

# PCR Starter Kit

## For 60-well blocks, 0.5 mL

Part. No. 4386261

Revision D



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.

©2019 Thermo Fisher Scientific Inc. All rights reserved.

# Contents

Product description .....	4
Kit contents .....	4
Limited product warranty .....	4
How to use the GeneAmp™ Thermal Isolation Frame .....	5
Prepare samples using GeneAmp™ reaction tubes .....	5
MicroAmp™ plastic compatibility chart.....	6



## 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
GeneAmp™ Thin-Walled Reaction Tube, with flat cap, 0.5 mL	N8010737	200 pc	
GeneAmp™ Thermal Isolation Frame	N/A	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.

## Limited product warranty

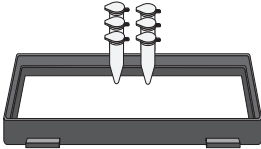
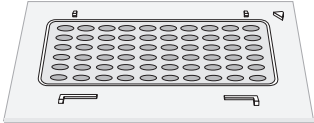
Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html).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

# How to use the GeneAmp™ Thermal Isolation Frame

The GeneAmp™ Thermal Isolation Frame is used to create a proper thermal and physical seal between the heated cover and reaction tubes to balance thermal uniformity across the block. The frame is also used to provide support and distribute pressure evenly across the sample block.

**IMPORTANT!** The retainer is essential for optimal performance, and its usage is recommended at all times for the Veriti™ 60-Well PCR System. The frame allows pressure to be distributed evenly over the tubes to ensure a proper seal, and prevent tubes from being crushed or deformed.

## Prepare samples using GeneAmp™ reaction tubes

Step		Action
1	Fill and cap tubes	<ul style="list-style-type: none"><li>a. Pipette the reaction mix into the tubes.</li><li>b. Cap tubes.</li></ul> 
2	Load tray	<ul style="list-style-type: none"><li>a. Ensure a frame is on the sample block.</li><li>b. Load the tubes into the sample block.</li></ul> 

# 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)

23 July 2019

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