

Product Information

TargeTron® Vector pNL9164

Catalog Number **T6701**

Storage Temperature $-20\text{ }^{\circ}\text{C}$

TECHNICAL BULLETIN

Product Description

TargeTron Vector pNL9164 is a 9,295 bp expression vector derived from pNL9162¹. pNL9164 is to be used in conjunction with the TargeTron Gene Knockout System, Catalog Number TA0100. This circularized vector can be used for targeted gene knockouts in gram-positive organisms such as *Staphylococcus aureus*¹. Expression of the group II intron RNA and protein is under the control of the cadmium inducible promoter, $P_{\text{cad-cadC}}$ ^{1,2}. To facilitate curing of the plasmid, pNL9164 has a temperature sensitive origin of replication, pT181 cop-634ts repC4^{1,2}. As supplied, the pNL9164 vector is re-targeted to the *S. aureus hsa* gene. The *hsa* gene encodes a member of the HU family of bacterial histone-like proteins³. For validation in a specific *S. aureus* strain, this vector can be used directly to knockout *hsa* without any further modifications using the enclosed protocol. In order to re-target this vector to knockout other genes, the *hsa* specific IBS-EBS fragment (350 bp) between the *Hind* III and *BsrG* I sites can be cut out and replaced with another gene specific fragment. For re-targeting protocols, substitute the restriction digest modification below when referring to the TargeTron Gene Knockout System User Guide, Catalog Number TA0100, at sigma-aldrich.com.

Digestion of 350 bp re-targeted PCR product with *Hind* III, *BsrG* I and *Dpn* I

1. Set up a restriction digestion as follows:

8 μl	Purified PCR product (~200 ng)
2 μl	10 \times Restriction Enzyme Buffer
1 μl	<i>Hind</i> III (20 U/ μl)
1 μl	<i>BsrG</i> I (10 U/ μl)
1 μl	<i>Dpn</i> I (20 U/ μl)
7 μl	Water (molecular biology reagent)
20 μl	Total volume

2. Incubate the reaction for:

40 minutes, $37\text{ }^{\circ}\text{C}$

20 minutes, $60\text{ }^{\circ}\text{C}$

10 minutes, $80\text{ }^{\circ}\text{C}$

Note: *BsrG*I is thermophilic and has increased activity at $60\text{ }^{\circ}\text{C}$.

Key Features of the TargeTron Vector pNL9164
pT181 cop-634ts repC4 <i>ori</i>
Erythromycin resistance
Ampicillin resistance
ColE1 <i>ori</i>
cadC inducible promoter
5' exon (IBS)
Intron RNA
EBS2
EBS1d
3' exon
LtrA ORF

Reagent

Supplied at a concentration of 100 ng/ μl in 10 mM Tris-HCl, pH 8.0, 1 mM EDTA.

Precautions/Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Storage

Store at $-20\text{ }^{\circ}\text{C}$.

Protocol for Validation of *hsa* Knockout in *S. aureus* RN4220

1. Electroporate pNL9164 into *S. aureus* RN4220.
Note: pNL9164 was propagated in *E. coli* and therefore might be sensitive to restriction-modification in certain *S. aureus* strains. RN4220 is a restriction-defective *S. aureus* strain that can be used to condition the pNL9164 vector before electroporation into other restriction-modification strains if necessary.
2. Grow cells at 32 °C in 1.0 ml BHI (brain heart infusion) medium for 1 hour.
3. Add 1.0 ml into 5.0 ml fresh BHI with erythromycin (10 µg/ml) and grow overnight at 32 °C.
4. **(Optional)** Make a 1:100 dilution of overnight culture into fresh BHI with erythromycin (10 µg/ml) and grow until early log phase (OD₅₉₅ = 0.5). Induce the culture with 10 µM CdCl₂ for 90 minutes.
Note: The *Pcad* promoter has low basal level transcription in the absence of cadmium. The *hsa-24s* Targetron is very efficient, routinely giving 100% knockouts even without cadmium induction in *S. aureus* RN4220¹. For other target genes, cadmium induction may be necessary to increase knockout efficiency. Therefore, the induction step is required.
5. Plate on BHI agar plates containing 10 µg/ml erythromycin.
6. Incubate plates at 32 °C.
7. Perform colony PCR to confirm TargeTron integration using the primers listed below.
8. The PCR results should yield a 0.4 kb amplicon for wild-type RN4220 *hsa* and a 1.3 kb amplicon for the mutant *hsa* disruptant containing the intron.

hsa forward primer

5'-CGGAATCAGGAGGTGAATGTCTAATGA

hsa reverse primer

5'-CGGAATCACTTAGAAGAACATTACAATTTAT

Hsa-24s target sequence

TGTCTAATGAACAAAACAGATTTAATCAAT - intron
GCAGTTGCAGAGCAA

In addition to strains RN4220 and RN8098¹, the *Hsa-24s* target sequence is present exactly as presented above in all sequenced strains of *S. aureus* (per NCBI microbial BLAST), namely:

Staphylococcus aureus RF122
Staphylococcus aureus COL
Staphylococcus aureus MRSA252
Staphylococcus aureus MSSA476

Staphylococcus aureus MW2
Staphylococcus aureus Mu50
Staphylococcus aureus N315
Staphylococcus aureus NCTC 8325
Staphylococcus aureus USA300

The *hsa-24s* sequence is also present in *Bacillus halodurans* C-125 and in *Bacillus clausii* KSM-K16 with three or fewer mutations.

Temperature sensitive plasmid curing

For intron insertions in the sense direction, it is possible for the intron to splice out of mRNA with the assistance of the LtrA protein encoded on pNL9164. Thus, for accurately observing phenotypic effects of a given disruption, a re-targeted pNL9164 plasmid should be cured from the host. Conversely, the intron cannot splice out when inserted in the antisense ORF orientation and curing is still recommended, but not absolutely required. In order to cure the pNL9164 donor plasmid after a knockout has been confirmed, grow the strain overnight at 43 °C in BHI medium (without antibiotic) and screen for erythromycin sensitive colonies^{1,2}. If the targeted gene is essential or partially essential for cell survival, curing of the donor plasmid could be more difficult or unattainable¹. For example, *hsa-24s* disruptants have been shown to be conditional and can grow at 32 °C but not at 43 °C¹.

Delivery of heterologous DNA

pNL9164 has an *Mlu* I restriction site located within group II intron RNA coding sequence. Digestion at the *Mlu* I site allows for insertion of other DNA such as promoters (for mitigating polar effects), reporter genes, *loxP* recombination sequences, other antibiotic RAM-type markers, etc. The *Mlu* I site has been used to successfully deliver *tetM* and *abiD* genes⁴, a trimethoprim –RAM⁵, a kanamycin-RAM (plasmid pACD4K-C TA0100 kit), a removable kanamycin-RAM flanked with *loxP* sites (plasmid T2826, pACD4K-C-*loxP*), and a *lacZα* gene⁶. The efficiency of the intron may be affected by insertions at the *Mlu* I site. A good starting point is to attempt to insert an intron containing heterologous DNA into an easily screenable or selectable gene. For instance, the *hsa-24s* targeting sequence in pNL9164 is very efficient and is a good control to test the effects of heterologous DNA inserted at the *Mlu* I site on intron mobility.

References:

1. Yao, J et al., *RNA*. (2006) (in press)
2. Charpentier, E., et al., *Appl Environ Microbiol.* **70**(10):6076-6085 (2004).

3. Viter S., et al., *Res. Microbiol.* **150**(4):287-290 (1999).
4. Frazier , C., et al., *Appl Environ Microbiol.* **69**(2):1121-1128 (2003).
5. Zhong, J., et al., *Nucleic Acids Res.* **31**(6):1656-1664 (2003).

6. Jones, J.P., et al., *Mol. Ther.* **11**(5):687-94 (2005).

TargeTron is a registered trademark of InGex LLC.

License Agreement

This product and its use are the subject of one or more of U.S. Patent Nos. 5,698,421, 5,804,418, 6,027,895, and 6,001,608. BEFORE OPENING OR USING THIS PRODUCT, PLEASE READ THE TERMS AND CONDITIONS SET FORTH BELOW. YOUR USE OF THIS PRODUCT SHALL CONSTITUTE ACKNOWLEDGMENT AND ACCEPTANCE OF THESE TERMS AND CONDITIONS. If you do not agree to use this product pursuant to the following terms and conditions, please contact Sigma Technical Services to return the unused and unopened product for a full refund.

The purchase of this product conveys to you, the buyer, the non-transferable right to use the purchased product in non-commercialized research conducted by you. Sigma does not have the right to grant you a license to use this product for any other purposes. If you wish to use this product for any non-research purposes, you must obtain a separate commercial license from InGex LLC. Commercial entities may use this product for research and evaluation purposes for one year from the date of purchase. IF YOU ARE A COMMERCIAL ENTITY, YOUR RIGHT TO USE THIS PRODUCT EXPIRES ONE YEAR FROM THE DATE OF PURCHASE. Any commercial entity that wishes to use this product beyond this one-year period, must obtain a commercial license from InGex, LLC.

You may not transfer the product, its components or any materials or information made through the use of this product, including, but not limited to cell lines, to any third party without prior written approval of Sigma and without the transferee entering into a use license with Sigma. Notwithstanding the foregoing, the product, its components or any materials or information made through use of the product may be transferred by you to your legal affiliates or bona fide third party contractors performing paid work on your behalf provided the use by such third party contractors is limited to performance of work for you and such third party contractors agree to abide by the terms of this license in writing. You may also (a) make such a transfer to a scientific collaborator for research purposes if such collaborator agrees to abide by the usage terms and conditions in writing,

(b) disclose in a peer reviewed scientific publication information made through the use of the product and (c) deposit materials made through use of the product as required in association with a peer reviewed scientific publication, provided the recipient is under an obligation not to use or distribute the materials. Your rights and the rights of any third party who obtains the product, its components or any materials or information made through use of the product from you, to use the product will terminate immediately if you or such third party fail to comply with these terms and conditions. You and any third party who obtains the product, its components or any materials or information made through use of the product from you shall, upon such termination of your rights, destroy all product and components thereof in your control, and notify Sigma of such in writing.

For information on purchasing a license to this product for purposes other than non-commercial research, contact Licensing Department, InGex, LLC, 3655 Vista Ave, St. Louis, MO 63110, 314-865-0113.

AH,PHC 10/10-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.