Technical Note: 51262

Automatic Power-Outage Recovery with an Antaris Process FT-NIR Analyzer and RESULT Software

Jeffrey Hirsch, Ph.D., Thermo Fisher Scientific, Madison, WI, USA

Key Words

- Antaris
- FT-NIR
- Power Outage
- Process
- Recovery
- Restart

One of the primary considerations in putting a Near-Infrared (NIR) analyzer online in a manufacturing process is how the analyzer will react in the event of a power outage on the factory grid. This document discusses the steps needed to set up an Antaris[™] FT-NIR process analyzer to recover automatically from a power outage. Once recovered, the instrument can be set to either begin collecting data automatically to monitor and record process information as the plant recovers, or it can be left in a "ready-to-run" state until other systems come online.

The first step configures the Binary Instrument Operating System (BIOS) settings of the computer attached to the instrument to auto-launch the computer after power loss. To set the BIOS options, turn on the computer and, when prompted to enter setup, hit the appropriate key. This will launch the BIOS setup utility. From there, navigate to the section titled "Power Management" and select the "AC Recovery" option (versions of BIOS differ in the location of this setting. If you cannot find power options contact your IT department). Set this option to "On". The computer will now automatically power on after power interrupt instead of requiring an operator to press a button. Figure 1 is a picture of how the "AC Recovery" option in the BIOS interface might look.



Figure 1: One version of the BIOS Setup utility for a PC

In highly-regulated environments, it is usually required that an operator log into a workstation using either a unique password or a pass key. If login security is required, then there are two main options to accomplish a recovery, remote desktop for the workstation or a remote computer. Remote desktop is the most efficient of the two, enabling a login by an authentic user over a network connection. To set up this feature, Microsoft[®] Windows[®] Remote Desktop must be enabled by going into the "System" control panel. The tab labeled "Remote" contains the activation for the Remote Desktop feature (Figure 2). This will allow a registered user to enter the required login over a network after power has been disrupted. Please check with your IT department to ensure compliance with company policy regarding this type of setup.



Figure 2: Microsoft Windows' Remote Desktop connection preferences

With the computer on and logged in, RESULT[™] software must be automatically started. To accomplish this task, simply place a copy of the RESULT Operation shortcut in the "Startup" folder. To get to this folder, you can either use Windows Explorer (typically found under C:\Documents and Settings\localuser\Start Menu\Programs) or you can go in through the Start Menu Properties. The latter method is explained in detail below.

Right click on the Windows Start Menu and select "Properties." The screen that comes up in Windows XP is shown in Figure 3a. Click "Customize" and then click "Advanced" on the next screen. You will get a Windows Explorer screen that is defaulted to a folder called "Start Menu" (Figure 3b). Open the "Start Menu" folder, expand the folder called "Programs", and click once on the folder called "Startup." The contents of this "Startup" folder will be launched automatically at computer power-up. Make a copy of the RESULT Operation icon on the desktop and place it in this folder. The next time the computer reboots, RESULT Operation will automatically launch. This same task can be accomplished using Windows Explorer to find the same "Startup" folder and copying in the icon.

At this point, both the computer and instrument are powered on, a legitimate user is logged in and RESULT Operation is open, all without operator interaction. The analyzer is now in a "ready-to-run" state.





There is another	Taskbar and Start Menu Properties	RESULTOPCServerDA and RESULTOPCServerAE - Thermo	o Electron Corp. R 🗖 🗖 🔀		
ten that can be taken	Taskber Stat Menu Ele OPC Yew Demo Help				
step that can be taken	6 3 M24				
o get the analyzer					
online and running a		Tag	Value		
finne and running a		RESULT.Mode	1		
pecific workflow		RESULT.RunningStatus	On		
pecific worknow		RESULT.Workflow.Name			
without operator		RESULT.Workflow.WorkflowPath	C: RESULT Data Workflows		
without operator	(Cana) State and the second se	RESULT.Workflow.ReadyToRun	On		
nteraction Using either		RESULT. Workflow.RunningState	U		
interaction. Using entiter	Start menu Customize	RESULT. Workflow CurrentEventStartTime			
he Industrial	Select this menu style for easy access to the Internet, e-mail, and your favorite programs	RESULT. Workflow. CurrentScanStatus			
ne maustriar		RESULT.Workflow.ErrorMessage			
Automation options in	Classic Start menu Customize	RESULT.Workflow.StartEvent.StartEventStatus	Off		
automation options in	Select this option to use the menu style from earlier versions of Windows	RESULT.Workflow.StopEvent.StopEventStatus	Off		
RESULT Operation or		RESULT.Workflow.MeasureEvent.NumberComponents	0		
CLOULI Operation of		RESULT.Workflow.RequestEvent.NumberRequests	0		
n OPC command a	OK Cancel Apply	RESULT.Workflow.CheckEvent.NumberPassFailResults	0		
in Of C command, a		RESULT. Workflow, ArchiveEvent, ArchiveEventName			
pecific workflow can	Figure 3a: Start Menu properties	RESULT. Workflow RequestIngutEvent IngutData Available	off		
pecific worknow can	rigure od. otdrt menu properties	RESULT. Workflow RequestingutEvent. InputDataAvailable	Off		
e loaded and launched		RESULT. Workflow.RequestInputEvent.NumberInputEields	0		
loaded and faunched.			-		
			2		
C:\Documents and Settings\Jeffrey	y.Hirsch\Start Menu\Programs\Startup 📘 🔲 🔀	Ready	1.		

Figure 4: RESULT OPC Test Client General Tab

menu. Figure 5 shows the "Automation Options" dialogue. To have RESULT initiate a workflow, set the "Start Workflow" line number under "Workflow Controls" to the desired number on our Process Communications Controller (this process controller works in tandem with any regular PLC but serves as a bridge to RESULT software). This line should be selected as digital. At the top of this dialogue, under workflow, select the workflow that you would like to run. This setting will allow RESULT, once opened, to automatically select a specified workflow and begin running.

So once the power comes back up after an outage, you can be assured that your Antaris FT-NIR process analyzer will be back online automatically to help with diagnostics or to simply continue normal operation quickly and smoothly.

😠 Industrial Automation Settings				l l	
Opto22 controller IP address:	10.0.0	0.0.1			
RESULT alive line frequency (seconds):	10	10			
Workflow selection:	Fixed	Fixed			
Workflow name:		C:\RESULT Data\Workflows\Fluid Bed Drier.wfl		Browse	
Workflow control:		Digital Lines			
RESULT and workflow running states:					
Item	[Point Number	Point Type		
RESULT Alive Line		12	Digital Point		
Workflow Running State		None Selected	None Selected		
StartStop Item Start State		None Selected	None Selected	None Selected	
StartStopItem Stop State		None Selected	None Selected		
Workflow controls:					
Item		Point Number	Point Type		
Start Workflow		1	Digital Point		
Stop Workflow		2	Digital Point		
StartStop Item Start Trigger		None Selected	None Selected		
StartStop Item Stop Trigger		None Selected	None Selected		
L			OK	Cancel	

Figure 5: RESULT Industrial Automation Options in RESULT Operation

In addition to these offices. Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

Australia Austria +43 1 333 50340 Belgium Canada 1 800 532 4752 China +86 10 5850 3588 Denmark +45 70 23 62 60 France +33 1 60 92 48 00 **Germany** +49 6103 408 1014 India +91 22 6742 9434 **Italy** +39 02 950 591 **Japan** +81 45 453 9100 Latin America +1 608 276 5659 **Netherlands** +31 76 587 98 88 South Africa +27 11 570 1840 Spain 91 657 4930 Sweden/Norway/ Finland +46 8 5<u>56 468 00</u> Switzerland UK +44 1442 233555 USA +1 800 532 4752 www.thermo.com 9001 Thermo Electron Scientific Instruments LLC, Madison, WI ©2006 Thermo Fisher

Scientific Inc. All rights reserved. Microsoft and Windows are registered trademarks of the Microsof Corporation, All other traḋemarks are the property of Thermo Fisher Scientific nc. and its subsidiaries

pricing are subject to change Not all products are availabl consult your local sa

TN51262_E 12/06M





Figure 3b: Copying the RESULT Operation icon into the Startup menu

To use the OPC option available in all configurations of RESULT software, edit the values for the tags labeled Name, WorkflowPath and RunningState (shown in Figure 4 using the RESULT OPC Test Client). This will allow a remote operator to load and initiate any RESULT workflow on the local computer automatically. The only workflows available to be called in this manner are the ones that already exist in RESULT Operation's "Select Workflow" menu.

After power interrupt, if a PLC line is still active, or if the PLC is already online, RESULT software can read its input and use it to do the same thing as above - load and initiate a RESULT workflow. The settings used in this case (as opposed to the OPC case) are found in RESULT's "Industrial Automation Options" in the "Administration"