

R&D systems™
by biotechne

TcBuster™ Non-Viral Gene Delivery System

Now Available Online in Research Grade and GMP

Derisk your genome engineering with TcBuster, a non-viral DNA transposon (plasmid) system that efficiently integrates large DNA cargos into various cell types, such as T, NK, and CHO.

Research grade TcBuster reagents are now available for cell therapy research and development without any licensing or MTA required. GMP TcBuster mRNA also available.

Experience TcBuster in your lab today!

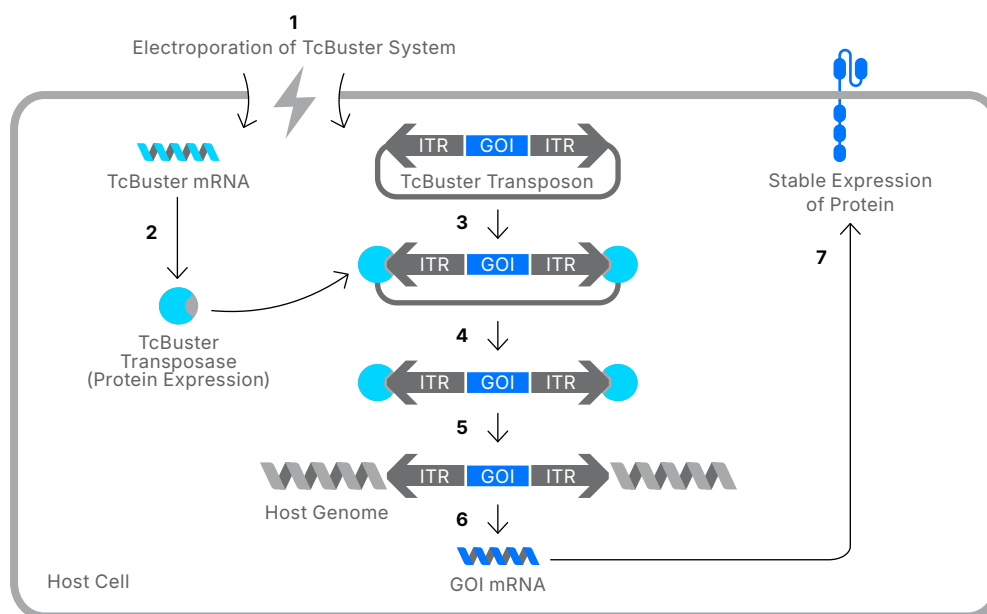


Buy It and Try It!

Scan the QR code or visit
www.rndsystems.com/tcbuster



TcBuster Mechanism Workflow:



1. TcBuster-M mRNA transposase and plasmid DNA transposon containing gene of interest (GOI) are delivered via electroporation.
2. mRNA is translated to TcBuster-M protein.
3. TcBuster-M transposase binds ITRs on plasmid DNA.
4. TcBuster-M transposase cuts plasmid DNA.
5. TcBuster-M randomly inserts DNA into host genome.
6. Inserted GOI is transcribed into mRNA.
7. GOI mRNA is translated into protein.

TcBuster Reagents and Use Cases

TcBuster mRNA Transposase

Catalog No.	Use Cases
TCB-001.1-100 (100 µg)	Useful for system evaluation.
TCB-001.1-500 (500 µg)	Useful for further development.
TCB-001.1-GMP	For clinical and commercial applications.

TcBuster Compatible Transposons

Catalog No.	Use Cases
TCBP001-100 (100 µg)	Multicistronic CAR-containing transposon. Useful in cell therapy applications.
TCBP002-100 (100 µg)	Whole plasmid insertion of eGFP with no untransposed plasmid expression. Useful in bioprocessing system evaluation.
Custom Transposons	For process development using your own therapeutic cargo. In partnership with Aldevron.

Anti-TcBuster-M Transposase Antibody

Catalog No.	Use Cases
MAB11511 (25 µg, 100 µg)	Detects TcBuster-M Transposase in Western blots.

Why Use TcBuster?

- ✓ Lower the cost and accelerate the timeline to cell-based therapy.
- ✓ Achieve greater gene delivery efficiencies compared to viral systems.
- ✓ Reduce risks associated with viral vectors, such as insertional mutagenesis.
- ✓ Streamline workflows with compatibility across multiple cell types, eliminating the need for additional tools or reagents.

TcBuster Common FAQs

Where should I go to purchase TcBuster online?

You can now purchase TcBuster online at:
www.rndsystems.com/tcbuster

What items do I need to start evaluating TcBuster in my lab?

You will need an access to an electroporator, TcBuster transposase (TCB-001.1-100), and an eGFP encoding experimental control transposon plasmid (TCBP001-100 or TCBP002-100) for system evaluation in your cells.

What if I need a custom transposon?

Bio-Techne does not provide custom transposons; instead we have partnered with Aldevron to generate your custom transposons. To discuss all your custom transposase needs, please get started by emailing:
nanoplasmids@aldevron.com.



Learn More About the TcBuster Non-Viral Gene Delivery System

Scan the QR code or visit
www.rndsystems.com/tcbuster