 a2="654">
    </T4>
  </T2>
  <![CDATA[ABCDEFGH]]>
</T1>
```

Path to the T3:

```
T1/T2/T3
```

Path to the text of the tag T3:

```
T1/T2/T3/text()
```

Path to the attribute "a1" of the tag T3:

```
T1/T2/T3/@a1
```

Path to the attribute "a2" of the third T5 tag with a value of "543"

```
T1/T2/T4/T5[3]@a2
```

Path to the CDATA of the tag T1

Note that XPath does not support addressing a CDATA element, we'll use this syntax:

```
T1[CDATA]
```

Thermo Scientific

KFModule.dll

Interface Specification

Contents

1	KFModuleDll Interface Specification	1
2	Using KFModuleDll	3
2.1	About	3
2.2	Exported functions	4
2.2.1	KFModule_OpenUsb	5
2.2.2	KFModule_OpenSerial	6
2.2.3	KFModule_OpenLan	7
2.2.4	KFModule_ListLanDevices	8
2.2.5	KFModule_OpenSimulator	9
2.2.6	KFModule_Close	10
2.2.7	KFModule_Connect	11
2.2.8	KFModule_Disconnect	12
2.2.9	KFModule_AttachEvent	13
2.2.10	KFModule_AttachMsg	14
2.2.11	KFModule_AttachCallback	15
2.2.12	KFModule_Send	16
2.2.13	KFModule_Abort	17
2.2.14	KFModule_ReadReceived	18
2.2.15	KFModule_ReadResponse	19
2.2.16	KFModule_ReadEvent	20
2.2.17	KFModule_UploadProtocol	21
2.2.18	KFModule_DownloadProtocol	22
2.2.19	KFModule_GetError	23
3	File Index	25
3.1	File List	25
4	File Documentation	27
4.1	api.h File Reference	27
4.2	KFModuleDll.c File Reference	27
4.2.1	Detailed Description	28
4.2.2	Function Documentation	28
4.2.2.1	KFModule_Abort	28
4.2.2.2	KFModule_AttachCallback	28
4.2.2.3	KFModule_AttachEvent	29
4.2.2.4	KFModule_AttachMsg	29
4.2.2.5	KFModule_Close	29
4.2.2.6	KFModule_Connect	30
4.2.2.7	KFModule_Disconnect	30
4.2.2.8	KFModule_DownloadProtocol	30
4.2.2.9	KFModule_GetError	31
4.2.2.10	KFModule_ListLanDevices	31
4.2.2.11	KFModule_OpenLan	32
4.2.2.12	KFModule_OpenSerial	32
4.2.2.13	KFModule_OpenSimulator	33

4.2.2.14	KFModule_OpenUsb	33
4.2.2.15	KFModule_ReadEvent	34
4.2.2.16	KFModule_ReadReceived	34
4.2.2.17	KFModule_ReadResponse	35
4.2.2.18	KFModule_Send	35
4.2.2.19	KFModule_UploadProtocol	36
4.3	KFModuleDll.h File Reference	36
4.3.1	Detailed Description	38
4.3.2	Macro Definition Documentation	38
4.3.2.1	KF_PRESTO_PID	38
4.3.2.2	KFMODULE_ERROR	38
4.3.2.3	KFMODULE_EVENT	38
4.3.2.4	KFMODULE_RECEIVE	38
4.3.2.5	KFMODULE_RESPONSE	38
4.3.2.6	KFMODULE_TRANSMIT	39
4.3.3	Enumeration Type Documentation	39
4.3.3.1	KFM_ERROR	39
4.3.3.2	KFM_SIMULATOR_PORT	39
4.3.4	Function Documentation	39
4.3.4.1	KFModule_Abort	39
4.3.4.2	KFModule_AttachCallback	40
4.3.4.3	KFModule_AttachEvent	40
4.3.4.4	KFModule_AttachMsg	40
4.3.4.5	KFModule_Close	41
4.3.4.6	KFModule_Connect	41
4.3.4.7	KFModule_Disconnect	41
4.3.4.8	KFModule_DownloadProtocol	42
4.3.4.9	KFModule_GetError	42
4.3.4.10	KFModule_ListLanDevices	42
4.3.4.11	KFModule_OpenLan	43
4.3.4.12	KFModule_OpenSerial	43
4.3.4.13	KFModule_OpenSimulator	44
4.3.4.14	KFModule_OpenUsb	44
4.3.4.15	KFModule_ReadEvent	45
4.3.4.16	KFModule_ReadReceived	45
4.3.4.17	KFModule_ReadResponse	46
4.3.4.18	KFModule_Send	47
4.3.4.19	KFModule_UploadProtocol	47
4.4	mainpage.h File Reference	48
Index		49

Chapter 1

KFModuleDll Interface Specification

About

The purpose of the KFModule.dll dynamic link library is to make interface to a KFModule instrument easy. For communication with the instrument the KFModule.dll uses another DLLs: ThermoUSB.dll, ThermoLAN.dll and ThermoCOM.dll. The user of the KFModule.dll does not have to know anything about the actual hardware port. Just use the functions provided by KFModule.dll.

Confidential

This document has been prepared by Thermo Fisher Scientific Oy to be used solely for the purposes defined by Thermo Fisher Scientific Oy. Use for other purposes is not authorized.

Please note that any and all information contained in this document is the property of Thermo Fisher Scientific Oy. This confidential information ("Confidential Information") shall not be reproduced in whole part or disclosed to any third party without the prior written approval of Thermo Fisher Scientific Oy. The receiving party shall ensure that its employees, officers, representatives and agents shall not disclose to third parties any Confidential Information.

Upon written request from Thermo Fisher Scientific Oy, the receiving party shall promptly return all Confidential Information or destroy all Confidential Information.

Chapter 2

Using KFModuleDll

2.1 About

Because it is not trivial to write PC software from scratch to communicate with a KFModule instrument, two Dynamic Link Libraries are provided which hide much of the complexity of the communication.

The first one is a DLL for the actual HW interface: ThermoUSB.dll, ThermoLAN.dll or ThermoCOM.dll. These DLLs are generic libraries for communicating with several different Thermo Scientific microplate instruments.

The second one is KFModule.dll, which uses DLLs mentioned above to communicate with the KFModule instrument.

To communicate with a KFModule instrument you use the exported functions of KFModule.dll. No knowledge of the other DLLs is required, they only need to be in a folder where Windows can find them (probably the same place where the KFModule.dll is loaded from).

Depending on your project setup, you may find useful a couple of other files which are also provided. The header file [KFModuleDll.h](#) contains the prototypes of the exported functions and definitions of constant values used by the dll. File KFModule.lib contains information about the dll the linker uses to add references to the library in the executable. This way the dll is automatically loaded and the exported functions of the library can be called as easy as the functions in the code using the dll.

2.2 Exported functions

- [KFModule_OpenUsb](#)
- [KFModule_OpenSerial](#)
- [KFModule_OpenLan](#)
- [KFModule_ListLanDevices](#)
- [KFModule_OpenSimulator](#)
- [KFModule_Close](#)
- [KFModule_Connect](#)
- [KFModule_Disconnect](#)
- [KFModule_AttachEvent](#)
- [KFModule_AttachMsg](#)
- [KFModule_AttachCallback](#)
- [KFModule_Send](#)
- [KFModule_Abort](#)
- [KFModule_ReadReceived](#)
- [KFModule_ReadResponse](#)
- [KFModule_ReadEvent](#)
- [KFModule_UploadProtocol](#)
- [KFModule_DownloadProtocol](#)
- [KFModule_GetError](#)

2.2.1 KFModule_OpenUsb

Open an USB communication channel to a KFModule instrument. Full declaration: [KFModule_OpenUsb\(\)](#).

Parameters

<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first device with matching VendorID and ProductID.
<i>productId</i>	Manufacturer product id number of the USB device.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the KFModule_OpenXxx() functions must be called first before using any other functions in the library. Only one connection per device is allowed.



Note. Function [KFModule_Connect\(\)](#) must be called after this function in order to use the XML interface to communicate with the instrument.

2.2.2 KFModule_OpenSerial

Open serial communication port to a KFModule instrument. Full declaration: [KFModule_OpenSerial\(\)](#).

Parameters

<i>port</i>	Serial port number.
<i>baud</i>	Serial port baudrate.
<i>DeviceName</i>	Pointer to the device name string of the device. The device must report an identical name to the VER command for the connection to succeed. This parameter may be NULL, in which case the device name is ignored.
<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number to the VER command for the connection to succeed. This parameter may be NULL, in which case the serial number is ignored.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the KFModule_OpenXxx() functions must be called first before using any other functions in the library. Only one connection per serial port is allowed. If both DeviceName and SerialNumber are NULL, no VER command is sent to the instrument.



Note. Function [KFModule_Connect\(\)](#) must be called after this function in order to use the XML interface to communicate with the instrument.

2.2.3 KFModule_OpenLan

Open a LAN communication channel to a KFModule instrument. Full declaration: [KFModule_OpenLan\(\)](#).

Parameters

<i>instrumentName</i>	Name of the instrument. This must match the device name the instrument uses for matching a WS-Discovery Probe and Resolve.
<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first found KFModule instrument.
<i>timeout</i>	Timeout which is used to search the instrument. If 0 is given then WS-Discovery will use default 4 sec timeout.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the KFModule_OpenXxx() functions must be called first before using any other functions in the library. Only one connection per device is allowed.



Note. Function [KFModule_Connect\(\)](#) must be called after this function in order to use the XML interface to communicate with the instrument.

2.2.4 KFModule_ListLanDevices

List instruments found on the LAN. Full declaration: [KFModule_ListLanDevices\(\)](#).

Parameters

<i>instrumentName</i>	Name of the instrument. Only instruments with a matching name are listed. May be NULL, in which case the found devices are not filtered by the name.
<i>SerialNumber</i>	The serial number of the instrument. Only instruments with a matching serial number are listed. May be NULL, in which case the found devices are not filtered by the serial number.
<i>buf</i>	The buffer to list the found devices to.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

The size in bytes of the complete list of found devices. If the returned value is equal or smaller than the given buffer size, the buffer contains the complete list of devices found on the LAN. If the returned value is higher than the given buffer size, no data is returned and the caller must call the function again with big enough buffer. Return value 0 means that no instruments were found.

On success, the caller's buffer contains zero terminated strings with a combined length of the return value. The string terminating zeros are included in the length.

For each found instrument, the first string is the instrument IPv4 address and the TCP port number it is listening, enclosed in square brackets, e.g. [10.32.196.210:49536].

The IP address string is followed by the WS-Discovery match strings, usually 3 of them. The first one is always 'Thermo-Device', the second one is the instrument name and the third one the instrument serial number string.

If the match strings are followed by a string in angle brackets, e.g. <10.32.196.154:57403>, it means that the instrument is currently connected that IP address and TCP port, and trying to connect to that instrument with function [KFModule_OpenLan\(\)](#) will fail. If there is no string in square brackets, the instrument will accept a connection.

2.2.5 KFModule_OpenSimulator

Open communication channel to KFModule simulator. Full declaration: [KFModule_OpenSimulator\(\)](#).

Parameters

<i>port</i>	The simulator port to connect to, one of KFM_SIMULATOR_USB , KFM_SIMULATOR_COM , KFM_SIMULATOR_DBG or KFM_SIMULATOR_LAN .
<i>productId</i>	(USB) product id of the instrument.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the [KFModule_OpenXxx\(\)](#) functions must be called first before using any other functions in the library. Only one connection per simulator communication channel is allowed.



Note. Function [KFModule_Connect\(\)](#) must be called after this function in order to use the XML interface to communicate with the instrument.

2.2.6 KFModule_Close

Close a communication channel. Full declaration: [KFModule_Close\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

This function should be called when the communication channel is no longer needed. Failing to close the channel prevents new connections to the channel as long as the dll stays loaded.



Note. When using the XML interface, function [KFModule_Disconnect\(\)](#) must be called before this function in order to disconnect the instrument from the open communication channel. Otherwise the instrument may refuse new connections.

2.2.7 KFModule_Connect

Connect to the instrument. Full declaration: [KFModule_Connect\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

This function sends the "Connect" command with a current date/time setting to the instrument. It is recommended to use this function to connect to the instrument so that the calendar of the instrument is always set to the local time automatically.



Note. This function returns before the actual connection is done, meaning that the return value indicates merely if the starting of the command is successful or not. The instrument will send a separate response after the connection is finished. See interface specification of the instrument in question for more details.

2.2.8 KFModule_Disconnect

Disconnect the instrument. Full declaration: [KFModule_Disconnect\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

This function sends the "disconnect" command to the instrument. This is only needed because of the symmetry in the list of API functions. Disconnect() is the pair to Connect() function.



Note. This function returns before the actual connection is closed, meaning that the return value indicates merely if the starting of the command is successful or not. The instrument will send a separate response just before the disconnection. See interface specification of the instrument in question for more details.

2.2.9 KFModule_AttachEvent

Attach an event object to a KFModule connection. Full declaration: [KFModule_AttachEvent\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>ev</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

Parameter 'ev' may be any combination of [KFMODULE_RECEIVE](#), [KFMODULE_TRANSMIT](#), [KFMODULE_RESPONSE](#), [KFMODULE_EVENT](#) and [KFMODULE_ERROR](#). The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event. If the same event object is used for more than one event, the application must check for all possible events after the event object is signalled.

2.2.10 KFModule_AttachMsg

Attach an event message to the KFModule connection. Full declaration: [KFModule_AttachMsg\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>hWnd</i>	Handle of the window to receive the message.
<i>msg</i>	Message id of the message to send.
<i>evCode</i>	Event(s) for which a message is sent.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

A message with message id 'msg' is sent to window 'hWnd' whenever any of the event(s) selected with parameter 'evCode' occurs. Parameter 'evCode' may be any combination of [KFMODULE_RECEIVE](#), [KFMODULE_TRANSMIT](#), [KFMODULE_RESPONSE](#), [KFMODULE_EVENT](#) and [KFMODULE_ERROR](#). The wParam of the sent message will be one of these event codes.

2.2.11 KFModule_AttachCallback

Attach an event callback function to the KFModule connection. Full declaration: [KFModule_AttachCallback\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>callback</i>	Address of the callback function.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

The callback function will be called whenever any of events [KFMODULE_RECEIVE](#), [KFMODULE_TRANSMIT](#), [KFMODULE_RESPONSE](#), [KFMODULE_EVENT](#) or [KFMODULE_ERROR](#) occurs. The connection handle and the event code are passed as parameters to the function.

2.2.12 KFModule_Send

Send a NUL terminated ASCII string to the instrument. Full declaration: [KFModule_Send\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>buf</i>	The NUL terminated string to send.

Returns

One of [KFM_ERROR](#) codes, [KFM_ERROR_SUCCESS](#) on success.

Use this function to send commands and acknowledges to the instrument. For uploading and downloading a protocol there are dedicated functions [KFModule_UploadProtocol\(\)](#) and [KFModule_DownloadProtocol\(\)](#). Also connecting/disconnecting is recommended to be done with separate [KFModule_Connect\(\)](#) and [KFModule_Disconnect\(\)](#) functions.

The data to be sent to the instrument is buffered in the dll. A [KFMODULE_TRANSMIT](#) event is sent to the application when all data is sent. You may call this function repeatedly without waiting for the event, but a [KFM_ERROR_OUT_OF_HEAP](#) error may be returned if you send a lot of data without waiting for the event.



Note. Remember to terminate commands with a new line character (ASCII 10). See interface specification of the instrument in question for more details.

2.2.13 KFModule_Abort

Send Abort command to the instrument. Full declaration: [KFModule_Abort\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

Aborts any command(s) being executed in the instrument.

2.2.14 KFModule_ReadReceived

Read a received data line from the received data chain. Full declaration: [KFModule_ReadReceived\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>buf</i>	Caller's buffer for the line.
<i>bufsize</i>	Size of the caller's buffer.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS if any data was copied to caller's buffer.

Call this function in response to the [KFMODULE_RECEIVE](#) event. The received data is neither a XML Response nor a XML Event. An empty string will be copied to caller's buffer when there is no more data to read. The returned string is always NUL terminated and does not contain the CR/LF characters sent by the instrument.



Note. Always read all data from the KFModuleDll when handling the event. Otherwise subsequent strings from the instrument may be missed. This can happen especially when the user application doesn't react to the events immediately. See example below:

```
void handler_KFMODULE_RECEIVE( HANDLE connection )
{
    KFM_ERROR    err;
    char         buf[100];

    while( (err = KFModule_ReadReceived( connection, buf, sizeof( buf ))) == KFM_ERROR_SUCCESS)
    {
        // Do something with the data
    }

    if( err > KFM_ERROR_NO_DATA)
    {
        // Handle the error
    }
}
```

2.2.15 KFModule_ReadResponse

Read a received XML response from the response chain. Full declaration: [KFModule_ReadResponse\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>buf</i>	Caller's buffer for the response.
<i>bufsize</i>	Size of the caller's buffer.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS if any data was copied to caller's buffer.

Call this function in response to the [KFMODULE_RESPONSE](#) event. The received data is a XML Response. The event is not sent until a whole XML Response is received. If the caller's buffer is not large enough to hold the whole response, a partial response is copied. An empty string is copied when there is no more data to read. The returned string is always NUL terminated and does not contain the CR/LF characters sent by the instrument.



Note. Always read all data from the KFModuleDll when handling the event. Otherwise subsequent responses may be missed. This can happen especially when the user application doesn't react to the events immediately. See example below:

```
void handler_KFMODULE_RESPONSE( HANDLE connection )
{
    KFM_ERROR    err;
    char         buf[100];

    while( (err = KFModule_ReadResponse( connection, buf, sizeof( buf ) )) == KFM_ERROR_SUCCESS)
    {
        // Do something with the data
    }

    if( err > KFM_ERROR_NO_DATA)
    {
        // Handle the error
    }
}
```

2.2.16 KFModule_ReadEvent

Read a received XML event from the event chain. Full declaration: [KFModule_ReadEvent\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>buf</i>	Caller's buffer for the event.
<i>bufsize</i>	Size of the caller's buffer.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS if any data was copied to caller's buffer.

Call this function in response to the [KFMODULE_EVENT](#) event. The received data is a XML Event. The event is not sent until a whole XML Event is received. If the caller's buffer is not large enough to hold the whole Event, a partial Event is copied. An empty string is copied when there is no more data to read. The returned string is always NUL terminated and does not contain the CR/LF characters sent by the instrument.



Note. Always read all data from the KFModuleDll when handling the event. Otherwise subsequent events from the instrument may be missed. This can happen especially when the user application doesn't react to the events immediately. See example below:

```
void handler_KFMODULE_EVENT( HANDLE connection )
{
    KFM_ERROR    err;
    char         buf[100];

    while( (err = KFModule_ReadEvent( connection, buf, sizeof( buf ) )) == KFM_ERROR_SUCCESS)
    {
        // Do something with the data
    }

    if( err > KFM_ERROR_NO_DATA)
    {
        // Handle the error
    }
}
```

2.2.17 KFModule_UploadProtocol

Export the requested protocol to the instrument. Full declaration: [KFModule_UploadProtocol\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>protocolName</i>	Name of the protocol to export.
<i>path</i>	Complete path of the .bdz file containing the protocol to export.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

The given protocol name must match exactly with a protocol name stored in the .bdz file. The .bdz file may contain several protocols, but only the requested protocol is sent to the instrument.



Note. This function returns before the actual transfer is completed, meaning that the return value indicates merely if the starting of the transfer is successful or not. The instrument will send a separate response after the transfer. See interface specification of the instrument in question for more details.

2.2.18 KFModule_DownloadProtocol

Import the requested protocol to a .bdz file. Full declaration: [KFModule_DownloadProtocol\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>protocolName</i>	Name of the protocol to export.
<i>path</i>	Complete path of the .bdz file to receive the protocol.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

The given protocol name must match exactly with a protocol name stored in the instrument. The dll must have write access to the given .bdz file. The file will be overwritten by the dll.



Note. This function returns before the actual transfer is completed, meaning that the return value indicates merely if the starting of the transfer is successful or not. The instrument will send a separate response after the transfer. See interface specification of the instrument in question for more details.

2.2.19 KFModule_GetError

Get the asynchronous error code. Full declaration: [KFModule_GetError\(\)](#).

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

Most dll functions return an error code, but an error may occur asynchronously in the receive or transmit thread of the dll. This kind of errors are reported with the KFMODULE_ERROR event. The application should then call this function to read the error code and take action. The error usually is KFM_ERROR_DISCONNECTED, in which case the application should close the current connection and try to open a new one.

Chapter 3

File Index

3.1 File List

Here is a list of all files with brief descriptions:

api.h	Part of KFModuleDll documentation	27
KFModuleDll.c	Interface to a Thermo Fisher Scientific KingFisher instrument that contains KingFisher module and uses ThermoUSB.dll, ThermoLAN.dll or ThermoCOM.dll	27
KFModuleDll.h	Functions exported from KFModule.dll	36
mainpage.h	KFModuleDll Interface Specification	48

Chapter 4

File Documentation

4.1 api.h File Reference

Part of KFModuleDll documentation.

4.2 KFModuleDll.c File Reference

Interface to a Thermo Fisher Scientific KingFisher instrument that contains KingFisher module and uses ThermoUSB.dll, ThermoLAN.dll or ThermoCOM.dll.

Functions

- HANDLE WINAPI [KFModule_OpenUsb](#) (LPCSTR SerialNumber, WORD productId)
Open an USB communication channel to a KFModule instrument.
- HANDLE WINAPI [KFModule_OpenSerial](#) (WORD port, DWORD baud, LPCSTR DeviceName, LPCSTR SerialNumber)
Open serial communication port to a KFModule instrument.
- HANDLE WINAPI [KFModule_OpenLan](#) (LPCSTR DeviceName, LPCSTR SerialNumber, UINT timeout)
Open a LAN communication channel to a KFModule instrument.
- DWORD WINAPI [KFModule_ListLanDevices](#) (LPCSTR DeviceName, LPCSTR SerialNumber, LPSTR buf, DWORD bufsize)
List instruments found on the LAN.
- HANDLE WINAPI [KFModule_OpenSimulator](#) (KFM_SIMULATOR_PORT port, WORD productId)
Open communication channel to KFModule simulator.
- [KFM_ERROR](#) WINAPI [KFModule_Close](#) (HANDLE hConn)
Close a communication channel.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_Connect](#) (HANDLE hConn)
Connect to the instrument.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_Disconnect](#) (HANDLE hConn)
Disconnect the instrument.
- [KFM_ERROR](#) WINAPI [KFModule_ReadReceived](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read a received data line from the received data chain.
- [KFM_ERROR](#) WINAPI [KFModule_ReadResponse](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read a received XML response from the response chain.
- [KFM_ERROR](#) WINAPI [KFModule_ReadEvent](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)

Read a received XML event from the event chain.

- **KFM_ERROR** WINAPI [KFModule_AttachEvent](#) (HANDLE hConn, UINT ev, HANDLE object)

Attach an event object to a KFModule connection.

- **KFM_ERROR** WINAPI [KFModule_AttachMsg](#) (HANDLE hConn, HANDLE hWnd, UINT msg, UINT evCode)

Attach an event message to the KFModule connection.

- **KFM_ERROR** WINAPI [KFModule_AttachCallback](#) (HANDLE hConn, void(*callback)(HANDLE conn, UINT ev))

Attach an event callback function to the KFModule connection.

- **KFM_ERROR** WINAPI [KFModule_Send](#) (HANDLE hConn, LPCSTR buf)

Send a NUL terminated ASCII string to the instrument.

- **KFM_ERROR** WINAPI [KFModule_Abort](#) (HANDLE hConn)

Send Abort command to the instrument.

- **KFM_ERROR** WINAPI [KFModule_UploadProtocol](#) (HANDLE hConn, LPCSTR protocolName, LPCSTR path)

Export the requested protocol to the instrument.

- **KFM_ERROR** WINAPI [KFModule_DownloadProtocol](#) (HANDLE hConn, LPCSTR protocolName, LPCSTR path)

Import the requested protocol to a .bdz file.

- **KFM_ERROR** WINAPI [KFModule_GetError](#) (HANDLE hConn)

Get the asynchronous error code.

4.2.1 Detailed Description

Note

Copyright by Thermo Fisher Scientific Oy 2015

Definition in file [KFModuleDll.c](#).

4.2.2 Function Documentation

4.2.2.1 KFM_ERROR WINAPI KFModule_Abort (HANDLE hConn)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, [KFM_ERROR_SUCCESS](#) on success.

Aborts any command(s) being executed in the instrument.

Definition at line 2553 of file [KFModuleDll.c](#).

4.2.2.2 KFM_ERROR WINAPI KFModule_AttachCallback (HANDLE hConn, void(*) (HANDLE conn, UINT ev) callback)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>callback</i>	Address of the callback function.

Returns

One of [KFM_ERROR](#) codes, [KFM_ERROR_SUCCESS](#) on success.

The callback function will be called whenever any of events [KFMODULE_RECEIVE](#), [KFMODULE_TRANSMIT](#), [KFMODULE_RESPONSE](#), [KFMODULE_EVENT](#) or [KFMODULE_ERROR](#) occurs. The connection handle and the event code are

passed as parameters to the function.

Definition at line 2451 of file KFModuleDll.c.

4.2.2.3 KFM_ERROR WINAPI KFModule.AttachEvent (HANDLE *hConn*, UINT *ev*, HANDLE *object*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>ev</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

Parameter 'ev' may be any combination of [KFMODULE_RECEIVE](#), [KFMODULE_TRANSMIT](#), [KFMODULE_RESPONSE](#), [KFMODULE_EVENT](#) and [KFMODULE_ERROR](#). The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event. If the same event object is used for more than one event, the application must check for all possible events after the event object is signaled.

Definition at line 2362 of file KFModuleDll.c.

4.2.2.4 KFM_ERROR WINAPI KFModule.AttachMsg (HANDLE *hConn*, HANDLE *hWnd*, UINT *msg*, UINT *evCode*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>hWnd</i>	Handle of the window to receive the message.
<i>msg</i>	Message id of the message to send.
<i>evCode</i>	Event(s) for which a message is sent.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

A message with message id 'msg' is sent to window 'hWnd' whenever any of the event(s) selected with parameter 'evCode' occurs. Parameter 'evCode' may be any combination of [KFMODULE_RECEIVE](#), [KFMODULE_TRANSMIT](#), [KFMODULE_RESPONSE](#), [KFMODULE_EVENT](#) and [KFMODULE_ERROR](#). The wParam of the sent message will be one of these event codes.

Definition at line 2417 of file KFModuleDll.c.

4.2.2.5 KFM_ERROR WINAPI KFModule.Close (HANDLE *hConn*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

This function should be called when the communication channel is no longer needed. Failing to close the channel prevents new connections to the channel as long as the dll stays loaded.



Note. When using the XML interface, function [KFModule_Disconnect\(\)](#) must be called before this function in order to disconnect the instrument from the open communication channel. Otherwise the instrument may refuse new connections.

Definition at line 2099 of file KFModuleDll.c.

4.2.2.6 DllExport KFM_ERROR WINAPI KFModule.Connect (HANDLE *hConn*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

This function sends the "Connect" command with a current date/time setting to the instrument. It is recommended to use this function to connect to the instrument so that the calendar of the instrument is always set to the local time automatically.



Note. This function returns before the actual connection is done, meaning that the return value indicates merely if the starting of the command is successful or not. The instrument will send a separate response after the connection is finished. See interface specification of the instrument in question for more details.

Definition at line 2133 of file KFModuleDll.c.

4.2.2.7 DllExport KFM_ERROR WINAPI KFModule.Disconnect (HANDLE *hConn*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

This function sends the "disconnect" command to the instrument. This is only needed because of the symmetry in the list of API functions. Disconnect() is the pair to Connect() function.



Note. This function returns before the actual connection is closed, meaning that the return value indicates merely if the starting of the command is successful or not. The instrument will send a separate response just before the disconnection. See interface specification of the instrument in question for more details.

Definition at line 2178 of file KFModuleDll.c.

4.2.2.8 KFM_ERROR WINAPI KFModule.DownloadProtocol (HANDLE *hConn*, LPCSTR *protocolName*, LPCSTR *path*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>protocolName</i>	Name of the protocol to export.
<i>path</i>	Complete path of the .bdz file to receive the protocol.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

The given protocol name must match exactly with a protocol name stored in the instrument. The dll must have write access to the given .bdz file. The file will be overwritten by the dll.



Note. This funtion returns before the actual transfer is completed, meaning that the return value indicates merely if the starting of the transfer is succesfull of not. The instrument will send a separate response after the transfer. See interface specification of the instrument in question for more details.

Definition at line 2627 of file KFMModuleDll.c.

4.2.2.9 KFM_ERROR WINAPI KFMModule_GetError (HANDLE *hConn*)

Parameters

<i>hConn</i>	A handle returned by one of the KFMModule_OpenXxx() functions.
--------------	--

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

Most dll functions return an error code, but an error may occur asynchronously in the receive or transmit thread of the dll. This kind of errors are reported with the KFMODULE_ERROR event. The application should then call this function to read the error code and take action. The error usually is KFM_ERROR_DISCONNECTED, in which case the application should close the current connection and try to open a new one.

Definition at line 2673 of file KFMModuleDll.c.

4.2.2.10 DWORD WINAPI KFMModule_ListLanDevices (LPCSTR *DeviceName*, LPCSTR *SerialNumber*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>instrumentName</i>	Name of the instrument. Only instruments with a matching name are listed. May be NULL, in which case the found devices are not filtered by the name.
<i>SerialNumber</i>	The serial number of the instrument. Only instruments with a matching serial number are listed. May be NULL, in which case the found devices are not filtered by the serial number.
<i>buf</i>	The buffer to list the found devices to.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

The size in bytes of the complete list of found devices. If the returned value is equal or smaller than the given buffer size, the buffer contains the complete list of devices found on the LAN. If the returned value is higher than the given buffer size, no data is returned and the caller must call the function again with big enough buffer. Return value 0 means that no instruments were found.

On success, the caller's buffer contains zero terminated strings with a combined length of the return value. The string terminating zeros are included in the length.

For each found instrument, the first string is the instrument IPv4 address and the TCP port number it is listening, enclosed in square brackets, e.g. [10.32.196.210:49536].

The IP address string is followed by the WS-Discovery match strings, usually 3 of them. The first one is always 'Thermo-Device', the second one is the instrument name and the third one the instrument serial number string.

If the match strings are followed by a string in angle brackets, e.g. <10.32.196.154:57403>, it means that the instrument is currently connected that IP address and TCP port, and trying to connect to that instrument with function [KFModule_OpenLan\(\)](#) will fail. If there is no string in square brackets, the instrument will accept a connection.

Definition at line 2000 of file KFModuleDll.c.

4.2.2.11 HANDLE WINAPI KFModule.OpenLan (LPCSTR DeviceName, LPCSTR SerialNumber, UINT timeout)

Parameters

<i>instrumentName</i>	Name of the instrument. This must match the device name the instrument uses for matching a WS-Discovery Probe and Resolve.
<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first found KFModule instrument.
<i>timeout</i>	Timeout which is used to search the instrument. If 0 is given then WS-Discovery will use default 4 sec timeout.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the KFModule_OpenXxx() functions must be called first before using any other functions in the library. Only one connection per device is allowed.



Note. Function [KFModule_Connect\(\)](#) must be called after this function in order to use the XML interface to communicate with the instrument.

Definition at line 1935 of file KFModuleDll.c.

4.2.2.12 HANDLE WINAPI KFModule.OpenSerial (WORD port, DWORD baud, LPCSTR DeviceName, LPCSTR SerialNumber)

Parameters

<i>port</i>	Serial port number.
<i>baud</i>	Serial port baudrate.
<i>DeviceName</i>	Pointer to the device name string of the device. The device must report an identical name to the VER command for the connection to succeed. This parameter may be NULL, in which case the device name is ignored.
<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number to the VER command for the connection to succeed. This parameter may be NULL, in which case the serial number is ignored.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the KFModule_OpenXxx() functions must be called first before using any other functions in the library. Only one connection per serial port is allowed. If both DeviceName and SerialNumber are NULL, no VER command is sent to the instrument.



Note. Function [KFModule_Connect\(\)](#) must be called after this function in order to use the XML interface to communicate with the instrument.

Definition at line 1880 of file KFModuleDll.c.

4.2.2.13 HANDLE WINAPI KFModule.OpenSimulator (KFM_SIMULATOR_PORT *port*, WORD *productId*)

Parameters

<i>port</i>	The simulator port to connect to, one of KFM_SIMULATOR_USB , KFM_SIMULATOR_COM , KFM_SIMULATOR_DBG or KFM_SIMULATOR_LAN .
<i>productId</i>	(USB) product id of the instrument.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the KFModule_OpenXxx() functions must be called first before using any other functions in the library. Only one connection per simulator communication channel is allowed.



Note. Function [KFModule_Connect\(\)](#) must be called after this function in order to use the XML interface to communicate with the instrument.

Definition at line 2030 of file KFModuleDll.c.

4.2.2.14 HANDLE WINAPI KFModule.OpenUsb (LPCSTR *SerialNumber*, WORD *productId*)

Parameters

<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first device with matching VendorID and ProductID.
<i>productId</i>	Manufacturer product id number of the USB device.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the KFModule_OpenXxx() functions must be called first before using any other functions in the library. Only one connection per device is allowed.



Note. Function [KFModule_Connect\(\)](#) must be called after this function in order to use the XML interface to communicate with the instrument.

Definition at line 1827 of file KFModuleDll.c.

4.2.2.15 KFM_ERROR WINAPI KFMModule_ReadEvent (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hConn</i>	A handle returned by one of the KFMModule_OpenXxx() functions.
<i>buf</i>	Caller's buffer for the event.
<i>bufsize</i>	Size of the caller's buffer.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS if any data was copied to caller's buffer.

Call this function in response to the [KFMODULE_EVENT](#) event. The received data is a XML Event. The event is not sent until a whole XML Event is received. If the caller's buffer is not large enough to hold the whole Event, a partial Event is copied. An empty string is copied when there is no more data to read. The returned string is always NUL terminated and does not contain the CR/LF characters sent by the instrument.



Note. Always read all data from the KFMModuleDll when handling the event. Otherwise subsequent events from the instrument may be missed. This can happen especially when the user application doesn't react to the events immediately. See example below:

```
void handler_KFMODULE_EVENT( HANDLE connection )
{
    KFM_ERROR    err;
    char         buf[100];

    while( (err = KFMModule_ReadEvent( connection, buf, sizeof( buf ))) == KFM_ERROR_SUCCESS)
    {
        // Do something with the data
    }

    if( err > KFM_ERROR_NO_DATA)
    {
        // Handle the error
    }
}
```

Definition at line 2333 of file KFMModuleDll.c.

4.2.2.16 KFM_ERROR WINAPI KFMModule_ReadReceived (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hConn</i>	A handle returned by one of the KFMModule_OpenXxx() functions.
<i>buf</i>	Caller's buffer for the line.
<i>bufsize</i>	Size of the caller's buffer.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS if any data was copied to caller's buffer.

Call this function in response to the [KFMODULE_RECEIVE](#) event. The received data is neither a XML Response nor a XML Event. An empty string will be copied to caller's buffer when there is no more data to read. The returned string is always NUL terminated and does not contain the CR/LF characters sent by the instrument.



Note. Always read all data from the KFMModuleDll when handling the event. Otherwise subsequent strings from the instrument may be missed. This can happen especially when the user application doesn't react to the events immediately. See example below:

```

void handler_KFMODULE_RECEIVE( HANDLE connection )
{
    KFM_ERROR    err;
    char         buf[100];

    while( (err = KFModule_ReadReceived( connection, buf, sizeof( buf ))) == KFM_ERROR_SUCCESS)
    {
        // Do something with the data
    }

    if( err > KFM_ERROR_NO_DATA)
    {
        // Handle the error
    }
}

```

Definition at line 2239 of file KFModuleDll.c.

4.2.2.17 KFM_ERROR WINAPI KFModule_ReadResponse (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>buf</i>	Caller's buffer for the response.
<i>bufsize</i>	Size of the caller's buffer.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS if any data was copied to caller's buffer.

Call this function in response to the [KFMODULE_RESPONSE](#) event. The received data is a XML Response. The event is not sent until a whole XML Response is received. If the caller's buffer is not large enough to hold the whole response, a partial response is copied. An empty string is copied when there is no more data to read. The returned string is always NUL terminated and does not contain the CR/LF characters sent by the instrument.



Note. Always read all data from the KFModuleDll when handling the event. Otherwise subsequent responses may be missed. This can happen especially when the user application doesn't react to the events immediately. See example below:

```

void handler_KFMODULE_RESPONSE( HANDLE connection )
{
    KFM_ERROR    err;
    char         buf[100];

    while( (err = KFModule_ReadResponse( connection, buf, sizeof( buf ))) == KFM_ERROR_SUCCESS)
    {
        // Do something with the data
    }

    if( err > KFM_ERROR_NO_DATA)
    {
        // Handle the error
    }
}

```

Definition at line 2286 of file KFModuleDll.c.

4.2.2.18 KFM_ERROR WINAPI KFModule_Send (HANDLE *hConn*, LPCSTR *buf*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>buf</i>	The NUL terminated string to send.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

Use this function to send commands and acknowledges to the instrument. For uploading and downloading a protocol there are dedicated functions [KFModule_UploadProtocol\(\)](#) and [KFModule_DownloadProtocol\(\)](#). Also connecting/disconnecting is recommended to be done with separate [KFModule_Connect\(\)](#) and [KFModule_Disconnect\(\)](#) functions.

The data to be sent to the instrument is buffered in the dll. A [KFMODULE_TRANSMIT](#) event is sent to the application when all data is sent. You may call this function repeatedly without waiting for the event, but a [KFM_ERROR_OUT_OF_HEAP](#) error may be returned if you send a lot of data without waiting for the event.



Note. Remember to terminate commands with a new line character (ASCII 10). See interface specification of the instrument in question for more details.

Definition at line 2495 of file KFModuleDll.c.

4.2.2.19 KFM_ERROR WINAPI KFModule_UploadProtocol (HANDLE *hConn*, LPCSTR *protocolName*, LPCSTR *path*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>protocolName</i>	Name of the protocol to export.
<i>path</i>	Complete path of the .bdz file containing the protocol to export.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

The given protocol name must match exactly with a protocol name stored in the .bdz file. The .bdz file may contain several protocols, but only the requested protocol is sent to the instrument.



Note. This function returns before the actual transfer is completed, meaning that the return value indicates merely if the starting of the transfer is successful or not. The instrument will send a separate response after the transfer. See interface specification of the instrument in question for more details.

Definition at line 2600 of file KFModuleDll.c.

4.3 KFModuleDll.h File Reference

Functions exported from KFModule.dll.

Macros

- #define [KFMODULE_RECEIVE](#) 1
Data line received.
- #define [KFMODULE_TRANSMIT](#) 2
All data transmitted.
- #define [KFMODULE_RESPONSE](#) 4
A response received from KFModule.

- #define [KFMODULE_EVENT](#) 8
An event received from KFModule.
- #define [KFMODULE_ERROR](#) 16
Fatal error.
- #define [KF_PRESTO_PID](#) 713
Thermo Fisher Scientific KingFisher Presto product id.

Enumerations

- enum [KFM_ERROR](#) {
[KFM_ERROR_SUCCESS](#), [KFM_ERROR_NO_DATA](#), [KFM_ERROR_INVALID_HANDLE](#), [KFM_ERROR_INVALID_ARGUMENT](#),
[KFM_ERROR_DISCONNECTED](#), [KFM_ERROR_FILE_NOT_FOUND](#), [KFM_ERROR_INVALID_BDZ_FILE](#), [KFM_ERROR_PROTOCOL_NOT_FOUND](#),
[KFM_ERROR_TEMP_FILE_CREATE](#), [KFM_ERROR_EXPORT_IMPORT_BUSY](#), [KFM_ERROR_OUT_OF_HEAP](#)
}

Error codes returned by KFModuleDll functions.
- enum [KFM_SIMULATOR_PORT](#) { [KFM_SIMULATOR_USB](#), [KFM_SIMULATOR_COM](#), [KFM_SIMULATOR_DBG](#), [KFM_SIMULATOR_LAN](#) }
KFModule simulator ports.

Functions

- DllExport HANDLE WINAPI [KFModule_OpenUsb](#) (LPCSTR SerialNumber, WORD productId)
Open an USB communication channel to a KFModule instrument.
- DllExport HANDLE WINAPI [KFModule_OpenSerial](#) (WORD port, DWORD baud, LPCSTR DeviceName, LPCSTR SerialNumber)
Open serial communication port to a KFModule instrument.
- DllExport HANDLE WINAPI [KFModule_OpenLan](#) (LPCSTR DeviceName, LPCSTR SerialNumber, UINT timeout)
Open a LAN communication channel to a KFModule instrument.
- DllExport DWORD WINAPI [KFModule_ListLanDevices](#) (LPCSTR DeviceName, LPCSTR SerialNumber, LPSTR buf, DWORD bufsize)
List instruments found on the LAN.
- DllExport HANDLE WINAPI [KFModule_OpenSimulator](#) ([KFM_SIMULATOR_PORT](#) port, WORD productId)
Open communication channel to KFModule simulator.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_Close](#) (HANDLE hConn)
Close a communication channel.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_Connect](#) (HANDLE hConn)
Connect to the instrument.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_Disconnect](#) (HANDLE hConn)
Disconnect the instrument.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_AttachEvent](#) (HANDLE hConn, UINT ev, HANDLE object)
Attach an event object to a KFModule connection.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_AttachMsg](#) (HANDLE hConn, HANDLE hWnd, UINT msg, UINT ev)
Attach an event message to the KFModule connection.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_AttachCallback](#) (HANDLE hConn, void(*callback)(HANDLE conn, UINT ev))
Attach an event callback function to the KFModule connection.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_Send](#) (HANDLE hConn, LPCSTR buf)
Send a NUL terminated ASCII string to the instrument.

- DllExport [KFM_ERROR](#) WINAPI [KFModule_Abort](#) (HANDLE hConn)
Send Abort command to the instrument.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_ReadReceived](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read a received data line from the received data chain.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_ReadResponse](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read a received XML response from the response chain.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_ReadEvent](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read a received XML event from the event chain.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_UploadProtocol](#) (HANDLE hConn, LPCSTR protocolName, LPCSTR path)
Export the requested protocol to the instrument.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_DownloadProtocol](#) (HANDLE hConn, LPCSTR protocolName, LPCSTR path)
Import the requested protocol to a .bdz file.
- DllExport [KFM_ERROR](#) WINAPI [KFModule_GetError](#) (HANDLE hConn)
Get the asynchronous error code.

4.3.1 Detailed Description

Note

Copyright by Thermo Fisher Scientific Oy 2016

Interface to a KingFisher module. Following instruments uses this interface:

- KingFisher Presto

Definition in file [KFModuleDll.h](#).

4.3.2 Macro Definition Documentation

4.3.2.1 #define KF_PRESTO_PID 713

Definition at line 65 of file [KFModuleDll.h](#).

4.3.2.2 #define KFMODULE_ERROR 16

Definition at line 28 of file [KFModuleDll.h](#).

4.3.2.3 #define KFMODULE_EVENT 8

Definition at line 27 of file [KFModuleDll.h](#).

4.3.2.4 #define KFMODULE_RECEIVE 1

Definition at line 24 of file [KFModuleDll.h](#).

4.3.2.5 #define KFMODULE_RESPONSE 4

Definition at line 26 of file [KFModuleDll.h](#).

4.3.2.6 #define KFMODULE_TRANSMIT 2

Definition at line 25 of file KFMModuleDII.h.

4.3.3 Enumeration Type Documentation

4.3.3.1 enum KFM_ERROR

These are the error codes returned by the API functions.

Enumerator

KFM_ERROR_SUCCESS No error.

KFM_ERROR_NO_DATA No more data to return.

KFM_ERROR_INVALID_HANDLE The connection handle is invalid.

KFM_ERROR_INVALID_ARGUMENT Invalid function argument.

KFM_ERROR_DISCONNECTED The instrument is disconnected.

KFM_ERROR_FILE_NOT_FOUND Protocol input or output file not found.

KFM_ERROR_INVALID_BDZ_FILE The given file is not a valid bdz file.

KFM_ERROR_PROTOCOL_NOT_FOUND Requested protocol not found in the given bdz file.

KFM_ERROR_TEMP_FILE_CREATE Failed to create a temporary file.

KFM_ERROR_EXPORT_IMPORT_BUSY Can only execute one Export or Import at a time.

KFM_ERROR_OUT_OF_HEAP Out of heap memory.

Definition at line 34 of file KFMModuleDII.h.

4.3.3.2 enum KFM_SIMULATOR_PORT

Enumerator

KFM_SIMULATOR_USB USB port.

KFM_SIMULATOR_COM Serial command port.

KFM_SIMULATOR_DBG Serial debug port.

KFM_SIMULATOR_LAN LAN port.

Definition at line 50 of file KFMModuleDII.h.

4.3.4 Function Documentation

4.3.4.1 DIExport KFM_ERROR WINAPI KFMModule_Abort (HANDLE hConn)

Parameters

<i>hConn</i>	A handle returned by one of the KFMModule_OpenXxx() functions.
--------------	--

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

Aborts any command(s) being executed in the instrument.

Definition at line 2553 of file KFMModuleDII.c.

4.3.4.2 DllExport KFM_ERROR WINAPI KModule_AttachCallback (HANDLE *hConn*, void (*)(HANDLE *conn*, UINT *ev*) *callback*)

Parameters

<i>hConn</i>	A handle returned by one of the KModule_OpenXxx() functions.
<i>callback</i>	Address of the callback function.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

The callback function will be called whenever any of events [KFMODULE_RECEIVE](#), [KFMODULE_TRANSMIT](#), [KFMODULE_RESPONSE](#), [KFMODULE_EVENT](#) or [KFMODULE_ERROR](#) occurs. The connection handle and the event code are passed as parameters to the function.

Definition at line 2451 of file KModuleDll.c.

4.3.4.3 DllExport KFM_ERROR WINAPI KModule_AttachEvent (HANDLE *hConn*, UINT *ev*, HANDLE *object*)

Parameters

<i>hConn</i>	A handle returned by one of the KModule_OpenXxx() functions.
<i>ev</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

Parameter 'ev' may be any combination of [KFMODULE_RECEIVE](#), [KFMODULE_TRANSMIT](#), [KFMODULE_RESPONSE](#), [KFMODULE_EVENT](#) and [KFMODULE_ERROR](#). The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event. If the same event object is used for more than one event, the application must check for all possible events after the event object is signalled.

Definition at line 2362 of file KModuleDll.c.

4.3.4.4 DllExport KFM_ERROR WINAPI KModule_AttachMsg (HANDLE *hConn*, HANDLE *hWnd*, UINT *msg*, UINT *evCode*)

Parameters

<i>hConn</i>	A handle returned by one of the KModule_OpenXxx() functions.
<i>hWnd</i>	Handle of the window to receive the message.
<i>msg</i>	Message id of the message to send.
<i>evCode</i>	Event(s) for which a message is sent.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

A message with message id 'msg' is sent to window 'hWnd' whenever any of the event(s) selected with parameter 'evCode' occurs. Parameter 'evCode' may be any combination of [KFMODULE_RECEIVE](#), [KFMODULE_TRANSMIT](#), [KFMODULE_RESPONSE](#), [KFMODULE_EVENT](#) and [KFMODULE_ERROR](#). The wParam of the sent message will be one of these event codes.

Definition at line 2417 of file KModuleDll.c.

4.3.4.5 DllExport KFM_ERROR WINAPI KFModule_Close (HANDLE *hConn*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

This function should be called when the communication channel is no longer needed. Failing to close the channel prevents new connections to the channel as long as the dll stays loaded.



Note. When using the XML interface, function [KFModule_Disconnect\(\)](#) must be called before this function in order to disconnect the instrument from the open communication channel. Otherwise the instrument may refuse new connections.

Definition at line 2099 of file KFModuleDll.c.

4.3.4.6 DllExport KFM_ERROR WINAPI KFModule_Connect (HANDLE *hConn*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

This function sends the "Connect" command with a current date/time setting to the instrument. It is recommended to use this function to connect to the instrument so that the calendar of the instrument is always set to the local time automatically.



Note. This function returns before the actual connection is done, meaning that the return value indicates merely if the starting of the command is successful or not. The instrument will send a separate response after the connection is finished. See interface specification of the instrument in question for more details.

Definition at line 2133 of file KFModuleDll.c.

4.3.4.7 DllExport KFM_ERROR WINAPI KFModule_Disconnect (HANDLE *hConn*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

This function sends the "disconnect" command to the instrument. This is only needed because of the symmetry in the list of API functions. Disconnect() is the pair to Connect() function.



Note. This function returns before the actual connection is closed, meaning that the return value indicates merely if

the starting of the command is succesfull of not. The instrument will send a separate response just before the disconnection. See interface specification of the instrument in question for more details.

Definition at line 2178 of file KFModuleDll.c.

4.3.4.8 DllExport KFM_ERROR WINAPI KFModule.DownloadProtocol (HANDLE *hConn*, LPCSTR *protocolName*, LPCSTR *path*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>protocolName</i>	Name of the protocol to export.
<i>path</i>	Complete path of the .bdz file to receive the protocol.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

The given protocol name must match exactly with a protocol name stored in the instrument. The dll must have write access to the given .bdz file. The file will be overwritten by the dll.



Note. This funtion returns before the actual transfer is completed, meaning that the return value indicates merely if the starting of the transfer is succesfull of not. The instrument will send a separate response after the transfer. See interface specification of the instrument in question for more details.

Definition at line 2627 of file KFModuleDll.c.

4.3.4.9 DllExport KFM_ERROR WINAPI KFModule.GetError (HANDLE *hConn*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
--------------	---

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

Most dll functions return an error code, but an error may occur asynchronously in the receive or transmit thread of the dll. This kind of errors are reported with the KFMODULE_ERROR event. The application should then call this function to read the error code and take action. The error usually is KFM_ERROR_DISCONNECTED, in which case the application should close the current connection and try to open a new one.

Definition at line 2673 of file KFModuleDll.c.

4.3.4.10 DllExport DWORD WINAPI KFModule.ListLanDevices (LPCSTR *DeviceName*, LPCSTR *SerialNumber*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>instrumentName</i>	Name of the instrument. Only instruments with a matching name are listed. May be NULL, in which case the found devices are not filtered by the name.
<i>SerialNumber</i>	The serial number of the instrument. Only instruments with a matching serial number are listed. May be NULL, in which case the found devices are not filtered by the serial number.
<i>buf</i>	The buffer to list the found devices to.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

The size in bytes of the complete list of found devices. If the returned value is equal or smaller than the given buffer size, the buffer contains the complete list of devices found on the LAN. If the returned value is higher than the given buffer size, no data is returned and the caller must call the function again with big enough buffer. Return value 0 means that no instruments were found.

On success, the caller's buffer contains zero terminated strings with a combined length of the return value. The string terminating zeros are included in the length.

For each found instrument, the first string is the instrument IPv4 address and the TCP port number it is listening, enclosed in square brackets, e.g. [10.32.196.210:49536].

The IP address string is followed by the WS-Discovery match strings, usually 3 of them. The first one is always 'Thermo-Device', the second one is the instrument name and the third one the instrument serial number string.

If the match strings are followed by a string in angle brackets, e.g. <10.32.196.154:57403>, it means that the instrument is currently connected that IP address and TCP port, and trying to connect to that instrument with function [KFModule_OpenLan\(\)](#) will fail. If there is no string in square brackets, the instrument will accept a connection.

Definition at line 2000 of file KFModuleDll.c.

4.3.4.11 DllExport HANDLE WINAPI KFModule_OpenLan (LPCSTR *DeviceName*, LPCSTR *SerialNumber*, UINT *timeout*)

Parameters

<i>instrumentName</i>	Name of the instrument. This must match the device name the instrument uses for matching a WS-Discovery Probe and Resolve.
<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first found KFModule instrument.
<i>timeout</i>	Timeout which is used to search the instrument. If 0 is given then WS-Discovery will use default 4 sec timeout.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the KFModule_OpenXxx() functions must be called first before using any other functions in the library. Only one connection per device is allowed.



Note. Function [KFModule_Connect\(\)](#) must be called after this function in order to use the XML interface to communicate with the instrument.

Definition at line 1935 of file KFModuleDll.c.

4.3.4.12 DllExport HANDLE WINAPI KFModule_OpenSerial (WORD *port*, DWORD *baud*, LPCSTR *DeviceName*, LPCSTR *SerialNumber*)

Parameters

<i>port</i>	Serial port number.
<i>baud</i>	Serial port baudrate.
<i>DeviceName</i>	Pointer to the device name string of the device. The device must report an identical name to the VER command for the connection to succeed. This parameter may be NULL, in which case the device name is ignored.

<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number to the VER command for the connection to succeed. This parameter may be NULL, in which case the serial number is ignored.
---------------------	--

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the `KFModule_OpenXxx()` functions must be called first before using any other functions in the library. Only one connection per serial port is allowed. If both `DeviceName` and `SerialNumber` are NULL, no VER command is sent to the instrument.



Note. Function `KFModule_Connect()` must be called after this function in order to use the XML interface to communicate with the instrument.

Definition at line 1880 of file `KFModuleDll.c`.

4.3.4.13 DllExport HANDLE WINAPI KFModule_OpenSimulator (KFM_SIMULATOR_PORT *port*, WORD *productId*)

Parameters

<i>port</i>	The simulator port to connect to, one of <code>KFM_SIMULATOR_USB</code> , <code>KFM_SIMULATOR_COM</code> , <code>KFM_SIMULATOR_DBG</code> or <code>KFM_SIMULATOR_LAN</code> .
<i>productId</i>	(USB) product id of the instrument.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the `KFModule_OpenXxx()` functions must be called first before using any other functions in the library. Only one connection per simulator communication channel is allowed.



Note. Function `KFModule_Connect()` must be called after this function in order to use the XML interface to communicate with the instrument.

Definition at line 2030 of file `KFModuleDll.c`.

4.3.4.14 DllExport HANDLE WINAPI KFModule_OpenUsb (LPCSTR *SerialNumber*, WORD *productId*)

Parameters

<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first device with matching VendorID and ProductID.
<i>productId</i>	Manufacturer product id number of the USB device.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

One of the `KFModule_OpenXxx()` functions must be called first before using any other functions in the library. Only one connection per device is allowed.



Note. Function `KFModule_Connect()` must be called after this function in order to use the XML interface to communicate with the instrument.

Definition at line 1827 of file `KFModuleDll.c`.

4.3.4.15 DllExport KFM_ERROR WINAPI KFModule_ReadEvent (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hConn</i>	A handle returned by one of the <code>KFModule_OpenXxx()</code> functions.
<i>buf</i>	Caller's buffer for the event.
<i>bufsize</i>	Size of the caller's buffer.

Returns

One of `KFM_ERROR` codes, `KFM_ERROR_SUCCESS` if any data was copied to caller's buffer.

Call this function in response to the `KFMODULE_EVENT` event. The received data is a XML Event. The event is not sent until a whole XML Event is received. If the caller's buffer is not large enough to hold the whole Event, a partial Event is copied. An empty string is copied when there is no more data to read. The returned string is always NUL terminated and does not contain the CR/LF characters sent by the instrument.



Note. Always read all data from the `KFModuleDll` when handling the event. Otherwise subsequent events from the instrument may be missed. This can happen especially when the user application doesn't react to the events immediately. See example below:

```
void handler_KFMODULE_EVENT( HANDLE connection )
{
    KFM_ERROR    err;
    char         buf[100];

    while( (err = KFModule_ReadEvent( connection, buf, sizeof( buf ) )) == KFM_ERROR_SUCCESS)
    {
        // Do something with the data
    }

    if( err > KFM_ERROR_NO_DATA)
    {
        // Handle the error
    }
}
```

Definition at line 2333 of file `KFModuleDll.c`.

4.3.4.16 DllExport KFM_ERROR WINAPI KFModule_ReadReceived (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hConn</i>	A handle returned by one of the <code>KFModule_OpenXxx()</code> functions.
<i>buf</i>	Caller's buffer for the line.
<i>bufsize</i>	Size of the caller's buffer.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS if any data was copied to caller's buffer.

Call this function in response to the [KFMODULE_RECEIVE](#) event. The received data is neither a XML Response nor a XML Event. An empty string will be copied to caller's buffer when there is no more data to read. The returned string is always NUL terminated and does not contain the CR/LF characters sent by the instrument.



Note. Always read all data from the KFMModuleDll when handling the event. Otherwise subsequent strings from the instrument may be missed. This can happen especially when the user application doesn't react to the events immediately. See example below:

```
void handler_KFMODULE_RECEIVE( HANDLE connection )
{
    KFM_ERROR    err;
    char         buf[100];

    while( (err = KFMModule_ReadReceived( connection, buf, sizeof( buf ))) == KFM_ERROR_SUCCESS)
    {
        // Do something with the data
    }

    if( err > KFM_ERROR_NO_DATA)
    {
        // Handle the error
    }
}
```

Definition at line 2239 of file KFMModuleDll.c.

4.3.4.17 DllExport KFM_ERROR WINAPI KFMModule_ReadResponse (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hConn</i>	A handle returned by one of the KFMModule_OpenXxx() functions.
<i>buf</i>	Caller's buffer for the response.
<i>bufsize</i>	Size of the caller's buffer.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS if any data was copied to caller's buffer.

Call this function in response to the [KFMODULE_RESPONSE](#) event. The received data is a XML Response. The event is not sent until a whole XML Response is received. If the caller's buffer is not large enough to hold the whole response, a partial response is copied. An empty string is copied when there is no more data to read. The returned string is always NUL terminated and does not contain the CR/LF characters sent by the instrument.



Note. Always read all data from the KFMModuleDll when handling the event. Otherwise subsequent responses may be missed. This can happen especially when the user application doesn't react to the events immediately. See example below:

```
void handler_KFMODULE_RESPONSE( HANDLE connection )
{
    KFM_ERROR    err;
    char         buf[100];

    while( (err = KFMModule_ReadResponse( connection, buf, sizeof( buf ))) == KFM_ERROR_SUCCESS)
    {
        // Do something with the data
    }
}
```

```

    }

    if( err > KFM_ERROR_NO_DATA)
    {
        // Handle the error
    }
}

```

Definition at line 2286 of file KFModuleDll.c.

4.3.4.18 DllExport KFM_ERROR WINAPI KFModule_Send (HANDLE *hConn*, LPCSTR *buf*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>buf</i>	The NUL terminated string to send.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

Use this function to send commands and acknowledges to the instrument. For uploading and downloading a protocol there are dedicated functions [KFModule_UploadProtocol\(\)](#) and [KFModule_DownloadProtocol\(\)](#). Also connecting/disconnecting is recommended to be done with separate [KFModule_Connect\(\)](#) and [KFModule_Disconnect\(\)](#) functions.

The data to be sent to the instrument is buffered in the dll. A [KFMODULE_TRANSMIT](#) event is sent to the application when all data is sent. You may call this function repeatedly without waiting for the event, but a [KFM_ERROR_OUT_OF_HEAP](#) error may be returned if you send a lot of data without waiting for the event.



Note. Remember to terminate commands with a new line character (ASCII 10). See interface specification of the instrument in question for more details.

Definition at line 2495 of file KFModuleDll.c.

4.3.4.19 DllExport KFM_ERROR WINAPI KFModule_UploadProtocol (HANDLE *hConn*, LPCSTR *protocolName*, LPCSTR *path*)

Parameters

<i>hConn</i>	A handle returned by one of the KFModule_OpenXxx() functions.
<i>protocolName</i>	Name of the protocol to export.
<i>path</i>	Complete path of the .bdz file containing the protocol to export.

Returns

One of [KFM_ERROR](#) codes, KFM_ERROR_SUCCESS on success.

The given protocol name must match exactly with a protocol name stored in the .bdz file. The .bdz file may contain several protocols, but only the requested protocol is sent to the instrument.



Note. This funtion returns before the actual transfer is completed, meaning that the return value indicates merely if the starting of the transfer is succesfull or not. The instrument will send a separate response after the transfer. See interface specification of the instrument in question for more details.

Definition at line 2600 of file KFModuleDll.c.

4.4 mainpage.h File Reference

KFModuleDll Interface Specification.

Index

api.h, [27](#)

KFM_ERROR_DISCONNECTED
 KFMModuleDll.h, [39](#)
KFM_ERROR_EXPORT_IMPORT_BUSY
 KFMModuleDll.h, [39](#)
KFM_ERROR_FILE_NOT_FOUND
 KFMModuleDll.h, [39](#)
KFM_ERROR_INVALID_ARGUMENT
 KFMModuleDll.h, [39](#)
KFM_ERROR_INVALID_BDZ_FILE
 KFMModuleDll.h, [39](#)
KFM_ERROR_INVALID_HANDLE
 KFMModuleDll.h, [39](#)
KFM_ERROR_NO_DATA
 KFMModuleDll.h, [39](#)
KFM_ERROR_OUT_OF_HEAP
 KFMModuleDll.h, [39](#)
KFM_ERROR_PROTOCOL_NOT_FOUND
 KFMModuleDll.h, [39](#)
KFM_ERROR_SUCCESS
 KFMModuleDll.h, [39](#)
KFM_ERROR_TEMP_FILE_CREATE
 KFMModuleDll.h, [39](#)
KFM_SIMULATOR_COM
 KFMModuleDll.h, [39](#)
KFM_SIMULATOR_DBG
 KFMModuleDll.h, [39](#)
KFM_SIMULATOR_LAN
 KFMModuleDll.h, [39](#)
KFM_SIMULATOR_USB
 KFMModuleDll.h, [39](#)
KFMModuleDll.h
 KFM_ERROR_DISCONNECTED, [39](#)
 KFM_ERROR_EXPORT_IMPORT_BUSY, [39](#)
 KFM_ERROR_FILE_NOT_FOUND, [39](#)
 KFM_ERROR_INVALID_ARGUMENT, [39](#)
 KFM_ERROR_INVALID_BDZ_FILE, [39](#)
 KFM_ERROR_INVALID_HANDLE, [39](#)
 KFM_ERROR_NO_DATA, [39](#)
 KFM_ERROR_OUT_OF_HEAP, [39](#)
 KFM_ERROR_PROTOCOL_NOT_FOUND, [39](#)
 KFM_ERROR_SUCCESS, [39](#)
 KFM_ERROR_TEMP_FILE_CREATE, [39](#)
 KFM_SIMULATOR_COM, [39](#)
 KFM_SIMULATOR_DBG, [39](#)
 KFM_SIMULATOR_LAN, [39](#)
 KFM_SIMULATOR_USB, [39](#)

KF_PRESTO_PID
 KFMModuleDll.h, [38](#)
KFM_ERROR
 KFMModuleDll.h, [39](#)
KFM_SIMULATOR_PORT
 KFMModuleDll.h, [39](#)
KFMODULE_ERROR
 KFMModuleDll.h, [38](#)
KFMODULE_EVENT
 KFMModuleDll.h, [38](#)
KFMODULE_RECEIVE
 KFMModuleDll.h, [38](#)
KFMODULE_RESPONSE
 KFMModuleDll.h, [38](#)
KFMODULE_TRANSMIT
 KFMModuleDll.h, [38](#)
KFMModule_Abort
 KFMModuleDll.c, [28](#)
 KFMModuleDll.h, [39](#)
KFMModule_AttachCallback
 KFMModuleDll.c, [28](#)
 KFMModuleDll.h, [39](#)
KFMModule_AttachEvent
 KFMModuleDll.c, [29](#)
 KFMModuleDll.h, [40](#)
KFMModule_AttachMsg
 KFMModuleDll.c, [29](#)
 KFMModuleDll.h, [40](#)
KFMModule_Close
 KFMModuleDll.c, [29](#)
 KFMModuleDll.h, [40](#)
KFMModule_Connect
 KFMModuleDll.c, [30](#)
 KFMModuleDll.h, [41](#)
KFMModule_Disconnect
 KFMModuleDll.c, [30](#)
 KFMModuleDll.h, [41](#)
KFMModule_DownloadProtocol
 KFMModuleDll.c, [30](#)
 KFMModuleDll.h, [42](#)
KFMModule_GetError
 KFMModuleDll.c, [31](#)
 KFMModuleDll.h, [42](#)
KFMModule_ListLanDevices
 KFMModuleDll.c, [31](#)
 KFMModuleDll.h, [42](#)
KFMModule_OpenLan

- KFModuleDll.c, [32](#)
- KFModuleDll.h, [43](#)
- KFModule_OpenSerial
 - KFModuleDll.c, [32](#)
 - KFModuleDll.h, [43](#)
- KFModule_OpenSimulator
 - KFModuleDll.c, [33](#)
 - KFModuleDll.h, [44](#)
- KFModule_OpenUsb
 - KFModuleDll.c, [33](#)
 - KFModuleDll.h, [44](#)
- KFModule_ReadEvent
 - KFModuleDll.c, [33](#)
 - KFModuleDll.h, [45](#)
- KFModule_ReadReceived
 - KFModuleDll.c, [34](#)
 - KFModuleDll.h, [45](#)
- KFModule_ReadResponse
 - KFModuleDll.c, [35](#)
 - KFModuleDll.h, [46](#)
- KFModule_Send
 - KFModuleDll.c, [35](#)
 - KFModuleDll.h, [47](#)
- KFModule_UploadProtocol
 - KFModuleDll.c, [36](#)
 - KFModuleDll.h, [47](#)
- KFModuleDll.c, [27](#)
 - KFModule_Abort, [28](#)
 - KFModule_AttachCallback, [28](#)
 - KFModule_AttachEvent, [29](#)
 - KFModule_AttachMsg, [29](#)
 - KFModule_Close, [29](#)
 - KFModule_Connect, [30](#)
 - KFModule_Disconnect, [30](#)
 - KFModule_DownloadProtocol, [30](#)
 - KFModule_GetError, [31](#)
 - KFModule_ListLanDevices, [31](#)
 - KFModule_OpenLan, [32](#)
 - KFModule_OpenSerial, [32](#)
 - KFModule_OpenSimulator, [33](#)
 - KFModule_OpenUsb, [33](#)
 - KFModule_ReadEvent, [33](#)
 - KFModule_ReadReceived, [34](#)
 - KFModule_ReadResponse, [35](#)
 - KFModule_Send, [35](#)
 - KFModule_UploadProtocol, [36](#)
- KFModuleDll.h, [36](#)
 - KF_PRESTO_PID, [38](#)
 - KFM_ERROR, [39](#)
 - KFMODULE_ERROR, [38](#)
 - KFMODULE_EVENT, [38](#)
 - KFMODULE_RECEIVE, [38](#)
 - KFMODULE_RESPONSE, [38](#)
 - KFMODULE_TRANSMIT, [38](#)
 - KFModule_Abort, [39](#)
 - KFModule_AttachCallback, [39](#)
 - KFModule_AttachEvent, [40](#)
 - KFModule_AttachMsg, [40](#)
 - KFModule_Close, [40](#)
 - KFModule_Connect, [41](#)
 - KFModule_Disconnect, [41](#)
 - KFModule_DownloadProtocol, [42](#)
 - KFModule_GetError, [42](#)
 - KFModule_ListLanDevices, [42](#)
 - KFModule_OpenLan, [43](#)
 - KFModule_OpenSerial, [43](#)
 - KFModule_OpenSimulator, [44](#)
 - KFModule_OpenUsb, [44](#)
 - KFModule_ReadEvent, [45](#)
 - KFModule_ReadReceived, [45](#)
 - KFModule_ReadResponse, [46](#)
 - KFModule_Send, [47](#)
 - KFModule_UploadProtocol, [47](#)
- mainpage.h, [48](#)

Thermo Scientific

ThermoUSB.dll

Interface Specification

Contents

1	ThermoUSB dll Interface Specification	1
2	Using ThermoUSB	3
2.1	About	3
2.2	Exported functions	4
2.2.1	ThermoUSBOpen	5
2.2.2	ThermoUSBClose	6
2.2.3	ThermoUSBAttachEvent	7
2.2.4	ThermoUSBAttachMsg	8
2.2.5	ThermoUSBRead	9
2.2.6	ThermoUSBReadBinary	10
2.2.7	ThermoUSBWrite	11
2.2.8	ThermoUSBWriteBinary	12
2.2.9	ThermoUSBGetError	13
2.2.10	ThermoUSBAbort	14
2.2.11	ThermoUSBGetThermoVendorId	15
3	File Index	17
3.1	File List	17
4	File Documentation	19
4.1	api.h File Reference	19
4.2	mainpage.h File Reference	19
4.3	ThermoUSB.c File Reference	19
4.3.1	Detailed Description	20
4.3.2	Function Documentation	20
4.3.2.1	ThermoUSBAbort	20
4.3.2.2	ThermoUSBAttachEvent	20
4.3.2.3	ThermoUSBAttachMsg	20
4.3.2.4	ThermoUSBClose	21
4.3.2.5	ThermoUSBGetError	21
4.3.2.6	ThermoUSBGetThermoVendorId	21
4.3.2.7	ThermoUSBOpen	21
4.3.2.8	ThermoUSBRead	22
4.3.2.9	ThermoUSBReadBinary	22
4.3.2.10	ThermoUSBWrite	23
4.3.2.11	ThermoUSBWriteBinary	23
4.4	ThermoUSB.h File Reference	23
4.4.1	Detailed Description	24
4.4.2	Macro Definition Documentation	24
4.4.2.1	THERMOUSB.ERROR	24
4.4.2.2	THERMOUSB.RECEIVE	24
4.4.2.3	THERMOUSB.TRANSMIT	24
4.4.3	Function Documentation	25
4.4.3.1	ThermoUSBAbort	25
4.4.3.2	ThermoUSBAttachEvent	25

4.4.3.3	ThermoUSBAttachMsg	25
4.4.3.4	ThermoUSBClose	26
4.4.3.5	ThermoUSBGetError	26
4.4.3.6	ThermoUSBGetThermoVendorId	26
4.4.3.7	ThermoUSBOpen	26
4.4.3.8	ThermoUSBRead	27
4.4.3.9	ThermoUSBReadBinary	27
4.4.3.10	ThermoUSBWrite	27
4.4.3.11	ThermoUSBWriteBinary	27

Chapter 1

ThermoUSB dll Interface Specification

About

The purpose of the ThermoUSB.dll dynamic link library is to make interface to the Thermo microplate instruments easy. The user of the ThermoUSB.dll does not have to know the details of USB communication. Knowing the USB Product Id of the instrument is enough.

Confidential

This document has been prepared by Thermo Fisher Scientific Oy to be used solely for the purposes defined by Thermo Fisher Scientific Oy. Use for other purposes is not authorized.

Please note that any and all information contained in this document is the property of Thermo Fisher Scientific Oy. This confidential information ("Confidential Information") shall not be reproduced in whole part or disclosed to any third party without the prior written approval of Thermo Fisher Scientific Oy. The receiving party shall ensure that its employees, officers, representatives and agents shall not disclose to third parties any Confidential Information.

Upon written request from Thermo Fisher Scientific Oy, the receiving party shall promptly return all Confidential Information or destroy all Confidential Information.

Chapter 2

Using ThermoUSB

2.1 About

Because it is not trivial to write PC software from scratch to communicate with the instrument, the ThermoUSB Dynamic Link Library is provided. It hides much of the complexity of the communication.

ThermoUSB.dll is a generic library for communicating with several different Thermo Scientific microplate instruments. To communicate with an instrument you use the exported functions of ThermoUSB.dll.

Depending on your project setup, you may find useful a couple of other files which are also provided. The header file [ThermoUSB.h](#) contains the prototypes of the exported functions and definitions of constant values used by the dll. File ThermoUSB.lib contains information about the dll the linker uses to add references to the library in the executable. This way the dll is automatically loaded and the exported functions of the library can be called as easy as the functions in the code using the dll.

2.2 Exported functions

- [ThermoUSBOpen](#)
- [ThermoUSBClose](#)
- [ThermoUSBAttachEvent](#)
- [ThermoUSBAttachMsg](#)
- [ThermoUSBRead](#)
- [ThermoUSBReadBinary](#)
- [ThermoUSBWrite](#)
- [ThermoUSBWriteBinary](#)
- [ThermoUSBGetError](#)
- [ThermoUSBAbort](#)
- [ThermoUSBGetThermoVendorId](#)

2.2.1 ThermoUSBOpen

Search for the requested USB device and open a communication channel to it. Full declaration: [ThermoUSBOpen\(\)](#).

Parameters

<i>VendorID</i>	The USB vendor id of the device manufacturer, 0x0AB6 for Thermo Fisher Scientific Oy.
<i>ProductID</i>	Product id number of the device.
<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first device with matching VendorID and ProductID.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

This function (or ThermoUSBOpenSimulator) must be called first before using any other functions in the library. Only one connection per device is allowed.

2.2.2 ThermoUSBClose

Close an USB connection. Full declaration: [ThermoUSBClose\(\)](#).

Parameters

<i>hConn</i>	Connection handle returned form a call to ThermoUSBOpen()
--------------	---

Closes the communication channel to the instrument. This function should be called when the application has finished communicating with the instrument. It is not possible to open a communication channel to the instrument while the previous channel is still open. When the dll is unloaded from the memory, all channels still open are automatically closed.

2.2.3 ThermoUSBAttachEvent

Attach an event object to an USB connection. Full declaration: [ThermoUSBAttachEvent\(\)](#).

Parameters

<i>hConn</i>	Connection handle returned from a call to ThermoUSBOpen() (pointer to a USB_CONNECTION structure).
<i>evCode</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Parameter 'evCode' may be any combination of THERMOUSB_RECEIVE, THERMOUSB_TRANSMIT and THERMOUSB_ERROR. The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event.

The THERMOUSB_RECEIVE event is signaled whenever something is written to the receive buffer. There is no guarantee that a whole response line is received or that only one response line is received. therefore upon receiving this event the application should call [ThermoUSBRead\(\)](#) or ThermoUSBReadBinary as long as they return data.

The THERMOUSB_TRANSMIT event is signaled when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOUSB_TRANSMIT event is received, functions [ThermoUSBWrite\(\)](#) and [ThermoUSBWriteBinary\(\)](#) will never fail.

2.2.4 ThermoUSBAttachMsg

Attach an event message to an USB connection. Full declaration: [ThermoUSBAttachMsg\(\)](#).

Parameters

<i>hConn</i>	Connection handle returned from a call to ThermoUSBOpen() (pointer to a USB_CONNECTION structure).
<i>hWnd</i>	Handle of the window to receive the message.
<i>msg</i>	Message id of the message to send.
<i>event</i>	Event(s) for which a message is sent.

A message with message id 'msg' is sent to window 'hWnd' whenever any of the event(s) selected with parameter 'evCode' occurs. Parameter 'evCode' may be any combination of THERMOUSB_RECEIVE, THERMOUSB_TRANSMIT and THERMOUSB_ERROR. The event code is passed in the wParam member of the sent Windows message.

The THERMOUSB_RECEIVE event is sent whenever something is written to the receive buffer. There is no guarantee that a whole response line is received or that only one response line is received. Therefore upon receiving this event the application should call [ThermoUSBRead\(\)](#) or [ThermoUSBReadBinary](#) as long as they return data.

The THERMOUSB_TRANSMIT event is sent when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOUSB_TRANSMIT event is received, functions [ThermoUSBWrite\(\)](#) and [ThermoUSBWriteBinary\(\)](#) will never fail.

2.2.5 ThermoUSBRead

Read a received response line. Full declaration: [ThermoUSBRead\(\)](#).

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
<i>buf</i>	Pointer to a buffer to receive the response.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

TRUE if a response line was retrieved, else FALSE.

Use this function to read full response lines returned from the instrument. The response line is returned without the terminating CRLF. Therefore TRUE may be returned even if the resulting line is empty. The returned line always ends to the NUL character even if FALSE is returned.

If the caller's buffer is too small to hold the whole response line, as much as fits to the buffer is returned and the rest is returned on subsequent call(s).

2.2.6 ThermoUSBReadBinary

Read data received from the USB device. Full declaration: [ThermoUSBReadBinary\(\)](#).

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
<i>buf</i>	Pointer to buffer receiving the data.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

Number of bytes copied to the user buffer.

Unlike function [ThermoUSBRead\(\)](#), this function returns all data received from the instrument without doing any interpretation on it. You can control the number of bytes returned with the bufsize parameter.

If the return value is less than the given bufsize parameter, there is no more data.

2.2.7 ThermoUSBWrite

Write a string to the transmit buffer. Full declaration: [ThermoUSBWrite\(\)](#).

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
<i>buf</i>	Buffer containing the ASCIIZ string to send.

Returns

TRUE if the string was written to the transmit buffer, else FALSE.

If the whole string does not fit in the transmit buffer, writes nothing and returns FALSE. The application may retry at a later time.

2.2.8 ThermoUSBWriteBinary

Write binary data to the transmit buffer. Full declaration: [ThermoUSBWriteBinary\(\)](#).

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
<i>buf</i>	Buffer containing the data to send.
<i>count</i>	Number of bytes to send.

Returns

TRUE if the data was written to the transmit buffer, else FALSE.

If all data does not fit in the transmit buffer, writes nothing and returns FALSE. The application may retry at a later time.

2.2.9 ThermoUSBGetError

Return the error code stored to the connection structure. Full declaration: [ThermoUSBGetError\(\)](#).

2.2.10 ThermoUSBAbort

Send Abort command to the device. Full declaration: [ThermoUSBAbort\(\)](#).

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
--------------	---

2.2.11 ThermoUSBGetThermoVendorId

Return Thermo Fisher Scientific Oy USB vendor id. Full declaration: [ThermoUSBGetThermoVendorId\(\)](#).

Returns

Thermo Fisher Scientific Oy USB vendor id, 0x0AB6.

Chapter 3

File Index

3.1 File List

Here is a list of all files with brief descriptions:

api.h	Part of ThermoUSB dll documentation	19
mainpage.h	Start page of ThermoUSB.dll Design Description	19
ThermoUSB.c	Implements USB interface to Thermo microplate instruments using either Windows HID driver or the libusb-win32 driver from http://libusb-win32.sourceforge.net	19
ThermoUSB.h	Functions exported from ThermoUSB.dll	23

Chapter 4

File Documentation

4.1 api.h File Reference

Part of ThermoUSB dll documentation.

4.2 mainpage.h File Reference

Start page of ThermoUSB.dll Design Description.

4.3 ThermoUSB.c File Reference

Implements USB interface to Thermo microplate instruments using either Windows HID driver or the libusb-win32 driver from <http://libusb-win32.sourceforge.net>.

Functions

- HANDLE WINAPI [ThermoUSBOpen](#) (WORD VendorID, WORD ProductID, LPCSTR SerialNumber)
Search for the requested USB device and open a communication channel to it.
- void WINAPI [ThermoUSBClose](#) (HANDLE hConn)
Close an USB connection.
- BOOL WINAPI [ThermoUSBAttachEvent](#) (HANDLE hConn, UINT evCode, HANDLE object)
Attach an event object to an USB connection.
- BOOL WINAPI [ThermoUSBAttachMsg](#) (HANDLE hConn, HANDLE hWnd, UINT msg, UINT evCode)
Attach an event message to an USB connection.
- BOOL WINAPI [ThermoUSBRead](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read a received response line.
- DWORD WINAPI [ThermoUSBReadBinary](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read data received from the USB device.
- BOOL WINAPI [ThermoUSBWrite](#) (HANDLE hConn, LPCSTR buf)
Write a string to the transmit buffer.
- BOOL WINAPI [ThermoUSBWriteBinary](#) (HANDLE hConn, LPCSTR buf, DWORD count)
Write binary data to the transmit buffer.
- void WINAPI [ThermoUSBAbort](#) (HANDLE hConn)
Send Abort command to the device.

- DWORD WINAPI [ThermoUSBGetError](#) (HANDLE hConn)
Return the error code stored to the connection structure.
- WORD WINAPI [ThermoUSBGetThermoVendorId](#) (void)
Return Thermo Fisher Scientific Oy USB vendor id.

4.3.1 Detailed Description

Note

Copyright by Thermo Fisher Scientific Oy 2015

Definition in file [ThermoUSB.c](#).

4.3.2 Function Documentation

4.3.2.1 void WINAPI ThermoUSBAbort (HANDLE hConn)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
--------------	---

Definition at line 2241 of file ThermoUSB.c.

4.3.2.2 BOOL WINAPI ThermoUSBAttachEvent (HANDLE hConn, UINT evCode, HANDLE object)

Parameters

<i>hConn</i>	Connection handle returned form a call to ThermoUSBOpen() (pointer to a USB_CONNECTION structure).
<i>evCode</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Parameter 'evCode' may be any combination of THERMOUSB_RECEIVE, THERMOUSB_TRANSMIT and THERMOUSB_ERROR. The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event.

The THERMOUSB_RECEIVE event is signaled whenever something is written to the receive buffer. There is no quarantee that a whole response line is received or that only one response line is received. therefore upon receiving this event the application should call [ThermoUSBRead\(\)](#) or ThermoUSBReadBinary as long as they return data.

The THERMOUSB_TRANSMIT event is signaled when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOUSB_TRANSMIT event is received, functions [ThermoUSBWrite\(\)](#) and [ThermoUSBWriteBinary\(\)](#) will never fail.

Definition at line 1853 of file ThermoUSB.c.

4.3.2.3 BOOL WINAPI ThermoUSBAttachMsg (HANDLE hConn, HANDLE hWnd, UINT msg, UINT evCode)

Parameters

<i>hConn</i>	Connection handle returned form a call to ThermoUSBOpen() (pointer to a USB_CONNECTION structure).
<i>hWnd</i>	Handle of the window to receive the message.
<i>msg</i>	Message id of the message to send.
<i>event</i>	Event(s) for which a message is sent.

A message with message id 'msg' is sent to window 'hWnd' whenever any of the event(s) selected with parameter 'ev-Code' occurs. Parameter 'evCode' may be any combination of THERMOUSB_RECEIVE, THERMOUSB_TRANSMIT and THERMOUSB_ERROR. The event code is passed in the wParam member of the sent Windows message.

The THERMOUSB_RECEIVE event is sent whenever something is written to the receive buffer. There is no guarantee that a whole response line is received or that only one response line is received. Therefore upon receiving this event the application should call [ThermoUSBRead\(\)](#) or ThermoUSBReadBinary as long as they return data.

The THERMOUSB_TRANSMIT event is sent when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOUSB_TRANSMIT event is received, functions [ThermoUSBWrite\(\)](#) and [ThermoUSBWriteBinary\(\)](#) will never fail.

Definition at line 1908 of file ThermoUSB.c.

4.3.2.4 void WINAPI ThermoUSBClose (HANDLE hConn)

Parameters

<i>hConn</i>	Connection handle returned from a call to ThermoUSBOpen()
--------------	---

Closes the communication channel to the instrument. This function should be called when the application has finished communicating with the instrument. It is not possible to open a communication channel to the instrument while the previous channel is still open. When the dll is unloaded from the memory, all channels still open are automatically closed.

Definition at line 1767 of file ThermoUSB.c.

4.3.2.5 DWORD WINAPI ThermoUSBGetError (HANDLE hConn)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
--------------	---

Returns

The current Windows error code stored for the connection.

Error conditions may be encountered asynchronously and not just as a direct result of a function call to the dll. Therefore the application should enable error reporting with [ThermoUSBAttachEvent\(\)](#) or [ThermoUSBAttachMsg\(\)](#). When an error is reported to the application, it may call this function to get the Windows error code, which may or may not be helpful in resolving the problem. In general, if an error is reported then some receive or transmit data is lost, and the best way of action is to close the port and then retry opening it again.

A call to [ThermoUSBGetError\(\)](#) resets the stored error code to ERROR_SUCCESS.

Definition at line 2308 of file ThermoUSB.c.

4.3.2.6 WORD WINAPI ThermoUSBGetThermoVendorId (void)

Returns

Thermo Fisher Scientific Oy USB vendor id, 0x0AB6.

Definition at line 2332 of file ThermoUSB.c.

4.3.2.7 HANDLE WINAPI ThermoUSBOpen (WORD VendorID, WORD ProductID, LPCSTR SerialNumber)

Parameters

<i>VendorID</i>	The USB vendor id of the device manufacturer, 0x0AB6 for Thermo Fisher Scientific Oy.
-----------------	---

<i>ProductID</i>	Product id number of the device.
<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first device with matching VendorID and ProductID.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

This function (or ThermoUSBOpenSimulator) must be called first before using any other functions in the library. Only one connection per device is allowed.

Definition at line 1738 of file ThermoUSB.c.

4.3.2.8 BOOL WINAPI ThermoUSBRead (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
<i>buf</i>	Pointer to a buffer to receive the response.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

TRUE if a response line was retrieved, else FALSE.

Use this function to read full response lines returned from the instrument. The response line is returned without the terminating CRLF. Therefore TRUE may be returned even if the resulting line is empty. The returned line always ends to the NUL character even if FALSE is returned.

If the caller's buffer is too small to hold the whole response line, as much as fits to the buffer is returned and the rest is returned on subsequent call(s).

Definition at line 1948 of file ThermoUSB.c.

4.3.2.9 DWORD WINAPI ThermoUSBReadBinary (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
<i>buf</i>	Pointer to buffer receiving the data.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

Number of bytes copied to the user buffer.

Unlike function [ThermoUSBRead\(\)](#), this function returns all data received from the instrument without doing any interpretation on it. You can control the number of bytes returned with the bufsize parameter.

If the return value is less than the given bufsize parameter, there is no more data.

Definition at line 2123 of file ThermoUSB.c.

4.3.2.10 BOOL WINAPI ThermoUSBWrite (HANDLE *hConn*, LPCSTR *buf*)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
<i>buf</i>	Buffer containing the ASCIIZ string to send.

Returns

TRUE if the string was written to the transmit buffer, else FALSE.

If the whole string does not fit in the transmit buffer, writes nothing and returns FALSE. The application may retry at a later time.

Definition at line 2169 of file ThermoUSB.c.

4.3.2.11 BOOL WINAPI ThermoUSBWriteBinary (HANDLE *hConn*, LPCSTR *buf*, DWORD *count*)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
<i>buf</i>	Buffer containing the data to send.
<i>count</i>	Number of bytes to send.

Returns

TRUE if the data was written to the transmit buffer, else FALSE.

If all data does not fit in the transmit buffer, writes nothing and returns FALSE. The application may retry at a later time.

Definition at line 2190 of file ThermoUSB.c.

4.4 ThermoUSB.h File Reference

Functions exported from ThermoUSB.dll.

Macros

- #define [THERMOUSB_RECEIVE](#) 1
Data received.
- #define [THERMOUSB_TRANSMIT](#) 2
All data transmitted.
- #define [THERMOUSB_ERROR](#) 4
Fatal error event.

Functions

- DllExport HANDLE WINAPI [ThermoUSBOpen](#) (WORD VendorID, WORD ProductID, LPCSTR SerialNumber)
Search for the requested USB device and open a communication channel to it.
- DllExport void WINAPI [ThermoUSBClose](#) (HANDLE *hConn*)
Close an USB connection.
- DllExport BOOL WINAPI [ThermoUSBAttachEvent](#) (HANDLE *hConn*, UINT *evCode*, HANDLE *object*)
Attach an event object to an USB connection.

- DllExport BOOL WINAPI [ThermoUSBAttachMsg](#) (HANDLE hConn, HANDLE hWnd, UINT msg, UINT event)
Attach an event message to an USB connection.
- DllExport BOOL WINAPI [ThermoUSBRead](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read a received response line.
- DllExport DWORD WINAPI [ThermoUSBReadBinary](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read data received from the USB device.
- DllExport BOOL WINAPI [ThermoUSBWrite](#) (HANDLE hConn, LPCSTR buf)
Write a string to the transmit buffer.
- DllExport BOOL WINAPI [ThermoUSBWriteBinary](#) (HANDLE hConn, LPCSTR buf, DWORD count)
Write binary data to the transmit buffer.
- DllExport DWORD WINAPI [ThermoUSBGetError](#) (HANDLE hConn)
Return the error code stored to the connection structure.
- DllExport void WINAPI [ThermoUSBAbort](#) (HANDLE hConn)
Send Abort command to the device.
- DllExport WORD WINAPI [ThermoUSBGetThermoVendorId](#) (void)
Return Thermo Fisher Scientific Oy USB vendor id.

4.4.1 Detailed Description

Note

Copyright by Thermo Fisher Scientific Oy 2015

USB interface to Thermo microplate instruments.

Definition in file [ThermoUSB.h](#).

4.4.2 Macro Definition Documentation

4.4.2.1 #define THERMOUSB_ERROR 4

In case of error the application receives an error event or an error message. The application may query the error code with function [ThermoUSBGetError\(\)](#). If the device is a HID device, the error code may be any of the Windows error codes. In case of a libusb-win32 device, the error code from libusb driver is translated to one of the following Windows error codes:

Windows error	Code	Description
ERROR_SUCCESS	0	No error.
ERROR_ACCESS_DENIED	5	I/O error.
ERROR_NOT_ENOUGH_MEMORY	12	Not enough memory.
ERROR_BAD_COMMAND	22	Invalid parameter.
WAIT_TIMEOUT	116	A transfer timed out.

What the application can do in case of an error is to first call [ThermoUSBClose\(\)](#) and then try to reopen with [ThermoUSBOpen\(\)](#).

Definition at line 47 of file ThermoUSB.h.

4.4.2.2 #define THERMOUSB_RECEIVE 1

Definition at line 18 of file ThermoUSB.h.

4.4.2.3 #define THERMOUSB_TRANSMIT 2

Definition at line 19 of file ThermoUSB.h.

4.4.3 Function Documentation

4.4.3.1 DllExport void WINAPI ThermoUSBAbort (HANDLE *hConn*)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
--------------	---

Definition at line 2241 of file ThermoUSB.c.

4.4.3.2 DllExport BOOL WINAPI ThermoUSBAttachEvent (HANDLE *hConn*, UINT *evCode*, HANDLE *object*)

Parameters

<i>hConn</i>	Connection handle returned form a call to ThermoUSBOpen() (pointer to a USB_CONNECTION structure).
<i>evCode</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Parameter 'evCode' may be any combination of THERMOUSB_RECEIVE, THERMOUSB_TRANSMIT and THERMOUSB_ERROR. The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event.

The THERMOUSB_RECEIVE event is signaled whenever something is written to the receive buffer. There is no quarantee that a whole response line is received or that only one response line is received. therefore upon receiving this event the application should call [ThermoUSBRead\(\)](#) or ThermoUSBReadBinary as long as they return data.

The THERMOUSB_TRANSMIT event is signaled when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOUSB_TRANSMIT event is received, functions [ThermoUSBWrite\(\)](#) and [ThermoUSBWriteBinary\(\)](#) will never fail.

Definition at line 1853 of file ThermoUSB.c.

4.4.3.3 DllExport BOOL WINAPI ThermoUSBAttachMsg (HANDLE *hConn*, HANDLE *hWnd*, UINT *msg*, UINT *evCode*)

Parameters

<i>hConn</i>	Connection handle returned form a call to ThermoUSBOpen() (pointer to a USB_CONNECTION structure).
<i>hWnd</i>	Handle of the window to receive the message.
<i>msg</i>	Message id of the message to send.
<i>event</i>	Event(s) for which a message is sent.

A message with message id 'msg' is sent to window 'hWnd' whenever any of the event(s) selected with parameter 'evCode' occurs. Parameter 'evCode' may be any combination of THERMOUSB_RECEIVE, THERMOUSB_TRANSMIT and THERMOUSB_ERROR. The event code is passed in the wParam member of the sent Windows message.

The THERMOUSB_RECEIVE event is sent whenever something is written to the receive buffer. There is no quarantee that a whole response line is received or that only one response line is received. Therefore upon receiving this event the application should call [ThermoUSBRead\(\)](#) or ThermoUSBReadBinary as long as they return data.

The THERMOUSB_TRANSMIT event is sent when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOUSB_TRANSMIT event is received, functions [ThermoUSBWrite\(\)](#) and [ThermoUSBWriteBinary\(\)](#) will never fail.

Definition at line 1908 of file ThermoUSB.c.

4.4.3.4 DllExport void WINAPI ThermoUSBClose (HANDLE *hConn*)

Parameters

<i>hConn</i>	Connection handle returned from a call to ThermoUSBOpen()
--------------	---

Closes the communication channel to the instrument. This function should be called when the application has finished communicating with the instrument. It is not possible to open a communication channel to the instrument while the previous channel is still open. When the dll is unloaded from the memory, all channels still open are automatically closed.

Definition at line 1767 of file ThermoUSB.c.

4.4.3.5 DllExport DWORD WINAPI ThermoUSBGetError (HANDLE *hConn*)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
--------------	---

Returns

The current Windows error code stored for the connection.

Error conditions may be encountered asynchronously and not just as a direct result of a function call to the dll. Therefore the application should enable error reporting with [ThermoUSBAttachEvent\(\)](#) or [ThermoUSBAttachMsg\(\)](#). When an error is reported to the application, it may call this function to get the Windows error code, which may or may not be helpful in resolving the problem. In general, if an error is reported then some receive or transmit data is lost, and the best way of action is to close the port and then retry opening it again.

A call to [ThermoUSBGetError\(\)](#) resets the stored error code to ERROR_SUCCESS.

Definition at line 2308 of file ThermoUSB.c.

4.4.3.6 DllExport WORD WINAPI ThermoUSBGetThermoVendorId (void)

Returns

Thermo Fisher Scientific Oy USB vendor id, 0x0AB6.

Definition at line 2332 of file ThermoUSB.c.

4.4.3.7 DllExport HANDLE WINAPI ThermoUSBOpen (WORD *VendorID*, WORD *ProductID*, LPCSTR *SerialNumber*)

Parameters

<i>VendorID</i>	The USB vendor id of the device manufacturer, 0x0AB6 for Thermo Fisher Scientific Oy.
<i>ProductID</i>	Product id number of the device.
<i>SerialNumber</i>	Pointer to the serial number string of the device. The device must report an identical serial number for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first device with matching VendorID and ProductID.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

This function (or ThermoUSBOpenSimulator) must be called first before using any other functions in the library. Only one connection per device is allowed.

Definition at line 1738 of file ThermoUSB.c.

4.4.3.8 DllExport BOOL WINAPI ThermoUSBRead (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
<i>buf</i>	Pointer to a buffer to receive the response.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

TRUE if a response line was retrieved, else FALSE.

Use this function to read full response lines returned from the instrument. The response line is returned without the terminating CRLF. Therefore TRUE may be returned even if the resulting line is empty. The returned line always ends to the NUL character even if FALSE is returned.

If the caller's buffer is too small to hold the whole response line, as much as fits to the buffer is returned and the rest is returned on subsequent call(s).

Definition at line 1948 of file ThermoUSB.c.

4.4.3.9 DllExport DWORD WINAPI ThermoUSBReadBinary (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
<i>buf</i>	Pointer to buffer receiving the data.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

Number of bytes copied to the user buffer.

Unlike function [ThermoUSBRead\(\)](#), this function returns all data received from the instrument without doing any interpretation on it. You can control the number of bytes returned with the bufsize parameter.

If the return value is less than the given bufsize parameter, there is no more data.

Definition at line 2123 of file ThermoUSB.c.

4.4.3.10 DllExport BOOL WINAPI ThermoUSBWrite (HANDLE *hConn*, LPCSTR *buf*)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
<i>buf</i>	Buffer containing the ASCII string to send.

Returns

TRUE if the string was written to the transmit buffer, else FALSE.

If the whole string does not fit in the transmit buffer, writes nothing and returns FALSE. The application may retry at a later time.

Definition at line 2169 of file ThermoUSB.c.

4.4.3.11 DllExport BOOL WINAPI ThermoUSBWriteBinary (HANDLE *hConn*, LPCSTR *buf*, DWORD *count*)

Parameters

<i>hConn</i>	Handle of the USB connection. This must be a handle returned from ThermoUSBOpen() .
--------------	---

<i>buf</i>	Buffer containing the data to send.
<i>count</i>	Number of bytes to send.

Returns

TRUE if the data was written to the transmit buffer, else FALSE.

If all data does not fit in the transmit buffer, writes nothing and returns FALSE. The application may retry at a later time.

Definition at line 2190 of file ThermoUSB.c.

Index

api.h, [19](#)

mainpage.h, [19](#)

THERMOUSB_ERROR
 ThermoUSB.h, [24](#)

THERMOUSB_RECEIVE
 ThermoUSB.h, [24](#)

THERMOUSB_TRANSMIT
 ThermoUSB.h, [24](#)

ThermoUSB.c, [19](#)
 ThermoUSBAbort, [20](#)
 ThermoUSBAttachEvent, [20](#)
 ThermoUSBAttachMsg, [20](#)
 ThermoUSBClose, [21](#)
 ThermoUSBGetError, [21](#)
 ThermoUSBGetThermoVendorId, [21](#)
 ThermoUSBOpen, [21](#)
 ThermoUSBRead, [22](#)
 ThermoUSBReadBinary, [22](#)
 ThermoUSBWrite, [22](#)
 ThermoUSBWriteBinary, [23](#)

ThermoUSB.h, [23](#)
 THERMOUSB_ERROR, [24](#)
 THERMOUSB_RECEIVE, [24](#)
 ThermoUSBAbort, [25](#)
 ThermoUSBAttachEvent, [25](#)
 ThermoUSBAttachMsg, [25](#)
 ThermoUSBClose, [25](#)
 ThermoUSBGetError, [26](#)
 ThermoUSBGetThermoVendorId, [26](#)
 ThermoUSBOpen, [26](#)
 ThermoUSBRead, [26](#)
 ThermoUSBReadBinary, [27](#)
 ThermoUSBWrite, [27](#)
 ThermoUSBWriteBinary, [27](#)

ThermoUSBAbort
 ThermoUSB.c, [20](#)
 ThermoUSB.h, [25](#)

ThermoUSBAttachEvent
 ThermoUSB.c, [20](#)
 ThermoUSB.h, [25](#)

ThermoUSBAttachMsg
 ThermoUSB.c, [20](#)
 ThermoUSB.h, [25](#)

ThermoUSBClose
 ThermoUSB.c, [21](#)
 ThermoUSB.h, [25](#)

ThermoUSBGetError
 ThermoUSB.c, [21](#)
 ThermoUSB.h, [26](#)

ThermoUSBGetThermoVendorId
 ThermoUSB.c, [21](#)
 ThermoUSB.h, [26](#)

ThermoUSBOpen
 ThermoUSB.c, [21](#)
 ThermoUSB.h, [26](#)

ThermoUSBRead
 ThermoUSB.c, [22](#)
 ThermoUSB.h, [26](#)

ThermoUSBReadBinary
 ThermoUSB.c, [22](#)
 ThermoUSB.h, [27](#)

ThermoUSBWrite
 ThermoUSB.c, [22](#)
 ThermoUSB.h, [27](#)

ThermoUSBWriteBinary
 ThermoUSB.c, [23](#)
 ThermoUSB.h, [27](#)

Thermo Scientific

ThermoLAN.dII

Interface Specification

Contents

1	ThermoLAN dll Interface Specification	1
2	Using ThermoLAN	3
2.1	About	3
2.2	Exported functions	4
2.2.1	ThermoLANListDevices	5
2.2.2	ThermoLANOpen	6
2.2.3	ThermoLANClose	7
2.2.4	ThermoLANWrite	8
2.2.5	ThermoLANRead	9
2.2.6	ThermoLANAttachEvent	10
2.2.7	ThermoLANAbort	11
2.2.8	ThermoLANGetError	12
3	File Index	13
3.1	File List	13
4	File Documentation	15
4.1	api.h File Reference	15
4.2	mainpage.h File Reference	15
4.3	ThermoLANWrapper.cpp File Reference	15
4.3.1	Detailed Description	16
4.3.2	Function Documentation	16
4.3.2.1	ThermoLANAbort	16
4.3.2.2	ThermoLANAttachEvent	16
4.3.2.3	ThermoLANClose	17
4.3.2.4	ThermoLANGetError	17
4.3.2.5	ThermoLANListDevices	17
4.3.2.6	ThermoLANOpen	18
4.3.2.7	ThermoLANRead	18
4.3.2.8	ThermoLANWrite	18
4.4	ThermoLANWrapper.h File Reference	19
4.4.1	Detailed Description	19
4.4.2	Function Documentation	19
4.4.2.1	ThermoLANAbort	19
4.4.2.2	ThermoLANAttachEvent	20
4.4.2.3	ThermoLANClose	20
4.4.2.4	ThermoLANGetError	20
4.4.2.5	ThermoLANListDevices	21
4.4.2.6	ThermoLANOpen	21
4.4.2.7	ThermoLANRead	22
4.4.2.8	ThermoLANWrite	22

Chapter 1

ThermoLAN dll Interface Specification

About

The purpose of the ThermoLAN.dll dynamic link library is to make interface to the Thermo microplate instruments easy and to offer a similar interface to the instrument as the ThermoUSB.dll library offers. The user of the ThermoLAN.dll does not have to know the details of serial port communication.

Confidential

This document has been prepared by Thermo Fisher Scientific Oy to be used solely for the purposes defined by Thermo Fisher Scientific Oy. Use for other purposes is not authorized.

Please note that any and all information contained in this document is the property of Thermo Fisher Scientific Oy. This confidential information ("Confidential Information") shall not be reproduced in whole part or disclosed to any third party without the prior written approval of Thermo Fisher Scientific Oy. The receiving party shall ensure that its employees, officers, representatives and agents shall not disclose to third parties any Confidential Information.

Upon written request from Thermo Fisher Scientific Oy, the receiving party shall promptly return all Confidential Information or destroy all Confidential Information.

Chapter 2

Using ThermoLAN

2.1 About

ThermoLAN.dll is a generic library for communicating with several different Thermo Scientific microplate instruments via a Ethernet port. To communicate with an instrument you use the exported functions of ThermoLAN.dll.

Depending on your project setup, you may find useful a couple of other files which are also provided. The header file ThermoLAN.h contains the prototypes of the exported functions and definitions of constant values used by the dll. File ThermoLAN.lib contains information about the dll the linker uses to add references to the library in the executable. This way the dll is automatically loaded and the exported functions of the library can be called as easy as the functions in the code using the dll.

2.2 Exported functions

- [ThermoLANListDevices](#)
- [ThermoLANOpen](#)
- [ThermoLANClose](#)
- [ThermoLANWrite](#)
- [ThermoLANRead](#)
- [ThermoLANAttachEvent](#)
- [ThermoLANAbort](#)
- [ThermoLANGetError](#)

2.2.1 ThermoLANListDevices

Search for existing instruments. Full declaration: [ThermoLANListDevices\(\)](#).

Parameters

<i>instrumentName</i>	Name of the instrument. Only instruments with a matching name are returned in the list. May be NULL, which matches all instruments.
<i>serialNumber</i>	Pointer to the serial number string of the instrument. Only instruments with a matching serial number are returned. May be NULL, which matches any serial number.
<i>buf</i>	Buffer to list found instruments to.
<i>bufSize</i>	Size of the buffer.

Returns

The size in bytes of the complete list of found devices. If the returned value is equal or smaller than the given buffer size, the buffer contains the complete list of devices found on the LAN. If the returned value is higher than the given buffer size, no data is returned and the caller must call the function again with big enough buffer. Return value 0 means that no instruments were found.

On success, the caller's buffer contains zero terminated strings with a combined length of the return value. The string terminating zeros are included in the length.

For each found instrument, the first string is the instrument IPv4 address and the TCP port number it is listening, enclosed in square brackets, e.g. [10.32.196.210:49536].

The IP address string is followed by the WS-Discovery match strings, usually 3 of them. The first one is always 'Thermo-Device', the second one is the instrument name and the third one the instrument serial number string.

If the match strings are followed by a string in angle brackets, e.g. <10.32.196.154:57403>, it means that the instrument is currently connected that IP address and TCP port, and trying to connect to that instrument with function [ThermoLANOpen\(\)](#) will fail. If there is no string in square brackets, the instrument will accept a connection.

2.2.2 ThermoLANOpen

Search for the requested instrument and open a communication channel to it. Full declaration: [ThermoLANOpen\(\)](#).

Parameters

<i>instrumentName</i>	Name of the instrument.
<i>serialNumber</i>	Pointer to the serial number string of the instrument. This must match with the actual serial number of the instrument for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first device with matching instrument name.
<i>timeout</i>	Timeout which is used to search the instrument. If 0 is given then WS-Discovery will use default 4 sec timeout.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

This function must be called first before using any other functions in the library. Only one connection per instrument is allowed.

To find the requested instrument on the LAN, the WS-Discovery protocol is used. The instrument responds with a Probe Match to a Probe message if all Target Service strings of the <Types> element of the Probe match the Target Service strings of the instrument. One of the strings, "ThermoDevice", is common to all instruments. Other strings are the instrument name string and the serial number string. Due to Windows limitations, a "SN_" prefix is added to the serial number string.

The <XAddr> element of the Probe Match response contains the IP address of the instrument and the TCP port number it is listening to.

2.2.3 ThermoLANClose

Close an LAN connection. Full declaration: [ThermoLANClose\(\)](#).

Parameters

<i>hConn</i>	Connection handle returned form a call to ThermoLANOpen()
--------------	---

Closes the communication channel to the instrument. This function should be called when the application has finished communicating with the instrument. It is not possible to open a communication channel to the instrument while the previous channel is still open. When the dll is unloaded from the memory, all channels still open are automatically closed.

2.2.4 ThermoLANWrite

Write a string to the transmit buffer. Full declaration: [ThermoLANWrite\(\)](#).

Parameters

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
<i>buf</i>	Buffer containing the ASCII string to send.

Returns

TRUE if the string was written to the transmit buffer, else FALSE.

If the whole string does not fit in the transmit buffer, writes nothing and returns FALSE. If FALSE is returned, the application can call function [ThermoLANGetError\(\)](#) to determine if the transmit buffer was full or if some other error occurred. In case of transmit buffer full, [ThermoLANGetError\(\)](#) returns ERROR_SUCCESS and the application may retry sending at a later time.

2.2.5 ThermoLANRead

Read a received response line. Full declaration: [ThermoLANRead\(\)](#).

Parameters

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
<i>buf</i>	Pointer to a buffer to receive the response.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

TRUE if a response line was retrieved, else FALSE.

Use this function to read full response lines returned from the instrument. The response line is returned without the terminating CRLF. Therefore TRUE may be returned even if the resulting line is empty. The returned line always ends to the NUL character even if FALSE is returned.

If the caller's buffer is too small to hold the whole response line, as much as fits to the buffer is returned and the rest is returned on subsequent call(s).

2.2.6 ThermoLANAttachEvent

Attach an event object to a LAN connection. Full declaration: [ThermoLANAttachEvent\(\)](#).

Parameters

<i>hConn</i>	Connection handle returned from a call to ThermoLANOpen() .
<i>evCode</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Returns

TRUE on success, else FALSE.

Parameter 'evCode' may be any combination of THERMOLAN_RECEIVE, THERMOLAN_TRANSMIT and THERMOLAN_ERROR. The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event.

The THERMOLAN_RECEIVE event is signaled whenever something is written to the receive buffer. There is no guarantee that a whole response line is received or that only one response line is received. Therefore upon receiving this event the application should repeatedly call [ThermoLANRead\(\)](#) as long it returns TRUE.

The THERMOLAN_TRANSMIT event is signaled when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOLAN_TRANSMIT event is received, function [ThermoLANWrite\(\)](#) will never fail.

The THERMOLAN_ERROR event is signaled when the dll detects an error, for example trying to send data when the TCP connection has been closed by the instrument side. If the application wants to continue communication with the instrument after an error, it should close and reopen the connection.

2.2.7 ThermoLANAbort

Send Abort command to the instrument. Full declaration: [ThermoLANAbort\(\)](#).

Parameters

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
--------------	---

The LAN Abort sequence is this:

- Send an UDP message containing text "Abort" (without quotes or newline) to the same UDP port number as is used for the TCP and wait a moment for an identical response from the instrument. The response is sent over the same UDP port.
- Retry two times more at half second interval if no response.
- If a response received or no response after retries, send Abort character (0x1B) to the TCP port.

2.2.8 ThermoLANGetError

Get the last error code recorded for the communication channel. Full declaration: [ThermoLANGetError\(\)](#).

Parameters

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
--------------	---

Returns

One of the Windows system error codes, ERROR_SUCCESS if no error has occurred.

The recorded error code is automatically cleared when this function is called. ERROR_INVALID_HANDLE is returned if the connection handle is invalid.

Chapter 3

File Index

3.1 File List

Here is a list of all files with brief descriptions:

api.h	Part of ThermoLAN dll documentation	15
mainpage.h	Start page of ThermoLAN.dll Design Description	15
ThermoLANWrapper.cpp	Implements ethernet interface to Thermo microplate instruments	15
ThermoLANWrapper.h	Functions exported from ThermoLAN.dll	19

Chapter 4

File Documentation

4.1 api.h File Reference

Part of ThermoLAN dll documentation.

4.2 mainpage.h File Reference

Start page of ThermoLAN.dll Design Description.

4.3 ThermoLANWrapper.cpp File Reference

Implements ethernet interface to Thermo microplate instruments.

Functions

- void * [ThermoLANOpen](#) (char *instrumentName, char *serialNumber, int timeout)
Search for the requested instrument and open a communication channel to it.
- DWORD [ThermoLANListDevices](#) (char *instrumentName, char *serialNumber, char *buf, unsigned long bufSize)
Search for existing instruments.
- void [ThermoLANClose](#) (void *hConn)
Close an LAN connection.
- BOOL __stdcall [ThermoLANWrite](#) (void *hConn, const char *str)
Write a string to the transmit buffer.
- BOOL __stdcall [ThermoLANRead](#) (void *hConn, char *str, unsigned long bufSize)
Read a received response line.
- BOOL __stdcall [ThermoLANAttachEvent](#) (void *hConn, unsigned int evCode, void *receiver)
Attach an event object to a LAN connection.
- void [ThermoLANAbort](#) (void *hConn)
Send Abort command to the instrument.
- DWORD [ThermoLANGetError](#) (void *hConn)
Get the last error code recorded for the communication channel.

4.3.1 Detailed Description

Note

Copyright by Thermo Fisher Scientific Oy 2015

Definition in file [ThermoLANWrapper.cpp](#).

4.3.2 Function Documentation

4.3.2.1 void ThermoLANAbort (void * *hConn*)

Parameters

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
--------------	---

The LAN Abort sequence is this:

- Send an UDP message containing text "Abort" (without quotes or newline) to the same UDP port number as is used for the TCP and wait a moment for an identical response from the instrument. The response is sent over the same UDP port.
- Retry two times more at half second interval if no response.
- If a response received or no response after retries, send Abort character (0x1B) to the TCP port.

Definition at line 379 of file ThermoLANWrapper.cpp.

4.3.2.2 BOOL __stdcall ThermoLANAttachEvent (void * *hConn*, unsigned int *evCode*, void * *receiver*)

Parameters

<i>hConn</i>	Connection handle returned form a call to ThermoLANOpen() .
<i>evCode</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Returns

TRUE on success, else FALSE.

Parameter 'evCode' may be any combination of THERMOLAN_RECEIVE, THERMOLAN_TRANSMIT and THERMOLAN_ERROR. The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event.

The THERMOLAN_RECEIVE event is signaled whenever something is written to the receive buffer. There is no quarantee that a whole response line is received or that only one response line is received. Therefore upon receiving this event the application should repeatedly call [ThermoLANRead\(\)](#) as long it returns TRUE.

The THERMOLAN_TRANSMIT event is signaled when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOLAN_TRANSMIT event is received, function [ThermoLANWrite\(\)](#) will never fail.

The THERMOLAN_ERROR event is signaled when the dll detects an error, for example trying to send data when the TCP connection has been closed by the instrument side. If the application wants to continue communication with the instrument after an error, it should close and reopen the connection.

Definition at line 354 of file ThermoLANWrapper.cpp.

4.3.2.3 void ThermoLANClose (void * *hConn*)**Parameters**

<i>hConn</i>	Connection handle returned from a call to ThermoLANOpen()
--------------	---

Closes the communication channel to the instrument. This function should be called when the application has finished communicating with the instrument. It is not possible to open a communication channel to the instrument while the previous channel is still open. When the dll is unloaded from the memory, all channels still open are automatically closed.

Definition at line 234 of file ThermoLANWrapper.cpp.

4.3.2.4 DWORD ThermoLANGetError (void * *hConn*)**Parameters**

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
--------------	---

Returns

One of the Windows system error codes, ERROR_SUCCESS if no error has occurred.

The recorded error code is automatically cleared when this function is called. ERROR_INVALID_HANDLE is returned if the connection handle is invalid.

Definition at line 400 of file ThermoLANWrapper.cpp.

4.3.2.5 DWORD ThermoLANListDevices (char * *instrumentName*, char * *serialNumber*, char * *buf*, unsigned long *bufSize*)**Parameters**

<i>instrumentName</i>	Name of the instrument. Only instruments with a matching name are returned in the list. May be NULL, which matches all instruments.
<i>serialNumber</i>	Pointer to the serial number string of the instrument. Only instruments with a matching serial number are returned. May be NULL, which matches any serial number.
<i>buf</i>	Buffer to list found instruments to.
<i>bufSize</i>	Size of the buffer.

Returns

The size in bytes of the complete list of found devices. If the returned value is equal or smaller than the given buffer size, the buffer contains the complete list of devices found on the LAN. If the returned value is higher than the given buffer size, no data is returned and the caller must call the function again with big enough buffer. Return value 0 means that no instruments were found.

On success, the caller's buffer contains zero terminated strings with a combined length of the return value. The string terminating zeros are included in the length.

For each found instrument, the first string is the instrument IPv4 address and the TCP port number it is listening, enclosed in square brackets, e.g. [10.32.196.210:49536].

The IP address string is followed by the WS-Discovery match strings, usually 3 of them. The first one is always 'Thermo-Device', the second one is the instrument name and the third one the instrument serial number string.

If the match strings are followed by a string in angle brackets, e.g. <10.32.196.154:57403>, it means that the instrument is currently connected that IP address and TCP port, and trying to connect to that instrument with function [ThermoLANOpen\(\)](#) will fail. If there is no string in square brackets, the instrument will accept a connection.

Definition at line 213 of file ThermoLANWrapper.cpp.

4.3.2.6 void* ThermoLANOpen (char * *instrumentName*, char * *serialNumber*, int *timeout*)

Parameters

<i>instrumentName</i>	Name of the instrument.
<i>serialNumber</i>	Pointer to the serial number string of the instrument. This must match with the actual serial number of the instrument for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first device with matching instrument name.
<i>timeout</i>	Timeout which is used to search the instrument. If 0 is given then WS-Discovery will use default 4 sec timeout.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

This function must be called first before using any other functions in the library. Only one connection per instrument is allowed.

To find the requested instrument on the LAN, the WS-Discovery protocol is used. The instrument responds with a Probe Match to a Probe message if all Target Service strings of the <Types> element of the Probe match the Target Service strings of the instrument. One of the strings, "ThermoDevice", is common to all instruments. Other strings are the instrument name string and the serial number string. Due to Windows limitations, a "SN_" prefix is added to the serial number string.

The <XAddr> element of the Probe Match response contains the IP address of the instrument and the TCP port number it is listening to.

Definition at line 161 of file ThermoLANWrapper.cpp.

4.3.2.7 BOOL __stdcall ThermoLANRead (void * *hConn*, char * *str*, unsigned long *bufSize*)

Parameters

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
<i>buf</i>	Pointer to a buffer to receive the response.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

TRUE if a response line was retrieved, else FALSE.

Use this function to read full response lines returned from the instrument. The response line is returned without the terminating CRLF. Therefore TRUE may be returned even if the resulting line is empty. The returned line always ends to the NUL character even if FALSE is returned.

If the caller's buffer is too small to hold the whole response line, as much as fits to the buffer is returned and the rest is returned on subsequent call(s).

Definition at line 299 of file ThermoLANWrapper.cpp.

4.3.2.8 BOOL __stdcall ThermoLANWrite (void * *hConn*, const char * *str*)

Parameters

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
<i>buf</i>	Buffer containing the ASCII string to send.

Returns

TRUE if the string was written to the transmit buffer, else FALSE.

If the whole string does not fit in the transmit buffer, writes nothing and returns FALSE. If FALSE is returned, the application can call function [ThermoLANGetError\(\)](#) to determine if the transmit buffer was full or is some other error occurred. In case of transmit buffer full, [ThermoLANGetError\(\)](#) returns ERROR_SUCCESS and the application may retry sending at a later time.

Definition at line 260 of file ThermoLANWrapper.cpp.

4.4 ThermoLANWrapper.h File Reference

Functions exported from ThermoLAN.dll.

Functions

- void * [ThermoLANOpen](#) (char *instrumentName, char *serialNumber, int timeout)
Search for the requested instrument and open a communication channel to it.
- DWORD [ThermoLANListDevices](#) (char *instrumentName, char *serialNumber, char *buf, unsigned long bufSize)
Search for existing instruments.
- void [ThermoLANClose](#) (void *hConn)
Close an LAN connection.
- BOOL __stdcall [ThermoLANWrite](#) (void *hConn, const char *str)
Write a string to the transmit buffer.
- BOOL __stdcall [ThermoLANRead](#) (void *hConn, char *str, unsigned long bufSize)
Read a received response line.
- BOOL __stdcall [ThermoLANAttachEvent](#) (void *hConn, unsigned int evCode, void *receiver)
Attach an event object to a LAN connection.
- void [ThermoLANAbort](#) (void *hConn)
Send Abort command to the instrument.
- DWORD [ThermoLANGetError](#) (void *hConn)
Get the last error code recorded for the communication channel.

4.4.1 Detailed Description

Note

Copyright by Thermo Fisher Scientific Oy 2015

Ethernet interface to Thermo microplate instruments.

Definition in file [ThermoLANWrapper.h](#).

4.4.2 Function Documentation

4.4.2.1 void ThermoLANAbort (void * hConn)

Parameters

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
--------------	---

The LAN Abort sequence is this:

- Send an UDP message containing text "Abort" (without quotes or newline) to the same UDP port number as is used for the TCP and wait a moment for an identical response from the instrument. The response is sent over the same

UDP port.

- Retry two times more at half second interval if no response.
- If a response received or no response after retries, send Abort character (0x1B) to the TCP port.

Definition at line 379 of file ThermoLANWrapper.cpp.

4.4.2.2 **BOOL __stdcall ThermoLANAttachEvent (void * *hConn*, unsigned int *evCode*, void * *receiver*)**

Parameters

<i>hConn</i>	Connection handle returned form a call to ThermoLANOpen() .
<i>evCode</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Returns

TRUE on success, else FALSE.

Parameter 'evCode' may be any combination of THERMOLAN_RECEIVE, THERMOLAN_TRANSMIT and THERMOLAN_ERROR. The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event.

The THERMOLAN_RECEIVE event is signaled whenever something is written to the receive buffer. There is no quarantee that a whole response line is received or that only one response line is received. Therefore upon receiving this event the application should repeatedly call [ThermoLANRead\(\)](#) as long it returns TRUE.

The THERMOLAN_TRANSMIT event is signaled when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOLAN_TRANSMIT event is received, function [ThermoLANWrite\(\)](#) will never fail.

The THERMOLAN_ERROR event is signaled when the dll detects an error, for example trying to send data when the TCP connection has been closed by the instrument side. If the application wants to continue communication with the instrument after an error, it should close and reopen the connection.

Definition at line 354 of file ThermoLANWrapper.cpp.

4.4.2.3 **void ThermoLANClose (void * *hConn*)**

Parameters

<i>hConn</i>	Connection handle returned form a call to ThermoLANOpen()
--------------	---

Closes the communication channel to the instrument. This function should be called when the application has finished communicating with the instrument. It is not possible to open a communication channel to the instrument while the previous channel is still open. When the dll is unloaded from the memory, all channels still open are automatically closed.

Definition at line 234 of file ThermoLANWrapper.cpp.

4.4.2.4 **DWORD ThermoLANGetError (void * *hConn*)**

Parameters

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
--------------	---

Returns

One of the Windows system error codes, ERROR_SUCCESS if no error has occurred.

The recorded error code is automatically cleared when this function is called. ERROR_INVALID_HANDLE is returned if the connection handle is invalid.

Definition at line 400 of file ThermoLANWrapper.cpp.

4.4.2.5 DWORD ThermoLANListDevices (char * *instrumentName*, char * *serialNumber*, char * *buf*, unsigned long *bufSize*)**Parameters**

<i>instrumentName</i>	Name of the instrument. Only instruments with a matching name are returned in the list. May be NULL, which matches all instruments.
<i>serialNumber</i>	Pointer to the serial number string of the instrument. Only instruments with a matching serial number are returned. May be NULL, which matches any serial number.
<i>buf</i>	Buffer to list found instruments to.
<i>bufSize</i>	Size of the buffer.

Returns

The size in bytes of the complete list of found devices. If the returned value is equal or smaller than the given buffer size, the buffer contains the complete list of devices found on the LAN. If the returned value is higher than the given buffer size, no data is returned and the caller must call the function again with big enough buffer. Return value 0 means that no instruments were found.

On success, the caller's buffer contains zero terminated strings with a combined length of the return value. The string terminating zeros are included in the length.

For each found instrument, the first string is the instrument IPv4 address and the TCP port number it is listening, enclosed in square brackets, e.g. [10.32.196.210:49536].

The IP address string is followed by the WS-Discovery match strings, usually 3 of them. The first one is always 'Thermo-Device', the second one is the instrument name and the third one the instrument serial number string.

If the match strings are followed by a string in angle brackets, e.g. <10.32.196.154:57403>, it means that the instrument is currently connected that IP address and TCP port, and trying to connect to that instrument with function [ThermoLANOpen\(\)](#) will fail. If there is no string in square brackets, the instrument will accept a connection.

Definition at line 213 of file ThermoLANWrapper.cpp.

4.4.2.6 void* ThermoLANOpen (char * *instrumentName*, char * *serialNumber*, int *timeout*)**Parameters**

<i>instrumentName</i>	Name of the instrument.
<i>serialNumber</i>	Pointer to the serial number string of the instrument. This must match with the actual serial number of the instrument for the connection to succeed. This parameter may be NULL, in which case the connection is made to the first devicee with matching instrument name.
<i>timeout</i>	Timeout which is used to search the instrument. If 0 is given then WS-Discovery will use default 4 sec timeout.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

This function must be called first before using any other functions in the library. Only one connection per instrument is allowed.

To find the requested instrument on the LAN, the WS-Discovery protocol is used. The instrument responds with a Probe Match to a Probe message if all Target Service strings of the <Types> element of the Probe match the Target Service strings of the instrument. One of the strings, "ThermoDevice", is common to all instruments. Other strings are the instrument name string and the serial number string. Due to Windows limitations, a "SN_" prefix is added to the serial number string.

The <XAddr> element of the Probe Match response contains the IP address of the instrument and the TCP port number it is listening to.

Definition at line 161 of file ThermoLANWrapper.cpp.

4.4.2.7 **BOOL** __stdcall ThermoLANRead (void * *hConn*, char * *str*, unsigned long *bufSize*)

Parameters

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
<i>buf</i>	Pointer to a buffer to receive the response.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

TRUE if a response line was retrieved, else FALSE.

Use this function to read full response lines returned from the instrument. The response line is returned without the terminating CRLF. Therefore TRUE may be returned even if the resulting line is empty. The returned line always ends to the NUL character even if FALSE is returned.

If the caller's buffer is too small to hold the whole response line, as much as fits to the buffer is returned and the rest is returned on subsequent call(s).

Definition at line 299 of file ThermoLANWrapper.cpp.

4.4.2.8 **BOOL** __stdcall ThermoLANWrite (void * *hConn*, const char * *str*)

Parameters

<i>hConn</i>	Handle of the LAN connection. This must be a handle returned from ThermoLANOpen() .
<i>buf</i>	Buffer containing the ASCIIZ string to send.

Returns

TRUE if the string was written to the transmit buffer, else FALSE.

If the whole string does not fit in the transmit buffer, writes nothing and returns FALSE. If FALSE is returned, the application can call function [ThermoLANGetError\(\)](#) to determine if the transmit buffer was full or is some other error occurred. In case of transmit buffer full, [ThermoLANGetError\(\)](#) returns ERROR_SUCCESS and the application may retry sending at a later time.

Definition at line 260 of file ThermoLANWrapper.cpp.

Index

[api.h](#), [15](#)

[mainpage.h](#), [15](#)

ThermoLANAbort

[ThermoLANWrapper.cpp](#), [16](#)

[ThermoLANWrapper.h](#), [19](#)

ThermoLANAttachEvent

[ThermoLANWrapper.cpp](#), [16](#)

[ThermoLANWrapper.h](#), [20](#)

ThermoLANClose

[ThermoLANWrapper.cpp](#), [16](#)

[ThermoLANWrapper.h](#), [20](#)

ThermoLANGetError

[ThermoLANWrapper.cpp](#), [17](#)

[ThermoLANWrapper.h](#), [20](#)

ThermoLANListDevices

[ThermoLANWrapper.cpp](#), [17](#)

[ThermoLANWrapper.h](#), [21](#)

ThermoLANOpen

[ThermoLANWrapper.cpp](#), [17](#)

[ThermoLANWrapper.h](#), [21](#)

ThermoLANRead

[ThermoLANWrapper.cpp](#), [18](#)

[ThermoLANWrapper.h](#), [22](#)

[ThermoLANWrapper.cpp](#), [15](#)

 ThermoLANAbort, [16](#)

 ThermoLANAttachEvent, [16](#)

 ThermoLANClose, [16](#)

 ThermoLANGetError, [17](#)

 ThermoLANListDevices, [17](#)

 ThermoLANOpen, [17](#)

 ThermoLANRead, [18](#)

 ThermoLANWrite, [18](#)

[ThermoLANWrapper.h](#), [19](#)

 ThermoLANAbort, [19](#)

 ThermoLANAttachEvent, [20](#)

 ThermoLANClose, [20](#)

 ThermoLANGetError, [20](#)

 ThermoLANListDevices, [21](#)

 ThermoLANOpen, [21](#)

 ThermoLANRead, [22](#)

 ThermoLANWrite, [22](#)

ThermoLANWrite

[ThermoLANWrapper.cpp](#), [18](#)

[ThermoLANWrapper.h](#), [22](#)

Thermo Scientific

ThermoCOM.dll

Interface Specification

Contents

1	ThermoCOM dll Interface Specification	1
2	Using ThermoCOM	3
2.1	About	3
2.2	Exported functions	4
2.2.1	ThermoCOMOpen	5
2.2.2	ThermoCOMOpenSimulator	6
2.2.3	ThermoCOMClose	7
2.2.4	ThermoCOMSetParam	8
2.2.5	ThermoCOMAttachEvent	9
2.2.6	ThermoCOMAttachMsg	10
2.2.7	ThermoCOMRead	11
2.2.8	ThermoCOMReadBinary	12
2.2.9	ThermoCOMWrite	13
2.2.10	ThermoCOMWriteBinary	14
2.2.11	ThermoCOMGetError	15
2.2.12	ThermoCOMAbort	16
3	File Index	17
3.1	File List	17
4	File Documentation	19
4.1	api.h File Reference	19
4.2	mainpage.h File Reference	19
4.3	ThermoCOM.c File Reference	19
4.3.1	Detailed Description	20
4.3.2	Function Documentation	20
4.3.2.1	ThermoCOMAbort	20
4.3.2.2	ThermoCOMAttachEvent	20
4.3.2.3	ThermoCOMAttachMsg	21
4.3.2.4	ThermoCOMClose	21
4.3.2.5	ThermoCOMGetError	21
4.3.2.6	ThermoCOMOpen	22
4.3.2.7	ThermoCOMOpenSimulator	22
4.3.2.8	ThermoCOMRead	22
4.3.2.9	ThermoCOMReadBinary	23
4.3.2.10	ThermoCOMSetParam	23
4.3.2.11	ThermoCOMWrite	23
4.3.2.12	ThermoCOMWriteBinary	24
4.4	ThermoCOM.h File Reference	24
4.4.1	Detailed Description	25
4.4.2	Macro Definition Documentation	25
4.4.2.1	HSK_NONE	25
4.4.2.2	HSK_RTSCTS	25
4.4.2.3	HSK_XONXOFF	25
4.4.2.4	THERMOCOM_ERROR	26

4.4.2.5	THERMOCOM_RECEIVE	26
4.4.2.6	THERMOCOM_TRANSMIT	26
4.4.3	Function Documentation	26
4.4.3.1	ThermoCOMAbort	26
4.4.3.2	ThermoCOMAttachEvent	26
4.4.3.3	ThermoCOMAttachMsg	27
4.4.3.4	ThermoCOMClose	27
4.4.3.5	ThermoCOMGetError	27
4.4.3.6	ThermoCOMOpen	28
4.4.3.7	ThermoCOMOpenSimulator	28
4.4.3.8	ThermoCOMRead	28
4.4.3.9	ThermoCOMReadBinary	29
4.4.3.10	ThermoCOMSetParam	29
4.4.3.11	ThermoCOMWrite	29
4.4.3.12	ThermoCOMWriteBinary	30

Chapter 1

ThermoCOM dll Interface Specification

About

The purpose of the ThermoCOM.dll dynamic link library is to make interface to the Thermo microplate instruments easy and to offer a similar interface to the instrument as the ThermoUSB.dll library offers. The user of the ThermoCOM.dll does not have to know the details of serial port communication.

Confidential

This document has been prepared by Thermo Fisher Scientific Oy to be used solely for the purposes defined by Thermo Fisher Scientific Oy. Use for other purposes is not authorized.

Please note that any and all information contained in this document is the property of Thermo Fisher Scientific Oy. This confidential information ("Confidential Information") shall not be reproduced in whole part or disclosed to any third party without the prior written approval of Thermo Fisher Scientific Oy. The receiving party shall ensure that its employees, officers, representatives and agents shall not disclose to third parties any Confidential Information.

Upon written request from Thermo Fisher Scientific Oy, the receiving party shall promptly return all Confidential Information or destroy all Confidential Information.

Chapter 2

Using ThermoCOM

2.1 About

ThermoCOM.dll is a generic library for communicating with several different Thermo Scientific microplate instruments via a PC serial port. To communicate with an instrument you use the exported functions of ThermoCOM.dll.

Depending on your project setup, you may find useful a couple of other files which are also provided. The header file [ThermoCOM.h](#) contains the prototypes of the exported functions and definitions of constant values used by the dll. File ThermoCOM.lib contains information about the dll the linker uses to add references to the library in the executable. This way the dll is automatically loaded and the exported functions of the library can be called as easy as the functions in the code using the dll.

2.2 Exported functions

- [ThermoCOMOpen](#)
- [ThermoCOMOpenSimulator](#)
- [ThermoCOMClose](#)
- [ThermoCOMSetParam](#)
- [ThermoCOMAttachEvent](#)
- [ThermoCOMAttachMsg](#)
- [ThermoCOMRead](#)
- [ThermoCOMReadBinary](#)
- [ThermoCOMWrite](#)
- [ThermoCOMWriteBinary](#)
- [ThermoCOMGetError](#)
- [ThermoCOMAbort](#)

2.2.1 ThermoCOMOpen

Open the requested serial port and check whether an instrument with the requested serial number is connected to the port.
Full declaration: [ThermoCOMOpen\(\)](#).

Parameters

<i>PortNumber</i>	The Windows serial port number to open.
<i>baud</i>	The baudrate to use.
<i>InstrumentName</i>	The name of the instrument. If not NULL, the name the instrument returns to a version query must match this string.
<i>SerialNumber</i>	The serial number of the instrument. If not NULL, the serial number the instrument returns to a version query must match this string.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

This function (or ThermoCOMOpenSimulator) must be called first before using any other functions in the library. Only one connection per communication port is allowed.

If either InstrumentName or SerialNumber, or both, are defined, a VER command is sent to the serial port, and these arguments are checked against the returned response. If there is a mismatch, the port is closed and NULL is returned.

If both InstrumentName and SerialNumber are NULL, the VER command is not sent. This makes it possible to open a channel to an instrument which does not support the VER command. In this case it is the responsibility of the application to check what if anything is connected to the serial port.

2.2.2 ThermoCOMOpenSimulator

Open a pipe to a communication port of an instrument simulator. Full declaration: [ThermoCOMOpenSimulator\(\)](#).

Parameters

<i>pipeName</i>	Name of the named pipe used for simulating the communication port. The pipe must have been created by the simulated port driver of the instrument simulator.
-----------------	--

Returns

Communication handle to the simulator, NULL if the function fails.

This function (or ThermoCOMOpen) must be called first before using any other functions in the library. Only one connection per device is allowed. Standard names of the pipes used for simulating communication ports are:

- Serial port: \\0xFFFFCOM
- Debug port: \\0xFFFFDBG
- USB port: \\0xFFFFUSB
- LAN port: \\0xFFFFLAN

The XXXX in all pipe names is the ProductID of the device in hex.

2.2.3 ThermoCOMClose

Close a serial port connection. Full declaration: [ThermoCOMClose\(\)](#).

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
--------------	---

Returns

None.

All open connections are automatically closed when the dll is unloaded, but it is good programming practise to close them explicitly when no longer used.

2.2.4 ThermoCOMSetParam

Set the baud rate of the serial port. Full declaration: [ThermoCOMSetParam\(\)](#).

Parameters

<i>hConn</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>baud</i>	The baud rate to set.
<i>handshake</i>	The handshake to set.

Returns

TRUE on success, else FALSE.

It is recommended to use one of the standard baudrates from 110 to 256000. Using a non standard baudrate may lead to too high bit time error for the communication to work.

The handshake must be one of HSK_NONE, HSK_XONXOFF and HSK_RTSCS. Thermo microplate instruments use the HSK_XONXOFF handshake with the exception of some very old models.

2.2.5 ThermoCOMAttachEvent

Attach an event object to a serial port connection. Full declaration: [ThermoCOMAttachEvent\(\)](#).

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>evCode</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Returns

None

Parameter 'evCode' may be any combination of THERMOCOM_RECEIVE, THERMOCOM_TRANSMIT and THERMOCOM_ERROR. The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event.

The THERMOCOM_RECEIVE event is signaled whenever something is written to the receive buffer. There is no guarantee that a whole response line is received or that only one response line is received. therefore upon receiving this event the application should call [ThermoCOMRead\(\)](#) or [ThermoCOMReadBinary](#) as long as they return data.

The THERMOCOM_TRANSMIT event is signaled when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOCOM_TRANSMIT event is received, functions [ThermoCOMWrite\(\)](#) and [ThermoCOMWriteBinary\(\)](#) will never fail.

2.2.6 ThermoCOMAttachMsg

Attach an event message to a serial port connection. Full declaration: [ThermoCOMAttachMsg\(\)](#).

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>hWnd</i>	Handle of the window to receive the message.
<i>msg</i>	Message id of the message to send.
<i>evCode</i>	Event(s) for which a message is sent.

Returns

None

A message with message id 'msg' is sent to window 'hWnd' whenever any of the event(s) selected with parameter 'evCode' occurs. Parameter 'evCode' may be any combination of THERMOCOM_RECEIVE, THERMOCOM_TRANSMIT and THERMOCOM_ERROR. The event code is passed to the application in the wParam member of the message structure.

The THERMOCOM_RECEIVE event is sent whenever something is written to the receive buffer. There is no guarantee that a whole response line is received or that only one response line is received. therefore upon receiving this event the application should call [ThermoCOMRead\(\)](#) or [ThermoCOMReadBinary](#) as long as they return data.

The THERMOCOM_TRANSMIT event is sent when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOCOM_TRANSMIT event is received, functions [ThermoCOMWrite\(\)](#) and [ThermoCOMWriteBinary\(\)](#) will never fail.

2.2.7 ThermoCOMRead

Read a received response line. Full declaration: [ThermoCOMRead\(\)](#).

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Pointer to buffer receiving the response.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

TRUE if a response line retrieved, else FALSE.

Use this function to read full response lines returned from the instrument. The response line is returned without the terminating CRLF. Therefore TRUE may be returned even if the resulting line is empty. The returned line always ends to the NUL character even if FALSE is returned.

If the caller's buffer is too small to hold the whole response line, as much as fits to the buffer is returned and the rest is returned on subsequent call(s).

2.2.8 ThermoCOMReadBinary

Read received data. Full declaration: [ThermoCOMReadBinary\(\)](#).

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Pointer to buffer receiving the data.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

Number of bytes copied to the user buffer.

Unlike function [ThermoCOMRead\(\)](#), this function returns all data received from the instrument without doing any interpretation on it. Note however that the XON and XOFF characters are filtered by the serial port driver when XON/XOFF flow control is used. You can control the number of bytes returned with the bufsize parameter.

If the return value is less than the given bufsize parameter, there is no more data.

2.2.9 ThermoCOMWrite

Write a string to the transmit buffer. Full declaration: [ThermoCOMWrite\(\)](#).

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Buffer containing the NUL terminated string to send.

Returns

TRUE if string written to the transmit buffer, else FALSE.

When this function returns, the string is not yet sent to the instrument but is queued for sending. Memory for transmit data is dynamically allocated and the only reason for this function returning FALSE is that no more memory is available. If FALSE is returned, the string is not even partially written to transmit buffer and the application should retry with the whole string at a later time.

2.2.10 ThermoCOMWriteBinary

Write binary data to the transmit buffer. Full declaration: [ThermoCOMWriteBinary\(\)](#).

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Buffer containing the data to send.
<i>count</i>	Number of bytes to send from the buffer.

Returns

TRUE if data written to the buffer, else FALSE.

When this function returns, the data is not yet sent to the instrument but is queued for sending. Memory for transmit data is dynamically allocated and the only reason for this function returning FALSE is that no more memory is available. If FALSE is returned, the data is not even partially written to transmit buffer and the application should retry with all data at a later time.

2.2.11 ThermoCOMGetError

Return the error code stored to the connection structure. Full declaration: [ThermoCOMGetError\(\)](#).

2.2.12 ThermoCOMAbort

Send Abort command to instrument. Full declaration: [ThermoCOMAbort\(\)](#).

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
--------------	---

Returns

None

This function first flushes the transmit buffer and also cancels transmission of any data already in the serial port driver. It then sends characters ESC (Abort) and XON to the serial port.

Chapter 3

File Index

3.1 File List

Here is a list of all files with brief descriptions:

api.h	Part of ThermoCOM dll documentation	19
mainpage.h	Start page of ThermoCOM.dll Design Description	19
ThermoCOM.c	Implements serial port interface to Thermo microplate instruments	19
ThermoCOM.h	Functions exported from ThermoCOM.dll	24

Chapter 4

File Documentation

4.1 api.h File Reference

Part of ThermoCOM dll documentation.

4.2 mainpage.h File Reference

Start page of ThermoCOM.dll Design Description.

4.3 ThermoCOM.c File Reference

Implements serial port interface to Thermo microplate instruments.

Functions

- HANDLE WINAPI [ThermoCOMOpen](#) (DWORD PortNumber, DWORD baud, LPCSTR InstrumentName, LPCSTR SerialNumber)
Open the requested serial port and check whether an instrument with the requested serial number is connected to the port.
- HANDLE WINAPI [ThermoCOMOpenSimulator](#) (LPCSTR pipeName)
Open a pipe to a communication port of an instrument simulator.
- BOOL WINAPI [ThermoCOMSetParam](#) (HANDLE hConn, DWORD baud, DWORD handshake)
Set the baud rate of the serial port.
- void WINAPI [ThermoCOMClose](#) (HANDLE hComm)
Close a serial port connection.
- BOOL WINAPI [ThermoCOMAttachEvent](#) (HANDLE hConn, UINT evCode, HANDLE object)
Attach an event object to a serial port connection.
- BOOL WINAPI [ThermoCOMAttachMsg](#) (HANDLE hConn, HWND hWnd, UINT msg, UINT evCode)
Attach an event message to a serial port connection.
- BOOL WINAPI [ThermoCOMRead](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read a received response line.
- DWORD WINAPI [ThermoCOMReadBinary](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read received data.
- BOOL WINAPI [ThermoCOMWrite](#) (HANDLE hConn, LPCSTR buf)
Write a string to the transmit buffer.

- BOOL WINAPI [ThermoCOMWriteBinary](#) (HANDLE hConn, LPCSTR buf, DWORD count)
Write binary data to the transmit buffer.
- void WINAPI [ThermoCOMAbort](#) (HANDLE hConn)
Send Abort command to instrument.
- DWORD WINAPI [ThermoCOMGetError](#) (HANDLE hConn)
Return the error code stored to the connection structure.

4.3.1 Detailed Description

Note

Copyright by Thermo Fisher Scientific Oy 2015

Definition in file [ThermoCOM.c](#).

4.3.2 Function Documentation

4.3.2.1 void WINAPI ThermoCOMAbort (HANDLE hConn)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
--------------	---

Returns

None

This function first flushes the transmit buffer and also cancels transmission of any data already in the serial port driver. It then sends characters ESC (Abort) and XON to the serial port.

Definition at line 1843 of file ThermoCOM.c.

4.3.2.2 BOOL WINAPI ThermoCOMAttachEvent (HANDLE hConn, UINT evCode, HANDLE object)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>evCode</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Returns

None

Parameter 'evCode' may be any combination of THERMOCOM_RECEIVE, THERMOCOM_TRANSMIT and THERMOCOM_ERROR. The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event.

The THERMOCOM_RECEIVE event is signaled whenever something is written to the receive buffer. There is no guarantee that a whole response line is received or that only one response line is received. therefore upon receiving this event the application should call [ThermoCOMRead\(\)](#) or [ThermoCOMReadBinary](#) as long as they return data.

The THERMOCOM_TRANSMIT event is signaled when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOCOM_TRANSMIT event is received, functions [ThermoCOMWrite\(\)](#) and [ThermoCOMWriteBinary\(\)](#) will never fail.

Definition at line 1464 of file ThermoCOM.c.

4.3.2.3 BOOL WINAPI ThermoCOMAttachMsg (HANDLE *hConn*, HWND *hWnd*, UINT *msg*, UINT *evCode*)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>hWnd</i>	Handle of the window to receive the message.
<i>msg</i>	Message id of the message to send.
<i>evCode</i>	Event(s) for which a message is sent.

Returns

None

A message with message id 'msg' is sent to window 'hWnd' whenever any of the event(s) selected with parameter 'evCode' occurs. Parameter 'evCode' may be any combination of THERMOCOM_RECEIVE, THERMOCOM_TRANSMIT and THERMOCOM_ERROR. The event code is passed to the application in the wParam member of the message structure.

The THERMOCOM_RECEIVE event is sent whenever something is written to the receive buffer. There is no guarantee that a whole response line is received or that only one response line is received. therefore upon receiving this event the application should call [ThermoCOMRead\(\)](#) or [ThermoCOMReadBinary](#) as long as they return data.

The THERMOCOM_TRANSMIT event is sent when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOCOM_TRANSMIT event is received, functions [ThermoCOMWrite\(\)](#) and [ThermoCOMWriteBinary\(\)](#) will never fail.

Definition at line 1518 of file ThermoCOM.c.

4.3.2.4 void WINAPI ThermoCOMClose (HANDLE *hComm*)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
--------------	---

Returns

None.

All open connections are automatically closed when the dll is unloaded, but it is good programming practise to close them explicitly when no longer used.

Definition at line 1377 of file ThermoCOM.c.

4.3.2.5 DWORD WINAPI ThermoCOMGetError (HANDLE *hConn*)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
--------------	---

Returns

Windows system error code.

The application may call this function after receiving the THERMOCOM_ERROR event. The returned error code may or may not give a clue about what the actual problem is. If an error is reported, data is already lost and the best action for the application to do is to close the connection and then try to reopen it.

A call to [ThermoCOMGetError\(\)](#) resets the stored error code to ERROR_SUCCESS.

Definition at line 1890 of file ThermoCOM.c.

4.3.2.6 HANDLE WINAPI ThermoCOMOpen (DWORD *PortNumber*, DWORD *baud*, LPCSTR *InstrumentName*, LPCSTR *SerialNumber*)

Parameters

<i>PortNumber</i>	The Windows serial port number to open.
<i>baud</i>	The baudrate to use.
<i>InstrumentName</i>	The name of the instrument. If not NULL, the name the instrument returns to a version query must match this string.
<i>SerialNumber</i>	The serial number of the instrument. If not NULL, the serial number the instrument returns to a version query must match this string.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

This function (or ThermoCOMOpenSimulator) must be called first before using any other functions in the library. Only one connection per communication port is allowed.

If either *InstrumentName* or *SerialNumber*, or both, are defined, a VER command is sent to the serial port, and these arguments are checked against the returned response. If there is a mismatch, the port is closed and NULL is returned.

If both *InstrumentName* and *SerialNumber* are NULL, the VER command is not sent. This makes it possible to open a channel to an instrument which does not support the VER command. In this case it is the responsibility of the application to check what if anything is connected to the serial port.

Definition at line 1237 of file ThermoCOM.c.

4.3.2.7 HANDLE WINAPI ThermoCOMOpenSimulator (LPCSTR *pipeName*)

Parameters

<i>pipeName</i>	Name of the named pipe used for simulating the communication port. The pipe must have been created by the simulated port driver of the instrument simulator.
-----------------	--

Returns

Communication handle to the simulator, NULL if the function fails.

This function (or ThermoCOMOpen) must be called first before using any other functions in the library. Only one connection per device is allowed. Standard names of the pipes used for simulating communication ports are:

- Serial port: \0xFFFFCOM
- Debug port: \0xFFFFDBG
- USB port: \0xFFFFUSB
- LAN port: \0xFFFFLAN

The XXXX in all pipe names is the ProductID of the device in hex.

Definition at line 1288 of file ThermoCOM.c.

4.3.2.8 BOOL WINAPI ThermoCOMRead (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Pointer to buffer receiving the response.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

TRUE if a response line retrieved, else FALSE.

Use this function to read full response lines returned from the instrument. The response line is returned without the terminating CRLF. Therefore TRUE may be returned even if the resulting line is empty. The returned line always ends to the NUL character even if FALSE is returned.

If the caller's buffer is too small to hold the whole response line, as much as fits to the buffer is returned and the rest is returned on subsequent call(s).

Definition at line 1557 of file ThermoCOM.c.

4.3.2.9 DWORD WINAPI ThermoCOMReadBinary (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)**Parameters**

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Pointer to buffer receiving the data.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

Number of bytes copied to the user buffer.

Unlike function [ThermoCOMRead\(\)](#), this function returns all data received from the instrument without doing any interpretation on it. Note however that the XON and XOFF characters are filtered by the serial port driver when XON/XOFF flow control is used. You can control the number of bytes returned with the bufsize parameter.

If the return value is less than the given bufsize parameter, there is no more data.

Definition at line 1733 of file ThermoCOM.c.

4.3.2.10 BOOL WINAPI ThermoCOMSetParam (HANDLE *hConn*, DWORD *baud*, DWORD *handshake*)**Parameters**

<i>hConn</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>baud</i>	The baud rate to set.
<i>handshake</i>	The handshake to set.

Returns

TRUE on success, else FALSE.

It is recommended to use one of the standard baudrates from 110 to 256000. Using a non standard baudrate may lead to too high bit time error for the communication to work.

The handshake must be one of HSK_NONE, HSK_XONXOFF and HSK_RTSCS. Thermo microplate instruments use the HSK_XONXOFF handshake with the exception of some very old models.

Definition at line 1347 of file ThermoCOM.c.

4.3.2.11 BOOL WINAPI ThermoCOMWrite (HANDLE *hConn*, LPCSTR *buf*)**Parameters**

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Buffer containing the NUL terminated string to send.

Returns

TRUE if string written to the transmit buffer, else FALSE.

When this function returns, the string is not yet sent to the instrument but is queued for sending. Memory for transmit data is dynamically allocated and the only reason for this function returning FALSE is that no more memory is available. If FALSE is returned, the string is not even partially written to transmit buffer and the application should retry with the whole string at a later time.

Definition at line 1781 of file ThermoCOM.c.

4.3.2.12 BOOL WINAPI ThermoCOMWriteBinary (HANDLE *hConn*, LPCSTR *buf*, DWORD *count*)**Parameters**

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Buffer containing the data to send.
<i>count</i>	Number of bytes to send from the buffer.

Returns

TRUE if data written to the buffer, else FALSE.

When this function returns, the data is not yet sent to the instrument but is queued for sending. Memory for transmit data is dynamically allocated and the only reason for this function returning FALSE is that no more memory is available. If FALSE is returned, the data is not even partially written to transmit buffer and the application should retry with all data at a later time.

Definition at line 1805 of file ThermoCOM.c.

4.4 ThermoCOM.h File Reference

Functions exported from ThermoCOM.dll.

Macros

- #define [THERMOCOM_RECEIVE](#) 1
Data received event.
- #define [THERMOCOM_TRANSMIT](#) 2
Data transmitted event.
- #define [THERMOCOM_ERROR](#) 4
Fatal error event.
- #define [HSK_NONE](#) 0
Do not use any handshake on serial port.
- #define [HSK_XONXOFF](#) 1
- #define [HSK_RTSC](#) 2
Use hardware handshake on serial port.

Functions

- DllExport HANDLE WINAPI [ThermoCOMOpen](#) (DWORD PortNumber, DWORD baud, LPCSTR InstrumentName, LPCSTR SerialNumber)
Open the requested serial port and check whether an instrument with the requested serial number is connected to the port.

- DllExport HANDLE WINAPI [ThermoCOMOpenSimulator](#) (LPCSTR pipeName)
Open a pipe to a communication port of an instrument simulator.
- DllExport void WINAPI [ThermoCOMClose](#) (HANDLE hComm)
Close a serial port connection.
- DllExport BOOL WINAPI [ThermoCOMSetParam](#) (HANDLE hConn, DWORD baud, DWORD handshake)
Set the baud rate of the serial port.
- DllExport BOOL WINAPI [ThermoCOMAttachEvent](#) (HANDLE hConn, UINT evCode, HANDLE object)
Attach an event object to a serial port connection.
- DllExport BOOL WINAPI [ThermoCOMAttachMsg](#) (HANDLE hConn, HWND hWnd, UINT msg, UINT evCode)
Attach an event message to a serial port connection.
- DllExport BOOL WINAPI [ThermoCOMRead](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read a received response line.
- DllExport DWORD WINAPI [ThermoCOMReadBinary](#) (HANDLE hConn, LPSTR buf, DWORD bufsize)
Read received data.
- DllExport BOOL WINAPI [ThermoCOMWrite](#) (HANDLE hConn, LPCSTR buf)
Write a string to the transmit buffer.
- DllExport BOOL WINAPI [ThermoCOMWriteBinary](#) (HANDLE hConn, LPCSTR buf, DWORD count)
Write binary data to the transmit buffer.
- DllExport DWORD WINAPI [ThermoCOMGetError](#) (HANDLE hConn)
Return the error code stored to the connection structure.
- DllExport void WINAPI [ThermoCOMAbort](#) (HANDLE hConn)
Send Abort command to instrument.

4.4.1 Detailed Description

Note

Copyright by Thermo Fisher Scientific Oy 2015

Serial port interface to Thermo microplate instruments.

Definition in file [ThermoCOM.h](#).

4.4.2 Macro Definition Documentation

4.4.2.1 #define HSK_NONE 0

Definition at line 47 of file ThermoCOM.h.

4.4.2.2 #define HSK_RTSCS 2

Definition at line 65 of file ThermoCOM.h.

4.4.2.3 #define HSK_XONXOFF 1

Definition at line 58 of file ThermoCOM.h.

4.4.2.4 #define THERMOCOM_ERROR 4

What the application can do in case of an error is to first call [ThermoCOMClose\(\)](#) and then try to reopen with [ThermoCOMOpen\(\)](#).

Definition at line 40 of file ThermoCOM.h.

4.4.2.5 #define THERMOCOM_RECEIVE 1

Definition at line 22 of file ThermoCOM.h.

4.4.2.6 #define THERMOCOM_TRANSMIT 2

Definition at line 29 of file ThermoCOM.h.

4.4.3 Function Documentation

4.4.3.1 DllExport void WINAPI ThermoCOMAbort (HANDLE *hConn*)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
--------------	---

Returns

None

This function first flushes the transmit buffer and also cancels transmission of any data already in the serial port driver. It then sends characters ESC (Abort) and XON to the serial port.

Definition at line 1843 of file ThermoCOM.c.

4.4.3.2 DllExport BOOL WINAPI ThermoCOMAttachEvent (HANDLE *hConn*, UINT *evCode*, HANDLE *object*)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>evCode</i>	Event(s) to signal.
<i>object</i>	Handle of an event object.

Returns

None

Parameter 'evCode' may be any combination of THERMOCOM_RECEIVE, THERMOCOM_TRANSMIT and THERMOCOM_ERROR. The event object is set to signaled state whenever any of the selected event(s) occurs.

Parameter 'object' is a handle of a Windows event object. If NULL, the selected events will not be signaled.

This function may be called repeatedly to set up a different event object for each event.

The THERMOCOM_RECEIVE event is signaled whenever something is written to the receive buffer. There is no guarantee that a whole response line is received or that only one response line is received. therefore upon receiving this event the application should call [ThermoCOMRead\(\)](#) or [ThermoCOMReadBinary](#) as long as they return data.

The THERMOCOM_TRANSMIT event is signaled when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOCOM_TRANSMIT event is received, functions [ThermoCOMWrite\(\)](#) and [ThermoCOMWriteBinary\(\)](#) will never fail.

Definition at line 1464 of file ThermoCOM.c.

4.4.3.3 DllExport BOOL WINAPI ThermoCOMAttachMsg (HANDLE *hConn*, HWND *hWnd*, UINT *msg*, UINT *evCode*)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>hWnd</i>	Handle of the window to receive the message.
<i>msg</i>	Message id of the message to send.
<i>evCode</i>	Event(s) for which a message is sent.

Returns

None

A message with message id 'msg' is sent to window 'hWnd' whenever any of the event(s) selected with parameter 'evCode' occurs. Parameter 'evCode' may be any combination of THERMOCOM_RECEIVE, THERMOCOM_TRANSMIT and THERMOCOM_ERROR. The event code is passed to the application in the wParam member of the message structure.

The THERMOCOM_RECEIVE event is sent whenever something is written to the receive buffer. There is no guarantee that a whole response line is received or that only one response line is received. therefore upon receiving this event the application should call [ThermoCOMRead\(\)](#) or [ThermoCOMReadBinary](#) as long as they return data.

The THERMOCOM_TRANSMIT event is sent when the last data from the transmit buffer is sent. It can be used as flow control: If new data is not sent until THERMOCOM_TRANSMIT event is received, functions [ThermoCOMWrite\(\)](#) and [ThermoCOMWriteBinary\(\)](#) will never fail.

Definition at line 1518 of file ThermoCOM.c.

4.4.3.4 DllExport void WINAPI ThermoCOMClose (HANDLE *hComm*)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
--------------	---

Returns

None.

All open connections are automatically closed when the dll is unloaded, but it is good programming practise to close them explicitly when no longer used.

Definition at line 1377 of file ThermoCOM.c.

4.4.3.5 DllExport DWORD WINAPI ThermoCOMGetError (HANDLE *hConn*)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
--------------	---

Returns

Windows system error code.

The application may call this function after receiving the THERMOCOM_ERROR event. The returned error code may or may not give a clue about what the actual problem is. If an error is reported, data is already lost and the best action for the application to do is to close the connection and then try to reopen it.

A call to [ThermoCOMGetError\(\)](#) resets the stored error code to ERROR_SUCCESS.

Definition at line 1890 of file ThermoCOM.c.

4.4.3.6 DllExport HANDLE WINAPI ThermoCOMOpen (DWORD *PortNumber*, DWORD *baud*, LPCSTR *InstrumentName*, LPCSTR *SerialNumber*)

Parameters

<i>PortNumber</i>	The Windows serial port number to open.
<i>baud</i>	The baudrate to use.
<i>InstrumentName</i>	The name of the instrument. If not NULL, the name the instrument returns to a version query must match this string.
<i>SerialNumber</i>	The serial number of the instrument. If not NULL, the serial number the instrument returns to a version query must match this string.

Returns

A handle to the opened communication channel. This handle must be passed to subsequent calls to the other functions in this library. If the channel could not be opened, NULL is returned.

This function (or ThermoCOMOpenSimulator) must be called first before using any other functions in the library. Only one connection per communication port is allowed.

If either *InstrumentName* or *SerialNumber*, or both, are defined, a VER command is sent to the serial port, and these arguments are checked against the returned response. If there is a mismatch, the port is closed and NULL is returned.

If both *InstrumentName* and *SerialNumber* are NULL, the VER command is not sent. This makes it possible to open a channel to an instrument which does not support the VER command. In this case it is the responsibility of the application to check what if anything is connected to the serial port.

Definition at line 1237 of file ThermoCOM.c.

4.4.3.7 DllExport HANDLE WINAPI ThermoCOMOpenSimulator (LPCSTR *pipeName*)

Parameters

<i>pipeName</i>	Name of the named pipe used for simulating the communication port. The pipe must have been created by the simulated port driver of the instrument simulator.
-----------------	--

Returns

Communication handle to the simulator, NULL if the function fails.

This function (or ThermoCOMOpen) must be called first before using any other functions in the library. Only one connection per device is allowed. Standard names of the pipes used for simulating communication ports are:

- Serial port: \0xFFFFCOM
- Debug port: \0xFFFFDBG
- USB port: \0xFFFFUSB
- LAN port: \0xFFFFLAN

The XXXX in all pipe names is the ProductID of the device in hex.

Definition at line 1288 of file ThermoCOM.c.

4.4.3.8 DllExport BOOL WINAPI ThermoCOMRead (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)

Parameters

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Pointer to buffer receiving the response.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

TRUE if a response line retrieved, else FALSE.

Use this function to read full response lines returned from the instrument. The response line is returned without the terminating CRLF. Therefore TRUE may be returned even if the resulting line is empty. The returned line always ends to the NUL character even if FALSE is returned.

If the caller's buffer is too small to hold the whole response line, as much as fits to the buffer is returned and the rest is returned on subsequent call(s).

Definition at line 1557 of file ThermoCOM.c.

4.4.3.9 DllExport DWORD WINAPI ThermoCOMReadBinary (HANDLE *hConn*, LPSTR *buf*, DWORD *bufsize*)**Parameters**

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Pointer to buffer receiving the data.
<i>bufsize</i>	Size of the buffer in bytes.

Returns

Number of bytes copied to the user buffer.

Unlike function [ThermoCOMRead\(\)](#), this function returns all data received from the instrument without doing any interpretation on it. Note however that the XON and XOFF characters are filtered by the serial port driver when XON/XOFF flow control is used. You can control the number of bytes returned with the bufsize parameter.

If the return value is less than the given bufsize parameter, there is no more data.

Definition at line 1733 of file ThermoCOM.c.

4.4.3.10 DllExport BOOL WINAPI ThermoCOMSetParam (HANDLE *hConn*, DWORD *baud*, DWORD *handshake*)**Parameters**

<i>hConn</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>baud</i>	The baud rate to set.
<i>handshake</i>	The handshake to set.

Returns

TRUE on success, else FALSE.

It is recommended to use one of the standard baudrates from 110 to 256000. Using a non standard baudrate may lead to too high bit time error for the communication to work.

The handshake must be one of HSK_NONE, HSK_XONXOFF and HSK_RTSCS. Thermo microplate instruments use the HSK_XONXOFF handshake with the exception of some very old models.

Definition at line 1347 of file ThermoCOM.c.

4.4.3.11 DllExport BOOL WINAPI ThermoCOMWrite (HANDLE *hConn*, LPCSTR *buf*)**Parameters**

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Buffer containing the NUL terminated string to send.

Returns

TRUE if string written to the transmit buffer, else FALSE.

When this function returns, the string is not yet sent to the instrument but is queued for sending. Memory for transmit data is dynamically allocated and the only reason for this function returning FALSE is that no more memory is available. If FALSE is returned, the string is not even partially written to transmit buffer and the application should retry with the whole string at a later time.

Definition at line 1781 of file ThermoCOM.c.

4.4.3.12 DllExport BOOL WINAPI ThermoCOMWriteBinary (HANDLE *hConn*, LPCSTR *buf*, DWORD *count*)**Parameters**

<i>hComm</i>	Connection handle returned from a call to ThermoCOMOpen() .
<i>buf</i>	Buffer containing the data to send.
<i>count</i>	Number of bytes to send from the buffer.

Returns

TRUE if data written to the buffer, else FALSE.

When this function returns, the data is not yet sent to the instrument but is queued for sending. Memory for transmit data is dynamically allocated and the only reason for this function returning FALSE is that no more memory is available. If FALSE is returned, the data is not even partially written to transmit buffer and the application should retry with all data at a later time.

Definition at line 1805 of file ThermoCOM.c.

Index

api.h, [19](#)

HSK_NONE

ThermoCOM.h, [25](#)

HSK_RTSCSTS

ThermoCOM.h, [25](#)

HSK_XONXOFF

ThermoCOM.h, [25](#)

mainpage.h, [19](#)

THERMOCOM_ERROR

ThermoCOM.h, [25](#)

THERMOCOM_RECEIVE

ThermoCOM.h, [26](#)

THERMOCOM_TRANSMIT

ThermoCOM.h, [26](#)

ThermoCOM.c, [19](#)

ThermoCOMAbort, [20](#)

ThermoCOMAttachEvent, [20](#)

ThermoCOMAttachMsg, [21](#)

ThermoCOMClose, [21](#)

ThermoCOMGetError, [21](#)

ThermoCOMOpen, [21](#)

ThermoCOMOpenSimulator, [22](#)

ThermoCOMRead, [22](#)

ThermoCOMReadBinary, [23](#)

ThermoCOMSetParam, [23](#)

ThermoCOMWrite, [23](#)

ThermoCOMWriteBinary, [24](#)

ThermoCOM.h, [24](#)

HSK_NONE, [25](#)

HSK_RTSCSTS, [25](#)

HSK_XONXOFF, [25](#)

THERMOCOM_ERROR, [25](#)

THERMOCOM_RECEIVE, [26](#)

ThermoCOMAbort, [26](#)

ThermoCOMAttachEvent, [26](#)

ThermoCOMAttachMsg, [27](#)

ThermoCOMClose, [27](#)

ThermoCOMGetError, [27](#)

ThermoCOMOpen, [27](#)

ThermoCOMOpenSimulator, [28](#)

ThermoCOMRead, [28](#)

ThermoCOMReadBinary, [29](#)

ThermoCOMSetParam, [29](#)

ThermoCOMWrite, [29](#)

ThermoCOMWriteBinary, [30](#)

ThermoCOMAbort

ThermoCOM.c, [20](#)

ThermoCOM.h, [26](#)

ThermoCOMAttachEvent

ThermoCOM.c, [20](#)

ThermoCOM.h, [26](#)

ThermoCOMAttachMsg

ThermoCOM.c, [21](#)

ThermoCOM.h, [27](#)

ThermoCOMClose

ThermoCOM.c, [21](#)

ThermoCOM.h, [27](#)

ThermoCOMGetError

ThermoCOM.c, [21](#)

ThermoCOM.h, [27](#)

ThermoCOMOpen

ThermoCOM.c, [21](#)

ThermoCOM.h, [27](#)

ThermoCOMOpenSimulator

ThermoCOM.c, [22](#)

ThermoCOM.h, [28](#)

ThermoCOMRead

ThermoCOM.c, [22](#)

ThermoCOM.h, [28](#)

ThermoCOMReadBinary

ThermoCOM.c, [23](#)

ThermoCOM.h, [29](#)

ThermoCOMSetParam

ThermoCOM.c, [23](#)

ThermoCOM.h, [29](#)

ThermoCOMWrite

ThermoCOM.c, [23](#)

ThermoCOM.h, [29](#)

ThermoCOMWriteBinary

ThermoCOM.c, [24](#)

ThermoCOM.h, [30](#)