



484PR

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

A Geno Technology, Inc. (USA) brand name

XIT™ Genomic DNA from Mouse Tail

For the Isolation of Genomic DNA from Mouse Tail

(Cat. # 786-350)



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

INTRODUCTION

The XIT™ Genomic DNA from Mouse Tail kit is designed for the isolation of genomic DNA from mouse tails. The XIT™ kit uses cell lysis, protein digestion and precipitation and finally DNA precipitation to isolate high quality genomic DNA.

XIT™ Genomic DNA from Mouse Tail kit is offered for the processing 5mm sections of mouse tail. The purified DNA has an A_{260}/A_{280} ratio between 1.7 and 1.9, and is up to 200kb in size.

ITEM(S) SUPPLIED

Description	Cat # 786-350 For 125mm Mouse Tail
XIT™ Lysis Buffer	10ml
LongLife™ Proteinase K	0.5ml
XIT™ Protein Precipitation Buffer	2.5ml
TE Buffer	1.5ml
LongLife™ RNase	0.5ml

STORAGE CONDITIONS

The kit is shipped at ambient temperature. Upon arrival, store the LongLife™ Proteinase K and LongLife™ RNase at -20°C, all other kit components can be stored at room temperature. The kit components are stable for 1 year, if stored properly.

ADDITIONAL ITEMS REQUIRED

Isopropanol, 70% ethanol

PREPARATION BEFORE USE

1. Preheat a waterbath or heating blocks to 55°C.
2. Equilibrate TE Buffer to 50-60°C.

PROTOCOL

1. Using a sharp razor or scalpel blade, chop 5mm (5-10mg) mouse tail into small pieces.
2. Transfer the ground or homogenized tissue to a 1.5ml microfuge tube and add 400µl XIT™ Lysis Buffer.
3. Add 10µl LongLife™ Proteinase K to the tube and mix by inverting the tube 10-20 times. Incubate at 55°C overnight and invert the tube periodically during the incubation.

NOTE: After incubation, the tube can be briefly centrifuged to remove the undigested vertebrae.

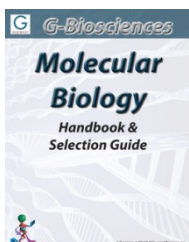
4. After incubation, incubate the sample on ice for 1 minute to quickly cool. Do not store on ice.
5. Add 90µl XIT™ Protein Precipitation Buffer to the sample and mix by inverting the tube 10-20 times.
6. Centrifuge at 14,000g for 3 minutes. Carefully, transfer the supernatant to a fresh tube.

NOTE: The precipitated protein should form a tight white pellet. If not, incubate the sample on ice for 5 minutes and repeat the centrifugation.

7. Add 400µl isopropanol to the supernatant and mix by gently inverting the sample 30-50 times.
8. Centrifuge at 14,000g for 5 minutes.
9. Discard the supernatant and use a pipette to carefully remove excess liquid.
10. Add 200µl 70% ethanol and invert the tube twice to wash the pellet.
11. Centrifuge at 14,000g for 2 minutes.
12. Discard the supernatant and drain the tube on a piece of clean absorbent paper. Allow to air dry for 15 minutes.
13. Add 50µl prewarmed TE buffer and 1µl LongLife™ RNase to remove the RNA (if required).
14. Rehydrate the genomic DNA by incubating at 55-65°C for one hour, followed by an overnight incubation at room temperature to ensure complete genomic DNA hydration.
15. Store DNA at 4°C, for long term storage store at -20 or -80°C.

RELATED PRODUCTS

Download our Molecular Biology Handbook.



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

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

Last saved: 8/29/2012 CMH



www.GBiosciences.com