

# PCR Starter Kit For 384-well blocks

Part. No. 100024353

Revision B



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.

**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.

**Trademarks:** All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2017 Thermo Fisher Scientific Inc. All rights reserved.

# Contents

Product description .....	4
Kit contents .....	4
How to use the MicroAmp™ Adhesive Film Applicator.....	5
Seal a PCR plate with adhesive film .....	5
How to use the MicroAmp™ Multi Removal Tool .....	6
Remove a PCR plate from a sample block with robotslots.....	6
Limited product warranty .....	6
MicroAmp™ plastic consumables for Applied Biosystems™ PCR systems .....	7
MicroAmp™ plastic compatibility chart.....	8

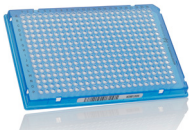



## Product description

The PCR Starter Kit contains sample consumables necessary to start using your new Applied Biosystems™ Thermal Cycler. These consumables are designed to promote efficient heat transfer to obtain fast, high-performance polymerase chain reaction (PCR).

The table below lists the samples included in the kit. It does not list the actual quantities of the starter kit components. To re-order any of these items refer to the catalog number.

**Note:** For safety and biohazard guidelines, refer to the “Safety” appendix in the user guide for your thermal cycler. Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

## Kit contents

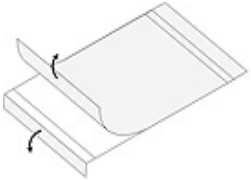


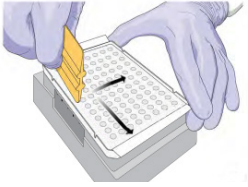

Component	Cat. No.	Quantity <sup>[1]</sup>	Image
MicroAmp™ EnduraPlate™ Optical 384-Well Clear Reaction Plates with Barcode	4483285	5 ea	
MicroAmp™ Clear Adhesive Film	4306311	5 pc	
MicroAmp™ Adhesive Film Applicator	4333183	1 ea	
MicroAmp™ Multi Removal Tool	4313950	1 ea	

[1] This number refers to the quantity included in the starter kit, and may differ from the quantity received when re-ordering components using the catalog numbers provided in the table.

# How to use the MicroAmp™ Adhesive Film Applicator

The MicroAmp™ Adhesive Film Applicator enables you to ergonomically push an adhesive film down upon a PCR plate to form a tight seal and minimize air bubbles. This tight seal will prevent well-to-well contamination and sample loss.

## Seal a PCR plate with adhesive film

Step		Action	
1	Apply film to plate	<ol style="list-style-type: none"><li>Place the PCR plate on a Splash-free 96-well base.</li><li>Remove the backing from the adhesive film.</li><li>Place the adhesive film on the PCR plate, making sure to cover all the wells.</li></ol>	
2	Seal plate	<ol style="list-style-type: none"><li>Rub the flat edge of the applicator back and forth along the long edge of the plate</li></ol>	
		<ol style="list-style-type: none"><li>Rub the flat edge of the applicator back and forth along the short edge (width) of the plate.</li></ol>	
		<ol style="list-style-type: none"><li>Rub the end of the applicator horizontally and vertically between all wells.</li></ol>	
		<ol style="list-style-type: none"><li>Rub the end of the applicator around all outside edges of the plate using small back and forth motions to form a complete seal around the outside wells.</li></ol>	

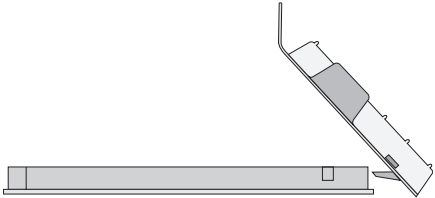
# How to use the MicroAmp™ Multi Removal Tool

The MicroAmp™ Multi Removal Tool makes it easy to open caps, and remove PCR plates that get stuck on the sample block due to heat expansion after performing PCR.

Use this multi-functional tool to open MicroAmp™ Tubes or GeneAmp™ Tubes with attached caps, remove PCR plates from a thermal cycler sample block, and remove cap strips from reaction tubes in a PCR plate and tray/retainer assembly.

## Remove a PCR plate from a sample block with robotslots

Use the tool to remove MicroAmp™ 384-Well Reaction Plate or 96-Well Fast Thermal Cycling Plate from a thermal cycler sample block.

Step	Action	
<b>1</b>	Insert the two hook protrusions into the short side of the PCR plate.	
<b>2</b>	Lift to remove the plate from the sample block.	

## 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](http://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](http://www.thermofisher.com/support).

## MicroAmp™ plastic consumables for Applied Biosystems™ PCR systems

These consumables are compatible with Veriti™ and ProFlex™ PCR systems with 384-well blocks. Go to [thermofisher.com/plastics](http://thermofisher.com/plastics) for additional details.

Product	Catalog No.
<b>384-well reaction plates</b>	
Optical 384-Well Plate	4343370
Optical 384-Well Plate with Barcode	4309849, 4326270, 4343814
EnduraPlate Optical 384-Well Clear Plate with Barcode <sup>[1]</sup>	4483285, 4483273
<b>Seals and covers</b>	
Clear Adhesive Film	4306311
Optical Adhesive Film	4360954, 4311971

[1] Multiple colors are available.

# MicroAmp™ plastic compatibility chart

Product	Catalog No.
<b>96-well 0.2 mL reaction plates</b>	
Optical 96-Well Plate	N8010560, 4316813
Optical 96-Well Plate with Barcode	4306737, 4326659
96-Well Plate with Barcode & Optical Caps	403012
Optical 96-Well Plate with Barcode & Optical Adhesive Films	4314320
EnduraPlate Optical 96-Well Clear Plate with Barcode <sup>[1]</sup>	4483354, 4483352
TriFlex 3 x 32-Well Reaction Plate <sup>[2]</sup>	A32810, A32811
<b>96-well 0.1 mL reaction plates</b>	
Fast Optical 96-Well Plate, 0.1 mL	4346907
Fast Optical 96-Well Plate with Barcode, 0.1 mL	4346906, 4366932
EnduraPlate Optical 96-Well Fast Clear Plate with Barcode <sup>[1]</sup>	4483485, 4483494
<b>384-well reaction plates</b>	
Optical 384-Well Plate	4343370
Optical 384-Well Plate with Barcode	4309849, 4326270, 4343814
EnduraPlate Optical 384-Well Clear Plate with Barcode <sup>[1]</sup>	4483285, 4483273
<b>Strip tubes and caps</b>	
Fast 8-Tube Strip, 0.1 mL	4358293
Optical 8-Tube Strip with Attached Optical Caps, 0.2 mL	A30588
8-Tube Strip with Attached Domed Caps, 0.2 mL	A30589
8-Tube Strip, 0.2 mL <sup>[1]</sup>	N8010580
Optical 8-Tube Strip, 0.2 mL	4316567
8-Cap Strip <sup>[1]</sup>	N8010535, N8011535
Optical 8-Cap Strip	4323032
12-Cap Strip <sup>[1]</sup>	N8010534, N8011534
<b>Single tubes</b>	
Fast Reaction Tube with Cap, 0.1 mL	4358297, 4358293
Reaction Tube with Cap, 0.2 mL <sup>[1]</sup>	N8010540, N8010612, N8011540
Reaction Tube without Cap, 0.2 mL <sup>[1]</sup>	N8010533, N8011533
Optical Tube without Cap, 0.2 mL	N8010933
<b>Seals and covers</b>	
Clear Adhesive Film	4306311
Optical Adhesive Film	4360954, 4311971
32-Well Clear Adhesive Film <sup>[2]</sup>	A32812
<b>Accessories</b>	
Splash-Free 96-Well Base	4312063
96-Well Support Base	4379590
96-Well Base	N8010531
96-Well Reaction Tube/Tray/Retainer Set, 0.2 mL	403083, 403086

[1] Multiple colors are available.

[2] Do not use MicroAmp™ 3 × 32-Well Retainer.



3 x 32-well	96-well, 0.2 mL	96-well, 0.1 mL (Fast)	384-well
ProFlex™	Veriti™, ProFlex™, SimpliAmp™	Veriti™	ProFlex™, Veriti™
96-well 0.2 mL reaction plates			
–	●	–	–
–	●	–	–
–	●	–	–
–	●	–	–
–	●	–	–
● <sup>[2]</sup>	●	–	–
96-well 0.1 mL reaction plates			
–	–	●	–
–	–	●	–
–	–	●	–
384-well reaction plates			
–	–	–	●
–	–	–	●
–	–	–	●
Strip tubes and caps			
–	–	●	–
●	●	–	–
●	●	–	–
●	●	–	–
●	●	–	–
●	●	●	–
●	●	●	–
–	●	●	–
Single tubes			
–	–	●	–
●	●	–	–
●	●	–	–
●	●	–	–
Seals and covers			
–	●	●	●
–	●	●	●
● <sup>[2]</sup>	●	–	–
Seals and covers			
–	●	●	–
–	●	●	–
–	●	–	–
–	–	–	–





[thermofisher.com/support](https://thermofisher.com/support) | [thermofisher.com/askaquestion](https://thermofisher.com/askaquestion)

[thermofisher.com](https://thermofisher.com)

4 December 2017

**ThermoFisher**  
S C I E N T I F I C