

## Epigenetic Services RIME Tissue Preparation and Best Practices

If you are planning on submitting tissue samples for RIME services, freeze tissue according to one of the protocols below. Tissue requirements for RIME services is a minimum of 800mg.

### Consumables

- Cryogenic vial(s) **or** 2 ml low-bind microcentrifuge tube(s)
- Liquid Nitrogen
- Dry ice

### Protocol A: Liquid Nitrogen

1. Excise the tissue and place in a cryogenic vial or microcentrifuge tube
2. Immediately submerge tube in liquid nitrogen for 2 minutes
3. Store at -80°C

### Protocol B: Dry Ice

1. Excise the tissue and place in cryogenic vial or microcentrifuge tube
2. Immediately place tube on dry ice for 15 minutes
3. Store at -80°C

### Best Practices for sending tissue samples to Active Motif

- Avoid overfilling tubes with tissue as this makes it very difficult to extract samples from tube
- Seal top of tube with parafilm to avoid tube from opening during transit
- Ensure that there is enough dry ice in package for transport
- Avoid shipping over a weekend or for Saturday delivery
- Ship samples Monday through Wednesday
- Ensure that a completed sample submission form is included in the shipment
- Ship samples to the following address

Active Motif, Inc.  
 Attention: Services  
 1914 Palomar Oaks Way, Suite 150  
 Carlsbad, CA 92008 USA  
 760-431-1263

Active Motif S.A. Waterloo Atrium  
 Drève Richelle 167 – boîte 4  
 BE-1410 Waterloo Belgium  
 +32 (0)2 653 0001