



025PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Arrest[™] Extraction Buffer

(Cat. #786-133)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

Arrest[™] Extraction Buffer is an optimized combination of various chaotropic agents and RNase inhibitors, which inhibits RNase in 5-10 minutes. *Arrest*[™] Extraction Buffer may be used in conjunction with any RNA extraction method, including extractions based on phenol, chloroform and other organic solvents and detergents. A quick 10 minutes single step protocol isolates high quality, total RNA from any tissue and is conveniently packed as two 50ml bottles. Enough reagents are provided to isolate RNA from a total of 10 grams of tissues.

ITEM(S) SUPPLIED (Cat# 786-133)

Description	Size
<i>Arrest</i> [™] Extraction Buffer	100ml

STORAGE CONDITION

It is shipped at ambient temperature. Upon arrival, store Refrigerated at 4⁰C. When stored and used properly the reagent is stable for 1 year.

PROTOCOL

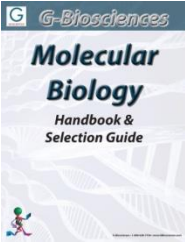
Preparation before Use:

1. Add 0.20ml beta-mercaptoethanol into the *Arrest*[™] Extraction Buffer. Prepare only one bottle at a time and place the date of use on the bottle label.
2. Use 0.5ml *Arrest*[™] Extraction Buffer for each 10-50mg tissue. For cell culture, add 0.5ml per 1-2 million cells.
3. Use any grinding method. After grinding, add 0.5ml chloroform and invert the tube several times to mix.
4. Centrifuge and collect aqueous phase supernatant for extraction of RNA. Take care not to disturb the debris in the inter-phase.
5. Add equal volume ice-cold isopropanol to precipitate RNA. Mix and incubate on ice for 5 minutes. Centrifuge 5 minutes at 4°C to collect the RNA pellet.
6. Wash the RNA pellet with 1ml ice-cold 70% ethanol, and centrifuge 3 minutes to collect the RNA pellet. Remove ethanol and invert the tube on a clean absorbent tissue for several minutes to allow the excess ethanol to drain away.
7. Rehydrate RNA pellet with a minimal volume (2-40ul) of water or other buffer.

NOTE: After grinding the tissue in *Arrest*[™] Extraction Buffer, the protocol can be modified for extraction of RNA, such as using different combination of chloroform and other organic solvents.

Related Products

Download our Molecular Biology Handbook.



<http://info2.gbiosciences.com/complete-molecular-biology-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 2/14/2014SM



www.GBiosciences.com