

Cat. # 9790A
9791A

For Research Use

TaKaRa

TaKaRa BioMasher Standard (Non-sterile)
TaKaRa BioMasher Standard (Sterile)

Product Manual

v201304_2_201803

Table of Contents

I. Description	3
II. Components	3
III. Storage	3
IV. Materials	3
V. Protocol.....	3
VI. Experimental Examples	4
VII. Related Products	4

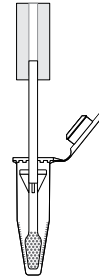
I. Description

TaKaRa BioMasher Standard is a disposable microtube homogenizer designed to efficiently crush small amounts of biological samples for nucleic acid and/or protein extraction. This product is a set of a 1.5-ml microtubes and micro stir bars. The inner wall of the tube and the tip of the stir bar are textured allowing effective disruption of the sample. In addition, there is a lid to prevent splashing of the sample and reagents during sample processing. After the sample has been homogenized, the tube can be centrifuged to minimize sample loss.

The TaKaRa BioMasher Standard is available either sterilized (Cat. #9791A) or nonsterilized (Cat.# 9790A).

II. Components

- | | |
|--|----|
| (1) TaKaRa BioMasher Standard (micro stir bar and microtube) | 50 |
| (2) Stir bar grip (PESTLE GRIP) | 1 |



- III. Storage** Room temperature
Keep out of sun- and UV light.

IV. Materials

TaKaRa BioMasher Standard	
Polypropylene microtubes	Autoclavable
Polyacetal stir bar	Not autoclavable*
Silicon rubber stir bar grip	Autoclavable

* : The stir bar can not be autoclaved. If you require sterilization, please use Cat. #9791A.

V. Protocol

NOTE: Wear a mask, gloves, lab coat, and safety goggles. If necessary, perform in a safety cabinet.

1. Insert the micro stir bar into the end of the PESTLE GRIP.
2. Add the sample to the microtube. Use less than 100 mg of sample. If necessary, add the extraction reagent (less than 250 μ l).
3. Insert the micro stir bar into the microtube. With the PESTLE GRIP, rotate and move the micro stir bar up and down, crushing the sample. If necessary, perform on ice.
4. After crushing, disconnect the PESTLE GRIP and discard the micro stir bar. The PESTLE GRIP can be used repeatedly; store for future use.
5. Close the lid of the microtube. The sample is ready for extraction.

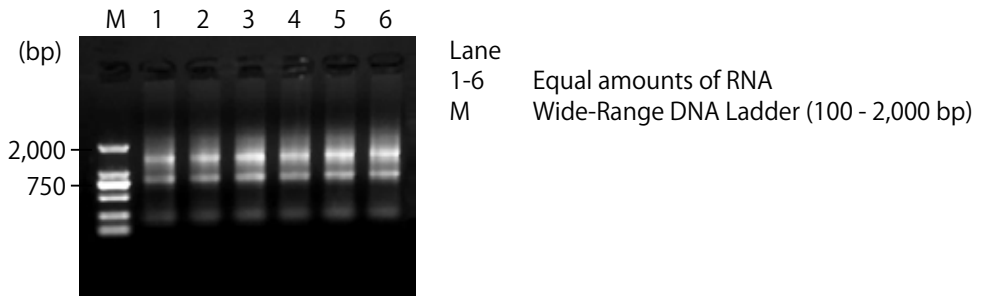
VI. Experimental Example

Total RNA extraction from mouse liver

100 mg of mouse liver was crushed using the TaKaRa BioMasher Standard (Sterile). As a comparison, 100 mg frozen mouse liver tissue was crushed with a mortar or with crusher beads. For all samples, total RNA was extracted using 1 ml RNAiso Plus (Cat. #9108/9109) according to the recommended protocol. Total RNA was quantified and analyzed on a 1% agarose gel.

Results: Processing the samples using the TaKaRa BioMasher Standard had similar extraction efficiency as the mortar and bead crushing protocols.

Crushing method	total RNA (μ g)	A ₂₆₀ /A ₂₈₀
1. Mortar	360	1.70
2. Mortar	424	1.80
3. Crusher beads	530	1.92
4. Crusher beads	494	1.93
5. TaKaRa BioMasher Standard	411	1.89
6. TaKaRa BioMasher Standard	431	1.90



VII. Related Products

RNAisoPlus	(Cat. #9108/9109) *
NucleoSpin® RNA II	(Cat. #740955.10/.50/.250)
NucleoSpin® RNA XS	(Cat. #740902.10/.50/.250)
NucleoSpin® Tissue	(Cat. #740952.10/.50/.250)
NucleoSpin® Tissue XS	(Cat. #740901.10/.50/.250)
Wide-Range DNA Ladder (100 - 2,000 bp)	(Cat. #3427A/B)

* : Not available in all geographic locations. Check for availability in your region.

NOTE: This product is for research use only. It is not intended for use in therapeutic or diagnostic procedures for humans or animals. Also, do not use this product as food, cosmetic, or household item, etc.

Takara products may not be resold or transferred, modified for resale or transfer, or used to manufacture commercial products without written approval from Takara Bio Inc.

If you require licenses for other use, please contact us by phone at +81 77 565 6973 or from our website at www.takara-bio.com.

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

All trademarks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions.
