

Terra™ PCR Direct Polymerase Mix Protocol-At-A-Glance

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

1. Terra PCR Direct is designed to amplify targets from whole tissues and tissue extracts. Use the sample amount—and/or preparation method—that is appropriate for your sample type.
2. On ice, combine the reagents shown in Table I in a PCR tube. Make sure you use enough of each reagent for all of your reactions plus an additional reaction to accommodate pipetting errors.

Table 1. Recommended Reagent Volumes

Reagent	Volume	Final concentration
2X Terra PCR Direct Buffer (with Mg ²⁺ , dNTP)	25 µl	1X
Primer 1	15 pmol	0.3 µM
Primer 2	15 pmol	0.3 µM
Tissue Sample/Extract	≤ 5 µl *	
Terra PCR Direct Polymerase Mix	1 µl	1.25 U
Sterile Water	to 50 µl	
Total volume per reaction	50 µl †	

* When using: 1) crude extracts from various animal tissues, add ≤ 5 µl; 2) blood treated with EDTA or heparin, add ≤ 5 µl; 3) mouse tail, add ≤ 1 mm; 4) mouse ear, add ≤ 1.5 mm²; or 5) plant leaves (e.g., tomato or spinach), add ≤ 1.2 mm diameter disk.

† For 25 µl reactions, be sure to add only half the amount indicated for each reagent.

3. Mix the contents of each tube by tapping the bottom of the tube, then centrifuge briefly.
4. Program your thermal cycler with the following cycling conditions:

- **3-Step PCR** (for amplification of standard targets < 2 kb):

98°C	2 min*	} 30–40 cycles
98°C	10 sec	
60°C	15 sec	
68°C	1 min/kb	

* The initial denaturation step must be performed at 98°C for 2 min in order to denature the hot start antibody.

- **2-Step PCR** (for amplification of targets that are GC-rich or ≥ 2 kb):

98°C	2 min*	} 30–40 cycles
98°C	10 sec	
68°C	1 min/kb	

* The initial denaturation step must be performed at 98°C for 2 min in order to denature the hot start antibody.

5. Post-PCR Considerations

- Use TAE running buffer when visualizing your PCR products by agarose gel electrophoresis.
- Add proteinase K to the gel loading buffer.
 - a. Add 5 µl of Proteinase K to 50 µl of 5–6X loading buffer.
 - b. Before loading your samples onto a gel, add 15 µl of the loading buffer-proteinase K mixture to the entire 50 µl PCR reaction, or add 1 µl of the loading buffer-proteinase K mixture to 4 µl of the PCR reaction.

Contact Us	
Customer Service/Ordering	Technical Support
tel: 800.662.2566 (toll-free)	tel: 800.662.2566 (toll-free)
fax: 800.424.1350 (toll-free)	fax: 800.424.1350 (toll-free)
web: www.clontech.com	web: www.clontech.com
e-mail: orders@clontech.com	e-mail: tech@clontech.com

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 prior written approval of Clontech Laboratories, Inc.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at <http://www.clontech.com>. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

Clontech, the Clontech logo, and Terra are trademarks of Clontech Laboratories, Inc. All other marks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions. Clontech is a Takara Bio Company. ©2013 Clontech Laboratories, Inc.

This document has been reviewed and approved by the Clontech Quality Assurance Department.