

# Advantage® GC Genomic LA Polymerase Mix Protocol-at-a-Glance

(PT3938-2)

Advantage GC Genomic LA Polymerase Mix is composed of a full-length *Taq* DNA polymerase, a small amount of proofreading enzyme, and a hot-start antibody. The neutralizing hot-start antibody binds and inhibits the *Taq* DNA polymerase at ambient temperatures, enabling reactions to be assembled at room temperature and preventing non-specific amplification due to mispriming and/or formation of primer-dimers before thermal cycling begins. The 2X Advantage GC-Melt Buffer is specifically designed to amplify DNA templates having high GC content or a significant level of secondary structure.

The protocol for using the Advantage GC Genomic LA Polymerase Mix is outlined below.

## PCR Protocol

1. Thaw all reagents on ice.
2. Assemble the PCR reactions by adding each of the components listed in Table I. This can be done at room temperature.

TABLE I. COMPONENTS OF PCR REACTIONS	
Reagents	Volume/Final Concentration per Reaction
Sterile deionized H <sub>2</sub> O	up to 25 $\mu$ l
2X Advantage GC-Melt LA Buffer <sup>1</sup>	12.5 $\mu$ l
dNTP Mixture (10 mM each)	1 $\mu$ l
Primer 1	0.1–0.5 $\mu$ M
Primer 2	0.1–0.5 $\mu$ M
Advantage GC Genomic LA Polymerase Mix (5 units/ $\mu$ l) <sup>2</sup>	0.25 $\mu$ l
Template DNA	10–100 ng

<sup>1</sup> 2 vials (1.25 ml each) of Advantage GC-Melt Buffer are supplied with the kit. This buffer contains 5 mM Mg<sup>2+</sup>.

<sup>2</sup> The enzyme is supplied in a storage buffer containing 20 mM Tris-HCl (pH 8.0), 100 mM KCl, 0.1 mM EDTA, 1 mM DTT, 0.5% Tween®20, 0.5% Nonidet P-40, and 50% glycerol.

3. Mix well and spin down briefly to collect all the liquid at the bottom of the wells.
4. Place the plates in a thermal cycler and begin thermal cycling immediately. The following parameters are useful starting points for amplifying a 2 kb fragment, but annealing temperatures will depend on the primers used.
  - 94°C for 1 min
  - 30 cycles
    - 94°C for 30 sec
    - 60°C for 30 sec
    - 72°C for 2 min\*
  - 72°C for 5 min

\*Note: The extension time should be determined by the size of the amplicon. As a guideline, use 1 min per 1 kb of amplicon.



**Clontech**

United States/Canada  
800.662.2566

Asia Pacific  
+1.650.919.7300

Europe  
+33.(0)1.3904.6880

Japan  
+81.(0)77.543.6116

Clontech Laboratories, Inc.  
A Takara Bio Company  
1290 Terra Bella Ave.  
Mountain View, CA 94043  
Technical Support (US)  
E-mail: tech@clontech.com  
www.clontech.com

(PR692088; published 14 November 2006)

**Notice to Purchaser**

Clontech products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, in vitro diagnostic purposes, therapeutics, or in humans. Clontech products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without written approval of Clontech Laboratories, Inc.

**NOTICE TO PURCHASER: LIMITED LICENSE**

Use of this product is covered

6,127,155. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim (such as method claims in U.S. Patents Nos. 5,994,056 and 6,171,785) and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is hereby conveyed by the purchase of this product expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses require a separate license from Roche. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

GC-Melt is licensed under U.S. Patent No. 5,545,539 and corresponding patents in other countries.

TaqStart® and other Hot Start Antibodies are licensed under U.S. Patent No. 5,338,671.

US Patent No. 5,436,149 for LA technology is owned by Takara Bio Inc.

Tween® is a registered trademark of ICI Americas, Inc.

Clontech, the Clontech logo and all other trademarks are the property of Clontech Laboratories, Inc.

Clontech is a Takara Bio Company. ©2006

Produced by Takara Biotechnology (Dalian) Co., Ltd.