

# GenCrispr T7 Endonuclease I

Cat. No.: Z03396

Version 2018-07-16

---

## Table of Contents

I Description	1
II Contents	1
III Applications	2
IV Quality Control Analysis	2
V Unit Definition	2
VI Storage	2
VII Activity test	3
VIII Protocol	3
IX Notes	4
X Ordering Information	5

---

## I Description

T7 Endonuclease I (T7E1) recognizes and cleaves non-perfectly matched DNA, cruciform DNA structures, Holliday structures or junctions, hetero duplex DNA and more slowly, nicked double-stranded DNA. The cleavage site is at the first, second or third phosphodiester bond that is 5' to the mismatch. The protein is the product of T7 gene 3. GenCrispr T7 Endonuclease I is a fusion protein produced from *E.coli*.

## II Contents

---

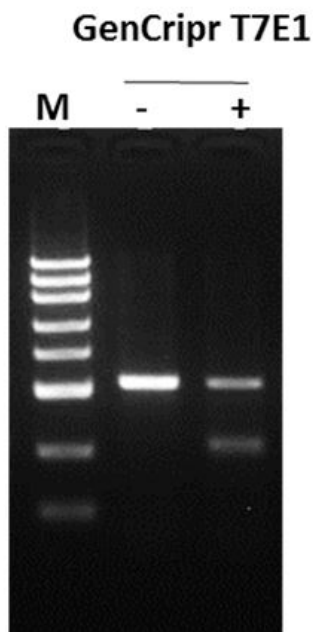
Contents	Quantity	Catalog No.	Components/Concentration
GenCrisprT7 Endonuclease I	250U	Z03396-250	10000 U/mL
	1250U	Z03396-1250	10000 U/mL
10X Reaction Buffer	1.0 mL		500 mM NaCl, 100mM Tris-HCl, 100 mM MgCl <sub>2</sub> , 10 mM DTT, pH 7.9 at 25°C

---



## VII Activity test

To test the activity of GenCrispr T7 Endonuclease I, GenScript used a control gRNA targeting HPRT, which is co-transfected with Cas9 protein into 293T cells. After 48 h, cells were lysed for genome PCR to amplify the specific target site. The PCR product ( 200 ng) was then annealed and incubated with 1  $\mu$ L T7E1 for 15 min at 37  $^{\circ}$ C. Add loading buffer to the reaction mixture directly, and detect the cleavage efficiency by agarose gel electrophoresis.



The T7E1 assay to detect the cleavage efficiency of control HPRT gRNA with Cas9 protein.

## VIII Protocol

### PCR

Set up a 50  $\mu$ l PCR reaction using 100 ng of genomic DNA as a template. For each amplicon set up 2 PCR reactions using the following templates:

- ◇ gDNA from targeted cells (e.g. Cas9, or TALEN transfected cells)
- ◇ gDNA from negative control cells (e.g. non-specific DNA transfected cells)

The amplification fragment is usually around 500 bp and the target site is better to avoid the middle of fragment.

## T7 Endonuclease I digestion:

Assemble reactions as follows:

PCR Products	200 ng
10 X reaction buffer	2 $\mu$ L
Nuclease-free Water	To 19 $\mu$ L

Anneal the above PCR products in a thermocycler using the following conditions:

Initial Denaturation	95°C	5 minutes
Annealing	95-85°C	-2°C/second
	85-25°C	-0.1°C/second
Hold	4°C	

Add 1  $\mu$ L T7 Endonuclease I to the annealed PCR products.  
Incubation at 37°C for 15 minutes.

## Detection:

Add loading buffer to the reaction mixture directly, and detect the cleavage efficiency by agarose gel electrophoresis.

## IX Notes

T7 Endonuclease I is a structure-selective enzyme. It acts on a variety of DNA substrates with different specific activities. It is important to control the amount of enzyme and the reaction time used for cleavage of a particular substrate. Temperatures above 42°C cause an increase in nonspecific nuclease activity and should be avoided.

## X Ordering Information

<b>Product Name</b>	<b>Cat. No.</b>
GenCrispr Cas9-C-NLS Nuclease	Z03385
GenCrispr Cas9 Nuclease	Z03386
GenCrispr Cas9-N-NLS Nuclease	Z03388
GenCrispr NLS-Cas9-NLS Nuclease	Z03389
GenCrispr NLS-Cas9-D10A Nickase	Z03390
GenCrispr NLS-Cas9-EGFP Nuclease	Z03393
GenCrispr T7 Endonuclease I	Z03396
GenCrispr Mutation Detection Kit	L00688
GenCrispr sgRNA Screening Kit	L00689
High-Efficiency gRNA-Cas9-GFP Plasmid (linear) Assembly Kit	L00690
High-Efficiency gRNA-Cas9-Puro Plasmid (linear) Assembly Kit	L00691
High-Efficiency gRNA-Cas9-GFP Plasmid Assembly Kit	L00692
High-Efficiency gRNA-Cas9-Puro Plasmid Assembly Kit	L00693

## Contact us

Web: <https://www.genscript.com>

Email: [product@genscript.com](mailto:product@genscript.com)

Fax: 1-732-518-5150