

I. Description

Prelude PreAmp Master Mix is a 2X PCR mix optimized for unbiased co-amplification of multiple (>100) targets in a single tube starting from low amounts of input sample (10 ng–100 pg cDNA). It contains an antibody-based hot start function for increased specificity, enabling room temperature reaction assembly. This premix works with intact gDNA and cDNA, as well as degraded DNA extracted from FFPE samples. Downstream applications include gene expression profiling and targeted resequencing; the formulation is compatible with both intercalating dye- and probe-based gene expression detection methods.

II. Preamplification of cDNA or gDNA targets

A. Pooling the assays

Combine equal volumes of each assay and dilute with low EDTA (0.1 mM) TE buffer to achieve a target assay concentration of 500 nM.

B. Reaction setup

A typical reaction volume of 50 µl is illustrated below; lower volumes can also be used:

25 µl	Prelude PreAmp Master Mix
5 µl	Primer pool (500 nM stock from Section II.A)
1 µl	Input cDNA (100 pg/µl)
19 µl	PCR-grade water
<hr/>	
50 µl	Total volume

C. Recommended cycling conditions

Cycle the reaction as needed for the sample input amount (10–14 cycles):

95°C	2 min	} 10–14 cycles
95°C	10 sec	
60°C	4 min	
4°C	Hold	

III. General Considerations

- For optimal performance, primers are pooled to target a final concentration of 25–50 nM in the amplification reaction; these can be preconfigured primer/probe sets or custom-designed assay primer sets. Lower concentrations (0.1 to 1 nM) have been tested and also work well.
- Depending on the sample input amount, the PCR program is cycled 10–14 times. The number of cycles can be extended further (up to 28 cycles), although minor loss of amplification uniformity may occur as primer concentrations become limiting (above 18 cycles).

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: takarabio.com	web: takarabio.com
e-mail: ordersUS@takarabio.com	e-mail: techUS@takarabio.com

Notice to Purchaser

Our 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. Our 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 Takara Bio USA, Inc.

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

© 2019 Takara Bio Inc. All Rights Reserved.

All trademarks are the property of Takara Bio Inc. or its affiliate(s) in the U.S. and/or other countries or their respective owners. Certain trademarks may not be registered in all jurisdictions. Additional product, intellectual property, and restricted use information is available at takarabio.com.

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