



480PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# XIT™ Genomic DNA from Yeast

For the Isolation of Genomic DNA  
From Yeast Overnight Cultures

(Cat. # 786-348, 786-349)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

The XIT™ Genomic DNA from Yeast kit is designed for the isolation of genomic DNA from yeast cultures. The XIT™ kit uses the principle of lytic digestion of cell walls, cell lysis, protein precipitation and finally DNA precipitation to isolate high quality genomic DNA. XIT™ Genomic DNA from Yeast kit is for the processing of a maximum of 25 or 250ml of culture. XIT™ Genomic DNA from Yeast Kit protocol is designed to use 1ml overnight culture, however the protocol can be easily adapted for larger tissue sample sizes. The purified DNA has an  $A_{260}/A_{280}$  ratio between 1.8-2.0 and has yields ranging between 1-6µg/ml depending on culture density.

## ITEM(S) SUPPLIED

Description	Cat # 786-348 For 25ml Culture	Cat # 786-349 For 250ml Culture
XIT™ Cell Suspension Buffer	10ml	100ml
XIT™ Lysis Buffer	10ml	100ml
LongLife™ Zymolyase®	0.5ml	3 x 0.5ml
XIT™ Protein Precipitation Buffer	2.5ml	25ml
TE Buffer	1.5ml	20ml
LongLife™ RNase	0.5ml	0.5ml

## STORAGE CONDITIONS

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

## ITEMS NEEDED BUT NOT SUPPLIED

Isopropanol, 70% ethanol

## PREPARATION BEFORE USE

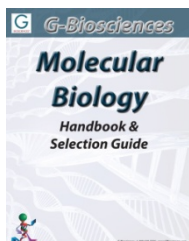
1. Preheat a waterbath or heating block to 37°C.
2. Equilibrate TE Buffer to 50-60°C.

## PROTOCOL

1. Remove 1ml of overnight yeast culture ( $\sim 1\text{-}2 \times 10^8$  cells) and transfer to a 1.5ml centrifuge tube.
2. Centrifuge tube at 5,000g for 5 minutes to pellet yeast.
3. Resuspend the yeast pellet in 400 $\mu$ l XIT™ Cell Suspension Buffer
4. Add 5 $\mu$ l LongLife™ Zymolyase® to the tube and mix by inverting the tube 10-20 times. Incubate at 37°C for 30 minutes with periodic inversion of the tube.
5. After incubation, centrifuge the tube at 5,000g for 5 minute to pellet the spheroplasts.
6. Add 400 $\mu$ l XIT™ Lysis Buffer and pipette up and down to lyze the yeast spheroplasts.
7. Add 90 $\mu$ l XIT™ Protein Precipitation Buffer to the sample and mix by inverting the tube 10-20 times.
8. Centrifuge at 14,000g for 5 minutes. Carefully, transfer supernatant to a new tube.  
**NOTE:** *The supernatant should be clear. If not, repeat the centrifugation.*
9. Add 400 $\mu$ l isopropanol to the supernatant and mix by gently inverting the sample at least 20-25 times.
10. Centrifuge at 14,000rpm for 5 minutes.
11. Discard the supernatant and use a pipette to carefully remove remaining liquid without disturbing the pellet.
12. Add 200 $\mu$ l 70% ethanol and invert the tube twice to wash the pellet.
13. Centrifuge at 14,000rpm for 5 minutes.
14. Discard the supernatant and drain the tube on a piece of clean absorbent paper. Allow to air dry for 15 minutes.
15. Add 50 $\mu$ l prewarmed TE buffer and 1 $\mu$ l LongLife™ RNase to remove the RNA (if required).
16. 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.
17. 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](http://www.GBiosciences.com) or contact us.

Last saved: 8/20/2012 CMH



**[www.GBiosciences.com](http://www.GBiosciences.com)**