



Tn5 Transposase User Manual

Tn5 Transposase

Cat. No: PC047 (20 µl) , PC048(100 µl)

User Manual

GeneCopoeia, Inc.
9620 Medical Center Drive, #101 Rockville, MD 20850
USA

301-762-0888

inquiry@genecopoeia.com

www.genecopoeia.com

Poly(A) Polymerase

- I. Description
- II. Contents and Storage
- III. Procedure
- IV. Limited Use License and Warranty

I. Description

Tn5 Transposase efficiently catalyzes the random insertion of Tn5 transposons into target sequences, exhibiting extremely high transposition insertion efficiency for both eukaryotic and prokaryotic DNA.

The enzyme specifically recognizes the inverted repeat Mosaic End (ME) sequences flanking the transposon. It recognizes DNA fragments bearing chimeric ME sequences at both ends (including primers containing ME sequences) and forms a Tn5 transposome. This transposome can randomly bind target DNA, cleave it, and insert the carried DNA fragment. Tn5 Transposase is widely used in in vitro transgenesis and next-generation sequencing (NGS) library preparation.

■ Source

Recombinant protein expressed in E. coli.

■ Quality Assurance

Free of host DNA contamination. Free of endonuclease and exonuclease contamination.

■ Applications

- ◇ Fragmentation and adapter addition for next-generation sequencing library preparation.
- ◇ Construction of random transposon insertion mutant libraries in vitro or in vivo.
- ◇ Rapid sequencing of large DNA molecules (e.g., BAC clones).
- ◇ Insertion of antibiotic resistance markers into target DNA.

II. Contents and Storage

Cat. No.	Contents	Part No.	Quantity
PC047	Tn5 Transposase	PC047-01	20 µl
	5× Tn5 Reaction Buffer	PC047-02	200 µl
	5×Stop Buffer	PC047-03	200 µl
PC048	Tn5 Transposase	PC047-01	20 µl×5
	5× Tn5 Reaction Buffer	PC047-02	200 µl×5
	5×Stop Buffer	PC047-03	200 µl×5

Store all components at -20°C (stable for at least 12 months). Avoid repeated freezing/ thawing.

III. Procedures

1. In Vitro Transposon Insertion Reaction

1. Tn5 transposon DNA is the inclusion of the 19-bp ME sequence flanking both ends of the DNA. This is appended to the transposon via PCR amplification with primers containing 5'-phosphorylated ME-sequence.

ME sequence: 5'-pCTGTCTCTTATACACATCT-3'

2. Prepare the transposon insertion mixture as follows:

Reagent	Volume
5× Reaction buffer	4 µl
Target DNA *	200-400ng
Tn5 Transposase	1 µl
Equimolar transposon	X µl
ddH2O	To 10 µl

Note: **Calculate the molar amount of target DNA in the reaction and add an equimolar amount of the Tn5 transposon fragment to avoid excess insertions.*

µmol target DNA = µg target DNA / [(base pairs in target DNA) × 660]

Example: 0.2 µg of a 1000-bp target DNA = 0.2µg / [1000bp×660]=0.3×10⁻⁶µmol = 0.3pmol.

Target DNA must be free of exogenous nucleic acid contamination.

3. Mix thoroughly and incubate at 37°C for 2 hours.
4. Add 2 µl of 5× Stop Buffer (thaw at room temperature), mix, and incubate at 55°C for 10 min to terminate the reaction..

2. Construction of Transposome Complex and Fragmentation for Next-Generation Sequencing

1. Construction of the Transposome Complex

- ① ME and adapter sequences:

Adapter 1: 5'-TCGTCGGCAGCGTCAGATGTGTATAAGAGACAG-3'

Adapter 2: 5'-GTCTCGTGGGCTCGGAGATGTGTATAAGAGACAG-3'

ME: 5'-pCTGTCTCTTATACACATCT-NH₂-3' (5'-phosphorylated, 3'-amino modifier)

- ② Preparation of the two adapters:

Mix equal volumes of 100 µM ME and 100 µM Adapter 1; separately, mix equal volumes of 100 µM ME and 100 µM Adapter 2. Anneal by slow cooling in a thermal cycler to obtain Adapter 1 and Adapter 2, each at 50 µM.

PCR program: 95°C for 3 min; ramp from 95°C to 25°C over 45 min; hold at 4°C.

- ③ Transposome assembly:

Mix equal volumes of Adapter 1 and Adapter 2 to obtain the adapter mix. Gently mix the enzyme and the adapter mix by pipetting and incubate at room temperature for 1 h.

A 1:1 molar ratio of enzyme to adapter is typically used; the adapter proportion may be increased if needed.

For a 1:1 molar ratio, prepare as follows:

Reagent	Volume
Double-stranded adapter mix (50 pmol/ μ l)	4 μ l
Tn5 Transposase (10pmol/ μ l)	20 μ l

The assembled transposome can be used directly for fragmentation assays or stored at -20°C .

2. Fragmentation Reaction

Set up a 20 μ l fragmentation reaction:

Reagent	Volume
5 \times Reaction buffer	4 μ l
DNA	50-100ng
Transposome complex	1 μ l
ddH ₂ O	To 20 μ l

- Mix by pipetting and incubate at 55°C for 10 min (use a heated lid).
- After the reaction, add 5 μ l of 5 \times Stop Buffer (thaw at room temperature), mix, and incubate at 55°C for 5 min to terminate the reaction.

The fragmented products can be analyzed directly or purified for library construction. If the resulting fragments are too large, increase the amount of transposome complex to reduce fragment size; conversely, decrease the amount of transposome complex if fragments are too small.

FAQ:

Q1: The fragment size distribution is unsatisfactory-fragments are too large or too small.

A: Test a gradient of enzyme amounts and incubation times to determine the optimal conditions for the desired fragment size.

Q2: Low transposition efficiency and insufficient library yield.

A: Improper storage or repeated freeze-thaw cycles of Tn5 may reduce activity; therefore, aliquot storage and avoidance of repeated freeze-thaw are recommended. Alternatively, inhibitors (e.g., SDS, EDTA, high salt concentrations) may be present in the sample-purify the sample before proceeding.s.

IV. Limited Use License and Warranty

Limited Use License

The following terms and conditions apply to the use of **Poly(A) Polymerase** (the Product). If the terms and conditions are not acceptable, the Product in its entirety must be returned to GeneCopoeia within 5 calendar days. A limited End-User license is granted to the purchaser of the Product. The Product shall be used by the purchaser for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged, or modified for resale, or used to manufacture commercial products without prior written consent from GeneCopoeia. This Product should be used in accordance with the NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

Limited Warranty

GeneCopoeia warrants that the Product meets the specifications described in the accompanying Product Datasheet. If it is proven to the satisfaction of GeneCopoeia that the Product fails to meet these specifications, GeneCopoeia will replace the Product. In the event a replacement cannot be provided, GeneCopoeia will provide the purchaser with a refund. This limited warranty shall not extend to anyone other than the original purchaser of the Product. Notice of nonconforming products must be made to GeneCopoeia within 30 days of receipt of the Product. GeneCopoeia's liability is expressly limited to replacement of Product, or a refund limited to the actual purchase price. GeneCopoeia's liability does not extend to any damages arising from use or improper use of the Product, or losses associated with the use of additional materials or reagents. This limited warranty is the sole and exclusive warranty. GeneCopoeia does not provide any other warranties of any kind, expressed or implied, including the merchantability or fitness of the Product for a particular purpose.

GeneCopoeia is committed to providing our customers with high-quality products. If you should have any questions or concerns about any GeneCopoeia products, please contact us at 301-762-0888.

© 2026, GeneCopoeia, Inc.

GeneCopoeia, Inc.

9620 Medical Center Drive, #101, Rockville, MD 20850

Tel: 301-762-0888 Fax: 301-762-3888,

Email: inquiry@genecopoeia.com

Web: www.genecopoeia.com