Automated Thermal Cycler (ATC) System

Installation and Use

Catalog Numbers A30207, A30208, A31486, A31487, A31488, A31489, A31490, A31491, A33976, A33977, A33978, A33979, A33980, A33981, and A33982

Pub. No. MAN0016422 Rev. A.0

Note: For safety and biohazard guidelines, see the "Safety" appendix in the *Automated Thermal Cycler (ATC) System User Guide* (Pub. no. MAN0014558). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Product information

Product description

The Automated Thermal Cycler (ATC) System is designed for integrating with laboratory automation products of other equipment manufacturers (OEM) for use in automation workflows. The ATC System can also be used as a stand-alone device in conjunction with desktop software.

Package contents

Item	Description	Quantity
Modules		
ATC block module	ATC block module You will receive one of the following: • ATC 96 well block module (Cat. No. A30207) • ATC 384 well block module (Cat. No.	
ATC control module (Cat. No. A30208)		1

ltem	Description	Quantity
Computer		
CMPTR,DELL E6540 RC (PN A25621)	8	1
Plates		
Full-skirted adaptor plate (Cat. No. A33045)		1
Semi-skirted plate adaptor (Cat. No. A33044)		1
Cables		
Connector cable	You will receive one of the following: • 10 cm cable (Cat. No. A31485) • 1 m cable (Cat. No. A31482) • 3 m cable (Cat. No. A31483)	1
Ethernet cable	_	1
AC power cable	_	1



Location of power points and ports



Figure 1 Back panel of the ATC control module

- 1 Power port
- Power button
- ③ Ethernet port
- (4) Automated Thermal Cycler (ATC) System DSUB
- ⑤ USB port



Figure 2 Back panel of the ATC block module

(1) Automated Thermal Cycler (ATC) System DSUB

Connect the hardware

 Connect the connector cable into the insertion points on the ATC block (Figure 1, item 4) and base module (Figure 2, item 1)

Note: Ensure the DSUB connector is securely fastened.

- 2. Connect the power cable into the back of the ATC control module (Figure 1, item 1)
- **3.** Connect the ethernet cable into the control module (Figure 1, item 3) and into the computer/network that you will be using to control the ATC.
- 4. Power on the ATC.

Use ATC plate adaptors

The plate adaptor helps to ensure that the supported consumables (full skirt and semi-skirt plates) are ejected properly, and that the temperature is kept uniform. The Automated Thermal Cycler comes factory configured with a full-skirted adaptor (amber color) already installed.

Note: ATC plate adaptors are only installed/required on the 96-well block. Adaptors are not required for the 384-well block.





• To uninstall the full-skirted adaptor, follow the instructions in Figure 4.



Figure 4 Uninstall the full-skirted plate adaptor

- ① Slightly push the tabs highlighted in red.
- 2 Grab the bottom of the adaptor, and tilt up.
- ③ Remove the adaptor.
- To install full-skirted plate adaptor follow the instructions in Figure 5.



Figure 5 Install full-skirted plate adaptor

- Remove half skirted adaptor, if any. place full skirted adaptor by tilting an angle allowing the front to insert under the frame.
- (2) Push the two tabs allowing it to be inserted into the frame.
- (3) Ensure the full-skirted plate adaptor sits freely under the frame.
- ④ Full-skirted plate adaptor is installed.
- To install the semi-skirted plate adaptor, follow the instructions in Figure 6.



Figure 6 Installation of semi-skirted plate adaptors and ejector buttons

- ① Slot the semi-skirted plate adaptor to the gap between the Alu top plate and drip pan.
- (2) Slot in the adaptor and center it.
- ③ Snap fit the ejector buttons.

④ There are four ejector buttons.

Operate the ATC as a stand alone instrument using a computer

Make sure all hardware connections are completed "Connect the hardware" on page 2.

1. Download the ATC desktop software at **thermofisher.com** and follow the installation instructions found in the software release notes.

Note: To fully install the software, the computer will need to be connected to the internet. Once the software is fully installed, disconnect the computer from the internet for best performance with the ATC desktop software.

2. Once the software is installed, click on the ATC Demo icon to launch the software.







Figure 8 ATC Demo software initial screen

1 Discover

- 2 A pop up window
- (3) Identified instrument name
- (4) Connect
- **3.** Make sure the ATC is powered on and connected via ethernet cable to your computer.

Note: It will take approximately 1.5 minutes for the ATC to initialize before it can connect to the software.

4. Once the ATC instrument has initialized, click **Discover** to identify all devices connected to the same local network.

A pop up window will appear.

5. Double click on the identified instrument name.

6. Click Connect.

nstrument Contr	ol	Real Time Monitor	ing	
Host Name or IP 2020 P Address 10.1	0816110076 Discover	Cover Temperature Sample Temperature	105.00 °C 25.00 °C	
Product Name ATC Serial Number 2021	: 384-well 0816110076	130.00 °C		
irmware Version 1.1.0 Connect Disconnect	Upgrade t Open Lid Close Lid	100.00 °C		
Run Progress	Stage - /Cycle - /Step -	80.00 °C 70.00 °C		
Remaining Time	C Elapsed Time	60.00 °C 50.00 °C		
:	:	40.00 °C 30.00 °C		
un Protocol: -		20.00 °C		

Figure 9 ATC Demo software connected to the ATC instrument

- **7.** Hover mouse over the real time monitoring temperature curves to see black window legend with graph related information.
- 8. Click Start to open Protocol Editor





Figure 10 Run Protocol

- 1) Add (+), Remove (-) stages
- Remove and arrange order
- ③ Run Mode drop down
- (4) Cover Temperature and Sample Volume
- 5 Temperature, Ramp Rate, Hold Time
- 6 Cancel
- ⑦ Start Run
- 8 Export SiLA
- 9 Save
- 10 Load
- 1. In **Stages**, add or remove protocol stages with the plus and minus signs.
- Click up and down arrows to arrange the sequence of stages. Specify the number of cycles required for the particular stage.

3. In the **Steps** field, specify the temperature set points, ramp rate, and duration.

The ramp rate is indicated as a % where 100% corresponds to a peak block ramp rate of 3.5C/s.

Type in −1 in the **Hold Time** field to achieve an infinity hold time.

- 4. Enter cover temperature and sample volume.
- 5. Click Start Run to start immediately, or Save for future use.
- 6. Press Load to display previously protocols.

Note: An empty protocol cannot be saved.

7. Click **Export SiLA** to export protocol for integration with SiLA based software or scheduler.

Note: Although SiLA exported protocol is an .xml file it should opened using a text editor in order to see the content, otherwise they might appear empty if opened with a default browser application. When exporting to SiLA format, no white space is allowed in stage label and protocol file name. Only alpha numeric characters and "-", "_" are accepted.

Note: A message will notify the user if the Protocol Editor is started while the instrument is already running (Edit Only mode). User can still edit and export a protocol while the software control and monitor the instrument running, without interfering with the current run.

Continue?		X
Instrument not ready for run. You may only edit and export pr Current active run protocol will	rotocol without starti not be affected.	ing a run.
	ОК	Cancel

Figure 11 Instrument not ready for run warning

Run Mode	Standard 👻	Cover Temperatur	ie 105	°C Sa	ample Volume	10	μL
Stages	Standard 9700 Simulation		Steps (Use	-1 as "Hold Ti	me" for Infinite Hol	ld step)	
Stage	Cycle		Temperatur	re (°C) Ran	np Rate (%)	Hold Time	e (s)

Figure 12 Run Protocol (Edit Only) mode

A run in progress will show the following display:



Figure 13 ATC Demo software -Run in progress monitoring

If the lid is opened while a run is in progress a warning message will open.

	X
Instrument is perfor	ming a run.
	ОК

Figure 14 Alert of run in progress

8. Click Abort to end the run.



Figure 15 Run ended notification

Connect to multiple instruments using the desktop software

 Connect multiple ATC instruments via ethernet cable to your ethernet network, or directly into your computer using a multiport ethernet/network switch.

Each ATC will be assigned its own IP address by the network.

- **2.** To contol multiple ATC instruments, open multiple instances of the ATC Demo software by repeatedly double clicking on the software shortcut. In each software instance, discover and connect to a different ATC.
- **3.** To launch a number of instances of the ATC Demo software at the same time, double click the ATC Demo desktop icon (FigureX) repeatedly to open multiple windows of the application. and then select from each specific application window (using Discovery button) the instrument you want to connect to.
- 4. Click **Discover** in each specific application window to identify the instrument that you want to connect to.



Figure 16 Parallel ATC instruments control and monitoring through multiple instances of ATC Demo software

Note: We recommend that you use one computer to control multiple ATC instruments, and we do not recommend that a single ATC is connected to multiple computers.

Connect Automated Thermal Cycler to a robotic system

Please contact your robotic system manufacturer for information on software compatibility, physical mounting of the ATC on your robot, and licensing information. For the most recent information regarding robot compatibility, driver information, and the full instrument user guide please visit **thermofisher.com/atc** or **thermofisher.com**.

Ordering information

Unless otherwise indicated, all materials are available through **thermofisher.com**.

ATC System

Item	Cat. No.
ATC 96 WELL SYSTEM, LAPTOP, 1m cable	A31486
ATC 96 WELL SYSTEM, LAPTOP, 3m cable	A31487
ATC 96 WELL SYSTEM, LAPTOP, 10 cm cable	A31488
ATC 96 WELL SYSTEM, 1m cable	A31489
ATC 96 WELL SYSTEM, 3m cable	A31490
ATC 96 WELL SYSTEM, 10 cm cable	A31491
ATC, 96 WELL BLOCK	A30207
ATC, 1m cable	A31482
ATC, 3m cable	A31483
ATC, 10 cm cable	A31485
ATC, CONTROL BOX	A30208
ATC, semi-skirted plate adaptor	A33044
ATC, full-skirted plate adaptor	A33045
ATC, 384 WELL BLOCK	A33976
ATC 384 WELL SYSTEM, LAPTOP,1m cable	A33977
ATC 384 WELL SYSTEM, LAPTOP, 3m cable	A33978
ATC 384 WELL SYSTEM, LAPTOP, 10 cm cable	A33979
ATC 384 WELL SYSTEM, 1m cable	A33980
ATC 384 WELL SYSTEM, 3m cable	A33981
ATC 384 WELL SYSTEM, 10 cm cable	A33982

Recommended plastics

Item	Cat. No.	Quantity
MicroAmp [™] EnduraPlate [™] Optical 96-	A31732	500
Well Full-Skirted Plates with Barcode, Clear	A31728	50
MicroAmp [™] EnduraPlate [™] Optical 96- Well Full-Skirted Plates with Barcode, Red	A31729	50
MicroAmp [™] EnduraPlate [™] Optical 96- Well Full-Skirted Plates with Barcode, Blue	A31727	50
MicroAmp [™] EnduraPlate [™] Optical 96- Well Full-Skirted Plates with Barcode, Green	A31731	50
MicroAmp™ EnduraPlate™ Optical 96- Well Blue Reaction Plates with Barcode	4483343	20
MicroAmp [™] EnduraPlate [™] Optical 96- Well Green Reaction Plates with Barcode	4483349	20
MicroAmp [™] EnduraPlate [™] Optical 96- Well Red Reaction Plates with Barcode	4483350	20
MicroAmp [™] EnduraPlate [™] Optical 96-	4483352	500
Barcode	4483354	20
MicroAmp [™] EnduraPlate [™] Optical 96-	4483355	5
Barcode	4483356	500
MicroAmp [™] EnduraPlate [™] Optical 96- Well Yellow Reaction Plates with Barcode	4483395	20

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at **www.thermofisher.com/us/en/home/global/terms-and-conditions.html**. If you have any questions, please contact Life Technologies at **www.thermofisher.com/support**.

Manufacturer's address: Life Technologies Holdings Pte Ltd | Block 33 | Marsiling Industrial Estate Road 3 | #07-06, Singapore 739256

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0016422

Revision	Date	Description
A.0	11 April 2017	New document

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.



thermofisher.com