

## Diversify® PCR Random Mutagenesis Kit Protocol-at-a-Glance (PT3393-2)

Please read the *User Manual* (PT3393-1) before using this abbreviated protocol. The Protocol-at-a-Glance is provided for your convenience, but is not intended for first-time users.

### Setting Up and Running Diversify® Mutagenesis Reactions

1. Consult the User Manual for guidelines on choosing the optimal buffer condition(s) for your experiment. Prepare reactions on ice; combine reagents in the order shown:

TABLE III: MUTAGENESIS REACTIONS

	Volumes by Buffer Condition (µl)									Std. <sup>a</sup>
	1	2	3	4	5	6	7	8	9	
Mutations per 1,000 bp	2.0	2.3	2.7	3.5	4.6	4.8	5.8	7.2	8.1	0.4
PCR Grade Water	40	39	38	37	36	35	34	33	32	41
10X TITANIUM <i>Taq</i> Buffer	5	5	5	5	5	5	5	5	5	5
MnSO <sub>4</sub> (8 mM)	0	1	2	3	4	4	4	4	4	0
dGTP (2 mM)	1	1	1	1	1	2	3	4	5	0
50X Diversify dNTP Mix	1	1	1	1	1	1	1	1	1	0
50X dNTP Mix	0	0	0	0	0	0	0	0	0	1
Primer mix <sup>b</sup>	1	1	1	1	1	1	1	1	1	1
Template DNA <sup>c</sup>	1	1	1	1	1	1	1	1	1	1
TITANIUM <i>Taq</i> Polym.	1	1	1	1	1	1	1	1	1	1
<b>Total volume</b>	<b>50</b>	<b>50</b>	<b>50</b>	<b>50</b>	<b>50</b>	<b>50</b>	<b>50</b>	<b>50</b>	<b>50</b>	<b>50</b>

<sup>a</sup> Standard PCR reaction using TITANIUM *Taq* DNA Polymerase

<sup>b</sup> Experimental or Control Primer Mix (10 µM each primer)

<sup>c</sup> Experimental or Control PCR Template (~1 ng/µl)

2. Mix well and spin briefly to collect all liquid at the bottom of the tubes.  
**Note:** If you are not using a hot-lid thermal cycler, overlay contents with mineral oil.
3. Commence thermal cycling using the following parameters for either hot-lid or non-hot-lid thermal cyclers:

- 94°C for 30 sec
- 25 cycles:
  - 94°C 30 sec
  - 68°C 1 min\*
- 68°C for 1 min
- 4°C soak

\* For experimental mutagenesis reactions with templates longer than 1 kb, add 1 min of extension time per additional kb.



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

(PR591061; published 7 December 2005)

**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.

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.

Clontech, the Clontech logo and all other trademarks are the property of Clontech Laboratories, Inc. Clontech is a Takara Bio Company. ©2006