

EnGen® sgRNA Synthesis Kit, *S. pyogenes*

NEB #E3322V/S

10/20 reactions

Version 6.0_8/26

Table of Contents

Section 1	
Protocol for Target-specific Oligo Design.....	4
Section 2	
Protocol for sgRNA Synthesis.....	5
Purification of sgRNAs.....	5
Evaluation of sgRNA Transcription Products	
Measuring the yield of sgRNA synthesis products.....	5
Analysis of sgRNAs by Gel Electrophoresis.....	6
<i>In vitro</i> Cas9 Nuclease Assay	6
Troubleshooting	7
Ordering Information	8
Revision History	8

The EnGen sgRNA Synthesis Kit, *S. pyogenes* Includes:

All kit components should be stored at –20°C. Each standard reaction yields between 4–25 µg of sgRNA.

EnGen sgRNA Enzyme Mix

EnGen 2X sgRNA Reaction Mix, *S. pyogenes*

DNase I (RNase-free)

EnGen sgRNA Control Oligo, *S. pyogenes*

Dithiothreitol (DTT)

Required Materials Not Included:

Target-specific DNA oligo(s)

Thermocycler/37°C heat block/incubator

Nuclease-free water

Equipment and reagents for RNA quantitation

Monarch® Spin RNA Cleanup Kit (50 µg) (NEB #T2040)

RNase-free tubes, aerosol tips

Microcentrifuge

Optional Materials:

Quick CIP

Cas9 Nuclease, *S. pyogenes*

EnGen Spy Cas9, NLS

Gels, running buffer, gel loading dye, RNA and DNA ladders, gel box

Introduction

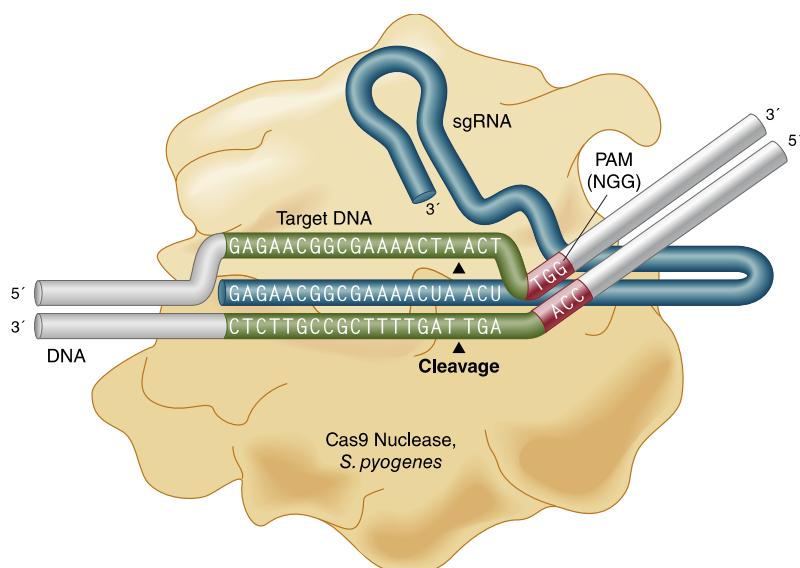
CRISPR is an acronym for Clustered Regularly Interspaced Short Palindromic Repeats which are genomic loci found in many bacteria and archaea. The CRISPR/Cas9 pathway, which naturally allows for the elimination of genomic material from invading sources in bacteria, has recently been adapted as a molecular biology tool to edit genomes in a target-specific manner. Cas9 (CRISPR-associated protein 9) is a double-stranded DNA endonuclease which forms an active ribonucleoprotein (RNP) when complexed with guide RNAs (gRNAs) encoded at the CRISPR loci. gRNAs provide sequence specificity to the RNP, directing the Cas9 nuclease to DNA targets resulting in double-strand breaks in the DNA.

In nature, *S. pyogenes* Cas9 is programmed with two separate RNAs, the crRNA and tracrRNA. The crRNA, or CRISPR RNA sequence contains approximately 20 nucleotides of homology complementary to the strand of DNA opposite and upstream of a PAM (Protospacer Adjacent Motif) (NGG) sequence. The tracrRNA, or transactivating crRNA, contains partial complementary sequence to the crRNA as well as the sequence and secondary structure that is recognized by Cas9. These sequences have been adapted for use in the lab by combining the tracrRNA and crRNA into one long single guide RNA (sgRNA) (1) species capable of complexing with Cas9 to recognize and cleave the target DNA (Figure 1).

When gRNA-programmed Cas9 RNPs are introduced into eukaryotic cells, dsDNA breaks are introduced at target sites. Activation of the cellular double-strand break machinery can lead to insertions and/or deletions (indels) through Non-Homologous End Joining (NHEJ), resulting in disruption of the gene at that specific locus. In the presence of a homologous repair template, the homology-directed repair (HDR) pathway can be activated leading to the introduction of specific changes to the DNA at the targeted site.

1. Jinek, M. et al. (2012) *Science* 816–821. PubMed ID: 22745249.

Figure 1. Cas9 Nuclease, *S. pyogenes*



*Cas9 nuclease, *S. pyogenes*, complexed with an sgRNA. Cleavage occurs three nucleotides upstream of the PAM sequence (shown in red). sgRNAs are complimentary to the strand of DNA opposite of the PAM.*

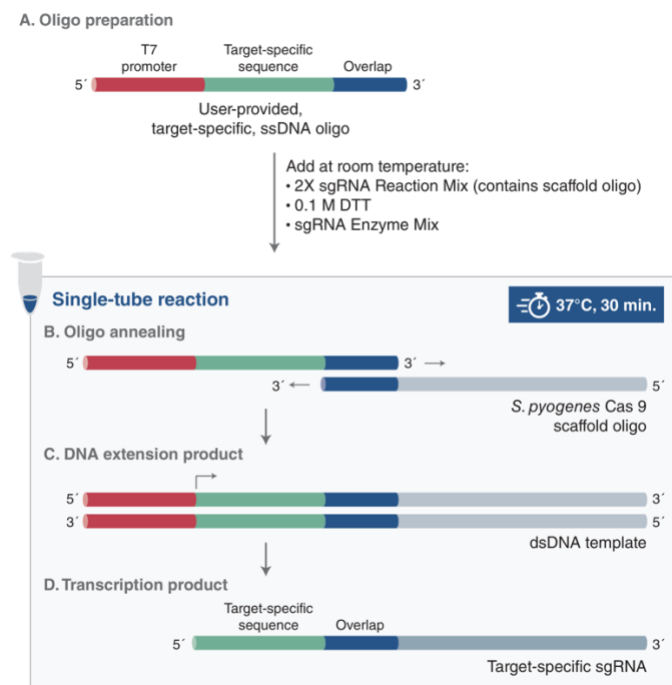
The EnGen sgRNA Synthesis Kit, *S. pyogenes* combines an *S. pyogenes* Cas9-specific Scaffold Oligo (included in the EnGen 2X sgRNA Reaction Mix) that is partially complementary to the target-specific oligos designed by the user. The two oligos anneal at the overlapping region and are filled in by the DNA polymerase, creating a double-stranded DNA (dsDNA) template for transcription. Synthesis of the dsDNA template and transcription of RNA occur in a single reaction, resulting in the generation of a functional sgRNA.

The EnGen sgRNA Synthesis Kit, *S. pyogenes* provides a simple and quick method for transcribing high yields of sgRNA in a single 30 minute reaction, using the supplied reagents and target-specific DNA oligos designed by the user.

Figure 2. General workflow for the EnGen sgRNA Synthesis Kit, *S. pyogenes*



Figure 3. EnGen sgRNA Synthesis Kit Overview



A. The target-specific oligo contains the T7 promoter sequence, ~20 nucleotides of target-specific sequence and a 14 nucleotide overlap sequence complementary to the *S. pyogenes* Cas9 Scaffold Oligo supplied in the reaction mix. Target-specific oligos are mixed with the EnGen 2X sgRNA Reaction Mix, 0.1 M DTT and the EnGen sgRNA Enzyme Mix (DNA and RNA polymerases) at room temperature. B. At 37°C the two oligos anneal at the 14-nucleotide overlap region of complementarity. C. The DNA polymerase contained in the EnGen sgRNA Enzyme Mix extends both oligos from their 3' ends creating a dsDNA template. D. The RNA polymerase contained in the EnGen sgRNA Enzyme Mix recognizes the dsDNA of the T7 promoter and initiates transcription, resulting in a target-specific sgRNA. All steps occur in a single reaction during a 30 minute incubation at 37°C.

Section 1

Protocol for Target-specific Oligo Design

This kit contains the *S. pyogenes* Cas9 Scaffold Oligo within the EnGen 2X sgRNA Reaction Mix, *S. pyogenes*. Target-specific oligos are designed by the user as follows:

1. Select a 20 nucleotide target sequence [not including the PAM (NGG) sequence]. We recommend using a target DNA selection program, such as [ChopChop](#).
2. Input the 20 nucleotide target sequence into our [EnGen® sgRNA Template Oligo Designer tool](#) (without the PAM site). An oligo sequence will be generated from the target sequence that you provide, which will include the T7 promoter sequence at the 5' end and a 14-nt overlap sequence at the 3' end. If there is no G at the 5' end of the target sequence (required for T7 RNA Polymerase transcription), the tool will add a G to the generated oligo.

Example using the provided EnGen sgRNA Control Oligo and *S. pyogenes* Cas9-specific Scaffold Oligo (Included in the EnGen 2X sgRNA).

Get started designing oligos.

EnGen sgRNA Template Oligo Designer can be used to design target-specific DNA oligos for use with the EnGen™ sgRNA Synthesis Kit, *S. pyogenes*. To determine your 20-nucleotide target sequence, we recommend using a target prediction program such as [ChopChop](#).

1. Enter your 20 nt target sequence.

Name/ID	Target DNA seq (5'→3') - NO PAM -	
EnGen sgRNA Control Oligo	TGCAACCTTCATTCCCTGC	[20 nt]
Note: G automatically added to 5' end if not present.		

2. Check generated oligo result.

Target-specific DNA oligo template (5'→3'):

TTCTAATACGACTCACTATAGTGCAACCTTCATTCCCTGC**GT**TTTAGAGCTAGA

Transcribed sgRNA sequence (5'→3'):

GUGCAACCUUCAUUUCCUGCGUUUUAGAGCUAGAAUAGCAAGUAAAAUAAGGCUAGUCCGUUAUCAACUUGAAAAAGUGGCACCGAGUCGGUGCUUUU

■ T7 Promoter ■ Added 5'G ■ Target DNA ■ Overlap/RNA Scaffold

1. Select target-specific DNA sequence (we recommend using a target DNA selection program):

TGCAACCTTCATTCCCTGCTGG

The PAM sequence (NGG, underlined) is required for Cas9 recognition of the target sequence and is **NOT** part of the sgRNA sequence. Remove these three nucleotides:

5' **TGCAACCTTCATTCCCTGC**

2. Check for "G" at the 5' end of the target sequence (at least one G is necessary for transcription). If no G is present, add one G:

5' **G**TGCAACCTTCATTCCCTGC 3'

3. To the 5' end, append the T7 promoter sequence (shown in red):

5' **TTCTAATACGACTCACTATAG**TGCAACCTTCATTCCCTGC 3'

4. To the 3' end, append the 14 nucleotide overlap sequence (shown in blue and underlined). This represents the oligo sequence to be ordered:

5' **TTCTAATACGACTCACTATAG**TGCAACCTTCATTCCCTGC**GT**TTTAGAGCTAGA 3'

5. The sequence of the *S. pyogenes* Cas9 Scaffold Oligo in the EnGen 2X sgRNA Reaction Mix (with reverse complimentary overlap shown in blue and underlined) is:

5' AAAAGCACCGACTCGGTGCCACTTTTTCAAGTTGATAACGGACTAGCCTTATTTAACTTGCTATT**TCTAGCTCTAAAC** 3'

In this example, the final sgRNA sequence of the provided control would be:

5' **GUGCAACCUUCAUUUCCUGCGUUUUAGAGCUAGAAUAGCAAGUAAAAUAAGGCUAGUCCGUUAUCAACUUGAAAAAGUGGCACCGAGUCGGUGCUUUU** 3'

Section 2

Protocol for sgRNA Synthesis

We strongly recommend wearing gloves and using nuclease-free tubes and reagents. Reactions should be assembled in microfuge tubes or PCR strip tubes.

1. Thaw the EnGen 2X sgRNA Reaction Mix, *S. pyogenes*, 0.1 M DTT, and the user-supplied target-specific oligo from Part I. Mix and pulse-spin each component in a microfuge prior to use. Keep the EnGen sgRNA Enzyme Mix on ice.
2. Assemble the reaction at **room temperature in the order listed**. Avoid master mixes, and add the enzyme last to each reaction:

COMPONENT	20 μ l REACTION	FINAL AMOUNT
Nuclease-free water	2 μ l	–
EnGen 2X sgRNA Reaction Mix, <i>S. pyogenes</i>	10 μ l	1X
Target-specific DNA Oligo (1 μ M)	5 μ l	250 nM
DTT	1 μ l	5 mM
EnGen sgRNA Enzyme Mix	2 μ l	–

3. Mix thoroughly and pulse-spin in a microfuge. Incubate at 37°C for 30 minutes.
4. Transfer the reaction to ice. To remove template DNA, add 30 μ l of Nuclease-free water and 2 μ l of the provided DNase I (RNase-free). Mix and incubate at 37°C for 15 minutes.
5. Optional – The EnGen sgRNA Kit reaction product is 5' triphosphorylated ssRNA, which can be converted to a 5' hydroxyl group by treating the sgRNA with Quick CIP (NEB #M0525). Phenol: chloroform extraction and/or spin-column cleanup is recommended to inactivate and remove the phosphatase.

Purification of sgRNAs

Spin columns will remove proteins, salts and most unincorporated nucleotides. We recommend following the standard protocol for the Monarch RNA Cleanup Kit (50 μ g) (NEB #T2040) to purify the reaction products. For columns from other manufacturers, please ensure that they are compatible with the size of sgRNAs (~100 nts) and expected RNA yields (4–25 μ g), and follow the manufacturer's instructions.

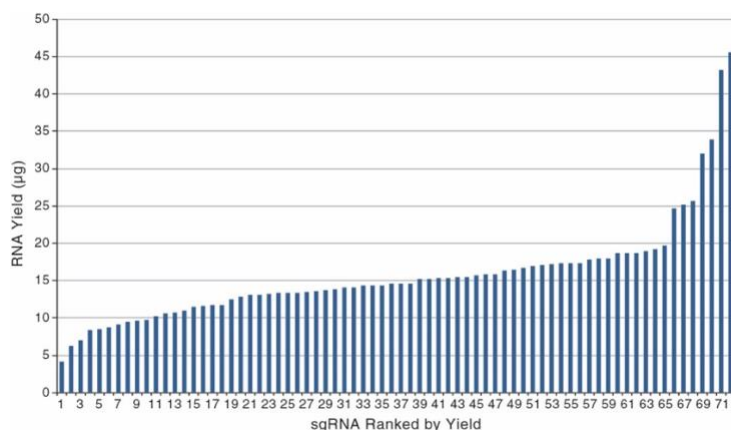
Evaluation of sgRNA Transcription Products

Measuring the yield of sgRNA transcription products

Following purification, we recommend determining sgRNA concentration directly on a NanoDrop®. The supplied EnGen *S. pyogenes* positive control performed using the protocol above should result in > 15 μ g sgRNA after purification. Alternatively, absorbance can be measured at 260 nm on a standard UV spectrophotometer following dilution of the sample 1 OD₂₆₀ unit for single-stranded RNA (ssRNA) is 40 μ g/ml. RNA concentration can be determined as follows:

$$A_{260} \times \text{dilution factor} \times 40 = "x" \mu\text{g/ml ssRNA.}$$

Figure 4. Range of sgRNA yield from the EnGen sgRNA Synthesis Kit, *S. pyogenes*.



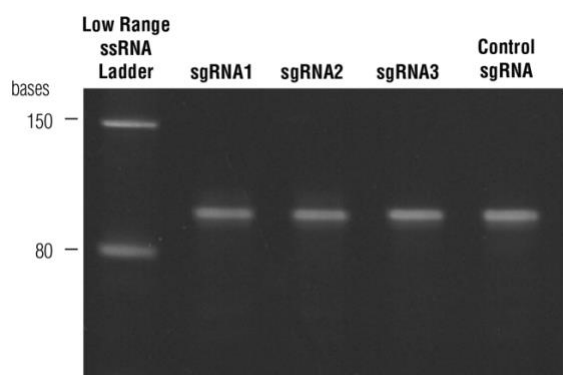
72 different target-specific oligos were designed and tested for RNA yield following DNase I treatment and column purification as determined by NanoDrop. RNA yields may be dependent on specific oligo sequence, with most falling between 4–25 μ g.

Analysis of sgRNAs by Gel Electrophoresis

Assessment of sgRNA quality and length can be evaluated by gel electrophoresis. Gels should be run under denaturing conditions to avoid formation of secondary structure.

1. Mix 20 ng of sgRNA with 20 μ l RNA Loading Dye, (2X) (NEB #B0363).
2. Denature at 70°C for 10 minutes, then load 5–10 μ l onto a TBE-urea gel, such as Novex® TBE-urea gels.
3. Visualize RNA by staining the gel with SYBR® Gold (preferred) or ethidium bromide. The supplied EnGen sgRNA Control Oligo, *S. pyogenes* positive control performed using the protocol above should result in a single band of ~100 nucleotides on a denaturing TBE-Urea gel under denaturing conditions.

Figure 5. Examples of sgRNAs synthesized using the EnGen sgRNA Synthesis Kit, *S. pyogenes* and multiple different target-specific oligos including the EnGen sgRNA Control Oligo, *S. pyogenes*.



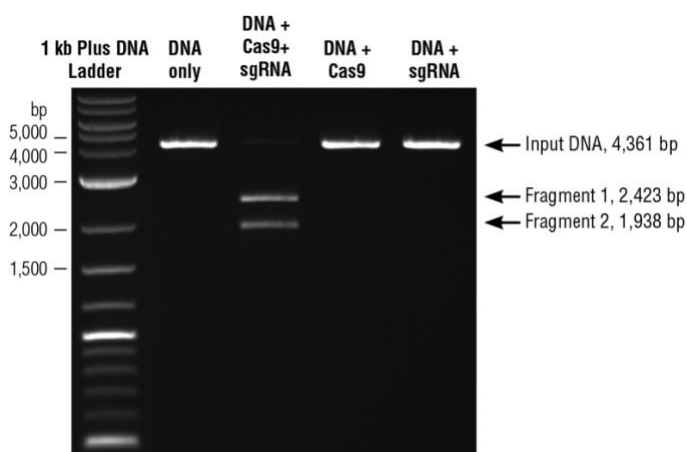
20 ng of RNA was run under denaturing conditions on a 10% Novex TBE-Urea gel and post-stained with SYBR Gold.

In vitro Cas9 Nuclease Assay

S. pyogenes Cas9 double-stranded DNA nuclease is guided to the target DNA based on sequence complementarity to the sgRNA that is loaded into the protein. Site-specific DNA cleavage occurs three nucleotides upstream of the PAM sequence (NGG), within the complementary region. sgRNAs synthesized with this kit can be used with Cas9 Nuclease, *S. pyogenes* (NEB #M0386) or EnGen Spy Cas9, NLS (NEB #M0646) for *in vitro* cleavage reactions.

Please see our detailed protocol, [In vitro digestion of DNA with Cas9 Nucleases \(NEB #M0386, #M0646, #M0650, #M0654, #M0667, #M0668\)](#).

Figure 6. Example of an *in vitro* Cas9 nuclease assay.



DNA target is PvuII-linearized pBR322, sgRNA was synthesized using the EnGen sgRNA Control Oligo, *S. pyogenes*. Reactions were set up following NEB® protocols with a ratio of 20:20:1 (Cas9:sgRNA:target site) in 30 μ l reactions. Aliquots of reactions were mixed with Gel Loading Dye, Purple (6X) (NEB #B7024). Cleavage products and 1 kb Plus DNA Ladder (NEB #N3200) were resolved on a 1% TBE agarose gel stained with ethidium bromide.

Troubleshooting

Positive Control Reaction

- The control oligo (EnGen sgRNA Control Oligo, *S. pyogenes*) provided with this kit can be used in place of the target-specific oligo to test the efficiency of the reaction. Following incubation for 30 minutes at 37°C, the control reaction should contain $\geq 15 \mu\text{g}$ sgRNA, post-purification, as determined by NanoDrop. Resulting sgRNA should run as a single band of ~ 100 nucleotides on a denaturing TBE-Urea gel under denaturing conditions.

Low Yield of RNA

- If the sgRNA yield using the control oligo is lower than expected, there may be issues during the reaction setup. It is very important to follow the directions exactly; components should be added in the exact order listed, making sure that the EnGen sgRNA Enzyme Mix is added last. Reactions should be set up at room temperature. Make sure that the heat block/incubator or thermocycler is at 37°C.

If the sgRNA yield using the control oligo is at the level expected, then there may be an issue with the design of the target-specific oligo(s). Make sure to follow the design protocol, and double-check the sequence before ordering. It is important to have at least one G following the T7 promoter sequence and to have a 14 nucleotide overlap with the *S. pyogenes* Cas9 Scaffold Oligo. If a higher yield is required, a longer incubation at 37°C is reasonable (1–2 hours). Use of a dry air incubator or thermocycler is recommended for incubations > 30 minutes.

As we have observed, transcription can be somewhat sequence dependent, though the yield of the majority of sgRNA synthesis reactions falls between 4–25 μg . Designing multiple sgRNAs per target is recommended.

sgRNA Smearing on Denaturing Gel

- We strongly recommend wearing gloves and using nuclease-free tubes and reagents. If the RNA appears degraded or smeared on a denaturing gel, it is likely that RNase contamination has been introduced at some step during the reaction. The sgRNAs synthesized from this kit should yield a single band of ~ 100 nucleotides on a TBE-Urea gel.

Multiple Bands are Detected on Denaturing Gel

- If more than one band appears on the gel it is likely that the RNA is not fully denatured due to the presence of strong secondary structure. It is important to run these reactions under denaturing conditions (using urea dye or formamide dye and heating the sample according to the manufacturer's directions prior to loading).

It is important to DNase treat the sample prior to gel analysis. DNA template present in the sgRNA synthesis reaction prior to DNase treatment may be detected on the gel as a slightly larger band compared to the RNA.

Ordering Information

NEB #	PRODUCT	SIZE
E3322V/S	EnGen sgRNA Synthesis Kit, <i>S. pyogenes</i>	10/20 reactions

COMPANION PRODUCTS

NEB #	PRODUCT	SIZE
M0386S/T/M	Cas9 Nuclease, <i>S. pyogenes</i>	90/500/2,500 pmol
M0646T/M	EnGen Spy Cas9, NLS	500/2,500 pmol
M0650S/T	EnGen Spy Cas9 Nickase	90/500 pmol
M0652S/T	EnGen Spy dCas9 (SNAP-tag®)	90/500 pmol
E3321S	EnGen Mutation Detection Kit	25 reactions
N0364S	Low Range ssRNA Ladder	100 gel lanes
B0363S	RNA Loading Dye, (2X)	4 ml
B7024S	Gel Loading Dye, Purple (6X)	4 ml
N3200S/L	1 kb Plus DNA Ladder	200/1,000 gel lanes
M0525S/L	Quick CIP	1,000/5,000 units
T2040S/L	Monarch® Spin RNA Cleanup Kit (50 µg)	10/100 preps

Revision History

REVISION #	DESCRIPTION	DATE
1.0	N/A	5/16
2.0		7/18
3.0		1/19
4.0	Apply new manual format and add a “V” size.	2/20
5.0	Apply new format and legal text and adjust an optional step to include Quick CIP	8/25
6.0	Cleaned up example sequence text and included the EnGen sgRNA Template Design Tool. Figure 3 updated. Minor protocol updates. Updated companion products list.	8/26

This product is intended for research purposes only. This product is not intended to be used for therapeutic or diagnostic purposes in humans or animals.

Products and content are covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc (NEB). The use of trademark symbols does not necessarily indicate that the name is trademarked in the country where it is being read; it indicates where the content was originally developed.

See www.neb.com/trademarks. The use of these products may require you to obtain additional third-party intellectual property rights for certain applications. For more information, please email busdev@neb.com.

B CORPORATION™ is a trademark of B Lab, Inc.

NANODROP® is a registered trademark of NanoDrop Technologies, Inc.

NOVEX® is a registered trademark of Life Technologies Corporation.

SYBR® is a registered trademark of Molecular Probes, Inc.

New England Biolabs is an ISO 9001, ISO 14001 and ISO 13485 certified facility

© Copyright 2026, New England Biolabs, Inc.; all rights reserved.

