
Product Manual

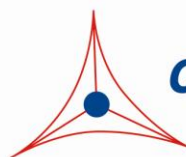
RAPAd[®] CMV Adenoviral Bicistronic Expression System (GFP)

Catalog Number

VPK-254

1 kit

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.

Creating Solutions for Life Science Research

Introduction

Recombinant adenoviruses have tremendous potential in both research and therapeutic applications. There are numerous advantages they provide when introducing genetic material into host cells. The permissive host cell range is very wide. The virus has been used to infect many mammalian cell types (both replicative and non-replicative) for high expression of the recombinant protein. Recombinant adenoviruses are especially useful for gene transfer and protein expression in cell lines that have low transfection efficiency with liposome. After entering cells, the virus remains epichromosomal (i.e. does not integrate into the host chromosome so does not activate or inactivate host genes). Recently, recombinant adenoviruses have been used to deliver RNAi into cells.

Two methods have traditionally been used to generate recombinant adenoviruses. The first involves homologous recombination of a shuttle vector containing gene of interest and an adenoviral backbone plasmid vector (restricted in E1/E3) in an adenovirus packaging cell line. The isolation of recombinant adenovirus by this method involves performing multiple plaque isolations to avoid wild-type virus and is extremely laborious and time consuming. The second method, pAdEasy system, employs the homologous recombination machinery in *E. coli*, a recombinant adenovirus is produced by a double-recombination event between cotransformed adenoviral backbone plasmid vector and a shuttle vector carrying the gene of interest. For the pAdEasy method, the system is high fidelity, but inefficient and requires the screening of many bacterial colonies. This results in a significant time commitment even before transfection of recombinant DNA into E1-expressing cells such as HEK293 cells.

Cell Biolabs' RAPAd® Adenoviral Expression System provides a much faster and safer method to generate RCA-free recombinant adenovirus at high titer (see Table 1). The RAPAd® system uses a novel Ad backbone devoid of the left-hand ITR, the packaging signal and E1 sequences. There is no need to perform the bacterial *in vitro* homologous recombination (pAdEasy method), and also the multiple plaque isolations (standard homologous recombination method in packaging cell line). The RAPAd® system allows for generation of a recombinant virus within 2 weeks and the virus produced contained virtually no contaminating E1a sequences or replication-competent virus (RCA).

Cell Biolabs' RAPAd® Adenoviral Expression System is simple to use. The method is straightforward and requires very limited 'hands on' time from shuttle/backbone cotransfection to the isolation of virus particles. It produces equivalent infectious titers as the standard viral genome/shuttle plasmid recombination method.

In Cell Biolabs' RAPAd® CMV Adenoviral Bicistronic Expression System (GFP), the shuttle vector contains a CMV ahead of the multiple cloning sites followed by IRES-GFP.

Standard Homologous Recombination	pAdEasy Expression System	RAPAd® Expression System
<p data-bbox="155 327 493 428">Cotransfect 293 cells with Shuttle Vector and Ad Backbone Vector</p> <p data-bbox="310 464 326 604">↓</p> <p data-bbox="261 621 380 722">Multiple Plaque Isolations</p> <p data-bbox="310 747 326 888">↓</p> <p data-bbox="233 911 412 978">Virus Amplification</p> <p data-bbox="310 989 326 1129">↓</p> <p data-bbox="250 1163 396 1192">Viral Stock</p>	<p data-bbox="610 327 799 428">Linearize Shuttle Vector using PmeI</p> <p data-bbox="691 464 708 604">↓</p> <p data-bbox="526 621 883 722">Cotransform <i>E.coli</i> BJ5183 cells with linearized Shuttle Vector and pAdEasy Vector</p> <p data-bbox="691 747 708 888">↓</p> <p data-bbox="529 911 880 978">Recombinant selection by restriction enzyme analysis</p> <p data-bbox="691 989 708 1129">↓</p> <p data-bbox="558 1163 850 1230">Linearize recombinant plasmid using PacI</p> <p data-bbox="691 1241 708 1381">↓</p> <p data-bbox="581 1415 828 1444">Transfect 293 cells</p> <p data-bbox="691 1472 708 1612">↓</p> <p data-bbox="630 1640 779 1669">Viral Stock</p>	<p data-bbox="915 327 1279 428">Linearize Shuttle Vector and RAPAd® Ad Backbone Vector using PacI</p> <p data-bbox="1097 464 1114 604">↓</p> <p data-bbox="1023 621 1175 680">Cotransfect 293 cells</p> <p data-bbox="1097 747 1114 888">↓</p> <p data-bbox="1023 911 1175 940">Viral Stock</p>
12-18 weeks	8-9 weeks	2-3 weeks

Table 1. Outline of Recombinant Adenovirus Systems

Related Products

1. AD-100: 293AD Cell Line
2. AD-200: ViraDuctin™ Adenovirus Transduction Reagent
3. VPK-099: ViraBind™ Adenovirus Miniprep Kit
4. VPK-110: QuickTiter™ Adenovirus Titer ELISA Kit
5. VPK-111: Rapid RCA Assay Kit
6. VPK-250: RAPAd® Universal Adenoviral Expression System
7. VPK-251: RAPAd® RSV Adenoviral Expression System
8. VPK-253: RAPAd® miRNA Adenoviral Expression System

Kit Components

1. pacAd5 CMV-IRES-GFP Shuttle Vector (Part No. 325401): One 40 µL vial at 0.25 mg/mL.
2. pacAd5 9.2-100 Vector (Part No. 325002): One 40 µL vial at 0.25 mg/mL.
3. pacAd5 CMV-GFP Control Vector (Part No. 325004): One 40 µL vial at 0.25 mg/mL.
4. pacAd5 CMV-ntLacZ Control Vector (Part No. 325202): One 40 µL vial at 0.25 mg/mL.

Materials Not Supplied

1. 293 cells: we recommend 293AD Cell Line (Cat.# AD-100) for high titer production of recombinant adenovirus.
2. 293 Cell Culture Medium
3. Transfection Reagents
4. PacI (New England Biolabs, Cat.# R0547L)

Storage

Upon receipt, store all kit components at -20°C.

Safety Considerations

Remember that you will be working with samples containing infectious virus. Follow the recommended NIH guidelines for all materials containing BSL-2 organisms.

Vector Features

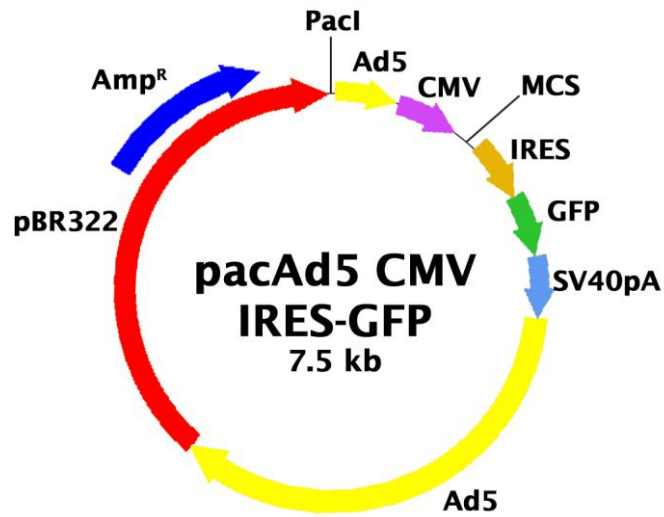


Figure 1. pacAd5 CMV-IRES-GFP shuttle Vector (7539 bp, Ampicillin-resistant).

pacAd5 CMV-IRES-GFP Features:

3-10:	PacI
16-368:	1-353 of Ad5
382-912	CMV Promoter
919-964:	MCS
983-1596:	IRES
1597-2316:	GFP
2317-2764:	SV40 pA
2759-5223:	3328-5792 of Ad5
6471-7331:	β Lactamase

Multiple Cloning Sites:

Pme I
EcoR I
 TTGGTACCGTTTAAACTCGAGGTCGACGGTATCGATAAGCTTGATATCGAATTCCTGCAGCCCGGGGATCCACTAGT
Cla I
EcoR V
BamH I

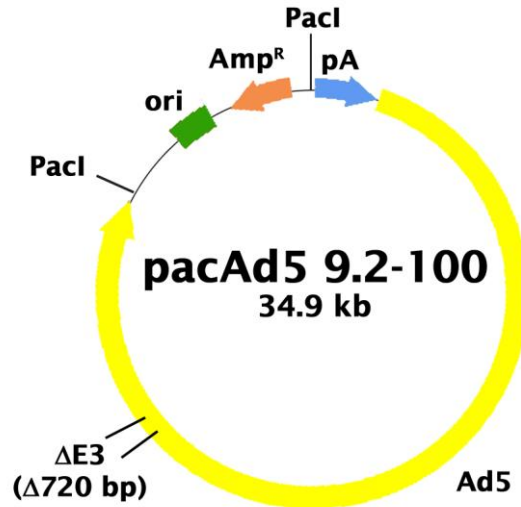


Figure 2. pacAd5 9.2-100 Vector (34947 bp, Ampicillin-resistant). The novel pacAd5 9.2-100 Ad backbone vector is devoid of the left-hand ITR, the packaging signal and E1 sequences.

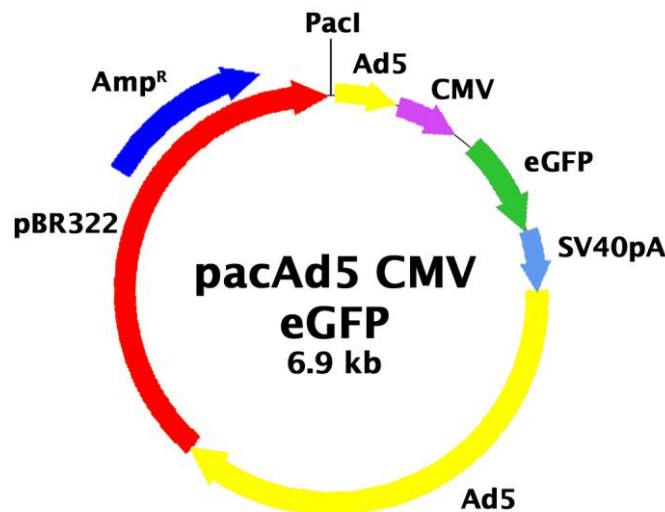


Figure 3. pacAd5 CMV-GFP Control Vector (6935 bp, Ampicillin-resistant).

pacAd5 CMV-GFP Features:

3-10:	PacI
16-368:	1-353 of Ad5
385-912:	CMV Promoter
992-1711:	GFP
1713-2160:	SV40 pA
2161-4615:	3328-5792 of Ad5
5867-6727:	β Lactamase

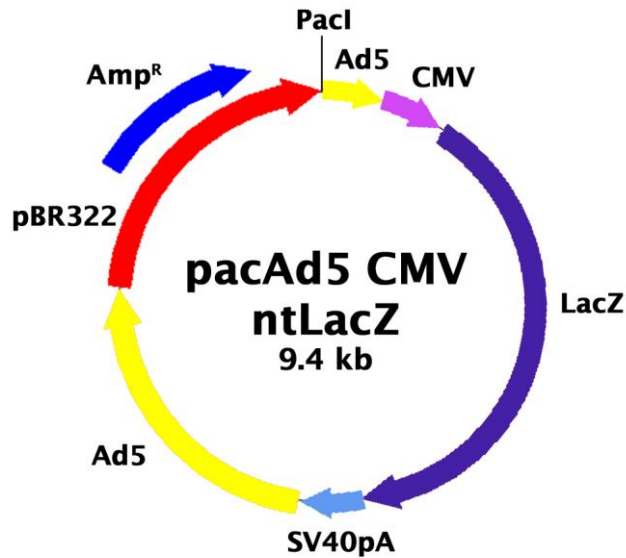


Figure 4. pacAd5 CMV-ntLacZ Control Vector (9278 bp, Ampicillin-resistant).

pacAd5 CMV-ntLacZ Features:

3-10:	PacI
16-368:	1-353 of Ad5
385-912:	CMV Promoter
1105-4148:	ntLacZ
4193-4640:	SV40 pA
4641-7095:	3328-5792 of Ad5
8347-9210:	β Lactamase

Preparation of Recombinant Adenovirus

I. Vector Linearization with PacI

1. Digest a sufficient amount of the pacAd5 CMV-IRES-GFP shuttle vector containing gene of interest and the pacAd5 9.2-100 Ad backbone vector with PacI.
2. Run 0.5 μ g of each digested DNA and undigested DNA on a 0.8% agarose gel to confirm the completion of PacI digestion (For pacAd5 9.2-100, one band of ~33 kb and a second band of 2.0 kb).
3. Remove buffer and enzyme from the remainder of the restriction reactions by phenol extraction/ethanol precipitation or using a similar DNA purification kit.
4. Resuspend the DNA in sterile dH₂O. Store the digested DNA at -20°C.

II. Transfection

1. Seed 2×10^6 cells in a 60 mm culture dish without antibiotics one day before transfection.

2. After 16 to 24 hours, start transfection when the culture becomes 70-80% confluence.
Note: We suggest transfecting cells with FuGENE® Transfection Reagent (Roche Applied Science) or Lipofectamine™ Plus (Invitrogen). For example, 4 µg of pacAd5 CMVK-NpA shuttle vector and 1 µg of pacAd5 9.2-100 Ad backbone vector are mixed with 9 µL FuGENE® Transfection Reagent according to the manufacturer's recommendation. The mixed DNA-FuGENE® complex is added by dropwise into the culture media.
3. Aspirate the media containing transfection reagent the next day and add 4 mL of complete culture medium.
4. After incubating for 7 days, check for the presence of plaques. If plate is ready for harvest, (>50% of cells lifted), then collect the Crude Viral Lysate. If not, feed the cells with 1 mL of complete culture medium, continue to incubate at 37°C with CO₂.
5. On day 10, check for the presence of plaques. If plate is ready for harvest, (>50% of cells lifted), then collect the Crude Viral Lysate. If not, feed the cells with 1 mL of complete culture medium, continue to incubate at 37°C with CO₂. Keep checking plate for the presence of plaques. Do not keep plate more than 15 days.

III. Harvesting the Crude Viral Lysate

1. Harvest adenovirus-containing cells by squirting cells off the plate with a 5 or 10 mL sterile serological pipette. Transfer cells and media to a sterile 15 ml tube. Scrape the cells into the medium with a cell lifter if necessary.
2. Release viruses from cells by three freeze/thaw cycles (10 minutes each in 37°C water bath and dry ice-methanol bath).
3. Centrifuge the cell lysate in a table-top centrifuge at 3000 rpm for 15 minutes at room temperature to pellet the cell debris.
4. Aliquot and store the Crude Viral Lysate (Initial Viral Stock) at -80°C.

IV. Amplification

Note: The following procedure is suggested for T75 flasks and may be optimized to suit individual needs.

1. Seed 3-5 x 10⁶ cells in a T75 flask one day before infection.
2. Add 50% of the above Crude Viral Lysate to the culture. We recommend using a multiplicity of ≥0.5 PFU (plaque forming units) or enough viruses that cells demonstrate cytopathic effects (CPEs) within 48 hrs.
3. During 24 - 48 hr infection, examine the monolayer twice per day under the microscope for CPE. When CPE is nearly complete (i.e. most cells rounded but not yet detached from the flask), harvest cells by pipetting media up and down to wash the infected cells from the flask into the media.
4. Pool infected cells and medium. Pellet cells by centrifugation at 1000 g for 5 minutes. Remove supernatant, resuspend cell pellet in medium or in 10 mM Tris, pH 8.0, 100 mM NaCl. (0.25-0.5 mL per T75 flask).
5. Release the adenoviruses from the cell suspension with three freeze/thaw cycles. Centrifuge at 3000 g for 10 minutes to pellet the cell debris. Discard the pellet and save supernatant as viral stock.
6. The viral supernatant can be stored at -80°C or immediately purified or titered.

Example of Results

The following figures demonstrate typical results of generating recombinant adenovirus. One should use the data below for reference only. This data should not be used to interpret actual results.

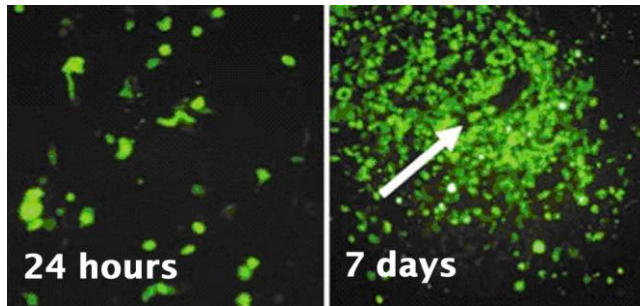


Figure 5. Generation of recombinant adenovirus using the RAPAd® Adenoviral Expression System. 293 cells were transfected with PacI linearized pacAd5 CMV-GFP vector and pacAd5 9.2-100 vector. Plates were examined for the presence of viral foci under inverted fluorescence microscope.

Appendix

pacAd5 CMV-IRES-GFP Plasmid Sequence

```
AATTAATTAAGCTAGCATCATCAATAATATACCTTATTTGGATTGAAGCCAATATGATAATGAGGGGGTGGAGTTTGTGACGTGGCGGGGGCGTGGGAACG
GGGCGGGTACGCTAGTAGTGTGGCGGAAGTGTGATGTTGCAAGTGTGGCGGAACACATGTAAGCGACGGATGTGGCAAAGTACGTTTGGTGTGCGCCGG
TGTACACAGGAAGTGACAATTTTCGCGCGGTTTGTAGCGGATGTTGTAGTAAATTTGGCGCTAACCGAGTAAGATTTGGCCATTTTCGCGGGAAAACCTGAATA
AGAGGAAGTGAATCTGAATAATTTGTGTTACTCATAGCGCGTAATATTTGTCTAGGGAGATCAGCCTGCAGGTCGTACATAACTTACGGTAAATGGCCCG
CTGGCTGACCGCCCAACGACCCCGCCCATTTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCCATTGACGTCAATGGGTGGAGT
ATTTACGGTAAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCCCGCTGGCATTATGC
CCAGTACATGACCTTATGGGACTTTCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGGTGTGCGGTTTTGGCAGTACATCAATGGGCGT
GGATAGCGGTTTACTCAGCGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGGTGGCACCAAAATCAACGGGACTTTCCAAAATGTGCGTA
ACAACCTCCGCCCATTTGACGCAAAATGGGCGGTAGGCGGTGTACGGTGGGAGGCTATATAAGCAGAGCTCGTTAGTGAACCGTCAGTTGGTACCGTTAAACT
CGAGGTCGACGGTATCGATAAGCTTGATATCGAATTCCTGCAGCCCGGGGGATCCTCGAGGGCCGGCGCCCGCGCCGCTACGTAATTTCCGCCCTCTCCC
TAACGTTACTGGCCGAAGCCGCTTGGAAATAAGGCCGGTGTGCGTTTTGTCTATATGTTATTTCCACCATAATGCGCGCTTTTGGCAATGTGAGGGCCCGGAAA
CCTGGCCCTGTCTTCTGACGAGCATTCCTAGGGGTCTTTCCCTCTCGCCAAAGGAATGCAAGGTCGTGTAATGTGCGTGAAGGAAGCAGTTCCTCTGGAAG
CTTCTTGAAGCAAAACACGCTGTGAGCGACCTTTGCAGGCGAGCGGAACCCCACTGGCGACAGGTGCCTCTGCGGCCAAAAGCCACGTATATAGATAC
ACCTGCAAAGGCGGCACAACCCAGTGCCACGTTGTGAGTTGGATAGTTGTGGAAGAGTCAAATGGCTCTCTCAAGCGTATTCAACAAGGGGTGAAGGAT
GCCAGAAGGTACCCATTTGATGGGATCTGATCTGGGGCCTCGGTGCACATGCTTTACATGTGTTTAGTCGAGGTTAAAAAACGCTAGGCCCCCGGAACC
ACGGGGACGTGGTTTTCTTTGAAAAACACGATGATAATATGGCCACAACCATGGTGTGCAAGGGCGAGGAGCTGTTACCGGGGTGGTGGCCATCTGGTGC
AGCTGGACGGGACGTAACGGCCACAAGTTCAGCGTGTCCGGCGAGGGCGAGGGCGATGCCACCTACGGCAAGCTGACCCGTAAGTTCATCTGCACCACCGG
CAAGCTGCCCGTGCCTTGGCCACCCCTCGTGACACCCCTGACCTACGGCGTGCAGTCTTACAGCCGTACCCCGACCCATGAAGCAGCAGCACTTCTTCAAG
TCCGCCATGCCGAAGGCTACGTCCAGGAGCGCACCATCTTCTTCAAGGACGACGGCAACTACAAGACCCCGCGCGAGGTGAAGTTCGAGGGCGACACCCCTGG
TGAACCGCATCGAGCTGAAGGGCATCGACTTCAAGGAGGACGGCAACATCTGGGGCACAAGCTGGAGTACAACATAACAGCCACACCTATATCATGGC
CGACAAGCAGAAGAAGCGCATCAAGGTGAACCTCAAGATCCGCCACAACATCGAGGACGGCAGCGTGCAGCTCGCCGACCACTACCAGCAGAACACCCCATC
GGCAGGCCCCGTGTGCTGCCGACAACCACTACCTGAGCACCCAGTCCGCCCTGAGCAAAGACCCCAACGAGAAGCGCGATCACATGGTCTGCTGGAGT
TCGTGACCGCCGCGCGGATCACTCTCGGCATGGACGAGCTGTACAAGTAAGCGGCCACCGCGGGGAGATCCAGACATGATAAGATACATTGATGAGTTTGA
GACAAACCAACTAGAAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATATAAGCTGCAATAAACAAGTTAA
CAACAACAATTCATTCATTTTATGTTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTTACAAAATGTGGTATGGCTGATATGAT
CCCGGTGCCTCGCGCGTTTTCGGTGATGACGGTGAACCTCTTGACACATGCAGCTCCCGGAGACGGTACAGCTTGTCTGTAAGCGGATGCCGGGAGCAGA
CAAGCCGTCAGGGCGGTGACGCGGTGTTGGCGGTGTCGGGGCGCAGCCATGAGGTCGACTCTAGTCCCGCGGTGGCAGATCTGGAAGGTGCTGAGGTAC
GATGAGACCCGACCAAGGTGACAGACCTGCGAGTGTGGCGTAAACATATTTAGGAACAGCCTGTGATGCTGGATGTGACCCAGGAGCTGAGGCCCGATCACT
TGGTGTGCGCTGCACCCGCTGAGTTTGGCTCTAGCGATGAAGATACAGATTGAGGTACGAAATGTGTGGCGTGGCTTAAAGGTGGGAAAGAATATATA
AGGTGGGGTCTTATGTAGTTTGTATCTGTTTTGACGAGCCCGCCCATGAGCACAACCTCGTTTGTGGAAGCATTTGAGCTCATATTTGACAACG
CGCATGCCCATGGCCGGGGTGGCTCAGAATGTGATGGGCTCCAGCATTTGATGGTGCACCCGCTCTGCCGCAAACCTACTACTACCTGACCTACGAGACCG
TGTCTGGAACCGGTTGGAGACTGCAGCCTCCGCCCGCGCTTACGGCGTGCAGCCACCGCCCGGGGATTTGACTGACTTTGCTTTCTGAGCCCGCTTGC
AAGCAGTGCAGCTTCCCGTTCATCCGCCCGGATGACAAGTTGACGGCTCTTTGGCACAATTTGGATTCTTTGACCCGGGAACCTAATGTGCTTTCTCAGCAG
CTGTTGGATCTGCGCCAGCAGGTTTCTGCCCTGAAGGCTTCTCCCTCCCAATGCGGTTTTAAACATAAATAAAAAACAGACTCTGTTGGATTTGGATCA
AGCAAGTGTCTGCTGCTTTATTTAGGGGTTTTGCGCGCGCGGTAGGCCCGGGACAGCGCTCTCGGTTCGTTGAGGGTCCGTGTATTTTTCCAGGACGTG
GTAAAGGTGACTCTGGATGTTTCAAGTACATGGGCATAAGCCGCTCTTGGGGTGGAGGTAGCACCCTGCAGAGCTTCATGCTGCGGGGTGGTGTGTAGATG
ATCCAGTCTGACAGGAGCGTGGGCGTGGTGCCTAAAATGTCTTTCAGTAGCAAGCTGATTGCCAGGGGACGGCCCTTGGTGAAGTGTTTACAAAGCGGT
```

TAAGCTGGGATGGGTGCATACGTGGGATATGAGATGCATCTTGGACTGTATTTTTAGGTTGGCTATGTGCCAGCCATATCCCTCCGGGATTCATGTTGTG
CAGAACCACCAGCACAGTGTATCCGGTGCACCTGGGAAATTTGTCATGTAGCTTAGAAGGAAATGCGTGGAAGAAGCTGGAGACGCGTTGTGACCTCCAAGA
TTTTCCATGCATTCGCATAATGATGGCAATGGGCCACGGGCGGCGGCTGGGCGAAGATATTTTGGGATCACTAACGTCATAGTTGTGTTCCAGGATGA
GATCGTCATAGGCCATTTTTACAAAAGCGCGGGCGGAGGTTGCCAGACTGCGGTATAATGGTTCATCCGGCCAGGGGCGTAGTTACCTCACAGATTTGCAT
TTCCACGCTTTGAGTTGAGTGGGGGATCATGTCTACCTGCGGGCGATGAAGAAAACGGTTTCCGGGGTAGGGGAGATCAGCTGGGAAGAAAGCAGGTTT
CTGAGCAGCTGCGACTTACCAGCAGCGGTGGGCCCTAAATCACACCTATTACCAGGTTGCAACTGGTAGTTAAGAGAGCTGCAGCTGCCGTATCCCTGAGCA
GGGGGGCCACTTCGTTAAGCATGTCCCTGACTCGCATGTTTTCCCTGACCAATCCGCCAGAAGGCGCTCGCCGCCAGCGATAGCAGTTCTTGAAGGAAGC
AAAGTTTTTCAACGGTTTGGACCGTCCGCGTAGGCATGCTTTTGGAGCGTTTGACCAAGCAGTTCCAGGCGGTCCCACAGCTCGGTACCTGCTCTACGGCA
TCTCGATCCAGCATATCTCCTCGTTTTGCGGGTTGGGGCGGCTTTCGCTGTACGGCAGTAGTCGGTCTCGTCCAGACGGGCCAGGGTCATGTCTTTCCACGG
GCGCAGGGTCTCGTACGCTAGTCTGGGTACGGTGAAGGGTGCCTCCGGGCTGCGCGTGGCCAGGGTGGCTTGGGCTGGTCTGCTGGTGTGAAG
CGCTGCCGGTCTCGCCCTGCGCGTGGCCAGGTAGCATTTGACCATGGTGTATAGTCCAGCCCTCCGCGGCGTGGCCCTTGGCGCGCAGCTTGCCTTGG
AGGAGGCGCCGACAGAGGGGCGAGTGCAGACTTTTGGGGCGTAGAGCTTGGGCGCGAGAAATACCGATTCCGGGGAGTAGGCATCCGCGCCGACGGCCCGCA
GACGGTCTCGCATCCACGACGAGGAGTTGAGCTTGGCGCTCCGGGTCAAAACAGGTTTCCCCCATGCTTTTTGATGCGTTTCTTACCTCGGTTTCCATG
AGCCGGTGTCCACGCTCGGTGACGAAAAGGCTGTCCGTGTCCCGTATACAGACTTGAGAGCCTGTCCCTGACCCGATGCGCTTGAAGCCTTCAACCCAGTC
AGCTCCTTCCGGTGGGCGGGGCGATGACTATCGTCCGCGCACTTATGACTGTCTTCTTTATCATGCAACTCGTAGGACAGGTGCCGGCAGCGCTCTGGGTCA
TTTTCCGGCAGGACCGCTTTCGCTGGAGCGGACGATGATCGCCCTGTGCGTGTGCGGTATTGGAATCTTGACGCGCTCGTCAAGCCTTGTGCTACTGGTCC
CGCCACCAACGTTTCCGGCAGAAAGCAGGCCATTATCGCCGGCATGGCGGGCAGCGCTGGGCTACGTTCTTGGTGGCTTCCGACGCGAGGCTGGATGGCC
TTCCTCATATGATTTCTCGCTTCCGGCGCATCGGGATGCCCGGTTGACGCGCATGCTGTCCAGGCGAGTAGATGACGACCATCAGGGACAGCTTCAAG
GCCAGCAAAAGGCCAGGAACCGTAAAAGGGCGGCTGTGCTGGCGTTTTTCCATAGGCTCCGCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAG
GTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTGGAAGCTCCCTCGTGCCTCTCCTGTCCGACCTGCCGCTTACCGGATACCTGTCC
GCCTTCTCCCTTCCGGAAGCGTGGCGCTTCTCATAGCTCACGCTGATAGGTATCTCAGTTCCGGTGTAGGTCGTTCCGCTCCAGCTGGGCTGTGTGACGAAC
CCCCGTTCCAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGGTCCAACCCGGTAAAGACAGCACTTATCGCCACTGGCAGCAGCCACTGGTAAACAG
GATTAGCAGAGCCGAGTATGAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCTTAACTACGGCTACACTAGAAGGACAGTATTTGGTATCTGCGCTGTGCTG
AAGCCAGTTTACCTTCCGAAAAGAGTTGGTAGCTCTTGTATCCGGCAAAACACCCGCTGGTAGCGGTTTGGTGGTTTGGTGGTGAAGCAGCAGATTACGCGCA
GAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTGGTTCATGAGATTATCAA
AAGGATCTTCCACTAGATCCTTTTAAATAAAAATGAAGTTTAAATCAATCTAAAGTATATATGAGTAAACTTGGTCTGACAGTTACCAATGCTTAAATCAGT
GAGGCACCTATCTCAGCGATCTGTCTATTTTCGTTTCCATAGTTGCTGACTCCCGTTCGTTAGATAACTACGATACGGGAGGGCTTACCATCTGGCCCA
GTGTCGAATGATACCGGAGACCCAGCTCACCGGCTCCAGATTTATCAGCAATAAACAGCCGCGGAGGGCGGAGCGCAGAAGTGGCTGCAACTTT
ATCCGCTCCATCCAGTCTATTAATTTGTTGCGGGGAGCTAGAGTAAAGTATGTTCCGCAAGTTAATAGTTTGGCAACGTTTGGCCATTGTGTCAGGCATCGTG
GTGTCAGCTCGTCTGTTGGTATGGCTTCAATTCAGTCCGGTTCACACGATCAAGGCGAGTTACATGATCCCCATGTTGTGCAAAAAAGCGGTTAGCTCCT
TCGGTCTCCGATCGTTGTGAGAAAGTGGCCCGAGTGTATCACTCATGTTTATGCGAGCAGTGCATAATCTCTTACTGTGATGCCATCCGTAAGATG
CTTTCTGTGACTGGTGAAGTACTCAACCAAGTCATTTCTGAGAATAGTGTATGCGGGCAGCGAGTTGCTTCTGCCGGCGTCAACACGGGATAATACCGCGCCA
CATAGCAGAACTTTAAAGTGCTCATCATTTGAAAACGTTCTTCCGGGCGAAAACCTCAAGGATCTTACCCTGTGAGATCCAGTTCCAGTGAATCAACCACTC
GTGCACCAACTGATCTTACGATCTTTTACTTTTCCACAGGCTTCTGGGTGAGCAAAAACAGGAAGGCAAAATGCCGCAAAAAGGGGAATAAGGGCGCACAG
GAAATGTTGAATACTCATACTCTTCTTTTCAATATATTTGAAGCATTTATCAGGTTTATGTCTCATGAGCGGATACATATTTGAATGTATTTAGAAAAAT
AAACAAATAGGGGTTCCGCGCACATTTCCCGAAAAGTGCCACCTGACGCTAAGAAACCATATTATCATGACATTAACCTATAAAAATAGCGGTATCACGA
GGCCCTTTCGTTCTCAAGAA

References

1. Feng, J. et al. (2015). SIRT6 suppresses glioma cell growth via induction of apoptosis, inhibition of oxidative stress and suppression of JAK2/STAT3 signaling pathway activation. *Oncol Rep.* doi:10.3892/or.2015.4477.
2. Bett AJ, Haddara W, Prevec L and Graham FL. (1994) *Proc Natl Acad Sci U S A.* **91**:8802-6 (Homologous recombination in packaging cell line)
3. He, T. C., Zhou, S., da Costa, L. T., Yu, J., Kinzler, K. W. et al. (1998) *Proc Natl Acad Sci U S A* **95**:2509-14. (pAdEasy System)
4. R D Anderson, R E Haskell, H Xia, B J Roessler and B L Davidson (2000) *Gene Ther.* **7**:1034-8. (RAPAd® System).

Recent Product Citations

1. Liu, Z. et al. (2020). Re-expression of DIRAS3 and p53 induces apoptosis and impaired autophagy in head and neck squamous cell carcinoma. *Mil Med Res.* **7**(1):48. doi: 10.1186/s40779-020-00275-3.
2. Park, J.E. et al. (2017). Hairy and enhancer of split 6 prevents hepatic lipid accumulation through inhibition of Pparg2 expression. *Hepatol Commun.* **1**(10):1085-1098. doi: 10.1002/hep4.1120.



3. Brunton, P. J. et al. (2015). 5α -reduced neurosteroids sex-dependently reverse central prenatal programming of neuroendocrine stress responses in rats. *J Neurosci.* **35**:666-677.
4. Lozić, M. et al. (2014). Overexpression of oxytocin receptors in the hypothalamic PVN increases baroreceptor reflex sensitivity and buffers BP variability in conscious rats. *Br J Pharmacol.* **171**:4385-4398.

Notice to Purchaser

This product is sold for research and development purposes only and is not to be incorporated into products for resale without written permission from Cell Biolabs. The patented RAPAd® technology is covered by a license from University of Iowa. By the use of this product you accept the terms and conditions of all applicable Limited Use Label Licenses. You may contact our Business Development department at busdev@cellbiolabs.com for information on sublicensing this technology.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS' sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2010-2020: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.