

Contents

1. Description
 - 1.1 Background information
 - 1.2 Reagent and instrument requirements
2. Protocol for homogenization of tissue for mRNA isolation

1. Description

1.1 Background information

The isolation of subcellular material such as mRNA from tissues or cells requires fast and thorough homogenization of the respective starting material. The gentleMACS™ Dissociators provide optimized programs that meet these requirements. In combination with M Tubes, the gentleMACS Dissociators allow the automated homogenization of tissues in a closed system, enabling sterile sample handling. A single sample or two samples can be processed in parallel.

1.2 Reagent and instrument requirements

- gentleMACS Dissociator (# 130-093-235)
- gentleMACS Octo Dissociator (# 130-095-937)
- gentleMACS M Tubes (# 130-093-236, # 130-096-335)
- μ MACS™ mRNA Isolation Kits: μ MACS mRNA Isolation Kit-Small Scale (# 130-075-201), μ MACS mRNA Isolation Kit-Small Scale (# 130-090-276), μ MACS mRNA Starting Kit (# 130-075-202), μ MACS mRNA Isolation Kit-Large Scale (# 130-075-101), μ MACS mRNA Isolation Kit-Large Scale (# 130-090-277)
- (Optional) Antifoam Y-30 emulsion (e.g. Sigma-Aldrich, A6457)

2. Protocol for homogenization of tissue for mRNA isolation

▲ The protocol has been tested successfully for a range of mouse tissues, such as liver, lung, brain, spleen, kidney, or heart.

▲ If working with fibrous and/or RNase-rich material, such as mouse tail, ear, skin, muscle, or pancreas it is required to prepare total RNA prior to mRNA isolation. Please refer to the gentleMACS Protocol “Homogenization of tissue for total RNA isolation”.

▲ **Note:** Very hard material such as bone should not be processed since it may damage the M Tubes.

▲ Sample volumes of 1 mL or 5 mL are processed per M Tube. For details, please refer to μ MACS mRNA Isolation Kit data sheets.

▲ Its molecular characteristics make RNA chemically unstable and inherently susceptible to ubiquitous RNases. It is therefore recommended to rapidly lyse samples in Lysis/Binding buffer without interruptions to minimize mRNA degradation. Avoid thawing of frozen samples before lysis.

▲ For details on the use of the gentleMACS Dissociators, refer to the gentleMACS Dissociator user manuals.

1. Choose one of the following gentleMACS Programs:
For fresh tissue: gentleMACS Program **RNA_01**
For frozen tissue: gentleMACS Program **RNA_02**
2. Adjust Lysis/Binding Buffer of μ MACS mRNA isolation kit to room temperature.
3. Add Antifoam A to a final concentration of 0.5% to the Lysis/Binding Buffer to prevent excessive foam formation during sample homogenization.
4. Pipette Lysis/Binding Buffer with 0.5 % Antifoam into the M Tube: 1 mL (small scale kit) or 5 mL (large scale kit). For details refer to the μ MACS mRNA isolation kit data sheets.
5. Transfer tissue sample into the Lysis/Binding Buffer in the M Tube.
▲ **Note:** Place sample directly into the buffer to avoid adherence of the tissue to the tube wall.
6. Tightly close M Tube and turn the tube upside down in one quick move ensuring that the sample material reaches the area of the rotor/stator.
7. Attach M Tube upside down onto the sleeve of the gentleMACS Dissociator.
8. Run one of the following gentleMACS Programs:
For fresh tissue: gentleMACS Program **RNA_01**
For frozen tissue: gentleMACS Program **RNA_02**
9. After termination of the program, detach M Tube from the gentleMACS Dissociator.
10. Centrifuge sample at 2000xg for one minute.
11. Remove the homogenized sample from the tube.
▲ **Note:** Homogenized tissue can be removed from the closed M Tube by pipetting through the septum-sealed opening in the center of the cap of the M Tube. Use ART 1000 REACH 1000 μ L pipette tips.
12. Proceed with mRNA isolation as described in the μ MACS mRNA Isolation Kit data sheets.

All gentleMACS Protocols are available at www.miltenyibiotec.com.

Warranty

The products sold hereunder are warranted only to be free from defects in workmanship and material at the time of delivery to the customer. Miltenyi Biotec GmbH makes no warranty or representation, either expressed or implied, with respect to the fitness of a product for a particular purpose. There are no warranties, expressed or implied, which extend beyond the technical specifications of the products. Miltenyi Biotec GmbH's liability is limited to either replacement of the products or refund of the purchase price. Miltenyi Biotec GmbH is not liable for any property damage, personal injury or economic loss caused by the product.

MACS is a registered trademark and gentleMACS and μ MACS are trademarks of Miltenyi Biotec GmbH.

Unless otherwise specifically indicated, Miltenyi Biotec products and services are for research use only and not for diagnostic or therapeutic use.

Copyright © 2011 Miltenyi Biotec GmbH. All rights reserved.