



PRODUCT INFORMATION & MANUAL

Retroviral Expression System

NBP2-29499

For research use only. Not for diagnostic or therapeutic procedures.

TABLE OF CONTENTS

I. Components	3
II. Introduction.....	4
III. Overview	6
<i>Outline of RetroMax Procedure</i>	
IV. Protocol	8
<i>Transfection of 293 cells with retroviral vector</i>	
<i>Determining the Viral Titer</i>	
<i>Selection for stable virus-producing cell lines</i>	
V. Maximizing Retrovirus Titers	13
VI. The Safe Use of Retrovirus Vectors.....	14
VII. References	16
VIII. Appendix	17
A. Expression Vectors	
B. Packaging Vectors	
C. LacZ Retroviral Reporter Vector	
D. Nucleotide Sequence.....	

I. COMPONENTS

Store at -20 °C		
Cat. No.	Description	Quantity
NBP2-29500	pCLXSN Expression Vector (0.5 mg/ml)	20 µl
NBP2-29502	pCLNCX Expression Vector (0.5 mg/ml)	20 µl
NBP2-29538	pCLNRX Expression Vector (0.5 mg/ml)	20 µl
NBP2-29539	pCLNDX Expression Vector (0.5 mg/ml)	20 µl
NBP2-29540	pCL-Eco Packaging Vector (0.5 mg/ml)	20 µl
NBP2-29541	pCL-Ampho Packaging Vector (0.5 mg/ml)	20 µl
NBP2-29542	pCL-10A1 Packaging Vector (0.5 mg/ml)	20 µl
NBP2-29543	pCLMFG-LacZ Retrovirus Reporter Vector (0.5 mg/ml)	20 µl

Store at 4 °C		
Description		Quantity
Retromax	Buffer (2x)	10 ml
2.0 M CaCl ₂		2 ml

Additional Materials Required:

- 293 Cells may be ordered from ATCC
- G418 selection antibiotic
- Cell culture reagents: DMEM, Fetal bovine serum, penicillin-streptomycin
- Trypsin-EDTA
- Phosphate buffered saline
- Cloning cylinders (for picking up stably transfected colonies)
- Polybrene for retroviral infection.

II. INTRODUCTION

Retrovirus vectors are very efficient tools for stably introducing genes into dividing cells. RetroMax™ retrovirus vector system is based on the pCL vector system developed by Naviaux et al. (1). The vectors used in this system have been designed to maximize recombinant-retrovirus titers in a simple, efficient, and flexible experimental system. All members of the RetroMax expression vector family (pCLXSN, pCLNCX, pCLNRX, and pCLNDX) have an extended packaging signal (ψ +) and are derived from safety-modified retrovirus vectors in which the gag open reading frame has been stopped by a point mutation, thereby minimizing the opportunity for replication-competent retrovirus production by recombination with packaging genome. The 5'-enhancer of Moloney murine sarcoma virus long terminal repeat (LTR) which is inhibited by E1A has been deleted and fused at the TATA box of the human CMV immediate early region (1). This results in initiation of viral RNA at or near the +1 position in the R region of the naturally programmed retrovirus. This results in transient-retrovirus titers in the range of $2-5 \times 10^6$ CFU/ml when 293 cells are used.

All three members of the RetroMax packaging vectors (pCL-Eco, pCL-Ampho, and pCL-10A1) have also been safety modified by deleting the packaging signal and the 3' LTR enhancer. This makes the RNAs of the helper genome virtually un-packageable. The advantage of these pCL packaging plasmids is a high level expression of gag, pol, and env proteins with a balanced stoichiometry that is not achieved with either transiently or stably expressed split-genome packaging constructs. Inclusion of these three packaging plasmids in the RetroMax kit allows the choice of expressing ectotropic, amphotropic, or 10A1 envelopes which leads to greater experimental flexibility.

The RetroMax system is designed for maximal virus titer in 293 cells. It takes advantages of two properties of 293 cells, i) high level of transfectibility, ii) strong E1A-mediated stimulation of CMV promoter controlled transcription. 293 cells are of nonmurine origin, hence the problem of selective packaging and transfer of VL30 genomes (present in all murine packaging cells) is avoided. Vector supernatants are free of helper virus and are of sufficiently high titer within 2 days of transient transfection in 293 cells to permit infection of more than 50% of dividing target cells in culture.

By introducing a retroviral vector into a cell expressing retroviral proteins, retroviral articles (virions) are shed into the culture medium at the rate of about one infectious particle/cell/day. Retrovirus tropism is determined at three levels. The first is simply a function of viral envelope protein, gp70. The envelope determines which cells the virus will enter. gp70 comes in three different flavors for gene therapy. A fourth one 10A1 is still experimental.

1. Ectotropic (usually (MoMuLV) mouse and rat cells only (not human)
2. Amphotropic (from 4070A MuLV) most mammalian cells (but not hamster)
3. Gibbon Ape Leukemia virus (GALV) many mammalian cells (including hamster)
4. 10A1 (MuLV) most mammalian cells (including hamster)

The second level of tropism is nuclear translocation and integration. This is defined by structural features of p30CA (but requires the full 160S nucleoprotein pre-integration complex, comprised of all the gag proteins and viral RNA and/or DNA). Naked DNA in the cytoplasm after retrovirus un-coating and reverse transcription is never seen.

The product of the FV-1 locus in murine cells interacts with p30CA, and can reduce the efficiency of translocation and integration (and thus apparent titer) 20-100 fold. Fortunately, the common Moloney-based packaging cells supply a p30CA form (NB tropic) that avoids this problem. The human equivalent of FV-1 has not yet been identified.

The third level of retrovirus tropism is determined by the transcriptional activity of the LTR (and/or internal promoter) in the transfected cell. In general, the Moloney (and MSV) LTR is active in most mammalian cell types, with the distinct exception of embryonic stem cells and teratocarcinoma cell lines (like F9), in which it is silenced. It is also potentially inhibited by E1A/p300 in 293 cells.

Ping-pong amplification is sometimes used to increase retrovirus vector titers, by co-culturing vector-producing ectotropic and amphotropic cell lines. This can increase vector titers 10 fold, but often at a heavy cost:

1. Frequent truncation, deletions, and point mutations may occur in the inserted cDNA.
2. You may generate helper virus if you are not using a safety-modified system.

III. OVERVIEW

Retrovirus particles are fragile. They are easily inactivated by 0.1% detergent, chloroform, phenol, 1% bleach, 70% ethanol, at 65°C for 30 min, pH <6.5 or >9.0, UV light, and autoclaving.

Simple high-speed centrifugation (100,000g x 90 min) produces enough hydrodynamic shear to strip many virions of their gp70, and thus infectivity (although reverse transcriptase activity is preserved).

Virus can be stored in culture medium (with 10% serum) at -70°C indefinitely. One freeze-thaw cycle reduces the titer about 2-3 fold compared to the fresh virus. The second freeze-thaw drops the titer another 5-10 fold. **Aliquot your virus for storage at -70°C.**

When filter sterilizing retrovirus, be sure to use non-detergent treated 0.22 or 0.45 um filters. Any trace of detergent will strip virus envelope and reduce your titers. Filter before freezing, and not after in order to avoid losses due to aggregation.

All murine retrovirus vectors produced in either mouse cells (like the NIH 3T3-based packaging cell lines) or primate cells (COS and 293) are rapidly inactivated by human serum complement, with kill kinetics of 2-3 logs in 5 min at 37°C. Human C1q initiates the cascade by binding p15ETM at the virion surface. This is an antibody-independent process.

Murine retroviruses are heat labile. They have an infectious half-life of only 6 h in culture medium at 4°C.

OUTLINE OF RETROMAX PROCEDURE

- Day 0** Seed 293 cells and grow overnight.
- Day 1** Transfect with retroviral vector containing gene of interest and an appropriate packaging vector.
- Day 2** Replace medium.
- Day 3** Harvest virus-containing supernatant. Virus may be stored at -70°C at this stage. Infect target cells, either for titer determination or for gene expression.
- Day 4** Split infected target cells and grow for selecting stable virus-producing cell lines. For transient expression experiments, you may harvest the cells at this stage.
- Day 5** Start selection by replacing the medium with G418 containing medium.
- Day 9** Change medium and continue selection.
- Day 14** Count antibiotic resistant colonies and calculate titer.

Note: If you are using retroviral expression system for the first time, we strongly recommend using the LacZ control plasmid included in the kit. The β -galactosidase expression can be monitored using β -gal staining kit (Cat.# NBP2-29546) or any other standard protocol.

IV. PROTOCOL

One Day Before Transfection

1. Seed 1×10^6 293 cells in 6 cm tissue culture plates. This should yield a cell density of about 30% confluence on the day of the experiment.
2. Incubate overnight at 37°C in DMEM supplemented with 10% fetal bovine serum, 1% pen-strep.

TRANSFECTION OF 293 CELLS WITH RETROVIRAL VECTOR

Day 1

1. Add 0.25 ml of RetroMax transfection buffer (previously tested for optimum transfection efficiency) to the required number of sterile 15 ml polypropylene tubes. Lipofectamine may have certain advantages in reproducibility, but this has not been tested extensively by us for overall virus titers.
2. Dilute 2 M CaCl_2 to final concentration of 0.25 M in sterile distilled water (30 μl of 2 M CaCl_2 and 220 μl of sterile water). Add the following to 1.5 ml sterile Eppendorf tube:

0.25 M CaCl_2 - 30 μl
pCL-ECO, pCL-Ampho, or pCL-10A1 - 10 μg
Retroviral vector containing your gene - 10 μg
Mix by vortexing.
3. Add the DNA/ CaCl_2 mix drop-wise to the transfection buffer tubes (step 1) while lightly vortexing. Incubate at room temperature for 20 min.
4. During this incubation aspirate the transfection media from plates to be transfected and add 2 ml DMEM containing 10% FBS/1% P/S and place back in the incubator until step 3 is completed.
5. Add the DNA/ CaPO_4 mix drop-wise to 293 cells on 6 cm T.C. plates. Place in humidified CO_2 incubator for 3-4 h. (Longer times may result in cells coming off the plates).
6. Carefully aspirate medium. Add 2 ml of warm PBS-15% Glycerol-no serum (glycerol shock medium) for 2 min. This step is optional. In some cases, it may increase the transfection efficiency by 2-fold.

- Aspirate glycerol shock medium. Carefully add 4 ml DMEM containing 10% FBS, 1% P/S along the side of the dish. Incubate for 12 h.

Culture vessel	Surface area/well (cm ²)	Volume of plating medium (ml)	Transfection buffer (ml)	Packaging vector	Retro-vector	2 M CaCl ₂ (μl) and dil. vol (μl)
6 well	10	2	0.125	5 μg	5 μg	15 μl + 110 μl
6 cm	20	4	0.25	10 μg	10 μg	30 μl + 220 μl
10 cm	60	10	0.5	20 μg	20 μg	60 μl + 440 μl

Day 2

- Aspirate medium and add fresh medium in the morning and incubate for 24 to 72 h.

Day 3 (24 h after addition of fresh medium).

- Filter sterilize (0.45 mm syringe filters are convenient) the virus-containing supernatant to remove any cells in suspension. The virus can now be used directly, or stored at -70°C until needed.
- Infect the desired target cells with 1 ml to 4 ml of 293 supernatant in 8 μg/ml Polybrene. The amount of supernatant you use depends on whether you are titrating virus or want to infect the maximum number of target cells possible. Do not forget the polybrene. Omission of polybrene will drop your apparent titers 100-1000 fold.

DETERMINING THE VIRAL TITER

Remember that for titrating, you must dilute the transfected supernatant at least 50 fold to stay in the linear part of the dilution curve. If you just want the maximum number of cells infected, then as little as a 2-fold dilution (equal volume mix) with the medium of the intended target cells is usually enough to prevent significant cell cycle inhibition.

- For titrating, prepare serial dilutions (four 10-fold dilutions) of vector supernatant in order to be sure that you are in the linear part of the titration curve (ie, out of the Poisson region). Infections for accurate titrating must be done at effective MOI is **0.1**. Target cells must be growing exponentially and only 30-50% confluent for maximum infection efficiencies.

2. Total virus-cell contact time should be a minimum of 12-24 h. This is because cycling cells are continuously entering and exiting the window of infectability. Even though the infective half life of the murine retrovirus particle is just 6-8 h at 37°C the rate of new cells entering the window is greater for the first 24 h, so longer contact times means more infected target cells.
3. Always test your titers on a standard control cell line (we use NIH 3T3) in parallel with infections of other desired target cells. Intrinsic infectability of many target cells can vary widely from 0.01-100% of the titers on NIH 3T3 cells.
 - Virus titers on NIH 3T3 cells for empty RetroMax vectors are typically $2-3 \times 10^6$ CFU/ml for ecotropic virus and 1×10^6 for amphotropic virus, assuming a typical 293 transfection efficiency of 30-50%.
 - When titrating virus on NIH 3T3 cells, infect 2×10^5 cells on a 6 cm plate (in 4 ml medium), overnight (16 h) with 1, 3, and 10 ml of pCL vector supernatant. You will need larger volumes for lower titer vectors, or cells that are more refractory to infection than NIH 3T3.
 - If virus stock is limiting: the most efficient use can be made by using 0.5-1 ml volumes to serially infect target cells in 6 cm plates (or 2-3 ml in 10 cm plates), and adding fresh virus every 4-6 h for 3-4 infection cycles. Continuous exposure to virus for about 24 h is necessary in order to ensure that all cells have cycled through their receptivity window (S-G2) for retroviral infection. Be sure to add polybrene to 8 μ g/ml.
4. Check your transfection efficiency by drawing a 1 cm square on the bottom of the plate of transfected 293 cells. Scrape harvest all the cells outside of this square (if desired) for RNA or protein analysis (CAT assays, ONPG-LacZ, Westerns, Northern, Hirts, etc.). Fix and stain the transfected cells remaining inside the 1 cm square with X-Gal to determine the transfection efficiency (TXE). Typical transfection efficiencies are 30-50% in this subline of 293 cells. The same DNA and reagents will give TXEs of 2-15% on COS cells.

SELECTION FOR STABLE CELL LINES**Day 4 (12-24 h after infection)**

1. If using a vector that confers G418 resistance, split the infected target cells at various dilutions (1:20 to 1:200) into 10 cm T.C. plates. A 1:20 dilution is about 10^5 NIH 3T3 cells. If 0.1% of the cells were infected, you will get about 100 colonies after 8-12 days of selection.

When infecting primary cells:

Accurate titers cannot be obtained when infecting primary fibroblasts, bone marrow or tumor cells because these cell types display density-dependent growth and typically have low plating efficiencies of 0.01%. This means that if 1000-10,000 cells are plated, only 1-100 colonies will actually clone out, *even if they are all infected and G418-resistant*. Therefore when infecting these cells, do not split them more than they will tolerate and only if they are >80% confluent (this is usually only a 1:2 to 1:4 dilution).

If you are selecting primary cells in G418, you will need to trypsinize and concentrate the cells by replating on sequentially smaller dishes until sufficient G418-resistant cells have grown out that you can begin expanding the infected pool of cells. This process can take 2 weeks. Effective titers for a particular primary cell type and vector will be a constant percentage of the titer observed on NIH 3T3 cells.

If using vectors that do not confer antibiotic resistance (like LacZ or GFP), simply change the medium today. Primary bone marrow cells should always be infected by co-cultivation of autochthonous stromal cells and virus producer cells in the presence of IL-3 (or WEHI-conditioned medium) and GM-CSF (a potent stromal cell growth factor). **Never select them in G418.**

Day 5 (2 days after infection)

1. Begin selection of cells infected with virus vectors conferring antibiotic resistance by adding 100 ml of a 100x stock to a 10 cm dish containing 10 ml of medium.

The correct concentration of G418 (or any antibiotic) varies widely for different cell types. You must determine the concentration empirically. For NIH 3T3 cells this is 400-1000 mg/ml (active) G418. For other cell types, the right concentration is that which results no observable death at Day 2 and about 30-50% on Day 4. Complete G418 selection is usually achieved in 7-10 days.

Retroviral Expression System Manual

2. If using a virus vector that does not contain a selectable marker (e.g., pCL-LacZ, MFG-GM-CSF, GFP), or if you would like a rapid assessment of gene expression in the infected target cells (for vectors expressing CAT, Luciferase, GFP, or LacZ), this can be tested today: b-gal staining of fixed cells in situ, (you can calculate the LacZ titer of your virus from this); CAT, Luciferase, or ONPG assays are done from cell lysates.
3. Because of the natural kinetics of retroviral infection, integration, and expression, no selection pressure (antibiotics) or assessment of gene expression should be made until 48 h after infection, i.e., if cells are infected on Day 3, gene expression cannot be accurately tested until Day 5.

Day 9 (4 days after starting selection).

1. Add fresh medium (and antibiotic) to cells under selection.
2. If infected cells were primary fibroblasts or primary tumor cells, you may need to increase the cell density (that has fallen due the death of uninfected cells under selection) by one of two methods, in order to avoid cell death due to densities falling below that tolerated by your particular primary cell type: concentrate the infected cells by trypsinization and plating on a smaller dish, or add uninfected primary cells (of the same type) to bring the density up to 50%, and continue selection. You must let the added (non-G418 resistant) cells attach to the plates for 3-4 h before adding G418 again.
3. Most primary cells will not grow as isolated clones because of density-dependent growth requirements. Attempts to pick clones frequently result in the loss of all infected cells.

Day 14 (10-13 days after infection)

1. Count the antibiotic resistant colonies, and calculate the titer (e.g., Neo titer) in your virus supernatants.

Example: Let us say you count 125 G418-resistant colonies on a 10 cm plate. If you infected (5×10^5) NIH 3T3 cells with 1 ml of virus supernatant, then split out the infected cells 1:20, your calculated titer is $125 \times 1000 \times 20 = 2.5 \times 10^6$ CFU/ml.

Note: Many cDNAs of interest may have either cytostatic or cytotoxic effects on infected cells, so that stable colony formation under G418 selection does not actually reflect the true number of cells initially infected. Only growing cells make colonies.

V. MAXIMIZING RETROVIRUS TITERS

1. The principal determinant of retrovirus titer is the abundance of packageable RNA, and not the abundance of viral proteins. Viral proteins are typically made in 20-fold stoichiometric excess. In fact, too much gp85 env can actually lower your titers because of impaired glycoprotein processing and assembly.
2. The RetroMax (pCL) system generates the highest abundance of packageable viral RNA of any known transient system by exploiting the power of the CMV IE enhancer-promoter in E1A-expressing 293 cells. The natural enhancer of the unmodified MuLV LTR is inhibited by E1A-p300 in 293 cells, so attempts to use non-pCL retroviral vectors in 293 cells will yield 20-50 fold lower titers, even with the same transfection efficiencies.
3. If you are studying cDNAs that do not have cytostatic or cytotoxic phenotypes, it may be possible to generate higher titer virus using traditional retrovirus packaging cells. This process takes 2 months (instead of 2 days for pCL). The highest titers are always obtained from stably infected (not transfected), cloned (not pooled) packaging cell lines. This is because transfected sequences are often inactivated by methylation, and because pro-virus integration position effects can influence gene expression from the same retrovirus vector in different clones of infected cells can vary over a 100-fold range (i.e., integration into heterochromatic regions of the genome gives poor expression, while integration into euchromatic regions gives high expression).
4. In deciding whether to go through the process of selecting and characterizing clones of packaging cells or simply preparing virus by the rapid pCL system, one must consider the intended applications. If you need a rapid test for the stable expression properties of a battery of mutant cDNAs that you have prepared, the pCL system is often adequate, or in the case of cytostatic and cytotoxic cDNAs, it is often the only way to produce usable amounts of virus.

Sometimes producing the virus (with a toxic or static cDNA) in cells from a different species can overcome the titer problems that result from cell growth inhibition.

If on the other hand, you plan to use the virus produced as a reagent that you can go back to many times over the next few years, then you need to pick clones of stable packaging cells.

5. pCL vectors reproducibly produce titers of $0.5-5 \times 10^6$ CFU/ml with good transfections, independent of phenotype and size (less than 4 kb) of the cDNA.
6. Typical retrovirus titers from cloned packaging cells are 10^4-10^6 CFU/ml (sometimes you can get 107), depending profoundly on the size and toxic properties of the cDNA expressed in these mouse fibroblast cell lines. cDNAs that are 2-4 kb long lead to modest reductions in titer because of packaging constraints. **cDNAs larger than 4 kb are subject to frequent spontaneous deletions and truncation during retroviral reverse transcription, and show large reductions in virus titers, and frequent non-expressing clones.**

Scaling Up:

1. Transfect 10 cm plates of 293 cells with 30-40 mg of pCL vector containing your gene of interest in 1 ml of CaCl_2 -HBS.
2. Replace the medium on Day 2.
3. Harvest and replace the medium every 24 h on Days 3, 4, and 5. This should give you 30 ml of virus supernatant from each transfected plate. The titers in supernatants harvested on Days 3 and 4 are equivalent. We suspect that Day 5 will be almost the same.

VI. THE SAFE USE OF MURINE RETROVIRUS VECTORS AND SAFETY PRECAUTIONS

Replication competent retroviruses (RCR) are called helper virus, or simply "Helper".

They require 3 trans- (gag, pol, and env), and 7 major, cis-active control elements (U3, R, U5, PBS, SD, ψ , and SA) in order to replicate.

The most common retrovirus vectors are based on the Moloney Murine Leukemia Virus (MoMuLV), encoding only the 7 cis elements.

These vectors are defective and cannot replicate without picking up 7.1 kb of sequence by homologous recombination with a helper genome (while simultaneously deleting your cDNA). Modern vectors are now "safety modified" by including a stop mutation early in "gag" (or a frame-shift) that pre-

vents gag translation and limits the sequence window available for productive recombination with helper genomes.

Packaging cells supply the 9 processed proteins encoded by gag, pol, and env (p15MA, p12 p30CA, p10NC, p14PRO, p85RT, p40IN, gp70SU, and p15ETM) necessary for virion assembly.

Modern packaging cells are safety-modified by dividing the gag-pol genes, and the env gene on two separate plasmids. These two plasmids are serially transfected (not co-transfected) into NIH 3T3 cells. The resulting safety modifications yield the modern split genome packaging cells.

Current evidence suggests that in order to initiate a pathogenic infection in primates with amphotropic murine retroviral vectors, three requirements must be met:

1. The infected host must be immunocompromised.
2. The vector preparation must contain helper virus.
3. Direct body fluid contact, e.g., intravenous inoculation is required for transfer.

However, for safe use of the RetroMax system, the user is strongly advised to follow the following guidelines:

1. According to NIH guidelines all retroviral production and transduction work must be done in a Biosafety Level 2 (BL2) facility.
2. Work in laminar flow, HEPA filtered hoods that receive annual maintenance and recertification.
3. Use sterile technique (flaming is not necessary and not recommended because of convection disturbances to airflow patterns).
4. Aspirate all liquid waste into flasks containing 5-10% (v/v) of a microbiocidal agent.
5. Discard spent plasticware in biohazard bags and autoclave before discarding.
6. Dispose spent glassware in detergent containers for cleaning and autoclaving.
7. Clean all surfaces with 70% ethanol at the end of the work.
8. Switch on the UV light immediately after work in laminar flow

Note: Retroviruses are not spread by aerosols.

VI. REFERENCES

1. Naviaux RK, Costanzi E, Haas M, and Verma I. The pCL vector system: Rapid production of helper-free, high titer, recombinant retroviruses. *J. Virol* 70: 5701-5705 (1996).
2. Cornetta K, Blaese RM, Anderson WF. Amphotropic leukemia retrovirus is not an acute pathogen for primates. *Human Gene Therapy*, 1: 15-30 (1990).
3. Naviaux RK and Verma IM. Retroviral vectors for persistent expression in vivo. *Curr. Opin. Biotechnol.* 3: 540-547 (1992).
4. Miller AD, Miller DG, Garcia JV, and Lynch C. Use of retroviral vectors for gene transfer and expression. *Methods in Enzymology* 217: 581-599 (1993).
5. Vanin EF, Kaloss M, Brosius C., and Nienhauis AW. Characterization of replication-competent retroviruses from nonhuman primates with virus-induced T-cell lymphomas and observations regarding the mechanism of oncogenesis. *J Virol.* 68: 4241-4250 (1994).
6. Ott D, Friedrich R and Rein A. Sequence analysis of amphotropic and 10A1 murine leukemia viruses: close relationship to mink cell focus-inducing viruses. *J. Virol.* 64: 757-766 (1990).

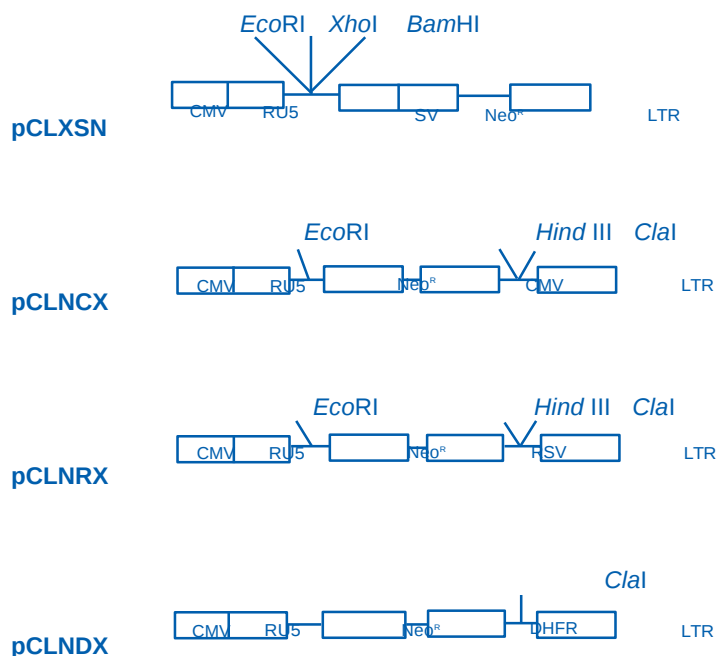
VII. APPENDIX

A. Expression Vectors

Choice of Vectors

The RetroMax expression vectors are designed to maximize recombinant-retrovirus titers in a simple, efficient, and flexible experimental system. All members of the RetroMax expression vector family (pCLXSN, pCLNCX, pCLNRX, and pCLNDX) have an extended packaging signal (ψ +) and are derived from safety-modified retrovirus vectors in which the gag open reading frame has been stopped by a point mutation (1), thereby minimizing the opportunity for replication competent retrovirus production by recombination with packaging genome. Four expression vectors are provided in the kit. Clone your gene of interest into one of these vectors:

Schematic Diagram of pCL Vectors



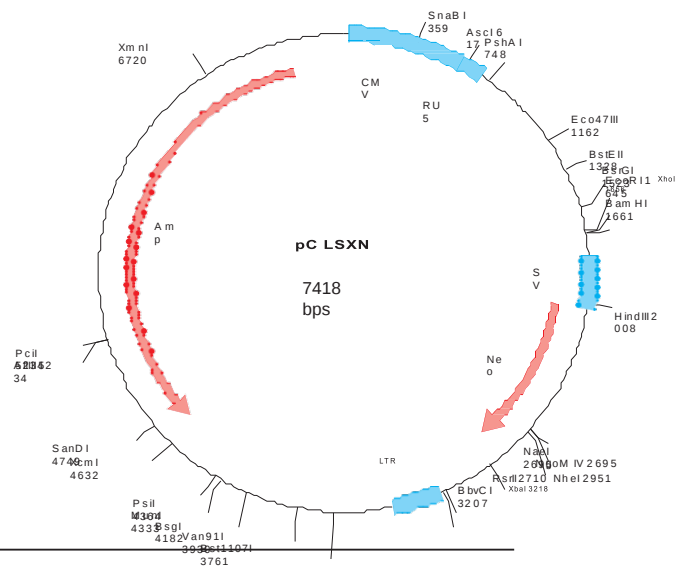
PCLXSN Retrovirus Expression Vector

Catalog No.: NBP2-29500

Quantity: 10 µg in 20 µl 1x TE (10 mM Tris, pH 7.5, 1 mM EDTA)

The inserted cDNA is under the control of LTR. The gene of interest can be cloned in *Eco* RI, *Xho* I and *Bam* HI cloning sites. In this case permanent cell lines can be selected. The complete vector sequence is available online at <http://www.novusbio.com/>.

Schematic presentation of pCLXSN vector



*CMV immediate early enhancer-promoter allows high-efficiency transcription in 293 cells. However, this is lost during viral replication. Hence, the gene of interest cloned into *Eco* RI, *Xho* I, or *Bam* HI cloning site is driven by Moloney MLV and murine sarcoma virus LTR (RU5).*

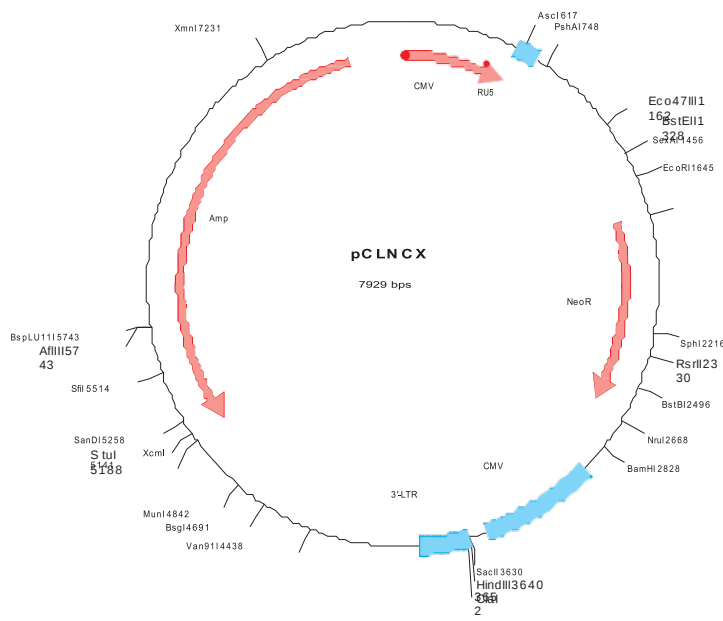
pCLNCX Retrovirus Expression Vector

Catalog No.: NBP2-29502

Quantity: 10 µg in 20 µl 1x TE (10 mM Tris, pH 7.5, 1 mM EDTA)

The gene of interest is cloned into *Hind* III and *Cla* I sites and is expressed under CMV promoter control. A second gene can be cloned into *Eco*RI site located upstream of the Neo^R gene. However, in this case permanent cell lines cannot be selected. Combined size of two inserts should not be more than 4 kb. The complete vector sequence is available online at <http://www.novusbio.com/>.

Schematic presentation of pCLNCX vector



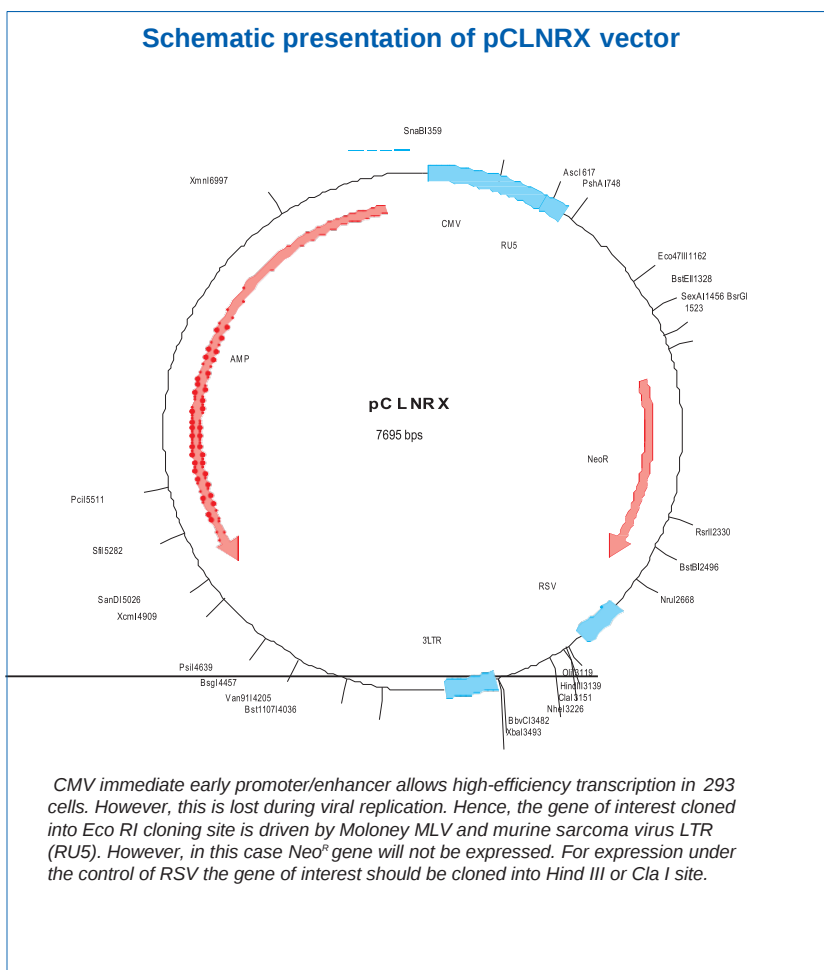
CMV immediate early promoter/enhancer allows high-efficiency transcription in 293 cells. However, this is lost during viral replication. Hence, the gene of interest cloned into *Eco*RI cloning site is driven by Moloney MLV and murine sarcoma virus LTR (RU5). However, in this case Neo^R gene will not be expressed. For expression under the control of CMV early promoter/enhancer, the gene of interest should be cloned into *Hind* III or *Cla* I site.

pCLNRX Retrovirus Expression Vector

Catalog No.: NBP2-29538

Quantity: 10 µg in 20 µl 1x TE (10 mM Tris, pH 7.5, 1 mM EDTA)

Similar to the pCLNCX vector, it differs only in that the foreign gene is cloned downstream of RSV promoter into *HindIII* or *ClaI*. A second gene can be cloned into *EcoRI* site located upstream of the Neo^R gene. However, in this case permanent cell lines cannot be selected. Combined size of two inserts should not be more than 4 kb. The complete vector sequence is available online at <http://www.novusbio.com/>.

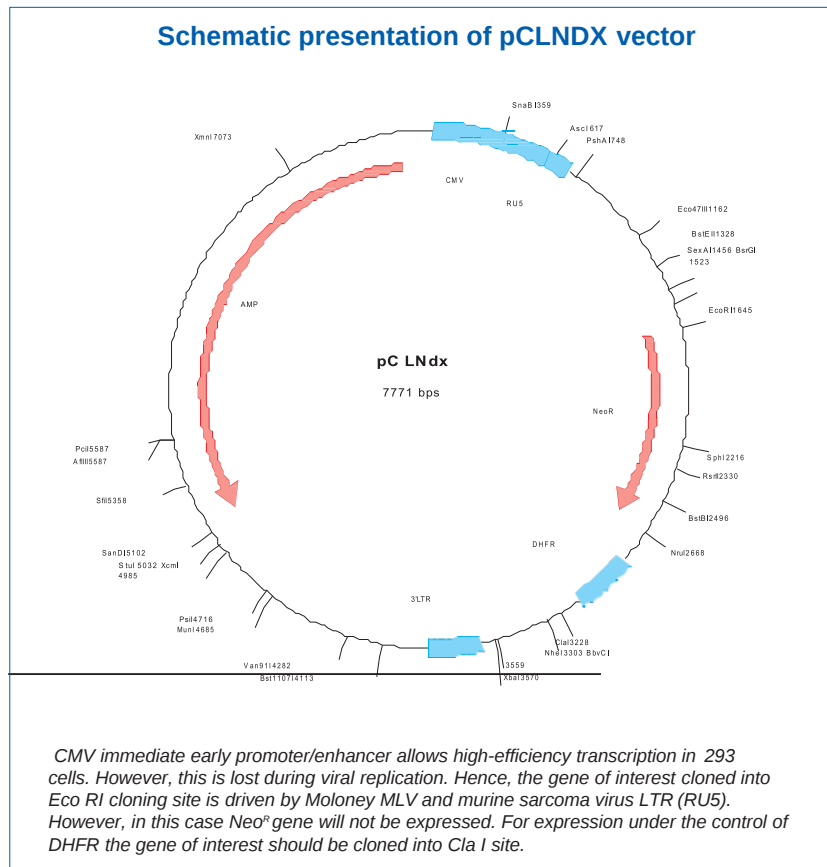


pCLNDX Retrovirus Expression Vector

Catalog No.: NBP2-29539

Quantity: 10 µg in 20 µl 1x TE (10 mM Tris, pH 7.5, 1 mM EDTA)

The gene of interest can be cloned into the single cloning site, *Cla*I, which puts it under the control of the DHFR promoter. However, in this case permanent cell lines cannot be selected. The complete vector sequence is available online at <http://www.novusbio.com/>.

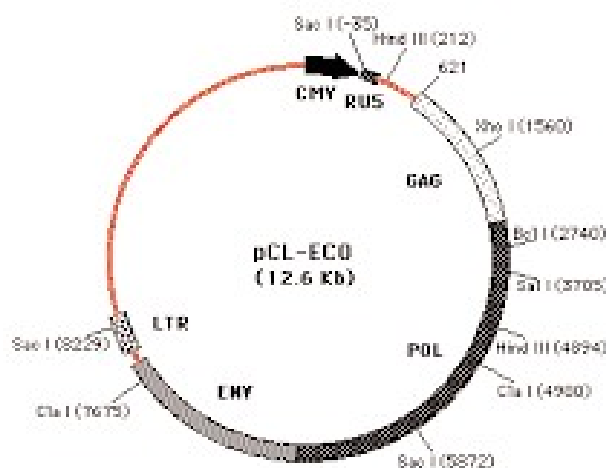


B. Packaging Vectors

The pCL packaging vectors is a part of the RetroMax Expression System (Cat# NBP2-29499) and has been designed to maximize recombinant retrovirus titers in a simple, efficient, and flexible experimental system. By introducing a retroviral vector into a cell expressing retroviral proteins, retroviral particles (virions) are shed into the culture medium at the rate of about 1 infectious particle/cell/day. Retrovirus tropism is determined at 3 levels. The first is simply a function of viral envelope protein, gp70. The envelope determines which cells the virus will enter. gp70 comes in three different flavors for gene therapy. Retroviruses obtained by co-transfection with pCL-Eco vector will infect mouse and rat cells, but not human cells.

Sequencing information is not available for any of our packaging vectors.

Schematic presentation of one of three packaging vectors: pCL-Eco



The gene coding for envelope protein was replaced with envelop gene from 4070A and 10A1 strain of MuLV to create pCL-Ampho and pCL-10A1 packaging vectors, respectively.

pCL-Eco Retrovirus Packaging Vector

Catalog No.: NBP2-29540

Quantity: 20 µg in 20 µl 1x TE (10 mM Tris, pH 7.5, 1 mM

Ectotropic (usually) (MoMuLV)	Mouse and rat cells only (not human)
-------------------------------	--------------------------------------

pCL-Ampho Retrovirus Packaging Vector

Catalog No.: NBP2-29541

Quantity: 20 µg in 20 µl (10 mM Tris, pH 7.5, 1 mM EDTA)

Amphotropic (from 4070A MuLV)	Most mammalian cells (but not hamster)
-------------------------------	--

pCL-10A1 Retrovirus Packaging Vector

Catalog No.: NBP2-29542

Quantity: 20 µg in 20 µl (10 mM Tris, pH 7.5, 1 mM EDTA)

10A1 (MuLV)	Most mammalian cells (including hamster)
-------------	--

Storage:

For long-term storage, store the Packaging at -20°C.

References:

1. Naviaux, RK, Costanzi, E, Haas, M and Verma, I. The pCL vector system: Rapid production of helper-free, high titer, recombinant retroviruses. J. Virol 70: 5701-5705 (1996).

Note: Packaging Vectors are for research use only. Not for use in humans. Use of this by commercial entities for any commercial purpose requires the user to obtain a commercial license.

C. LacZ Retroviral Reporter Vector

pCL-MFG-LacZ Retrovirus Reporter Vector

Catalog No.: NBP2-29543

Quantity: 20 µg in 20 µl (10 mM Tris, pH 7.5, 1 mM EDTA)

Sequencing information is not available for any of our packaging vectors.

Background:

The pCL-MFG-LacZ reporter vector is a part of the RetroMax expression system (Cat# NBP2-29499) and has been designed to assay beta-galactosidase activity. This plasmid has an ampicillin resistance gene and should be grown in LB-ampicillin media before use.

Novus also provides Beta-Galactosidase Staining Kit (Catalog no: NBP2-29546) as well as Beta-Galactosidase Quantitation Kit (Catalog no: NBP2-29547).

Storage:

For long-term storage, store at -20°C.

References:

1. Naviaux, RK, Costanzi, E, Haas, M and Verma, I. The pCL vector system: Rapid production of helper-free, high titer, recombinant retroviruses. J. Virol 70: 5701-5705 (1996).

Note: For research use only. Not for use in humans. Use of this by commercial entities for any commercial purpose requires the user to obtain a commercial license.

D. Nucleotide Sequence

Nucleotide Sequence pCLXSN

CGCGTTGACATTGATTATTGACTAGTTATTAATAGTAATCAATTACGGGGTCATTAGTTC
ATAGCCCATATATGGAGTTCGCGTTACATAACTTACGGTAAATGGCCCCGCTGGCTGAC
CGCCCAACGACCCCCGCCCATTGACGTCAATAATGACGTATGTTCCATAGTAACGCCAAT
AGGGACTTTCCATTGACGTCAATGGGTGGACTATTTACGGTAAACTGCCCACTTGGCAGT
ACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC
CGCCTGGCATTATGCCCAGTACATGACCTTATGGGACTTTCTACTTGGCAGTACATCTA
CGTATTAGTCATCGCTATTACCATGTGATGCGGTTTTTGGCAGTACATCAATGGGCGTGGA
TAGCGGTTTGACTACGCGGGATTTCCAAGTCTCCACCCCATTGACGTCAATGGGAGTTTG
TTTTGGCACCAAAATCAACGGGACTTTCCAAAATGTCGTAACAACCTCCGCCCCATTGACG
CAAATGGGCGGTAGGCGTGTACGGTGGGAGGTCTATATAAGCAGAGCTCAATAAAAGAGC
CCACAACCCCTCACTCGGCGCGCCAGTCTTCCGATAGACTGCGTCGCCCCGGGTACCCGTA
TTCCCAATAAAAGCCTCTTGCTGTTTGCATCCGAATCGTGGTCTCGCTGTTCTTGGGAGG
GTCTCCTCTGAGTGATTGACTACCCACGACGGGGGTCTTTCATTTGGGGGCTCGTCCGGG
ATTTGGAGACCCCTGCCCAGGACCAACCGACCCACCGGGAGGTAAGCTGGCCAGCAA
CTTATCTGTGTCTGTCCGATTGTCTAGTGTCTATGTTTGATGTTATGCGCCTGCGTCTGT
ACTAGTTAGCTAACTAGCTCTGTATCTGGCGGACCCGTGGTGGAAGTACGAGTTCTGAA
CACCCGGCCGCAACCCTGGGAGACGTCCCAGGGACTTTGGGGGCCGTTTTTGTGGCCCGA
CCTGAGGAAGGGAGTCGATGTGGAATCCGACCCCGTCAGGATATGTGGTTCTGGTAGGAG
ACGAGAACCTAAAACAGTTCCCGCCTCCGTCTGAATTTTTGCTTTCGGTTTGAACCGAA
GCCGCGCGTCTTGTCTGCTGCAGCGCTGCAGCATCGTTCTGTGTTGTCTCTGTCTGACTG
TGTTTCTGTATTTGTCTGAAAATTAGGGCCAGACTGTTACCACTCCCTTAAGTTTGACCT
TAGGTCACTGGAAAGATGTGAGCGGATCGCTCACAACCAAGTCGGTAGATGTCAAGAAGA
GACGTTGGGTTACCTTCTGCTCTGCAGAATGGCCAACCTTTAACGTCGGATGGCCGCGAG
ACGGCACCTTTAACCGAGACCTCATCACCCAGGTAAAGATCAAGGTCTTTTACCTGGCC
CGCATGGACACCCAGACCAGGTCCCCTACATCGTGACCTGGGAAGCCTTGGCTTTTGACC
CCCCTCCCTGGGTCAAGCCCTTTGTACACCCTAAGCCTCCGCCTCCTCTTCTCCATCCG
CCCCGTCTCTCCCCCTTGAACCTCCTCGTTCGACCCCGCCTCGATCCTCCCTTTATCCAG
CCCTCACTCCTTCTCTAGGCGCCGGAATTCGTTAACTCGAGGATCCGGCTGTGGAATGTG
TGTCAGTTAGGGTGTGGAAGTCCCCAGGCTCCCCAGCAGGCAGAAGTATGCAAAGCATG
CATCTCAATTAGTCAGCAACCAGGTGTGGAAGTCCCCAGGCTCCCCAGCAGGCAGAAGT
ATGCAAAGCATGCATCTCAATTAGTCAGCAACCATAGTCCCGCCCCTAACTCCGCCCATC
CCGCCCTAACTCCGCCAGTTCGGCCCATTTCTCCGCCCATGGCTGACTAATTTTTTTT
ATTTATGCAGAGGCCGAGGCCGCTCGCCCTCTGAGCTATTCCAGAAGTAGTGAGGAGGC
TTTTTTGGAGGCTTAGGCTTTTGCAAAAAGCTTGGGCTGCAGGTCGAGGCGGATCTGAT
AAGAGACAGGATGAGGATCGTTTCGCATGATTGAACAAGATGGATTGCACGCAGGTTCTC
CGGCCGCTTGGGTGGAGAGGCTATTCCGGCTATGACTGGGCACAACAGACAATCGGCTGCT
CTGATGCCGCCGTGTTCCGGCTGTCAGCGCAGGGGCGCCCGGTTCTTTTTGTCAAGACCG
ACCTGTCCGGTGCCCTGAATGAACTGCAGGACGAGGCAGCGCGGCTATCGTGGCTGGCCA
CGACGGGCGTTCTTGGCAGCTGTGCTCGACGTTGTCACTGAAGCGGGAAGGGACTGGC
TGCTATTGGGCGAAGTGCCGGGGCAGGATCTCCTGTCATCTCACCTTGCTCCTGCCGAGA
AAGTATCCATCATGGCTGATGCAATGCGGCGGCTGCATACGTTGATCCGGCTACCTGCC
CATTCGACCACCAAGCGAAACATCGCATCGAGCGAGCACGTACTCGGATGGAAGCCGGTC
TTGTCGATCAGGATGATCTGGACGAAGAGCATCAGGGGCTCGCGCCAGCCGAAGTGTTCG
CCAGGCTCAAGGCGCGCATGCCCGACGGCGAGGATCTCGTCGTGACCCATGGCGATGCCT
GCTTGCCGAATATCATGGTGGAATGGCCGCTTTTCTGGATTCATCGACTGTGGCCGGC
TGGGTGTGGCGGACCGCTATCAGGACATAGCGTTGGCTACCCGTGATATTGCTGAAGAGC
TTGGCGGCGAATGGGCTGACCGCTTCTCGTGCTTTACGGTATCGCCGCTCCCGATTTCG
AGCGCATCGCCTTCTATCGCCTTCTTGACGAGTTCTTCTGAGCGGGACTCTGGGGTTTCA
TAAAATAAAAGATTTTATTTAGTCTCCAGAAAAAGGGGGGAATGAAAGACCCACCTGTA
GGTTTGGCAAGCTAGCTTAAGTAACGCCATTTTGAAGGCATGAAAAATACATAACTGA

GAATAGAGAAGTTCAGATCAAGGTCAGGAACAGATGGAACAGCTGAATATGGGCCAAACA
GGATATCTGTGGTAAGCAGTTCCTGCCCCGGCTCAGGGCCAAGAACAGATGGAACAGCTG
AATATGGGCCAAACAGGATATCTGTGGTAAGCAGTTCCTGCCCCGGCTCAGGGCCAAGAA
CAGATGGTCCCCAGATGCGGTCCAGCCCTCAGCAGTTTCTAGAGAACCATCAGATGTTTC
CAGGGTGCCCCAAGGACCTGAAATGACCCTGTGCCTTATTTGAACTAACCAATCAGTTCG
CTTCTCGCTTCTGTTTCGCGCGCTTCTGCTCCCCGAGCTCAATAAAAGAGCCACAACCCC
TCACTCGGGGCGCCAGTCTCCGATTGACTGAGTCGCCCCGGGTACCCGTGTATCCAATAA
ACCCTCTTGACAGTTGCATCCGACTTGTGGTCTCGCTGTTCTTGGGAGGGTCTCCTCTGA
GTGATTGACTACCCGTCAGCGGGGGTCTTTCATTTGGGGGCTCGTCCGGGATCGGGAGAC
CCCTGCCCAGGACCACCGACCCACCACCGGGAGGTAAGCTGGCTGCCTCGCGCGTTTCG
GTGATGACGGTGAAAACCTCTGACACATGCAGCTCCCGGAGACGGTCACAGCTTGTCTGT
AAGCGGATGCCGGGAGCAGACAAGCCCGTCAGGGCGCGTCAGCGGGTGTGGCGGGTGTG
GGGGCGCAGCCATGACCCAGTCACGTAGCGATAGCGGAGTGTATACCTAGCTAGGTAGCT
AGAGGATCTTTGTGAAGGAACCTTACTTCTGTGGTGTGACATAATTGGACAACTACCTA
CAGAGATTTAAAGCTCTAAGGTAAATATAAAATTTTTAAGTGTATAATGTGTTAACTAC
TGATTCTAATTGTTTGTGATTTTAGATTCCAACCTATGGAAGTGAATGGGAGCAGT
GGTCCAATGCCTTTAATGAGGAAAACCTGTTTTGCTCAGAAGAAATGCCTCTAGTGATGA
TGAGGCTACTGCTGACTCTCAACATTCTACTCCTCCAAAAAAGAAGAGAAAAGGTAGAAGA
CCCCAAGGACTTTCCCTTCAGAATTGCTAAGTTTTTTGAGTCATGCTGTGTTTAGTAATAG
AACTCTTGCTTGCTTTGCTATTTACACCACAAAGGAAAAAGCTGCACTGCTATACAAGAA
AATTATGGAAAAATATTTGATGTATAGTGCCTTGACTAGAGATCATAATCAGCCATACCA
CATTTGTAGAGGTTTTACTTGCTTTAAAAAACCTCCCACACCTCCCCCTGAACCTGAAAC
ATAAAATGAATGCAATTGTTGTTGTTAACTTGTTTATTGCAGCTTATAATGGTTACAAAT
AAAGCAATAGCATCACAAATTTACAAATAAAGCATTTTTTTTCACTGCATTCTAGTTGTG
GTTTGTCCAAACTCATCAATGTATCTTATCATGTCTGATCAGCTTCAGAAGATGGCGGAG
GGCCTCCAACACAGTAATTTTCCCTCCCGACTCTTAAAATAGAAAATGTCAAGTCAGTTAA
GCAGGAAGTGGACTAACTGACGCAGCTGGCCGTGCGACATCCTCTTTTAATTAGTTGCTA
GGCAACGCCCTCCAGAGGGCGTGTGGTTTTGCAAGAGGAAGCAAAAGCCTCTCCACCCAG
GCCTAGAATGTTTCCACCCAATCATTACTATGACAACAGCTGTTTTTTTTAGTATTAAGC
AGAGGCCGGGGACCCCTGGCCCGCTTACTCTGGAGAAAAAAAACATTGTAGAGGCTTCCA
GAGGCAACTTGTCAAAACAGGACTGCTTCTATTTCTGTACACTGTCTGGCCCTGTCA
AGGTCCAGCACCTCCATACCCCTTTAATAAGCAGTTTGGGAACGGGTGCGGGTCTTACT
CCGCCATCCGCCCCTAACTCCGCCCAGTTCGCCCATTTCTCCGCCCCATGCTGACTAAT
TTTTTTTATTTATGCAGAGGCCGAGGCCGCTCGGCCTCTGAGCTATTCCAGAAGTAGTG
AGGAGGCTTTTTTGGAGGCTGCATTAATGAATCGGCCAACGCGCGGGGAGAGGCGGTTG
CGTATTGGGCGCTCTTCCGCTTCTCGCTCACTGACTCGCTGCGCTCGGTCTGTTCCGCTG
CGGCGAGCGGTATCAGCTCACTCAAAGGCGGTAAATACGGTTATCCACAGAATCAGGGGAT
AACGCAGGAAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCC
GCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAACTCGACG
TCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTCGGA
AGCTCCCTCGTGCGCTCTCCTGTTCGGACCTGCCGCTTACCGGATACCTGTCCGCTTT
CTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTG
TAGGTGCTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCCGTTCAGCCCGACCGCTGC
GCCTTATCCGGTAACATATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTG
GCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTC
TTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGGACAGTATTTGGTATCTGCGCTCTG
CTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACC
GCTGGTAGCGGTGGTTTTTTTTGTTTGCAAGCAGCAGATTACGCGCAGAAAAAAAGGATCT
CAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGGAACGAAAACCTCACGT
TAAGGGATTTTGGTCATGAGATTATCAAAAAGGATCTTCACCTAGATCCTTTTAAATTAA
AAATGAAGTTTTAAATCAATCTAAAGTATATATGAGTAAACTTGGTCTGACAGTTACCAA
TGCTTAATCAGTGAGGCACCTATCTCAGCGATCTGTCTATTTTCGTTTCATCCATAGTTGCC
TGACTCCCCGTCGTGTAGATAACTACGATACGGGAGGGCTTACCATCTGGCCCCAGTGCT
GCAATGATACCGCGAGACCCACGCTCACCGGCTCCAGATTTATCAGCAATAAACAGCCA
GCCGGAAGGGCCGAGCGCAGAAGTGGTCTGCAACTTTATCCGCTCCATCCAGTCTATT
AATTGTTGCCGGAAGCTAGAGTAAGTAGTTCGCCAGTTAATAGTTTGCGCAACGTTGTT

GCCATTGCTGCAGGCATCGTGGTGTCACGCTCGTCGTTTGGTATGGCTTCATTCAGCTCC
GGTTCCCAACGATCAAGGCGAGTTACATGATCCCCCATGTTGTGCAAAAAAGCGGTTAGC
TCCTTCGGTCCTCCGATCGTTGTCAGAAGTAAGTTGGCCGCAGTGTTATCACTCATGGTT
ATGGCAGCACTGCATAATTCTCTTACTGTCATGCCATCCGTAAGATGCTTTTCTGTGACT
GGTGAACCTCAACCAAGTCATTCTGAGAATAGTGTATGCGGCGACCGAGTTGCTCTTGCCC
GGCGTCAATACGGGATAATACCGCGCCACATAGCAGAACTTTAAAAGTGCTCATCATTGG
AAAACGTTCTTCGGGGCGAAAACTCTCAAGGATCTTACCGCTGTTGAGATCCAGTTCGAT
GTAACCCACTCGTGCACCCAACTGATCTTCAGCATCTTTTACTTTTACCAGCGTTTCTGG
GTGAGCAAAAACAGGAAGGCAAAATGCCGCAAAAAAGGGAATAAGGGCGACACGGAAATG
TTGAATACTCATACTCTTCCTTTTTCAATATTATTGAAGCATTTATCAGGGTTATTGTCT
CATGAGCGGATACATATTTGAATGTATTTAGAAAAATAAACAAATAGGGGTTCGCGCAC
ATTTCCCGAAAAAGTGCCACCTGACGTCTAAGAAACCATTATTATCATGACATTAACCTA
TAAAAATAGGCGTATCACGAGGCCCTTTTCGTCTCGCGCGTTTCGGTGATGACGGTGAAAA
CCTCTGACACATGCAGCTCCCGGAGACGGTCACAGCTTGTCTGTAAGCGGATGCCGGGAG
CAGACAAGCCCGTCAGGGCGCGTCAGCGGGTGTTGGCGGGTGTCGGGGCTGGCTTAATA
TGCGGCATCAGAGCAGATTGTACTGAGAGTGCACCATATGCGGTGTGAAATACCGCACAG
ATGCGTAAGGAGAAAAATACCGCATCAGGCGCCATTTCGCCATTTCAGGCTGCGCAACTGTTG
GGAAGGGCGATCGGTGCGGGCCTCTTCGCTATTACGCC

CMV - 21-584

RU5 – 584 -700

SV- 1770-2030

NeoR – 2030- 2880

3'LTR- 3230 – 3471

AMP- 7135-4666C

Nucleotide Sequence pCLNCX

```

1  cgcggttgaca ttgattattg actagttatt aatagtaatc aattacgggg tcattagttc
61  atagcccata tatggagttc cgcggttaca aacttacggt aaatggcccc cctggctgac
121 cgcccaacga ccccgccca ttgacgtcaa taatgacgta tgttccatag taacgccaat
181 agggactttc cattgacgtc aatgggtgga ctatttacgg taaactgccc acttggcagt
241 acatcaagtg tatcatatgc caagtacgcc ccctattgac gtcaatgacg gtaaattggc
301 cgcttgcat tatgcccagt acatgacctt atgggacttt cctacttggc agtacatcta
361 cgtattagtc atcgctatta ccatgtgatg cggttttggc agtacatcaa tgggcgtgga
421 tagcggtttg actcacgggg atttccaagt ctccaccca ttgacgtcaa tgggagtttg
481 ttttggcacc aaaatcaacg ggactttcca aaatgtcgta acaactccgc cccattgacg
541 caaatgggcg gtaggcgtgt acgggtgggag gtctatataa gcagagctca ataaaagagc
601 ccacaacccc tactcggcg cgccagtctt ccgatagact gcgtcgcccc ggtacccgta
661 ttcccaataa agcctcttgc tgtttgcata cgaatcgtgg tctcgtgtgt ccttgggagg
721 gtctcctctg agtgattgac taccacgacg gggggtcttt catttggggg ctcgtccggg
781 atttggagac ccctgcccag ggaccaccga cccaccaccg ggaggttaagc tggccagcaa
841 cttatctgtg tctgtccgat tgtctagtgt ctatgtttga tgttatgcgc ctgctctgtg
901 actagtttag taactagctc tgtatctggc ggaccctgtg tggaaactgac gatttctgaa
961 caccgggccg caaccctggg agacgtccca gggacttttg gggcggtttt tgtggcccga
1021 cctgaggaag ggagtcgatg tggaaatcca cccggtcagg atatgtggtt ctggtaggag
1081 acgagaacct aaacagttc ccgcctccgt ctgaattttt gctttcgggt tggaaaccgaa
1141 gccgcgcgtc ttgtctgtcg cagcgtgcca gcatcgttct gtgttgtctc tgtctgactg
1201 tgtttctgta tttgtctgaa aattaggggc agactgttac cactccctta agtttgacct
1261 taggtcactg gaaagatgtc gagcggatcg ctcaacaacc gtcggtagat gtcaagaaga
1321 gacgttgggt taccttctgc tctgcagaat ggccaacctt taacgtcgga tggccgcgag
1381 acggcacctt taaccgagac ctcatcacc aggttaagat caaggctttt tcacctggcc
1441 cgcattggaca ccagaccag gtcccctaca tcgtgacctg ggaagccttg gcttttgacc
1501 cccctccctg ggtcaagccc tttgtacacc ctaagcctcc gcctcctctt cctccatccg
1561 ccccgctctc ccccttgaa cctcctcgtt cgaccccgcc tcgacccctc ctttatccag
1621 ccctcactcc ttctctaggc gccggaattc cgatctgacg aagagacagg atgaggatcg
1681 tttcgcatga ttgaacaaga tggattgcac gcaggttctc cggccgcttg ggtggagagg
1741 ctattcgggt atgactgggc acaacagaca atcgggtgct ctgatgccgc cgtgttccgg
1801 ctgtcagcgc aggggcgccc ggttcttttt gtcaagaccg acctgtccgg tggcctgaat
1861 gaactgcagg acgaggcagc gcggctatcg tggctggcca cgacgggctg tccttgcgca
1921 gctgtgctcg acgttgtcac tgaagcggga agggactggc tgctattggg cgaagtgcg
1981 gggcaggatc tcctgtcatc tcaccttgct cctgccgaga aagtatccat catggctgat
2041 gcaatgcggc ggctgcatac gcttgatccg gctacctgcc cattcgacca ccaagcga
2101 catcgcatcg agcagcacg tactcggatg gaagccggtc ttgtcgatca ggaatgatctg
2161 gccgaagagc atcaggggct cgcgccagcc gaactgttcg ccaggctcaa ggcgcgcatg
2221 cccgacggcg aggatctcgt cgtgacccat ggcgatgcct gcttgccgaa tatcatgggtg
2281 gaaaatggcc gcttttctgg attcatcgac tgtggccggc tgggtgtggc ggaccgctat
2341 caggacatag cgttggctac ccgtgatatt gctgaagagc ttggcggcga atgggctgac
2401 cgcttccctg tgctttacgg tatcgccgct cccgattcgc agcgcacgc cttctatcgc
2461 cttcttgacg agttcttctg agcgggactc tggggttcga aatgaccgac caagcgacgc
2521 ccaacctgcc atcacgagat ttcgattcca ccgccgctt ctatgaaagg ttgggcttcg
2581 gaatcgtttt ccgggacgcc ggctggatga tcctccagcg cggggatctc atgctggagt
2641 tcttcgcca ccccgggctc gatccctcgt cgagttggtt cagctgctgc ctgaggctgg
2701 acgacctcgc ggagttctac cggcagtgca aatccgtcgg catccaggaa accagcagcg
2761 gctatccgcg catccatgcc cccgaactgc aggagtgagg aggcacgatg gccgctttgg
2821 tcgaggcgga tccggccatt agccatatta ttcatgggtt atatagcata aatcaatatt
2881 ggctattggc cattgcatac gttgtatcca tatcataata tgtacattta tattggctca
2941 tgtccaacat taccgcatg ttgacattga ttattgacta gttattaata gtaatcaatt
3001 acgggggtcat tagttcatag cccatatatg gagttccgcg ttacataact tacggtaaat
3061 ggcccgccgt gctgaccgcc caacgacccc cgccattga cgtcaataat gacgtatggt

```

```

3121 cccatagtaa cgccaatagg gactttccat tgacgtcaat ggggtggagta tttacggtaa
3181 actgcccact tggcagtaca tcaagtgtat catatgccaa gtacgcccc tattgacgtc
3241 aatgacggta aatggcccg ctagcattat gcccagtaca tgaccttatg ggactttcct
3301 acttggcagt acatctacgt attagtcatc gctattacca tggatgatgcg gttttggcag
3361 tacatcaatg ggcgtggata gcggtttgac tcacggggat ttccaagtct ccacccatt
3421 gacgtcaatg ggagtttggt ttggcaccaa aatcaacggg actttccaaa atgtcgtaac
3481 aactccgccc cattgacgca aatgggcggg aggcattgtac ggtgggaggt ctatataagc
3541 agagctcggt tagtgaaccg tcagatcgcc tggagacgcc atccacgctg ttttgacctc
3601 catagaagac accgggaccg atccagcctc cgcgcccca agcttggttaa catcgataaa
3661 ataaaagatt ttatttagtc tccagaaaaa ggggggaatg aaagacccca cctgtaggtt
3721 tggcaaaagt agagaacct cagatgtttc caggttgccc caaggacctg aaatgacctt
3781 gtgccttatt tgaactaacc aatcagttcg ctctcgctt ctgttcgcgc gcttctgctc
3841 cccgagctca ataaaagagc ccacaacccc tctactgggg cgccagtcct ccgattgact
3901 gagtcgcccg ggtacccgtg tatccaataa accctcttgc agttgcatcc gacttgtggt
3961 ctgctgtttc cttgggaggg tctcctctga gtgattgact acccgtcagc gggggtcttt
4021 cttttggggg ctgctccggg atcgggagac ccctgccag ggaccaccga cccaccaccg
4081 ggaggtaaag tggctgcctc gcgctttcg gtgatgacgg tgaaaacctc tgacacatgc
4141 agctcccga gacggtcaca gcttgtctgt aagcggatgc cgggagcaga caagcccgtc
4201 agggcgctgc agcgggtgtt ggcgggtgtc gggcgcgagc catgaccag tcacgtagcg
4261 atagcggagt gtatctagct aggtagctag aggatctttg tgaaggaacc ttacttctgt
4321 ggtgtgacat aattggacaa actacctaca gagatttaa gctctaaggt aaatataaaa
4381 tttttaagtg tataatgtgt taaactactg attctaattg tttgtgtatt ttagattcca
4441 acctatggaa ctgatgaatg ggagcagtg tggaatgcct ttaatgagga aaacctgttt
4501 tgctcagaag aaatgccatc tagtgatgat gaggctactg ctgactctca acattctact
4561 cctccaaaaa agaagagaaa ggtagaagac cccaaggact ttccttcaga attgctaagt
4621 tttttgagtc atgctgtgtt tagtaataga actcttgctt gctttgctat ttacaccaca
4681 aaggaaaaag ctgactgct atacaagaaa attatggaaa aatatttgat gtatagtgcc
4741 ttgactagag atcataatca gccataccac atttgtagag gttttacttg ctttaaaaaa
4801 cctcccacac ctccccctga acctgaaaca taaaatgaat gcaattgttg ttgttaactt
4861 gtttatttga gcttataatg gttacaaata aagcaatagc atcacaaata tcacaaataa
4921 agcatttttt tctactgcatt ctagtgttgg tttgtccaaa ctcataatg tatcttatca
4981 tgtctgatca gcttcagaag atggcggagg gcctccaaca cagtaatttt cctccgact
5041 cttaaaatag aaaatgtcaa gtcagttaag caggaagtgg actaactgac gcagctggcc
5101 gtgcgacatc ctcttttaat tagttgctag gcaacgccct ccagaggcg tgtggttttg
5161 caagaggaag caaaagcctc tccaccagg cctagaatgt ttccacccaa tcattactat
5221 gacaacagct gtttttttta gtattaagca gaggccggg acccctggcc cgcttactct
5281 ggagaaaaaa aacattgtag aggcctccag aggcaacttg tcaaaacagg actgcttcta
5341 tttctgtcac actgtctggc cctgtcacia ggtccagcac ctccataccc cctttaataa
5401 gcagtttggg aacgggtgcg ggtcttactc cgcccatccg cccctaactc cgccagttc
5461 cgccattct ccgccccatg ctgactaatt ttttttatt atgcagaggc cgaggccgcc
5521 tcggcctctg agctattcca gaagtagtga ggaggctttt ttggaggctg cattaatgaa
5581 tcggccaacg cgcggggaga ggcggtttgc gtattgggag ctcttcgctc tcctcgctca
5641 ctgactcgct gcgctcggtc gttcggctgc ggcgagcgg atcagctcac tcaaaggcgg
5701 taatacgggt atccacagaa tcaggggata acgcaggaaa gaacatgtga gcaaaaggcc
5761 agcaaaaggc caggaaaccgt aaaaaggccg cgttgctggc gtttttccat aggcctcgcc
5821 cccctgacga gcatcacaaa aatcgacgct caagtcagag gtggcgaaac ccgacaggac
5881 tataaagata ccaggcggtt cccctggaa gctccctcgt gcgctctcct gttccgacct
5941 tgccgcttac cggatacctg tccgccttcc tcccttcggg aagcgtggcg ctttctcata
6001 gctcacgctg taggtatctc agttcgggtg aggtcggtcg ctccaagctg ggctgtgtgc
6061 acgaaccccc cgttcagccc gaccgctgcg cttatccgg taactatcgt cttgagtcca
6121 acccggtaa gacacgactta tcgccactgg cagcagccac tggtaacagg attagcagag
6181 cgagggtatg aggcgggtgt acagagttct tgaagtgggt gcctaactac ggctacacta
6241 gaaggacagt atttgggtat tgcgctctgc tgaagccagt taccttcgga aaaagagttg
6301 gtagctcttg atccggcaaa caaaccaccg ctggtagcgg tggttttttt gtttgcaagc
6361 agcagattac gcgcagaaaa aaaggatctc aagaagatcc tttgatcttt tctacggggt
6421 ctgacgctca gtggaacgaa aactcacgtt aagggatttt ggtcatgaga ttatcaaaaa
6481 ggatcttcac ctagatcctt ttaaatataa aatgaagttt taaatcaatc taaagtatat

```

```

6541 atgagtaaac ttggtctgac agttaccaat gcttaatcag tgaggcacct atctcagcga
6601 tctgtctatt tcgttcatcc atagttgcct gactccccgt cgtgtagata actacgatac
6661 gggagggcctt accatctggc cccagtgtctg caatgatacc gcgagaccca cgctcaccgg
6721 ctccagatctt atcagcaata aaccagccag ccggaagggc cgagcgcaga agtggctctg
6781 caactttatc cgcctccatc cagtctatta attgttgccg ggaagctaga gtaagtagtt
6841 cgccagttaa tagtttgcg cagctgtgtg ccattgctgc aggcacgtg gtgtcacgct
6901 cgtcgttttg tatggcttca ttcagctccg gttcccaacg atcaaggcga gttacatgat
6961 ccccatgtt gtgcaaaaaa gcggttagct ccttcggtcc tccgatcgtt gtcagaagta
7021 agttggccgc agtgttatca ctcatgggta tggcagcact gcataattct cttactgtca
7081 tgccatccgt aagatgcttt tctgtgactg gtgagtactc aaccaagtca ttctgagaat
7141 agtgtatgcy gcgaccgagt tgctcttgcc cggcgtcaat acgggataat accgcgccac
7201 atagcagaac tttaaaagtg ctcatcattg gaaaacgttc ttcggggcga aaactctcaa
7261 ggatcttacc gctgttgaga tccagttcga tgtaaccac tcgtgcaccc aactgatctt
7321 cagcatcttt tactttcacc agcgtttctg ggtgagcaaa aacaggaagg caaaatgccg
7381 caaaaaaggg aataaggggc acacggaaat gttgaatact catactcttc ctttttcaat
7441 attattgaag catttatcag ggttattgtc tcatgagcgg atacatattt gaatgtattt
7501 agaaaaataa acaaataggg gttccgcgca catttccccg aaaagtgcc cctgacgtct
7561 aagaaaccat tattatcatg acattaacct ataaaaatag gcgtatcacg agggcccttc
7621 gtctcgcgcg tttcgggtgat gacggtgaaa acctctgaca catgcagctc ccggagacgg
7681 tcacagcttg tctgtaagcg gatgccggga gcagacaagc ccgtcagggc gcgtcagcgg
7741 gtgttgggcg gtgtcggggc tggcttaact atgcggcatc agagcagatt gtactgagag
7801 tgcaccatat gcggtgtgaa ataccgcaca gatgcgtaag gagaaaatac cgcacagggc
7861 gccattcgcc attcaggctg cgcaactgtt ggaagggcg atcgggtcgg gcctcttcgc
7921 tattacgcc

```

Molecule Features:

Name	Start	End
CMV prom	21	584
RU5 and packaging signal	585	700
NeoR	1645	2645
3'-LTR	3637	3878
AmpR	7624	5155

Nucleotide Sequence pCLNRX

CGCGTTGACATTGATTATTGACTAGTTATTAATAGTAATCAATTACGGGGTCATTAGTTC
 ATAGCCCATATATGGAGTTCCGCGTTACATAACTTACGGTAAATGGCCCGCTGGCTGAC
 CGCCCAACGACCCCGCCCATTTGACGTCAATAATGACGTATGTTCCATAGTAACGCCAAT
 AGGGACTTTCCATTGACGTCAATGGGTGGACTATTTACGGTAAACTGCCCACTTGGCAGT
 ACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC
 CGCCTGGCATTATGCCCAGTACATGACCTTATGGGACTTTCTACTTGGCAGTACATCTA
 CGTATTAGTCATCGCTATTACCATGTGATGCGGTTTTGGCAGTACATCAATGGGCGTGGA
 TAGCGGTTTGACTCACGGGGATTTCCAAGTCTCCACCCCATTTGACGTCAATGGGAGTTTG
 TTTTGGCACCAAAATCAACGGGACTTTCCAAAATGTCGTAACAACTCCGCCCCATTGACG
 CAAATGGGCGGTAGGCGTGTACGGTGGGAGGTCTATATAAGCAGAGCTCAATAAAAGAGC
 CCACAACCCCTCACTCGGCGCGCCAGTCTTCCGATAGACTGCGTCGCCCCGGGTACCCGTA
 TTCCCAATAAAGCCTCTTGCTGTTTGCATCCGAATCGTGGTCTCGCTGTTCTTGGGAGG
 GTCTCCTCTGAGTGATTGACTACCCACGACGGGGGTCTTTCATTTGGGGGCTCGTCCGGG
 ATTTGGAGACCCCTGCCAGGGACCAACCGACCCACCACGGGAGGTAAGCTGGCCAGCAA
 CTTATCTGTGTCTGTCCGATTGTCTAGTGTCTATGTTTGATGTTATGCGCCTGCGTCTGT
 ACTAGTTAGCTAACTAGCTCTGTATCTGGCGGACCCGTGGTGGAAGTACGAGTTCTGAA
 CACCCGGCCGCAACCCTGGGAGACGTCCCAGGGACTTTGGGGGCCGTTTTTGTGGCCGA
 CCTGAGGAAGGGAGTCGATGTGGAATCCGACCCCGTCAGGATATGTGGTTCTGGTAGGAG
 ACGAGAACCTAAAACAGTTCCCGCCTCCGTCTGAATTTTTGCTTTCGGTTTGAACCGAA
 GCCGCGCGTCTTGTCTGCTGCAGCGCTGCAGCATCGTTCTGTGTTGCTCTGTCTGACTG
 TGTTTCTGTATTTGTCTGAAAATTAGGGCCAGACTGTTACCACTCCCTTAAGTTTGACCT
 TAGGTCACCTGGAAAGATGTCGAGCGGATCGTCCACAACCAGTCGGTAGATGTCAAGAAGA
 GACGTTGGGTTACCTTCTGCTCTGCAGAATGGCCAACCTTTAACGTCGGATGGCCGCGAG
 ACGGCACCTTTAACCGAGACCTCATCACCCAGGTAAAGATCAAGGTCTTTTACCTGGCC
 CGCATGGACACCCAGACAGGTCCCTACATCGTGACCTGGGAAGCCTTGGCTTTTGACC
 CCCCTCCCTGGGTCAAGCCCTTTGTACACCTAAGCCTCCGCTCCTCTTCCCTCCATCCG
 CCCGTCTCTCCCCCTTGAACCTCCTCGTTGACCCCGCCTCGATCCTCCCTTTATCCAG
 CCCTCACTCCTTCTCTAGGCGCCGGAATTCGATCTGATCAAGAGACAGGATGAGGATCG
 TTTTCGATGATTGAACAAGATGGATTGCACGCAGGTTCTCCGGCCGCTTGGGTGGAGAGG
 CTATTCGGCTATGACTGGGCACAACAGACAATCGGCTGCTCTGATGCCGCCGTGTTCCGG
 CTGTCAGCGCAGGGGCGCCCGGTTCTTTTTGTCAAGACCGACCTGTCCGGTGCCCTGAAT
 GAACTGCAGGACGAGGCAGCGCGGCTATCGTGGCTGGCCACGACGGGCGTTCCTTGCGCA
 GCTGTGCTCGACGTTGTCACTGAAGCGGGAAGGGACTGGCTGCTATTGGGCGAAGTGCCG
 GGGCAGGATCTCCTGTCTATCTACCTTGCTCCTGCCGAGAAAGTATCCATCATGGCTGAT
 GCAATGCGGCGGCTGCATACGCTTGATCCGGCTACCTGCCCATTCGACCACCAAGCGAAA
 CATCGCATCGAGCGAGCACGTACTCGGATGGAAGCCGGTCTTGTGATCAGGATGATCTG
 GACGAAGAGCATCAGGGGCTCGCGCCAGCCGAAGTTCGCCAGGCTCAAGGCGCGCATG
 CCCGACGGCGAGGATCTCGTCGTGACCCATGGCGATGCCTGCTTGCCGAATATCATGGTG
 GAAAATGGCCGCTTTTCTGGATTATCGACTGTGGCCGGCTGGGTGTGGCGGACCGCTAT
 CAGGACATAGCGTTGGCTACCCGTGATATTGCTGAAGAGCTTGGCGGCGAATGGGCTGAC
 CGCTTCCTCGTGCTTTACGGTATCGCCGCTCCCGATTTCGACGCGCATCGCCTTCTATCGC
 CTTCTTGACGAGTTCTTCTGAGCGGGACTCTGGGGTTCGAAATGACCGACCAAGCGACGC
 CCAACCTGCCATCACGAGATTCGATTCCACCGCCGCTTCTATGAAAGGTTGGGCTTCG
 GAATCGTTTTCCGGGACGCCGGCTGGATGATCTCCAGCGCGGGGATCTCATGCTGGAGT
 TCTTCGCCCACCCCGGGCTCGATCCCTCGCGAGTTGGTTTCAGCTGCTGCCTGAGGCTGG
 ACGACCTCGCGGAGTTCTACCGGCAGTGCAAATCCGTGCGCATCCAGGAACAGCAGCG
 GCTATCCGCGCATCCATGCCCCGAACTGCAGGAGTGGGGAGGCACGATGGCCGCTTTGG
 TCGAGGCTCCCTCAGGATATAGTAGTTTCGCTTTTGCATAGGGAGGGGAAATGTAGTC
 TTATGCAATACACTTGTAGTCTTGCAACATGGTAACGATGAGTTAGCAACATGCCTTACA
 AGGAGAGAAAAAGCACCGTGCATGCCGATTGGTGGAAGTAAGGTGGTACGATCGTGCTT
 ATTAGGAAGGTAACAGACAGGTCTGACATGGATTGGACGAACCACTGAATTCGCTATGC
 AGAGATAATTGTATTTAAGTGCCTAGCTCGATACAATAAACGCCATTTGACCATTACCA

CATTGGTGTGCACCTCCAAAGCTTGTTAACATCGATAAAATAAAAGATTTTATTTAGTCT
 CCAGAAAAAGGGGGGAATGAAAGACCCACCTGTAGGTTTGGCAAGCTAGCTTAAGTAAC
 GCCATTTTGAAGGCATGGAAAAATACATAACTGAGAATAGAGAAGTTCAGATCAAGGTC
 AGGAACAGATGGAACAGCTGAATATGGGCCAAACAGGATATCTGTGGTAAGCAGTTCCTG
 CCCC GGCTCAGGGCCAAGAACAGATGGAACAGCTGAATATGGGCCAAACAGGATATCTGT
 GGTAAGCAGTTCCTGCCCCGGCTCAGGGCCAAGAACAGATGGTCCCCAGATGCGGTCCAG
 CCCTCAGCAGTTTCTAGAGAACCATCAGATGTTTCCAGGGTGCCCCAAGGACCTGAAATG
 ACCCTGTGCCTTATTTGAACTAACCAATCAGTTCGCTTCTCGCTTCTGTTGCGCGCTTC
 TGCTCCCCGAGCTCAATAAAAGAGCCACAACCCCTCACTCGGGGCGCCAGTCTCCGAT
 TGACTGAGTCGCCCCGGTACCCGTGTATCCAATAAACCCCTCTTGCAGTTGCATCCGACTT
 GTGGTCTCGCTGTTCTTGGGAGGGTCTCTCTGAGTGATTGACTACCCGTGAGCGGGG
 TCTTTTCAATTGGGGGCTCGTCCGGGATCGGGAGACCCCTGCCAGGGACCACCGACCCAC
 CACCGGGAGGTAAGCTGGCTGCCTCGCGCGTTTCGGTGATGACGGTGAAAACCTCTGACA
 CATGCAGTCCCCGAGACGGTCACAGCTTGTCTGTAAGCGGATGCCGGGAGCAGCAAGC
 CCGTCAGGGCGCTCAGCGGGTGTGGCGGGTGTGCGGGCGCAGCCATGACCCAGTCAG
 TAGCGATAGCGGAGTGATACCTAGCTAGGTAGCTAGAGGATCTTTGTGAAGGAACCTTA
 CTTCTGTGGTGTGACATAATTGGACAAACTACCTACAGAGATTTAAAGCTCTAAGGTA
 TATAAAATTTTAAAGTGATAATGTGTTAACTACTGATTCTAATTGTTGTGATTTTA
 GATTCCAACCTATGGAAGTATGAATGGGAGCAGTGGTCCAATGCCTTTAATGAGGAAAA
 CCTGTTTTGCTCAGAAGAAATGCCTCTAGTGATGATGAGGCTACTGCTGACTCTCAACAT
 TCTACTCCTCCAAAAAAGAAGAGAAAGGTAGAAGACCCCAAGGACTTTCCTTCAGAATTG
 CTAAGTTTTTTGAGTCATGCTGTGTTTAGTAATAGAACTCTTGCTTGCTTTGCTATTTAC
 ACCACAAAGGAAAAAGCTGCACTGCTATACAAGAAAATTATGAAAAATATTTGATGTAT
 AGTGCCTTGACTAGAGATCATAATCAGCCATACCACATTTGTAGAGGTTTTACTTGCTTT
 AAAAAACCTCCACACCTCCCCCTGAACCTGAAACATAAAATGAATGCAATTGTTGTTGT
 TAACTTGTTTTATTGCAGCTTATAATGGTTACAAATAAAGCAATAGCATCACAAATTTTAC
 AAATAAAGCATTTTTTTTCACTGCATTCTAGTTGTGGTTTGTCCAACTCATCAATGTATC
 TTATCATGTCTGTGATCAGCTTCAGAAGATGGCGGAGGGCCTCCAACACAGTAATTTCC
 TCCCGACTCTTAAAAATAGAAAATGTCAAGTCAGTTAAGCAGGAAGTGGACTAACTGACGC
 AGCTGGCCGTGCGACATCCTCTTTTAATTAGTTGCTAGGCAACGCCCTCCAGAGGGCGTG
 TGGTTTTGCAAGAGGAAGCAAAAGCCTCTCCACCCAGGCCTAGAATGTTTCCACCCAATC
 ATTACTATGACAACAGCTGTTTTTTTTAGTATTAAGCAGAGGCCGGGGACCCCTGGCCCG
 CTTACTCTGGAGAAAAAAAACATTGTAGAGGCTTCCAGAGGCAACTTGTCAAAACAGGAC
 TGCTTCTATTTCTGTACACTGTCTGGCCCTGTCACAAGGTCCAGCACCTCCATACCCCC
 TTTAATAAGCAGTTTGGGAACGGGTGCGGGTCTTACTCCGCCCATCCGCCCTAACTCCG
 CCCAGTTCCGCCCATCTCCGCCCATGCTGACTAATTTTTTTTATTTATGCAGAGGCCG
 AGGCCGCTCGGCCTCTGAGCTATTCCAGAAGTAGTGAGGAGGCTTTTTTGGAGGCTGCA
 TTAATGAATCGGCCAACGCGCGGGGAGAGGCGGTTTGCCTATTGGGCGCTCTTCCGCTTC
 CTCGCTCACTGACTCGCTGCGCTCGGTCGCTCGGTCGCGCGAGCGGTATCAGCTCACT
 AAAGGCGGTATACGGTTATCCACAGAACTCAGGGGATAACGCAGGAAAGACATGTGAGC
 AAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCGTTGCTGGCGTTTTTCCATAG
 GCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCC
 GACAGGACTATAAAGATACCAGGCGTTTCCCCCTGGAAGCTCCCTCGTGCGCTCTCCTGT
 TCCGACCCTGCCGCTTACCGGATACCTGTCCGCCTTTCTCCCTTCGGGAAGCGTGCGCT
 TTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGAGGTCGTTGCTCCAAGCTGGG
 CTGTGTGCACGAACCCCCCGTTACGCCCCACCGCTGCGCCTTATCCGTAACCTATCGTCT
 TGAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGAT
 TAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGG
 CTACACTAGAAGGACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGA
 AAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGT
 TTGCAAGCAGCAGATTACGCGCAGAAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTC
 TACGGGGTCTGACGCTCAGTGGAACGAAACTCACGTAAAGGGATTTTGGTCATGAGATT
 ATCAAAAAGGATCTTACCTAGATCCTTTTAAATTAATAAAGTAAAAATCAATCTA
 AAGTATATATGAGTAACTTGGTCTGACAGTTACCAATGCTTAATCAGTGAGGCACCTAT
 CTCAGCGATCTGTCTATTTTCGTTTCATCCATAGTTGCCTGACTCCCCGTGCTGTAGATAAC
 TACGATACGGGAGGGCTTACCATCTGGCCCCAGTGCTGCAATGATACCGCGAGACCCACG

CTCACCGGCTCCAGATTTATCAGCAATAAACCAGCCAGCCGGAAGGGCCGAGCGCAGAAG
TGGTCCTGCAACTTTATCCGCCTCCATCCAGTCTATTAATTGTTGCCGGGAAGCTAGAGT
AAGTAGTTCGCCAGTTAATAGTTTTCGCAACGTTGTTGCCATTGCTGCAGGCATCGTGGT
GTCACGCTCGTCGTTTGGTATGGCTTCATTCAGCTCCGGTCCCAACGATCAAGGCGAGT
TACATGATCCCCCATGTTGTGCAAAAAAGCGGTAGCTCCTTCGGTCTCCGATCGTTGT
CAGAAGTAAGTTGGCCGCAAGTGTATCACTCATGGTTATGGCAGCACTGCATAATTCTCT
TACTGTCATGCCATCCGTAAGATGCTTTTCTGTGACTGGTGAACCTCAACCAAGTCATTCT
GAGAATAGTGTATGCGGCGACCGAGTTGCTCTTGCCCGGCGTCAATACGGGATAATACCG
CGCCACATAGCAGAACTTTAAAAGTGCTCATCATTGGAAAACGTTCTTCGGGGCGAAAAC
TCTCAAGGATCTTACCGCTGTTGAGATCCAGTTCGATGTAACCCACTCGTGCACCCAACT
GATCTTCAGCATCTTTTACTTTTACCAGCGTTTCTGGGTGAGCAAAAACAGGAAGGCAAA
ATGCCGCAAAAAAGGGAATAAGGGCGACACGGAAATGTTGAATACTCATACTCTTCCTTT
TTCAATATTATTGAAGCATTTATCAGGGTATTGTCTCATGAGCGGATACATATTTGAAT
GTATTTAGAAAAATAACAAATAGGGGTTCGCGCACATTTCCCGAAAAGTGCCACCTG
ACGTCTAAGAAACCATATTATCATGACATTAACTATAAAAAATAGGCGTATCACGAGGC
CCTTTTCGTCTCGCGCTTTCGGTGATGACGGTGAAAACCTCTGACACATGCAGCTCCCGG
AGACGGTCACAGCTTGTCTGTAAGCGGATGCCGGGAGCAGACAAGCCCGTCAGGGCGCGT
CAGCGGGTGTGCGGGGTGTCGGGGCTGGCTTAACTATGCGGCATCAGAGCAGATTGTAC
TGAGAGTGACCATATGCGGTGTGAAATACCGCACAGATGCGTAAGGAGAAAAATACCGCA
TCAGGCGCCATTTCGCCATTCAGGCTGCGCAACTGTTGGGAAGGGCGATCGGTGCGGGCCT
CTTCGCTATTACGCC

CGCGTTGACATTGATTATTGACTAGTTATTAATAGTAATCAATTACGGGGTCATTAGTTC
ATAGCCCATATATGGAGTTCGCGTTACATAACTTACGGTAAATGGCCCGCCTGGCTGAC
CGCCCAACGACCCCCGCCCATTGACGTCAATAATGACGTATGTTCCATAGTAACGCCAAT
AGGGACTTTCCATTGACGTCAATGGGTGGACTATTTACGGTAAACTGCCCACTTGGCAGT
ACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC
CGCCTGGCATTATGCCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTA
CGTATTAGTCATCGCTATTACCATGTGATGCGGTTTTGGCAGTACATCAATGGGCGTGGA
TAGCGGTTTGACTCACGGGGATTTCCAAGTCTCCACCCCATTTGACGTCAATGGGAGTTTG
TTTTGGCACCAAAATCAACGGGACTTTCCAAAATGTCGTAACAACTCCGCCCCATTGACG
CAAATGGGCGGTAGGCGTGTACGGTGGGAGGTCTATATAAGCAGAGCTCAATAAAAGAGC
CCACAACCCCTCACTCGGCGCGCCAGTCTTCCGATAGACTGCGTCGCCCCGGGTACCCGTA
TTCCCAATAAAGCCTCTTGCTGTTTGCATCCGAATCGTGGTCTCGCTGTTCTTGGGAGG
GTCTCCTCTGAGTGATTGACTACCCACGACGGGGGTCTTTCAATTTGGGGGCTCGTCCGGG
ATTTGGAGACCCCTGCCCAGGGACCAACCGACCCACCGGGAGGTAAGCTGGCCAGCAA
CTTATCTGTGTCTGTCCGATTGTCTAGTGTCTATGTTTGATGTTATGCGCCTGCGTCTGT
ACTAGTTAGCTAACTAGCTCTGTATCTGCGGACCCGTGGTGGAAGTACGAGATTCTGAA
CACCCGGCGCAACCCCTGGGAGACGTCGCCAGGACTTTGGGGGCGGTTTTTGTGGCCGA
CCTGAGGAAGGGAGTCGATGTGGAATCCGACCCCGTCAGGATATGTGTTTCTGTGTAGGAG
ACGAGAACCTAAAACAGTTCCCGCCTCCGTCTGAATTTTTGCTTTCGGTTTTGGAACCGAA
GCCGCGCGTCTTGTCTGCTGCAGCGCTGCAGCATCGTTCGTGTGTGTCTGTCTGACTG
TGTTTTCTGTATTTGTCTGAAAATTAGGGCCAGACTGTTACCACTCCCTTAAAGTTTGACCT
TAGGTCACCTGGAAAGATGTCGAGCGGATCGCTCACAACCACTCGGTAGATGTCAAGAAGA
GACGTTGGGTACCTTCTGCTCTGCAGAATGGCCAACCTTTAACGTCGGATGGCCGCGAG
ACGGCACCTTTAACCGAGACCTCATCACCCAGGTTAAGATCAAGGTCTTTTACCTGGCC
CGCATGGACACCCAGACAGGTCCCCTACATCGTGACCTGGGAAGCCTTGGCTTTTGACC
CCCCTCCCTGGGTCAAGCCCTTTGTACACCCTAAGCCTCCGCCTCCTCTTCTCCATCCG
CCCCGTCTCTCCCCCTTGAACCTCCTCGTTCGACCCCGCCTCGATCCTCCCTTTATCCAG
CCCTCACTCCTTCTCTAGGCGCCGGAATTCCGATCTGATCAAGAGACAGGATGAGGATCG
TTTCGCATGATTGAACAAGATGGATTGCACGCAGGTTCTCCGGCCGCTTGGGTGGAGAGG
CTATTCGGCTATGACTGGGCACAACAGACAATCGGCTGCTCTGATGCCGCCGTGTTCCGG
CTGTCAGCGCAGGGGCGCCCGGTTCTTTTTGTCAAGACCGACCTGTCCGGTGCCCTGAAT
GAACTGCAGGACGAGGCAGCGCGGCTATCGTGGCTGGCCACGACGGGCGTTCCTTGCGCA
GCTGTGCTCGACGTTGTCACTGAAGCGGGAAGGGACTGGCTGCTATTGGGCGAAGTGCCG
GGGCAGGATCTCCTGTCATCTCACCTTGCTCCTGCCGAGAAAGTATCCATCATGGCTGAT

GCAATGCGGCGGCTGCATACGCTTGATCCGGCTACCTGCCCATTCGACCACCAAGCGAAA
CATCGCATCGAGCGAGCACGTACTCGGATGGAAGCCGGTCTTGTGATCAGGATGATCTG
GACGAAGAGCATCAGGGGCTCGCGCCAGCCGAAGTTCGCCAGGCTCAAGGCGCGCATG
CCCGACGGCGAGGATCTCGTCGTGACCCATGGCGATGCCTGCTTGCCGAATATCATGGTG
GAAAATGGCCGCTTTTCTGGATTATCGACTGTGGCCGGCTGGGTGTGGCGGACCGCTAT
CAGGACATAGCGTTGGCTACCCGTGATATTGCTGAAGAGCTTGGCGGCGAATGGGCTGAC
CGCTTCCTCGTGCTTTACGGTATCGCCGCTCCCGATTTCGCAGCGCATCGCCTTCTATCGC
CTTCTTGACGAGTTCTTCTGAGCGGGACTCTGGGGTTCGAAATGACCGACCAAGCGACGC
CCAACCTGCCATCACGAGATTCGATTCCACCGCCGCCTTCTATGAAAGGTTGGGCTTCG
GAATCGTTTTCCGGGACGCCGGCTGGATGATCCTCCAGCGCGGGGATCTCATGCTGGAGT
TCTTCGCCCACCCCGGGCTCGATCCCCTCGCGAGTTGGTTCAGCTGCTGCCTGAGGCTGG
ACGACCTCGCGGAGTTCTACCGGCAGTGCAAATCCGTCGGCATCCAGGAAACCAGCAGCG
GCTATCCGCGCATCCATGCCCCGAACTGCAGGAGTGGGGAGGCACGATGGCCGCTTTGG
TCGAGGCTCCCCTCAGGATATAGTAGTTTCGCTTTTGCATAGGGAGGGGAAATGTAGTC
TTATGCAATACACTTGTAGTCTTGCAACATGGTAACGATGAGTTAGCAACATGCCTTACA
AGGAGAGAAAAAGCACCGTGCCATGCCGATTGGTGGAAGTAAGGTGGTACGATCGTGCCCT
ATTAGGAAGGCAACAGACAGGTCTGACATGGATTGGACGAACCACTGAATTCCGCATTGC
AGAGATAATTGTATTTAAGTGCCTAGCTCGATACAATAAACGCCATTTGACCATTACCA
CATTGGTGTGCACCTCCAAAGCTTGTTAACATCGATAAAAATAAAAGATTTTATTTAGTCT
CCAGAAAAAGGGGGGAATGAAAGACCCACCTGTAGGTTTGGCAAGCTAGCTTAAGTAAC
GCCATTTTGAAGGCATGGAAAAATACATAACTGAGAATAGAGAAGTTCAGATCAAGGTC
AGGAACAGATGGAACAGCTGAATATGGGCCAAACAGGATATCTGTGGTAAGCAGTTCCTG
CCCCGGCTCAGGGCCAAGAACAGATGGAACAGCTGAATATGGGCCAAACAGGATATCTGT
GGTAAGCAGTTCCTGCCCGGCTCAGGGCCAAGAACAGATGGTCCCCAGATGCGGTCCAG
CCCTCAGCAGTTTCTAGAGAACCATCAGATGTTTCCAGGGTGCCCCAAGGACCTGAAATG
ACCCTGTGCCTTATTTGAACTAACCAATCAGTTCGCTTCTCGCTTCTGTTTCGCGCGCTTC
TGCTCCCCGAGCTCAATAAAAGAGCCCAACCCCTCACTCGGGGCGCCAGTCCTCCGAT
TGACTGAGTCGCCCCGGTACCCGTGTATCCAATAAACCCCTCTGCAGTTGCATCCGACTT
GTGGTCTCGCTGTTCTTGGGAGGGTCTCCTCTGAGTGATTGACTACCCGTCAGCGGGG
TCTTTTCAATTGGGGGCTCGTCCGGGATCGGGAGACCCCTGCCCAGGGACCACCGACCCAC
CACCGGGAGGTAAGCTGGCTGCCTCGCGCGTTCGGTGATGACGGTGAAAACCTCTGACA
CATGCAGCTCCCGGAGACGGTCACAGCTTGTCTGTAAGCGGATGCCGGGAGCAGACAAGC
CCGTCAGGGCGCGTCAGCGGGTGTGGCGGGTGTGCGGGCGCAGCCATGACCCAGTCACG
TAGCGATAGCGGAGTGTATACCTAGCTAGGTAGCTAGAGGATCTTTGTGAAGGAACCTTA
CTTCTGTGGTGTGACATAATTGGACAACTACCTACAGAGATTTAAAGCTCTAAGGTA
TATAAAATTTTAAAGTGTATAATGTGTTAACTACTGATTCTAATTGTTTGTGTATTTTA
GATTCCAACCTATGGAAGTGTGAATGGGAGCAGTGGTCCAATGCCTTTAATGAGGAAAA
CCTGTTTTGTCTCAGAAGAAATGCCTCTAGTGATGATGAGGCTACTGCTGACTCTCAACAT
TCTACTCCTCCAAAAAAGAGAAAGTAGAAGACCCCAAGGACTTTCCTTCAGGAATTG
CTAAGTTTTTTGAGTCATGCTGTGTTTGTAGTAATAAGAACTCTTGCTTGCTTTGCTTATC
ACCACAAAGGAAAAAGCTGCACTGCTATACAAGAAAATTATGAAAAAATATTTGATGTAT
AGTGCCTTGACTAGAGATCATAATCAGCCATACCACATTTGTAGAGGTTTTACTTGCTTT
AAAAAACCTCCACACCTCCCCCTGAACCTGAAACATAAAATGAATGCAATTGTTGTTGT
TAACTTGTTTATTGCAGCTTATAATGGTTACAAATAAAGCAATAGCATCACAAATTTTAC
AAATAAAGCATTTTTTTTCACTGCATTCTAGTTGTGGTTTGTCCAACTCATCAATGTATC
TTATCATGTCTGTGATCAGCTTCAGAAGATGGCGGAGGGCCTCCAACACAGTAATTTTCC
TCCGACTCTTAAAATAGAAAATGTCAAGTCAGTTAAGCAGGAAGTGGACTAACTGACGC
AGCTGGCCGTGCGACATCCTCTTTAATTAGTTGCTAGGCAACGCCCTCCAGAGGGCGTG
TGGTTTTGCAAGAGGAAGCAAAAGCCTCTCCACCCAGGCCTAGAATGTTTCCACCCAATC
ATTACTATGACAACAGCTGTTTTTTTTAGTATTAAGCAGAGGCCGGGGACCCCTGGCCCG
CTTACTCTGGAGAAAAAAAACATTGTAGAGGCTTCCAGAGGCAACTTGTCAAAACAGGAC
TGCTTCTATTTCTGTACACTGTCTGGCCCTGTCACAAGGTCCAGCACCTCCATACCCCC
TTTAATAAGCAGTTTGGGAACGGGTGCGGGTCTTACTCCGCCCATCCGCCCTAACTCCG
CCCAGTTCCGCCCATTTCTCCGCCCATGCTGACTAATTTTTTTTATTTATGCAGAGGCCG
AGGCCGCTCGGCCTCTGAGCTATTCCAGAAGTAGTGAGGAGGCTTTTTTGGAGGCTGCA
TTAATGAATCGGCCAACGCGCGGGGAGAGGCGGTTTGCCTATTGGGCGCTCTTCCGCTTC

CTCGCTCACTGACTCGCTGCGCTCGGTTCGGCTGCGGCGAGCGGTATCAGCTCACTC
AAAGGCGGTAATACGGTTATCCACAGAATCAGGGGATAACGCAGGAAAGAACATGTGAGC
AAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCGTTGCTGGCGTTTTTCCATAG
GCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCC
GACAGGACTATAAAGATACCAGGCGTTTTCCCCCTGGAAGCTCCCTCGTGCGCTCTCCTGT
TCCGACCCTGCCGCTTACCGGATACCTGTCCGCCTTTCTCCCTTCGGGAAGCGTGCGCT
TTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTTCGCTCCAAGCTGGG
CTGTGTGCACGAACCCCCCGTTCAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCT
TGAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGAT
TAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGG
CTACACTAGAAGGACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAA
AAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGT
TTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTC
TACGGGTCTGACGCTCAGTGGAACGAAAAACTCAGTTAAGGGATTTTGGTCATGAGATT
ATCAAAAAGGATCTTACCTAGATCCTTTTAAATTAATAAATGAAGTTTTAAATCAATCTA
AAGTATATATGAGTAAACTTGGTCTGACAGTTACCAATGCTTAATCAGTGAGGCACCTAT
CTCAGCGATCTGTCTATTTTCGTTTCATCCATAGTTGCCTGACTCCCCGTCGTGTAGATAAC
TACGATACGGGAGGGCTTACCATCTGGCCCCAGTGCTGCAATGATACCGCGAGACCCACG
CTCACCGGCTCCAGATTTATCAGCAATAAACCAGCCAGCCGGAAGGGCCGAGCGCAGAAG
TGGTCTGCAACTTTATCCGCCTCCATCCAGTCTATTAATTGTTGCCGGAAGCTAGAGT
AAGTAGTTCGCCAGTTAATAGTTTGCACAACGTTGTTGCCATTGCTGCAGGCATCGTGGT
GTCACGCTCGTCGTTTGGTATGGCTTCATTCAGTCCGGTTCCCAACGATCAAGGCGAGT
TACATGATCCCCATGTTGTGCAAAAAAGCGGTAGCTCCTTCGGTCTCCTCCGATCGTTGT
CAGAAGTAAGTTGGCCGAGTGTTATCACTCATGGTTATGGCAGCACTGCATAATTCTCT
TACTGTGTCATGCCATCCGTAAGATGCTTTTCTGTGACTGGTGAACCAACCAAGTCATTCT
GAGAATAGTGTATGCGGCGACCGAGTTGCTCTTGCCCGGCGTCAATACGGGATAATACCG
CGCCACATAGCAGAACTTTAAAAGTGCTCATCATTGGAAAACGTTCTTCGGGGCGAAAAC
TCTCAAGGATCTTACCGCTGTTGAGATCCAGTTCGATGTAACCCACTCGTGCACCCAACT
GATCTTCAGCATCTTTTACTTTACCAAGCGTTTCTGGGTGAGCAAAAACAGGAAGGCAAA
ATGCCGCAAAAAAGGGAATAAGGGCGACACGGAAATGTTGAATACTCATACTCTTCCTTT
TTCAATATTATTGAAGCATTTATCAGGGTTATTGTCTCATGAGCGGATACATATTTGAAT
GTATTTAGAAAAATAAACAATAGGGGTTCCGCGCACATTTCCCCGAAAAGTGCCACCTG
ACGCTAAGAAACCATATTATCATGACATTAACCTATAAAAAATAGGCGTATCACGAGGC
CCTTTCGTCTCGCGCGTTTCGGTGATGACGGTGAAAACCTCTGACACATGCAGCTCCCGG
AGACGGTCACAGCTTGTCTGTAAGCGGATGCCGGGAGCAGACAAGCCCGTCAGGGCGCGT
CAGCGGGTGTGGCGGGTGTGGGGCTGGCTTAACCTATGCGGCATCAGAGCAGATTGTAC
TGAGAGTGCACCATATGCGGTGTGAAATACCGCACAGATGCGTAAGGAGAAAATACCGCA
TCAGGCGCCATTCGCCATTCAGGCTGCGCAACTGTTGGGAAGGGCGATCGGTGCGGGCCT
CTTCGCTATTACGCC

CMV - 21-584

RU5 – 584 -700

NeoR – 1645 – 2645

RSV- 2827 – 3039

3'LTR- 3502 – 3743

AMP- 7495-5026C

Nucleotide Sequence pCLNDX

CGCGTTGACATTGATTATTGACTAGTTATTAATAGTAATCAATTACGGGGTCATTAGTTCATAG
CCCATATATGGAGTTCCGCGTTACATAACTTACGGTAAATGGCCCGCCTGGCTGACCGCCCAA
CGACCCCGCCCATTTGACGTCAATAATGACGTATGTTCCATAGTAACGCCAATAGGGACTTTC
CATTGACGTCAATGGGTGGACTATTTACGGTAAACTGCCCACTTGGCAGTACATCAAGTGTAT
CATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCCCGCCTGGCATTATGCC
CAGTACATGACCTTATGGGACTTTCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTA
CCATGTGATGCGGTTTTTGGCAGTACATCAATGGGCGTGGATAGCGGTTTGACTCACGGGGATT
TCCAAGTCTCCACCCCATTTGACGTCAATGGGAGTTTGTGTTTGGCACCAAAATCAACGGGACTT
TCCAAAATGTCGTAACAACCTCCGCCCCATTGACGCAAATGGGCGGTAGGCGTGTACGGTGGG
AGGTCTATATAAGCAGAGCTCAATAAAAGAGCCCAACCCCTCACTCGGCGCGCCAGTCTTC
CGATAGACTGCGTCGCCCCGGTACCCGTATTCCCAATAAAGCCTCTTGCTGTTTGCATCCGAA
TCGTGGTCTCGCTGTTCTTGGGAGGGTCTCCTCTGAGTGATTGACTACCCACGACGGGGGTCT
TTCATTTGGGGGCTCGTCCGGGATTTGGAGACCCCTGCCCAGGGACCACCGACCCACCACCGG
GAGGTAAGCTGGCCAGCAACTTATCTGTGTCTGTCCGATTGTCTAGTGTCTATGTTTGATGTTA
TGCGCCTGCGTCTGTACTAGTTAGCTAACTAGCTCTGTATCTGGCGGACCCGTGGTGGAACTG
ACGAGTTCTGAACACCCGCGCCGAACCTGGGAGACGTCCAGGGACTTTGGGGGCCGTTTTT
GTGGCCCGACCTGAGGAAGGGAGTGCATGTGGAATCCGACCCCGTCAGGATATGTGGTTCTG
GTAGGAGACGAGAACCTAAAACAGTTCCCGCCTCCGTCTGAATTTTTGCTTTCGGTTTTGGAAC
CGAAGCCGCGCGTCTTGCTGCTGCAGCGCTGCAGCATCGTTCTGTGTTGTCTCTGTCTGACTG
TGTTTTCTGTATTTGTCTGAAAATTAGGGCCAGACTGTTACCACTCCCTTAAGTTTGACCTTAGG
TCACTGGAAAGATGTCGAGCGGATCGCTCACAACCAGTCGGTAGATGTCAAGAAGAGACGTT
GGGTTACCTTCTGCTCTGCAGAATGGCCAACTTTAACGTTCGGATGGCCGCGAGACGGCACCT
TTAACCGAGACCTCATCACCCAGGTAAAGATCAAGGTCTTTTACCTGGCCCGCATGGACACC
CAGACCAGGTCCCCTACATCGTGACCTGGGAAGCCTTGCTTTTGACCCCTCCCTGGGTCA
AGCCCTTTGTACACCCTAAGCCTCCGCCTCCTCTTCTCCATCCGCCCCGTCTCTCCCTTGA
ACCTCTCGTTTCGACCCCGCCTCGATCCTCCCTTTATCCAGCCCTCACTCCTTCTCTAGGCGCC
GGAATTCCGATCTGATCAAGAGACAGGATGAGGATCGTTTCGCATGATTGAACAAGATGGAT
TGCACGCAGGTTCTCCGGCCGCTTGGGTGGAGAGGCTATTCCGGCTATGACTGGGCACAACAGA
CAATCGGCTGCTCTGATGCCGCCGTGTTCCGGCTGTCAGCGCAGGGGCGCCCGGTTCTTTTTGT
CAAGACCGACCTGTCCGGTGCCCTGAATGAACTGCAGGACGAGGCAGCGCGGCTATCGTGGC
TGGCCACGACGGGCGTTCTTGCGCAGCTGTGCTCGACGTTGTCACTGAAGCGGGAAGGGACT
GGCTGCTATTGGGCGAAGTGCCGGGGCAGGATCTCCTGTCTCTCACCTTGCTCCTGCCGAGA
AAGTATCCATCATGGCTGATGCAATGCGGCGGCTGCATACGCTTGATCCGGCTACCTGCCCAT
TCGACCACCAAGCGAAACATCGCATCGAGCGAGCACGTAAGTTCGGATGGAAGCCGGTCTTGTC
GATCAGGATGATCTGGACGAAGAGCATCAGGGGCTCGCGCCAGCCGAACTGTTCCGCCAGGCT
CAAGGCGCGCATGCCCGACGGCGAGGATCTCGTCTGACCCATGGCGATGCCTGCTTGCCGA
ATATCATGGTGGAAAATGGCCGCTTTTCTGGATTTCATCGACTGTGGCCGGCTGGGTGTGGCGG
ACCGCTATCAGGACATAGCGTTGGCTACCCGTGATATTGCTGAAGAGCTTGGCGGCGAATGGG
CTGACCGCTTCTCGTGCTTTACGGTATCGCCGCTCCCGATTTCGCAGCGCATCGCCTTCTATCG
CCTTCTTGACGAGTTCTTCTGAGCGGGACTCTGGGGTTCGAAATGACCGACCAAGCGACGCCC
AACCTGCCATCACGAGATTTGATTCCACCGCCGCTTCTATGAAAGGTTGGGCTTCGGAATC
GTTTTCCGGGACGCCGGCTGGATGATCCTCCAGCGCGGGGATCTCATGCTGGAGTTCTTCGCC
CACCCCGGGCTCGATCCCCTCGCGAGTTGGTTCAGCTGCTGCCTGAGGCTGGACGACCTCGCG
GAGTTCTACCGGCAGTGCAAATCCGTCGGCATCCAGGAAACCAGCAGCGGCTATCCGCGCAT
CCATGCCCCCGAACTGCAGGAGTGGGGAGGCACGATGGCCGCTTTGGTTCGAGGCGGATCAAT
TCGATAGCTTGGGCTGCTAGGAGCGCGAGCGCGCGGCCGCACTTTCTCGCGCCTGCGCGCGC
GCACGCCTCAACCTGTGCGGGACCGGCCTTGGGGGCGGAGCCTTAGCTACACAAATAGAATG
CGCGGCGGGCCTTGGTGGGGGCGGGGCCTTAGCTGCACAAATAGGATGCGCGGCGGGCCTTG
TAGGGGCGGAGCCTTAGCTGCACAAATAGGATGCGCGGCGGGCCTTGGTGGGGGCGGGGCC
TAAGCTGCGCAAGTGGTACACAGCTCAGGGCTGCGATTTCGCGCCAAACTTGACGGCAAATCC
TAGCGTGAAGGCTGGTAGGATTTTATCCCCGTGCCATCATGGTTTCGACCTGCAGCCAAGCTA
TCAAGAGCTTGTTTAACATCGATAAAATAAAAGATTTTATTTAGTCTCCAGAAAAGGGGGGA

ATGAAAGACCCACCTGTAGGTTTGGCAAGCTAGCTTAAGTAACGCCATTTTGAAGGCATGG
 AAAAATACATAACTGAGAATAGAGAAGTTCAGATCAAGGTCAGGAACAGATGGAACAGCTG
 AATATGGGCCAAACAGGATATCTGTGGTAAGCAGTTCCTGCCCGGCTCAGGGCCAAGAACA
 GATGGAACAGCTGAATATGGGCCAAACAGGATATCTGTGGTAAGCAGTTCCTGCCCGGCTC
 AGGGCCAAGAACAGATGGTCCCCAGATGCGGTCCAGCCCTCAGCAGTTTCTAGAGAACCATC
 AGATGTTTCCAGGGTGCCCCAAGGACCTGAAATGACCTGTGCCTTATTTGAACTAACCAATC
 AGTTCGCTTCTCGCTTCTGTTCGCGCGCTTCTGCTCCCCGAGCTCAATAAAAGAGCCCACAACC
 CCTCACTCGGGGCGCCAGTCCCTCCGATTGACTGAGTCGCCCCGGGTACCCGTGTATCCAATAAA
 CCCTCTTGCAAGTTGCATCCGACTTGTGGTCTCGCTGTTTCTTGGGAGGGTCTCCTCTGAGTGAT
 TGACTACCCGTCAGCGGGGGTCTTTCATTTGGGGGCTCGTCCGGGATCGGGAGACCCCTGCCC
 AGGGACCACCGACCCACCACCGGGAGGTAAGCTGGCTGCCTCGCGCGTTTCGGTGATGACGG
 TGAAAACCTCTGACACATGCAGCTCCCGGAGACGGTCACAGCTTGTCTGTAAGCGGATGCCGG
 GAGCAGACAAGCCCGTCAGGGCGCGTCAGCGGGTGTGGCGGGTGTGCGGGGCGCAGCCATGA
 CCCAGTACGTAGCGATAGCGGAGTGTATACCTAGCTAGGTAGCTAGAGGATCTTTGTGAAGG
 AACCTTACTTCTGTGGTGTGACATAAATTGGACAACTACCTACAGAGATTTAAAGCTCTAAGG
 TAAATATAAAATTTTTAAGTGTATAATGTGTTAACTACTGATTCTAATTGTTTGTGATTTTTA
 GATTCCAACCTATGGAAGTATGAATGGGAGCAGTGGTCCAATGCCTTTAATGAGGAAAACCT
 GTTTTGTCTCAGAAGAAATGCCTCTAGTGATGATGAGGCTACTGCTGACTCTCAACATTCTACTC
 CTCCAAAAAAGAAGAGAAAGGTAGAAGACCCCAAGGACTTTCCTTCAGAATTGCTAAGTTTTT
 TGAGTCATGCTGTGTTTAGTAATAGAAGTCTTGCTTGCTTTGCTATTTACACCACAAAGGAAAA
 AGCTGCACTGCTATACAAGAAAATTATGGAAGAAATATTTGATGTATAGTGCCTTGACTAGAGA
 TCATAATCAGCCATACCACATTTGTAGAGGTTTTACTTGCTTTAAAAAACCTCCCACACCTCCC
 CCTGAACCTGAAACATAAAATGAATGCAATTGTTGTTGTTAACTTGTTTATTGCAGCTTATAAT
 GGTACAAATAAAGCAATAGCATCACAAATTCACAAATAAAGCATTTTTTTTTACTGCATTCT
 AGTTGTGGTTTTGTCCAACTCATCAATGTATCTTATCATGTCTGGATCAGCTTCAGAAGATGGC
 GGAGGGCCTCCAACACAGTAATTTTCTCCCGACTCTTAAATAGAAAATGTCAAGTCAGTTA
 AGCAGGAAGTGGACTAACTGACGCAGCTGGCCGTGCGACATCCTCTTTTAATTAGTTGCTAGG
 CAACGCCCTCCAGAGGGCGTGTGGTTTTGCAAGAGGAAGCAAAAGCCTCTCCACCCAGGCCT
 AGAATGTTTCCACCCAATCATTACTATGACAACAGCTGTTTTTTTTTAGTATTAAGCAGAGGCCG
 GGGACCCCTGGCCCGCTTACTCTGGAGAAAAAAAACATTGTAGAGGCTTCAGAGGCAACTT
 GTCAAAACAGGACTGCTTCTATTTCTGTACACTGTCTGGCCCTGTCAAGGTCCAGCACCTC
 CATACCCCTTTAATAAGCAGTTTGGGAACGGGTGCGGGTCTTACTCCGCCCATCCGCCCTA
 ACTCCGCCCAGTTCGCCCATTTCTCCGCCCATGCTGACTAATTTTTTTTATTATGCAGAGGC
 CGAGGCCGCTCGGCCTCTGAGCTATTCCAGAAGTAGTGAGGAGGC
 TTTTTTGGAGGCTGCATTAATGAATCGGCCAACGCGCGGGGAGAGGCGGTTTGCATTTGGGC
 GCTCTTCGCTTCTCGCTCACTGACTCGCTGCGCTCGGTGCTTCGGCTGCGGCGAGCGGTATC
 AGCTCACTCAAAGGCGGTAAATACGGTTATCCACAGAATCAGGGGATAACGCAGGAAAGAACA
 TGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCGTTGCTGGCGTTTTTC
 CATAGGCTCCGCCCTGACGAGCATCAAAAAATCGACGCTCAAGTCAGAGGTGGCGAAA
 CCCGACAGGATAAAAGATACACAGGCGTTTCCCCCTGGAAGCTCCCTCGTGCCTCTCTCTGT
 TCCGACCCTGCCGCTTACCGGATACCTGTCCGCTTTTCTCCCTTCGGAAGCGTGCGCTTTCT
 CATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTGCTTCGCTCCAAGCTGGGCTGTGTG
 CACGAACCCCCCGTTACGCGGACCGCTGCGCCTTATCCGGTAAGTATCGTCTTGAGTCCAAC
 CCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAG
 GTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGGAC
 AGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAAGAGTTGGTAGCTCTTG
 ATCCGGCAAACAACACCGCTGGTAGCGGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCG
 CAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGA
 CGAAAACCTACGTTAAGGGATTTTGGTCATGAGATTATCAAAAAGGATCTTCACCTAGATCCT
 TTTAAATTAATAAATGAAGTTTTAAATCAATCTAAAGTATATATGAGTAACTTGGTCTGACAG
 TTACCAATGCTTAATCAGTGAGGCACCTATCTCAGCGATCTGTCTATTTTCGTTTCATCCATAGTT
 GCCTGACTCCCGTCTGTGTAGATAACTACGATACGGGAGGGCTTACCATCTGGCCCCAGTGCT
 GCAATGATACCGCGAGACCCACGCTCACCGGCTCCAGATTTATCAGCAATAAACCAGCCAGC
 CGGAAGGGCCGAGCGCAGAAAGTGGTCTGCAACTTTATCCGCCTCCATCCAGTCTATTAATTG
 TTGCCGGGAAGCTAGAGTAAGTAGTTCCGCAAGTTAATAGTTTGCAGCAACGTTGTTGCCATTGC
 TGCAGGCATCGTGGTGTACGCTCGTCTTGGTATGGCTTCATTACGCTCCGGTTCCCAACGA

TCAAGGCGAGTTACATGATCCCCCATGTTGTGCAAAAAAGCGGTTAGCTCCTTCGGTCCTCCG
ATCGTTGTCAGAAGTAAGTTGGCCGCAGTGTTATCACTCATGGTTATGGCAGCACTGCATAAT
TCTCTTACTGTCATGCCATCCGTAAGATGCTTTTCTGTGACTGGTGAACCTCAACCAAGTCATTC
TGAGAATAGTGTATGCGGCGACCGAGTTGCTCTTGCCCGGCGTCAATACGGGATAATACCGCG
CCACATAGCAGAACTTTAAAAGTGCTCATCATTGGAAAACGT
TCTTCGGGGCGAAAACTCTCAAGGATCTTACCGCTGTTGAGATCCAGTTCGATGTAACCCACT
CGTGCACCCAACTGATCTTCAGCATCTTTTACTTTACCAGCGTTTCTGGGTGAGCAAAAACA
GGAAGGCAAAATGCCGCAAAAAAGGGAATAAGGGCGACACGGAAATGTTGAATACTCATAC
TCTTCCTTTTTCAATATTATTGAAGCATTTATCAGGGTTATTGTCTCATGAGCGGATACATATTT
GAATGTATTTAGAAAAATAAACAAATAGGGGTTCGCGCACATTTCCCGAAAAAGTGCCACCT
GACGTCTAAGAAACCATTATTATCATGACATTAACCTATAAAAAATAGGCGTATCACGAGGCCC
TTTCGTCTCGCGCGTTTCGGTGATGACGGTGAAAACCTCTGACACATGCAGCTCCCGGAGACG
GTCACAGCTTGTCTGTAAGCGGATGCCGGGAGCAGACAAGCCCGTCAGGGCGCGTCAGCGGG
TGTTGGCGGGTGTCGGGGCTGGCTTAACCTATGCGGCATCAGAGCAGATTGTAAGTACTGAGAGTGCA
CCATATGCGGTGTGAAATACCGCACAGATGCGTAAGGAGAAAAATACCGCATCAGGCGCCATT
CGCCATTCAGGCTGCGCAACTGTTGGGAAGGGCGATCGGTGCGGGCCTCTTCGCTATTACGCC

CMV – 21- 584 bp
RU5 – 584- 700 bp
Neo – 1645 – 2643
DHFR – 2832 – 3615
3'LTR – 3637 – 3878
AMP – 7624 – 5155C