

GeneHero™ Cas9 NLS, *S. pyogenes* protein

Catalog number:	GE001
Product:	Cas9 Nuclease, <i>Streptococcus pyogenes</i>
Expressed in:	<i>E. coli</i>
Description:	GeneHero™ Cas9 Nuclease is a recombinant protein, purified for genome engineering experiments. It contains the nuclear localization sequence (NLS), which aids in transport of Cas9 protein to the nucleus. Cas9 protein forms a stable complex with sgRNA and can be used for both <i>in vitro</i> and <i>in vivo</i> applications.
Size:	25 µg
Concentration:	1 mg/ml
Storage buffer:	10 mM Tris-HCl , 0.3 M NaCl , 1 mM DTT , 0.1 mM EDTA , 50 % Glycerol, (pH 7.4)
Storage:	-20 °C
Protein tag/fusion:	Untagged
Shelf life:	24 months from the date of manufacture
Quality control:	Each lot is tested by SDS-PAGE for purity, <i>in vitro</i> and <i>in vivo</i> cleavage assays, endonuclease activity, DNase/RNase activity and bacteria/fungi/mycoplasma contamination assessment. Please refer to the CoA for details.

Citation of product: If use of this item results in a publication, please use this information: Cas9 NLS, *S. pyogenes* (GE001, GeneCopoeia, Inc., Rockville, MD).

Limited Use License

A limited use license is granted to the Buyer of the Product. The Product shall be used by the Buyer for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale, or used to manufacture commercial products without prior written consent from GeneCopoeia. This Product should be used in accordance with NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

Copyright ©2018 GeneCopoeia, Inc.

C9SCL-DS-060618

GeneCopoeia, Inc.

9620 Medical Center Drive, #101

Rockville, MD 20850 USA

Tel: 301-762-0888; Fax: 301-762-3888

Email: support@genecopoeia.com

Web: www.genecopoeia.com