

GenCrispr NLS-Cas9-EGFP Nuclease Cat. No. Z03467



I	Description.....	1
II	Kit Contents.....	2
III	Key Features.....	2
IV	Quality Control Analysis.....	2
V	Utilities of Product.....	2
VI	Storage.....	2
VII	Diluent Compatibility.....	3
VIII	Activity Test.....	3
IX	References.....	4



I Description

GenCrispr NLS-Cas9-EGFP is a fusion protein developed by GenScript. It contains a nuclear localization sequence (NLS) on its N terminal and EGFP on the C terminal. Cas9 nuclease is an RNA-guided endonuclease that can catalyze cleavage of double stranded DNA. This kind of targeted nuclease is a powerful tool for genome editing with high precision. Cas9 protein forms a very stable ribonucleoprotein (RNP) complex with the guide RNA (gRNA) component of the CRISPR/Cas9 system. The Cas9 RNP complex can localize to the nucleus immediately upon entering the cell with the addition of a nuclear localization signal (NLS). There is no requirement for transcription and translation compared with mRNA or plasmid systems. Additionally, the Cas9 RNP complex is rapidly cleared from the cells minimizing the chance of off-target cleavage when compared to other systems (Kim, et al. 2014). The EGFP can be used as a reporter for tracking or sorting transfected cells, which enables the possibility of enriching cell populations for desired genome edits via fluorescence activated cell sorting (FACS). It significantly reduces the labor and cost associated with single cell cloning and genotyping in genome editing applications.

Product Source: GenCrispr NLS-Cas9-EGFP is produced by expression from an *E. coli* strain.

II Kit Contents

Kit Contents	Quantity	Catalog No.	Components/Concentration
GenCrispr NLS- Cas9-EGFP Nuclease	100 µg 500 µg 1 mg	Z03467-100 Z03467-500 Z03467-1	10 mg/mL
10X Reaction Buffer	1.5 mL		Reagent composition confidentiality, 1000µg/mL BSA pH 7.9 at 25°C

III Key Features

- **DNA-free: no external DNA added to system**
- **High cleavage efficiency: NLS ensures the entry of Cas9 protein into nuclei**
- **Low off target: transient expression of Cas9 nuclease**
- **Time-saving: no need for transcription and translation**

IV Quality Control Analysis

- **High Protein purity:** GenCrispr Cas9 is >95% pure as determined by SDS-PAGE using Coomassie Blue detection and analysis by Image Lab.
- **Low Endotoxin:** Endotoxin level is <0.1EU/µg test by Tachypleus Amebocyte Lysate (TAL) gel-clot method: limit test.
- **Non-specific DNase activity:** A 20 µL reaction in Cas9 reaction buffer containing 100 ng linearized pUC57 plasmid and 0.1 µg GenCrispr Cas9 incubated for 16h at 37°C. No DNA degradation is determined by agarose gel electrophoresis.
- **Non-specific RNase activity:** A 10 µL reaction in Cas9 reaction buffer containing 1800 ng total RNA and 0.1 µg of GenCrispr Cas9 incubated for 2 h at 37°C. No RNA degradation as determined by Agarose gel electrophoresis.

High Bioactivity: 50 ng GenCrispr Cas9 incubated for 2 hours at 37°C result in 90% digestion of the substrate DNA as determined by agarose gel electrophoresis and compared with competing products.

V Utilities of Product

1. Screening for highly efficient and specific targeting gRNAs by *in vitro* DNA cleavage.
2. *In vivo* gene editing when combined with a specific gRNA by electroporation or injection.

VI Storage

GenCrispr NLS-Cas9-EGFP Nuclease is supplied with 1X storage buffer (10 mM Tris, 300 mM NaCl, 0.1 mM EDTA, 1 mM DTT, 50% Glycerol PH7.4, at 25°C) and recommended to be stored at -20°C. Guaranteed stable for 18 months when properly stored.

VII Diluent Compatibility

In *in vitro* assays, it is recommended to dilute Cas9 nuclease with Diluent Buffer B (validated in our *in vitro* tests).

Diluent Buffer B: 300 mM NaCl, 10 mM Tris-HCl, 0.1 mM EDTA, 1 mM DTT, 50% glycerol. (pH 7.4, at 25°C).

VIII Activity Test

Cas9 site-specific digestion:

GenScript used *in vitro* digestion of a linearized plasmid to determine the activity of the Cas9 nuclease. It is a sensitive assay for GenCrispr Cas9 quality control. The linearized plasmid containing the target site:

(CATCATTGGAAAACGTTCTT)

can be digested with gRNA:

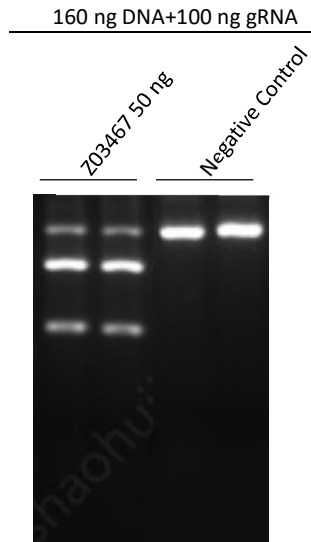
(CAUCAUUGGAAAACGUUCUUGUUUUAGAGCUAGAAAUAGCAAGUAAAAUAA
GGCUAGUCCGUUAUCAACUUGAAAAGUGGCACCGAGUCGGUGCUUUUUUU)

and GenCrispr Cas9. Two cleavage DNA fragments (526 bp and 1053 bp) are determined by agarose gel electrophoresis. A 20 µL reaction in 1xCas9 Nuclease Reaction Buffer containing 160 ng linearized pUC57, 100 ng gRNA and 50 ng GenCrispr Cas9 for 2 h at 37°C results in over 90% digestion of linearized pUC57, as determined by agarose gel electrophoresis and compared with competing products.

Reagent	Volume
10xCas9 Nuclease Reaction Buffer	2 µL
gRNA	100 ng
GenCrispr Cas9 Nuclease	50 ng
DEPC water	Up to 19 µL

Procedure:

1. The above components should be gently mixed, centrifuged, placed in a PCR machine and incubated at 37°C for 10 minutes.
2. 1 µL (160 ng) of linearized pUC57 was added, mixed and centrifuged, incubated at 37°C for 2h.
3. 5 µL of cleavage products are resolved on a 1.5% agarose gel.



In vitro DNA cleavage assay with GenCrispr NLS-Cas9-EGFP Nuclease

Reactions were set up according to recommended conditions, and cleavage products were resolved on a 1.5% agarose gel.

IX References

1. Jinek et al. A Programmable Dual-RNA–Guided DNA Endonuclease in Adaptive Bacterial Immunity. (2012) *Science* 337 (6096) 816-821 (2012).
2. Larson, M. H., et al. CRISPR interference (CRISPRi) for sequence-specific control of gene expression. *Nature Protocols*. 8, (11), 2180-2196 (2013).
3. Ran, F. A., et al. Genome engineering using the CRISPR-Cas9 system. *Nature Protocols*. 8, (11), 2281-2308 (2013).
4. Kim, S., Kim, D., Cho, S.W., Kim, J., Kim, J.S, (2014) Highly efficient RNA-guided genome editing in human cells via delivery of purified Cas9 ribonucleoprotein. *Genome Res*. 24(6), 1012-1019.

Note:

1. This is a basic protocol. The reagent concentrations, conditions, and parameters may need to be optimized.
2. 1000 nM is equal to 190 ng/ μ L.

GenScript US
860 Centennial Ave., Piscataway, NJ 08854
Tel: 732-885-9188, 732-885-9688
Fax: 732-210-0262, 732-885-5878
Email: product@genscript.com
Web: <http://www.genscript.com>
For Research Use Only.

For laboratory research use only. Direct human use, including taking orally and injection and clinical use are forbidden.