

SureGuide Cas9 Programmable Nuclease Kit

Data Sheet

Key Features

- CRISPR/Cas *in vitro* system in a complete and ready-to-use kit
 - Perform unrestricted *in* vitro cloning of large DNA fragments without the limitation imposed by common Restriction Enzymes or PCR
 - Clone complex DNA libraries with high specificity
 - Blunt ends ready for the next molecular step
- Fast & Easy synthesis and purification of guide RNAs with a single kit



The SureGuide Cas9 Programmable Nuclease Kit is the basis of the next-generation genetic engineering tool.

The full kit contains a purified, recombinant Cas9 *S. pyogenes* nuclease, the SureGuide gRNA in vitro expression and purification system, control target and sgRNA templates, and all the necessary reagents to get you started with the CRISPR/Cas9 system. Agilent's SureGuide product line provides a set of tools for genetic engineering that are extensively validated by our scientists, allowing you to obtain results you can be sure of.

The only complete, in vitro CRISPR/Cas9 system contains everything you need to immediately explore and benefit from this new technology.

Enjoy faster start-up times, and better results.



Cas9 programmable nuclease experimental workflow



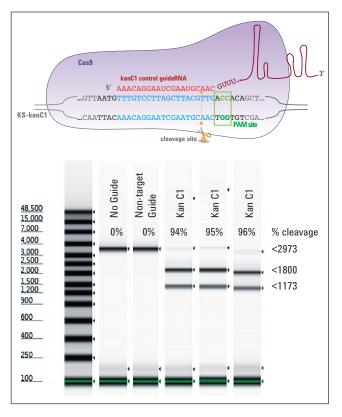


Figure 2. Cas9 digestion of KanC1 linear plasmid with three KanC1 Guide RNA controls (5190-7718) analyzed on the Agilent TapeStation. Cleavage is typically greater than 90%.

Ordering Instructions

PN	Description	Enzyme & Buffer	sgRNA Control	IVT
5190-7714	SureGuide CRISPR/Cas Complete Kit, 40rx	~	~	~
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, September 1, 2014 Publication Number 5991-5149EN



Agilent Technologies