

SureGuide gRNA Synthesis Kit

Data Sheet

Key Features

- Complete, easy-to-use gRNA expression kit; obtain ready-to-use guide RNAs in less than 2.5 hours
- Validated synthetic system to provide ready to use guide RNAs
- Enzymes and reagents optimized for maximum gRNA yields and consistent performance.



Complete, reliable, and validated kit to generate high concentrations of gRNA for CRISPR/Cas applications.

The SureGuide gRNA synthesis kit is designed for the rapid and simple preparation of gRNAs to be used in conjunction with the SureGuide Cas9 system (part no 5190-7714). It has been optimized for in vitro expression and contains all necessary components to prepare purified gRNA from a user-provided DNA template. gRNA expression using the SureGuide system can take as little as an hour to obtain micro-molar yields, enough for around 400 in vitro Cas9 digests.



High gRNA yields are obtained even with low DNA template concentrations



Figure 1. The effect of template input on yield of gRNA. Decreasing concentrations of template were subjected to 1 hour of vitro transcription, showing that even low concentrations of template DNA are sufficient to obtain high yields of gRNA.



1 hour is sufficient to obtain gRNA for CRISPR/Cas experiment

Figure 2. Three gene fragments were subjected to in vitro transcription for 1, 2, or 20 hours. gRNA yields after only 1 hour provided sufficient material for completing nearly 400 in vitro Cas9 digests. Additionally, yield continued to increase linearly over time indicating that PCR amplification would be unnecessary even for low concentration templates.

Agilent SureGuide Product Portfolio

PN	Description	Enzyme & Buffer	sgRNA Control	Ιντ
5190-7714	SureGuide CRISPR/Cas Complete Kit, 40rx	\checkmark	\checkmark	\checkmark
5190-7719	SureGuide gRNA Synthesis Kit, 50rx			~
5190-7717	SureGuide Cas9 Programmable Nuclease, 100rx	✓		
5190-7718	SureGuide gRNA Control Kit, 20rx		~	
5190-7715	SureGuide Cas9 Programmable Nuclease Kit, 20rx	✓	~	
5190-7716	SureGuide Cas9 Programmable Nuclease Kit, 100rx	✓	✓	

www.agilent.com/genomics/sureguide

For Research Use only. Not for use with diagnostics procedures. © Agilent Technologies, Inc., 2014 Published in USA, August 29, 2014 Publication Number 5991-5148EN

