

Product Information

pACD4 TargetTron® Vector Set

No selectable marker

Catalog Number **TV0010**

Storage Temperature -20°C

TECHNICAL BULLETIN

Product Description

The pACD4 vector set (no selectable marker) is a set of four linearized vectors intended for use with the TargetTron Gene Knockout System when the use of a selectable marker is not desired. Each vector in the pACD4 vector set is similar to the pACD4K-C TargetTron vector, but without a kanamycin RAM marker to select for chromosomal insertion. When these non-RAM TargetTron vectors express the ribonucleoprotein complex (RNP) and intron RNA is inserted into the specific sequence of interest, there is no selectable marker present in the group II intron insertion sequence. The pACD4 vector set also differs from the pACD4K-C TargetTron vector in that each vector in this set contains either an A, C, G, or T base in the $\delta+1$ position.¹ This base variation facilitates proper base pairing of group II intron precursor RNA for different target site designs and results in more efficient precursor RNA splicing to form active TargetTron RNPs.¹

Reagents

Set contains 2 μg of each vector supplied at 20 ng/ μl in 10 mM Tris-HCl, pH 8.0, with 1 mM EDTA. The vectors have been linearized with *Hind* III and *Bsr*GI.

Precautions and Disclaimer

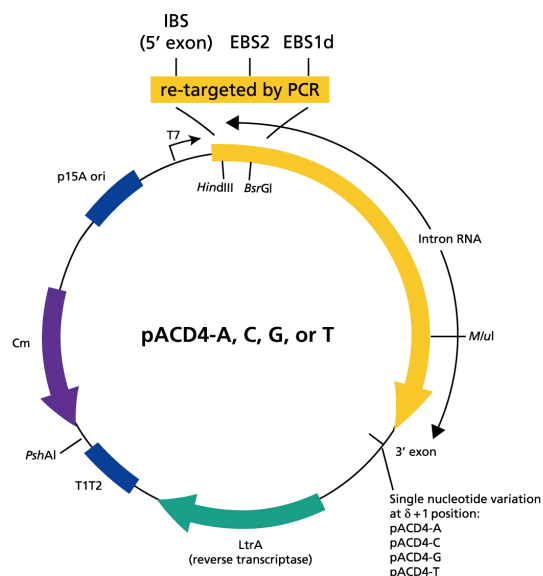
This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Storage

Store at -20°C . Shipped on wet ice.

Vector Features

Location (bp)	Feature
585-1497	p15A ori
1741-1762	T7 promoter
1802-1826	5' exon (IBS)
1827-2741	intron RNA
2049-2053	EBS2 (exon binding sequence 2)
2102-2111	EBS1d
2742	$\delta+1$ position
2742-2751	3' exon
2988-4787	LtrA ORF
4920-5190	T1/T2 transcriptional terminator
5861-219	chloramphenicol (Cm) resistance



Procedure

A. Vector selection

In general, the pACD4-C vector can be used. In some cases, upon performing colony PCR to locate a mutant, the efficiency of insertion may be just below what is reasonable to screen (for example, >100 colonies must be screened). If a mutant cannot be found, cloning into an alternative vector (pACD4-A, T, or G) may increase the insertion efficiency so that <100 colonies could be screened by colony PCR.

To choose the appropriate vector, match the vector name with the nucleotide at the $\delta+1$ position in the target site. For example, if the algorithm suggests the target site shown below, vector pACD4-A would be the appropriate vector to use.

CCATCTTGCGATAGACCGGATCTTTAGCGC –
intron – **A**GTTTTTACGTCCAC

↑
 $\delta + 1$ position

The $\delta+1$ position is the nucleotide immediately after the intron insertion in the 3' exon.

B. Additional plasmid information

Vectors are provided pre-linearized with *Hind* III and *Bsr*G I. Chloramphenicol (25 μ g/ml) is used to propagate each pACD4 vector. These vectors require the use of T7 RNA polymerase for intron expression. When using DE3 strains or a T7 RNA polymerase under the control of the *lac* UV5 promoter, glucose is typically included to provide additional suppression of the *lac* UV5 promoter prior to IPTG-induction. T7 RNA polymerase may also be supplied by co-transforming plasmid pAR1219 that expresses T7 RNA polymerase, Catalog Number T2076.

C. Sub-cloning into host specific shuttle vectors

Prior to sub-cloning, the pACD4 vectors need to be circularized. This may be accomplished using the *lacZ* control reaction in the TargeTron Kit (Product No. TA0100). For adaptation to another host, it is advised to sub-clone intron components first from pACD4-C and target an easily screenable gene. Sub-cloning the region bound by *Hind* III and *Psh*A I into an alternative host shuttle vector downstream of a host specific promoter should result in a functional TargeTron expression vector. The *Hind* III-*Psh*A I region should contain the intron RNA and LtrA reverse transcriptase coding regions, as well as the transcriptional terminator. In the final TargeTron shuttle vector, the *Hind* III and *Bsr*G I sites should be unique since these are used to routinely re-target the intron to knockout specific genes.

D. Delivery of heterologous DNA

Since these vectors do not contain the kan-RAM marker, heterologous DNA may be cloned into the *Mlu* I site in the intron for delivery to specific DNA sites. The *Mlu* I site has been used to successfully deliver a trimethoprim-RAM, a kanamycin-RAM (plasmid pACD4K-C in the TA0100 kit), and a *lacZ* α gene.³ 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, such as *lacZ*.

For detailed TargeTron gene disruption protocols, see the latest User Guide for the TargeTron Gene Knockout System, Catalog Number TA0100, at www.sigma-aldrich.com

References

1. Perutka, J., *et al.*, Use of computer-designed group II introns to disrupt *Escherichia coli* DExH/D-box protein and DNA helicase genes. *J. Mol. Biol.*, **336** (2), 421-39 (2004)
2. Zhong, J., *et al.*, Targeted and random bacterial gene disruption using a group II intron (targetron) vector containing a retrotransposition-activated selectable marker. *Nucleic Acids Res.*, **31**(6), 1656-64 (2003).
3. Jones, J.P., *et al.*, Retargeting mobile group II introns to repair mutant genes. *Mol. Ther.*, **11**(5), 687-94 (2005).

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.

Academic and Non-Profit Laboratory Assurance Letter

The T7 system is based on technology developed at Brookhaven National Laboratory under contract with the U.S. Department of Energy and is the subject of U.S. Patent No. 5,693,489 (expiration date, December 2, 2014) assigned to Brookhaven Science Associates, LLC. (BSA). BSA will grant a nonexclusive license for the use of this technology, including the enclosed material, based upon the following assurances:

1. These materials are to be used for noncommercial research purposes only. A separate license is required for any commercial use, including use of these materials for research purposes or production purposes by any commercial entity. Information about commercial licenses may be obtained from the Office of Technology Commercialization and Partnerships, Brookhaven National Laboratory, Bldg. 490-C, P.O. Box 5000, Upton, New York 11973-5000, telephone (631)-344-7134.
2. No materials that contain the cloned copy of T7 gene 1, the gene for T7 RNA polymerase, may be distributed further to third parties outside of your laboratory, unless the recipient receives a copy of this license and agrees to be bound by its terms. This limitation applies to strains of BL21(DE3), BL21(DE3)pLysS, and BL21(DE3)pLysE, and any derivatives.
3. You may refuse this license by returning the enclosed materials unused. By keeping or using the enclosed materials, you agree to be bound to the terms of this license.

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.