

Services ChIP-Seq Sample Preparation for Tissue

If submitting tissues for ChIP-Seq services, freeze according to one of the protocols below. Tissue samples will be fixed by Active Motif Services. You will have the option of selecting “Standard Fixation” or “Double Fixation” in the Sample Submission Portal. Tissue requirements for ChIP-Seq services are 100 - 500 mg. However, for fatty tissue such as mammary gland, Active Motif recommends 1 g per sample.

Required Reagents and Materials:

- Cryogenic vial(s) or 2 ml low-bind microcentrifuge tubes
- Liquid nitrogen or dry ice

Protocol A: Liquid Nitrogen

1. Excise the tissue from the animal and place it 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 from the animal 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 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 complete sample submission form is included in the shipment