



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

A Geno Technology, Inc. (USA) brand name

GET™ DNA Template

Genomic Efficient Technology for DNA Template

(Cat. # 786-353, 786-354)



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

INTRODUCTION 3

ITEM(S) SUPPLIED 4

STORAGE CONDITION 4

ADDITIONAL ITEM(S) REQUIRED 4

IMPORTANT INFORMATION 4

PREPARATION BEFORE USE 4

PROTOCOL 5

 FROM CULTURED CELLS 5

 FROM PARAFFIN EMBEDDED TISSUE 6

 FROM ETHANOL OR FORMALIN FIXED TISSUE 7

 FROM NUCLEATED BLOOD CELLS FROM BIRD, FISH & FROG 8

 FROM GRAM NEGATIVE BACTERIA 9

 FROM GRAM POSITIVE BACTERIA 9

 FROM PLANT TISSUE (FRESH OR FROZEN) 10

 FROM MOUSE TAIL TISSUE 11

 FROM YEAST 12

 FROM FUNGAL TISSUE 13

 FROM BLOOD (≤ 0.2 ML) 14

 FROM BLOOD STAINED & BODY FLUID STAINED MATERIAL 14

 FROM BODY FLUIDS 15

TROUBLESHOOTING 16

RELATED PRODUCTS 17

INTRODUCTION

Genomic Efficient Technology (GET) for purification of DNA templates from diverse sample types, including cells, tissues, mouse tails, plant, blood, nucleated blood cells, body fluid, fixed and embedded tissues, stained specimen, bacteria, yeast, fungal tissues, and so forth.

GET™ is based on a highly efficient Genomic lysis buffer that liberates nucleic acid from cellular protein complexes, making nucleic acids free and available for purification in pure form. Free nucleic acids, DNA templates, are immobilized, in the presence of high concentration of chaotropic agents, on silica solid phase membrane. Following the capture of DNA template on the silica membrane, a series of washing steps removes interfering impurities. In the final step, pure DNA template is eluted in concentrated form with elution buffer (Fig.1).

The eluted DNA template is highly pure and does not require any further processing, making it suitable for a wide variety of applications including PCR, library construction, southern blotting, SNP analysis and molecular diagnostic assays.

GET™ DNA Template kit is supplied in two sizes, 50 (Cat. # 786-353) and 100 (Cat. # 786-354) preps, suitable for processing 50 and 100 samples respectively

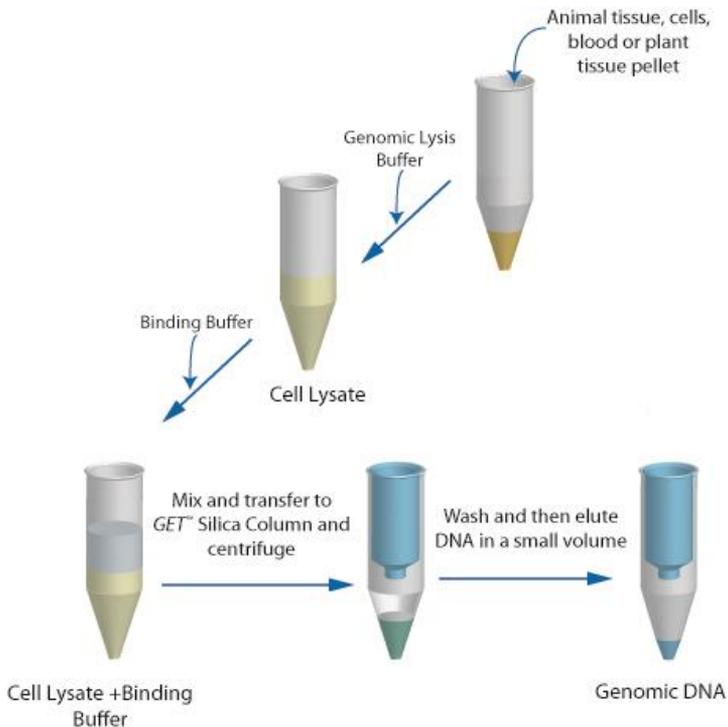


Fig.1

ITEM(S) SUPPLIED

Description	Cat. # 786-353 50 preps	Cat. # 786-354 100 preps
Genomic Lysis Buffer	30 ml	2 x 30 ml
GET Binding Buffer	50 ml	2 x 50 ml
Longlife™ Proteinase K	0.5 ml	0.5 ml
GET Silica Columns	50	2 x 50
GET Wash I	30 ml	2 x 30ml
GET Wash II	20ml	2 x 20ml
GET Elution Buffer	10ml	10ml

STORAGE CONDITION

The kit is shipped at ambient temperature. Upon arrival, store the kit components as recommended on the label. The kit components are stable for 1 year, if stored as directed.

ADDITIONAL ITEM(S) REQUIRED

- Absolute ethanol
- Nuclease-free microfuge tube, 1.5ml
- LongLife™ RNase (Cat. # 786-040) for removal of RNA
- *Longlife™* Lysozyme (Cat. # 786-037) for Gram Positive bacterial culture
- LongLife™ Zymolyase® (Cat # 786-036) for yeast culture

IMPORTANT INFORMATION

- Bring the buffers to room temperature if stored at 4°C. Check the GET Binding Buffer for any precipitate. Warm the buffer at 37°C to dissolve any precipitate.
- Thaw and keep the enzymes Longlife™ Proteinase K on ice when performing DNA isolation. Avoid repeated thawing of enzymes. Invert the vial a few to mix the enzyme suspension. Make small aliquots and store at -20°C for long term storage. Before use, invert the enzyme tube 3-4 times to create uniform suspension, then remove an aliquot for use.

PREPARATION BEFORE USE

1. Add 18 ml of molecular grade ethanol to 30 ml GET Wash I bottle and check the box on the bottle label to indicate ethanol has been added.
2. Add 80ml molecular grade ethanol to the GET Wash II bottle (20 ml) and check the box on the bottle label to indicate ethanol has been added.
3. Equilibrate Elution Buffer to 70°C.

PROTOCOL

From Animal Tissue

1. Transfer 10-50mg tissue to a microcentrifuge tube. Add 200µl Genomic Lysis Buffer. Homogenize the tissue with a microfuge pestle.
2. Add 5 µl Longlife™ Proteinase K suspension into the sample, mix and incubate at 55°C-60°C for 1 hr.
NOTE: Before use, Invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.
3. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 400 µl of GET Binding Buffer and vortex to mix.
4. Transfer the sample to a GET Silica column into the column, positioned in a microfuge tube.
5. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
6. Apply 0.6ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
7. Repeat the wash step as above with 0.6ml Wash II. Centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
8. Repeat the wash step 7, one more time.
9. After discarding the flow through from the last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual Wash II buffer.
10. Transfer the column to a clean nuclease-free 1.5ml microfuge tube for elution of DNA template.
11. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column membrane and incubate at room temperature for 15 minutes.
12. Centrifuge the spin column at 12,000xg for 1 minute to collect the eluted DNA. Store the eluted DNA at 4°C or -20°C until use.
13. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From Cultured Cells

1. Transfer 1-5x10⁶ cells, suspension or trypsinized cells into a 1.5ml microfuge tube.
2. Centrifuge the cells at 1000 x g for 5 minutes. Discard the supernatant.
3. Add 200µl Genomic Lysis Buffer to the pellet. Gently pipette up and down several times to release nuclei from the cells. Transfer lysate to a clean tube.
4. Add 5 µl Longlife™ Proteinase K suspension into the sample and incubate at 55°C-60°C for 1 hr.

NOTE: Before use, Invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.

5. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 400 µl of GET Binding Buffer and vortex to mix.
6. Transfer the sample to a GET Silica Column, positioned in a microfuge tube.
7. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
8. Apply 0.6ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
9. Apply 0.6ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
10. Repeat step 9 for a second wash with GET Wash II.
11. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.
12. Transfer the column to a clean nuclease-free 1.5ml microfuge tube for elution of DNA template.
13. Add 50µl prewarmed (70°) Elution Buffer on the top of the spin column matrix and incubate at room temperature for 15 minutes.
14. Centrifuge the spin column at 12000xg for 1 minute to collect the eluted DNA. Store DNA at 4°C or -20°C for later use or treat it with RNase.
15. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From paraffin embedded tissue

1. Finely chop 0.5-2mg paraffin-embedded tissue and place in 1.5ml microfuge tube with 100µl xylene or safe xylene substitute and incubate at room temperature for 5 minutes with constant mixing.
2. Centrifuge at 14,000xg for 2 minutes and discard xylene or xylene substitute. Repeat steps 1-2 twice to achieve a total of three washes.
3. Add 100µl 100% ethanol and incubate for 5 minutes at room temperature with constant mixing.
4. Centrifuge at 14,000xg for 2 minutes and discard ethanol. Repeat steps 3-4 to achieve a total of two washes.
5. Add 100µl Genomic Lysis Buffer and homogenize the sample with a microfuge pestle until a homogenous suspension is acquired, approximately 30-60 strokes.
6. Add 2.5 µl Longlife™ Proteinase K suspension into the sample and incubate at 55°C-60°C for 1 hr.
NOTE: *Before use, Invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.*
7. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 200 µl of GET Binding Buffer and vortex to mix.
8. Transfer the sample to a GET Silica Column, positioned in a microfuge tube.

9. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
10. Apply 0.6ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
11. Apply 0.6ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
12. Repeat step 11.
13. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.
14. Transfer the column to a clean nuclease-free 1.5ml microfuge tube for elution of DNA template.
15. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate for 15 minutes as room temperature.
16. Centrifuge the spin column at 12,000xg for 1 minute at 25°C to collect the eluted DNA. Store DNA at 4°C or -20°C for later use or treat it with RNase.
17. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From ethanol or formalin fixed tissue

1. Blot excess fixative from tissue with clean absorbent paper.
2. Transfer 10-50mg tissue to a microcentrifuge tube containing 200µl Genomic Lysis Buffer and incubate for 15 minutes at 55-65°C to soften tissue.
3. Homogenize the sample with a microfuge pestle until a homogenous suspension is acquired, approximately 30-60 strokes.
4. Add 5µl Longlife™ Proteinase K suspension into the sample and incubate at 55°C-60°C for 1 hr.
NOTE: *Before use, Invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.*
5. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 400 µl of GET Binding Buffer and vortex to mix.
6. Transfer the sample to a GET Silica Column, positioned in a microfuge tube.
7. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
8. Apply 0.6ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute. Discard the flow through.
9. Apply 0.6ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute. Discard the flow through.
10. Repeat step 9.
11. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.

12. Discard the collection tube and place the column on a clean nuclease-free 1.5ml microfuge tube.
13. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate at room temperature for 15 minutes.
14. Centrifuge the spin column at 12,000xg for 1 minute to collect the eluted DNA. Store DNA at 4°C or -20°C for later use or treat it with RNase.
15. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From nucleated blood cells from bird, fish & frog

1. Add 10µl nucleated blood to a 1.5ml microfuge tube containing 200µl Genomic Lysis Buffer.
2. Add 5 µl Longlife™ Proteinase K suspension into the sample and incubate at 55°C-60°C for 1 hr.
NOTE: *Before use, invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.*
3. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 400 µl of GET Binding Buffer to the sample and vortex to mix.
4. Transfer the sample to a GET Silica Column, positioned in a microfuge tube.
5. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
6. Apply 0.6ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
7. Apply 0.6ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
8. Repeat step 7.
9. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.
10. Discard the collection tube and place the column on a clean nuclease-free 1.5ml microfuge tube.
11. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate at room temperature for 15 minutes.
12. Centrifuge the spin column at 12,000xg for 1 minute to collect the eluted DNA. Store DNA at 4°C or -20°C for later use or treat it with RNase.
13. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From gram negative bacteria

1. Add 0.5 ml of an overnight culture to a 1.5ml microfuge tube.
2. Centrifuge at 16,000xg for 2-3 minutes to pellet the cells. Remove and discard the supernatant. Vortex the tube to re-suspend the cells in residual supernatant.
3. Add 200µl Genomic Lysis Buffer and mix by inverting a few times.
4. Add 5 µl Longlife™ Proteinase K suspension into the sample and incubate at 55°C-60°C for 1 hr.

NOTE: *Before use, Invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.*

5. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 400µl of GET Binding Buffer to the sample and vortex to mix.
6. Transfer the sample to a GET Silica Column, positioned in a microfuge tube.
7. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
8. Apply 0.6ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
9. Apply 0.6ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
10. Repeat step 9.
11. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.
12. Discard the collection tube and place the column on a clean nuclease-free 1.5ml microfuge tube.
13. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate at room temperature for 15 minutes.
14. Centrifuge the spin column at 12,000xg for 1 minute to collect the eluted DNA. Store DNA at 4°C or -20°C for later use.
15. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From gram positive bacteria

1. Aliquot 0.5ml Gram positive bacteria overnight culture into a 1.5ml microfuge tube and centrifuge at 14,000xg for 30 seconds. Discard the supernatant.
2. Add 450µl sterile water and 50µl EDTA to the pellet and gently vortex to resuspend.
3. Add 50µl Longlife™ Lysozyme, invert to mix and incubate at 37°C for 45 minutes with periodic mixing.

NOTE: *Before use, Invert the Longlife™ Lysozyme tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.*

4. Centrifuge the sample tube for 5 minutes at 14,000xg and discard the supernatant. Gently vortex the tube to resuspend the pellet in the residual liquid.
5. Add 200µl Genomic Lysis Buffer and mix by inverting the tube several times.
6. Add 5 µl Longlife™ Proteinase K suspension into the sample and incubate at 55°C-60°C for 1 hr.
NOTE: *Before use, Invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.*
7. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 400 µl of GET Binding Buffer to the sample and vortex to mix.
8. Transfer the sample to a GET Silica Column, positioned in a microfuge tube.
9. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
10. Apply 0.6ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
11. Apply 0.6 ml GET Wash II to the column and centrifuge at 8,000xg for 1 minute at 25°C. Discard the flow through.
12. Repeat step 11.
13. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.
14. Discard the collection tube and place the column on a clean nuclease-free 1.5ml microfuge tube.
15. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate at room temperature for 15 minutes.
16. Centrifuge the spin column at 12,000xg for 1 minute at 25°C to collect the eluted DNA. Store DNA at 4°C or -20°C for later use.
17. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From plant tissue (fresh or frozen)

1. Transfer 10-50mg finely ground dried tissue, frozen tissue or fresh leave tissue to a microcentrifuge tube.
NOTE: *Most plant tissues are best prepared by freezing in liquid nitrogen. Grinding samples in liquid nitrogen to a fine powder and quickly add to an appropriate volume of Genomic Lysis Buffer.*
2. Add 200µl Genomic Lysis Buffer.
3. If ground, vortex for 5 seconds; if unground, homogenize the sample with a microfuge pestle until a homogenous suspension is acquired, approximately 15-30 strokes.
4. Add 5µl Longlife™ Proteinase K suspension into the sample and incubate at 55°C-60°C for 1 hr.

NOTE: Before use, Invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.

5. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 400 µl of GET Binding Buffer and vortex to mix.
6. Transfer the sample to a GET Silica Column, positioned in a microfuge tube.
7. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
8. Apply 0.6 ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute. Discard the flow through.
9. Apply 0.6 ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute. Discard the flow through.
10. Repeat step 9.
11. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.
12. Discard the collection tube and place the column on a clean nuclease-free 1.5ml microfuge tube.
13. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate at room temperature for 15 minutes.
14. Centrifuge the spin column at 12,000xg for 1 minute to collect the eluted DNA. Store DNA at 4°C or -20°C for later use.
15. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From mouse tail tissue

1. Add 0.25-0.5cm, approximately 20-50mg, mouse tail in to a 1.5ml microcentrifuge tube with 200µl Genomic Lysis Buffer.
2. Add 5µl Proteinase K solution and incubate at 60°C for 3-4 hours to overnight. Invert the tube periodically if possible.

NOTE: Before use, Invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.
3. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 400 µl of GET Binding Buffer and vortex to mix.
4. Transfer the sample to a GET Silica Column, positioned in a microfuge tube.
5. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
6. Apply 0.6 ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
7. Apply 0.6 ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute. Discard the flow through.
8. Repeat step 7.

9. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.
10. Discard the collection tube and place the column on a clean nuclease-free 1.5ml microfuge tube.
11. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate at room temperature for 15 minutes.
12. Centrifuge the spin column at 12,000xg for 1 minute to collect the eluted DNA. Store DNA at 4°C or -20°C for later use.
13. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From Yeast

1. Aliquot 1.5ml yeast overnight culture into a 1.5ml microfuge tube and centrifuge at 14,000xg for 30 seconds. Discard the supernatant.
2. Add 150µl PBS, 5µl LongLife™ Zymolyase® and 1µl β-mercaptoethanol to the pellet and gently vortex to resuspend.
NOTE: *Before use, invert the Longlife™ Zymolyase® tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.*
3. Incubate at 37°C for 30 minutes with periodic mixing.
4. Centrifuge the sample tube for 5 minutes at 14,000xg and pour off the supernatant. Gently vortex the tube to resuspend the pellet in the residual liquid.
5. Add 200µl Genomic Lysis Buffer to the tube and mix by inverting the tube several times.
6. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 400 µl of GET Binding Buffer to the sample and vortex to mix.
7. Transfer the sample to a GET Silica Spin Column, positioned in a microfuge tube.
8. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
9. Apply 0.6 ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute. Discard the flow through.
10. Apply 0.6 ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute. Discard the flow through.
11. Repeat step 10.
12. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.
13. Discard the collection tube and place the column on a clean nuclease-free 1.5ml microfuge tube.
14. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate at room temperature for 15 minutes.

15. Centrifuge the spin column at 12,000xg for 1 minute to collect the eluted DNA. Store DNA at 4°C or -20°C for later use or treat it with RNase.
16. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From fungal tissue

1. Collect fungal tissue from liquid culture and wash 2-3 times in sterile water.
2. Fungal mycelia are best prepared by grinding samples using Molecular Grinding Resin™ in Genomic Lysis Buffer. For fungal teliospores, grinding samples in liquid nitrogen to a fine powder and quickly add to an appropriate volume of Genomic Lysis Buffer is recommended.
3. Transfer 10-20mg fungal mycelia to a microcentrifuge tube. Add 200µl Genomic Lysis Buffer. Add 30µl Molecular Grinding Resin™ using a wide bore pipette tips and grind with a microcentrifuge pestle. For teliospores, add ground powder to 200µl Genomic Lysis Buffer and vortex to wet sample.
4. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 400 µl of GET Binding Buffer to the sample and vortex to mix.
5. Transfer the sample to a GET Silica Column, positioned in a microfuge tube.
6. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
7. Apply 0.6 ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute. Discard the flow through.
8. Apply 0.6 ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute. Discard the flow through.
9. Repeat step 8.
10. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.
11. Discard the collection tube and place the column on a clean nuclease-free 1.5ml microfuge tube.
12. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate at room temperature for 15 minutes.
13. Centrifuge the spin column at 12,000xg for 1 minute to collect the eluted DNA. Store DNA at 4°C or -20°C for later use.
17. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From blood (≤0.2ml)

1. To ≤200µl whole blood, buffy coat, bone marrow or packed cells in a 2ml microfuge add 200 µl Genomic Lysis Buffer.
2. Add 10 µl Longlife™ Proteinase K suspension into the sample and incubate at 55°C-60°C for 1 hr.
3. Add 800 µl of GET Binding Buffer to the sample and vortex to mix.
4. Transfer 0.6 ml to a GET Silica Column, positioned in a microfuge tube.
5. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
6. Repeat step 4 and 5 once, with residual sample, to allow all sample to pass through GET Silica Column.
7. Apply 0.6 ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
8. Apply 0.6 ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
9. Repeat step 8.
10. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.
11. Discard the collection tube and place the column on a clean nuclease-free 1.5ml microfuge tube.
12. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate the column at room temperature for 15 minutes.
13. Centrifuge the spin column at 12,000xg for 1 minute to collect the eluted DNA. Store DNA at 4°C or -20°C for later use.
18. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From blood stained & body fluid stained material

1. To 10-30mm² section of stained material in a 2ml microfuge tube, add 500µl Genomic Lysis Buffer and 10µl Proteinase K. Invert to mix.
NOTE: *Before use, Invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.*
2. Incubate the sample at 65°C for 4 hours with periodic inversions.
3. Allow to cool to room temperature, remove the stained material and remove excess buffer from the material with a pipette and return to the tube.
4. Centrifuge the sample tube for 5 minutes at 5000 x g and transfer the clear supernatant to a clean tube. Add 1ml of GET Binding Buffer and vortex to mix.
5. Transfer 0.5 ml to a GET Silica Column, positioned in a microfuge tube.
6. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
7. Repeat step 5 and 6 twice, till all sample is passed through GET Silica Column.

8. Apply 0.6 ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
9. Apply 0.6 ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
10. Repeat step 9.
11. After discarding the flow through from last wash, replace the spin column on the microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.
12. Discard the collection tube and place the column on a clean nuclease-free 1.5ml microfuge tube.
13. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate the column at room temperature for 15 minutes.
14. Centrifuge the spin column at 12,000xg for 1 minute to collect the eluted DNA. Store DNA at 4°C or -20°C for later use.
15. **Optional:** Removal of RNA. Invert the LongLife™ RNase vial 3-4 times to uniformly suspend the enzyme mix. Add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C until use.

From body fluids

This includes CSF, plasma, saliva, serum, sputum, synovial fluid, urine and whole blood

1. Transfer 50µl body fluid to a 1.5ml microfuge tube.
NOTE: For body fluids with a low cell number, concentrate the cells by centrifuging 5-40ml sample at 2,000xg for 10 minutes.
2. For samples with a normal protein concentration, add 200µl Genomic Lysis Buffer and mix by pipetting up and down.
3. Add 5 µl Longlife™ Proteinase K suspension into the sample and incubate at 55°C-60°C for 1 hr.
NOTE: Before use, Invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.
4. Add 500µl of GET Binding Buffer to the sample and vortex to mix.
5. Transfer 0.5 ml to a GET Silica Column, positioned in a microfuge tube.
6. Centrifuge the column at 12,000 x g for 1 minute at 25°C. Discard the flow through.
7. Transfer the remaining sample to the column and repeat step 6 once.
8. Apply 0.6ml GET Wash I to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
9. Apply 0.6ml GET Wash II to the column and centrifuge at 12,000xg for 1 minute at 25°C. Discard the flow through.
10. Repeat step 9.
11. After discarding the flow through from last wash, replace the spin column on microfuge tube and spin at 14,000xg for 3 minutes to remove residual GET Wash II buffer.

12. Discard the collection tube and place the column on a clean nuclease-free 1.5ml microfuge tube.
13. Add 50µl prewarmed (70°C) Elution Buffer on the top of the spin column matrix and incubate the column at room temperature for 15 minutes.
14. Centrifuge the spin column at 12,000xg for 1 minute to collect the eluted DNA. Store DNA at 4°C or -20°C for later use.
15. **Optional:** Briefly centrifuge the LongLife™ RNase and add 1µl LongLife™ RNase per 50 µl of eluted DNA. Incubate at room temperature for 15 minutes. Store the DNA at 4°C or -20°C for later use.

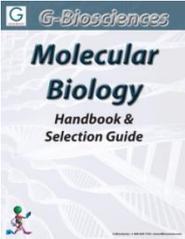
TROUBLESHOOTING

Issue	Suggested reason	Possible solution
Low nucleic acid yield or purity	Kit components are not stored properly	Store kit components as indicated in the label.
	Ethanol not added to the Wash Buffers	Add absolute ethanol to the Wash buffers before using
	Reagent and samples not properly mixed	Mix the sample tube well after addition of each reagent.
Low yield of Tissue DNA	Incomplete Proteinase K digestion	Homogenize tissue sample in Genomic Lysis buffers completely before adding Proteinase K
		Incubate tissue with proteinase K for longer period 2-4 hours
Low yield from bacteria or yeast	Bacteria and Yeast not lysed efficiently with lysozyme and zymolyase respectively	Lyse bacteria and Yeast as mentioned in the protocol
Eluted DNA is degraded	Nuclease activity in unlysed tissue	Tissue should be frozen from the time of extraction to lysis
		Use small pieces of tissue or homogenize in Genomic Lysis buffer
Eluted DNA from blood is slightly colored	Incomplete wash step	Wash GET Silica Column until flow through is colorless
		Add 200 µl of GET Binding Buffer to the eluted DNA and reload on a fresh GET Silica Column and perform

		wash steps and elution as mentioned in protocol
--	--	---

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.

Upgraded: 5/28/20



www.GBiosciences.com