

STOP

Before using this product, please read the Limited Use License statement below:

Important Limited Use License information for pCpGfree-basic-Lucia

The purchase of the pCpGfree-basic-Lucia vector conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes.

The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes.

Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic, or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research.

If the purchaser is unwilling to accept the limitations of this limited use statement, InvivoGen is willing to accept return of the product with a full refund. The product must be returned in resaleable condition. For information on purchasing a license to this product for purposes other than research, contact InvivoGen, 10515 Vista Sorrento Parkway San Diego, CA 92121 USA. Tel: 858-457-5873 Fax: 858-457-5843.

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3-622-34-80

E-mail: info@invivogen.com



pCpGfree-basic-Lucia

A Lucia® reporter plasmid without a promoter and devoid of CpG dinucleotides

Catalog # pcpgf-baslC

For research use only

Version 21F04-MMv02

PRODUCT INFORMATION

Content:

- 20 µg of pCpGfree-basic-Lucia plasmid provided as lyophilized DNA
- *E. coli* GT115 strain provided lyophilized on a paper disk
- 1 ml of Zeocin™ (100 mg/ml)

Storage and Stability:

- Products are shipped at room temperature.
- Store lyophilized DNA at -20°C.
- Resuspended DNA is stable 6 months when stored at -20°C.
- Bacteria should be stored at -20°C and are stable up to 1 year.
- Store Zeocin™ at 4 °C or at -20 °C. The expiry date is specified on the product label.

Quality control:

Plasmid construct has been confirmed by restriction analysis and sequencing. Plasmid DNA was purified by ion exchange chromatography and lyophilized. Viability of the lyophilized bacteria upon resuspension has been verified.

GENERAL PRODUCT USE

Methylation of CpG dinucleotides within the promoter/enhancer region of genes is often associated with transcriptional silencing. This epigenetic event plays an important role in the regulation of gene activity in normal and cancer cells. Recently, it has been confirmed that the activity of enhancers is correlated with DNA methylation¹.

InvivoGen provides pCpGfree-basic-Lucia, a secreted luciferase reporter plasmid that is completely devoid of CpG dinucleotides and lacks the entire promoter region. It contains a multiple cloning site upstream of the Lucia luciferase reporter gene. Expression of Lucia luciferase in cells transfected with this plasmid depends on the insertion of a functional promoter or enhancer/promoter cassette upstream from the Lucia luciferase gene. Thus, pCpGfree-basic-Lucia allows to study the effect of CpG methylation on a promoter, alone or combined with enhancer elements.

PLASMID FEATURES

All the elements required for replication and selection of the plasmid in *E. coli* and gene expression in mammalian cells are completely devoid of CpG dinucleotides. Furthermore, all Dam methylation sites (GATC) have been removed to prevent prokaryotic methylation.

Elements for expression in *E. coli*

- Origin of replication: The *E. coli* R6K gamma ori has been modified to remove all CpGs. This origin is activated by the R6K specific initiator protein π , encoded by the *pir* gene².
- Bacterial promoter: EM2K is a CpG-free version of the bacterial EM7 promoter.
- Selectable marker: The Zeocin™ resistance gene is a small gene (<400 bp) that contains numerous CpG dinucleotides. A synthetic new allele was created that contains no CpGs.

Elements for expression in mammalian cells

- Lucia luciferase is a synthetic CpG-free gene that codes for a secreted coelenterazine-utilizing luciferase.
ORF size (from the ATG to the stop codon): 634 bp
- Polyadenylation signal: The polyadenylation signal is a CpG-free form of the late SV40 polyadenylation signal.

- MAR: Matrix attached regions (MARs) are sequences typically AT-rich that are able to form barriers between independently regulated domains³. pCpGfree plasmids contains two MARs, from the 5' region of the human IFN- β gene or β -globin gene that were chosen because they are naturally CpG-free. The MARs are placed between the bacterial and mammalian transcription units.
- MCS: The multiple cloning site contains several commonly used restriction sites for convenient cloning of a gene of interest.
5' Sda I, Bsp 120I, Avr II, Nsi I, Ppu 10I, Sca I, Bam HI, Spe I, Hind III 3'

Due to the presence of the R6K γ origin of replication, pCpG plasmids can only be amplified in *E. coli* mutant strain expressing a *pir* mutant gene. They will not replicate in standard *E. coli* strains. Therefore, pCpG plasmids are provided with the *E. coli* GT115 strain, a *pir* mutant also deficient in *Dcm* methylation.

1. Hoivik EA. et al., 2011. DNA Methylation of Intronic Enhancers Directs Tissue-Specific Expression of Steroidogenic Factor 1/Adrenal 4 Binding Protein (SF-1/Ad4BP). *Endocrinology*. 152(5):2100-12. 22. 2. Wu F. et al. 1995. A DNA segment conferring stable maintenance on R6K gamma-origin core replicons. *J Bacteriol*. 177(22):6338-45. 3. Bode J. et al., 1996. Scaffold/matrix-attached regions: topological switches with multiple regulatory functions. *Crit Rev Eukaryot Gene Expr*. 6(2-3):115-38.

METHODS

Plasmid resuspension

Quickly spin the tube containing the lyophilized plasmid to pellet the DNA. To obtain a plasmid solution at 1 µg/µl, resuspend the DNA in 20 µl of sterile H₂O. Store resuspended plasmid at -20 °C.

Reconstitution of *E. coli* GT115 strain

Use sterile conditions to do the following:

1. Reconstitute *E. coli* GT115 by adding 1 ml of Luria-Bertani (LB) medium in the tube containing the paper disk. Let sit for 5 minutes.
2. Mix gently by vortexing for 1-2 minutes.
3. Streak bacteria taken from this suspension on a LB agar plate.
4. Place the plate in an incubator at 37°C overnight.
5. Isolate a single colony and grow the bacteria in LB or terrific broth (TB) medium.
6. Prepare competent cells utilizing protocol of choice.

Plasmid amplification and cloning

Plasmid amplification and cloning can be performed in *E. coli* GT115.

Zeocin™ usage

This antibiotic can be used for *E. coli* at 25 µg/ml in liquid or solid media.

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Hong Kong: +852 3-622-34-80
E-mail: info@invivogen.com



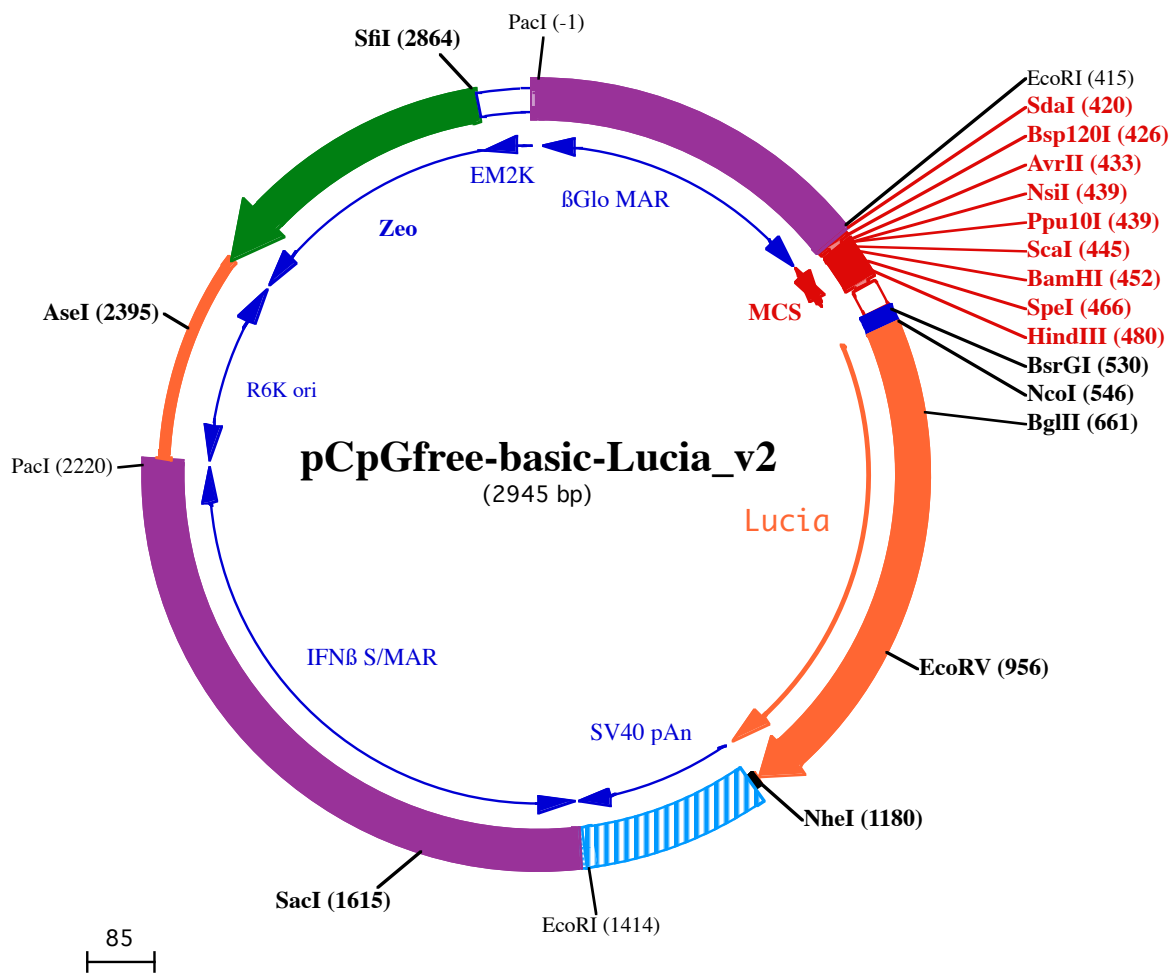
RELATED PRODUCTS

Product	Catalog Code
QUANTI-Luc™	rep-qlc1
ChemiComp GT115	gt115-11
pCpGfree-promoter-Lucia	pcpgf-promlc
pCpGfree-promoter (mSEAP)	pcpgf-prom
pCpGfree-basic (mSEAP)	pcpgf-bas

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Hong Kong: +852 3-622-34-80
E-mail: info@invivogen.com





PacI (-1)

1 TTAATTA~~AAAA~~TATCTCTAAGGCATGTGAACGGCTGTCTGGTTTTATCTGACTTCATCTGCTACCTCTGTGACCTGAAACATATTTATAATCCAT

101 TAAGCTGTGCATATGATAGATTTATCATATGATATTTTCTTAAAGGATTTTTGAAGAACTAATTGAATTGATACCTGTAAAGCTTTATCACACTACCC

201 AATAAATAATAAATCTCTTTGTTGAGCTCTCTGTTTCTATAAATATGTACCAGTTTTATTGTTTTAGTGGTAGTGATTTTATTCTCTTTCTATATATAT

301 ACACACACATGTGTGCATTATAAATATATACAATTTTTATGAATAAAAAATTATTAGCAATCAATATTGAAAACCACTGATTTTTGTTTATGTGAGCAA

Bsp120I (426) Ppu10I (439)
SdaI (420) NsiI (439) BamHI (452)
EcoRI (415) AvrII (433) ScaI (445) SpeI (466) HindIII (480)

401 ACAGCAGATTA~~AAAGGAAT~~CCTGCAGGGCCACCTAGGATGCATAGTACTAGGATCCAACATGTAAGCTAGTAGCATGCAAAGCTTAGAAAttgtactaac

501 cttcttctctttctctcctgacagGTTGGTGTACAGTAGCTCCACCATGGAATCAAGGTGCTGTTTGCCCTCATCTGTATTGCTGTTGCTGAGGCAA
 1 M E I K V L F A L I C I A V A E A

BsrGI (530) NcoI (546)
BglII (661)

601 AACCCACTGAAATCAATGAAGACCTCAATATAGCTGCTGTGGCCTCCAACCTTGGCCACCACAGATCTTGAGACTGACCTGTTCCACCACTGGGAGACCAT
 18 K P T E I N E D L N I A A V A S N F A T T D L E T D L F T N W E T M
 701 GAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACTCCCCAGATGTCCTGAGGGAGCTGGAG
 51 N V I S T D T E Q V N T D A D R G K L P G K K L P P D V L R E L E
 801 GCCAATGCCAGAAGGGCTGGTTGCACAAGAGGCTGCTCATTTGCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCC
 85 A N A R R A G C T R G C L I C L S H I K C T P K M K K F I P G R C

EcoRV (956)

901 ACACCTATGAAGGTGAAAGGAGTCTGCTCAGGGAGGATTGGAGAGCAATTTGATATCCCAGAGATTCCTGGCTCAAGGATAAGGAGCCACTGGA
 118 H T Y E G E K E S A Q G G I G E A I V D I P E I P G F K D K E P L D
 1001 CCAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCAACACTGGCTGTCTGAAGGGCCTTCCAATGTCCAGTGCCTGACCTCCTGAAGAAATTTATCCCTGGCAGGTGCC
 151 Q F I A Q V D L C A D C T T G C L K G L A N V Q C S D L L K K W L

NheI (1180)

1101 CCCAGAGGTGTACCACCTTTTGGCAGCAAGATTCAGGGTAGGGTGACAAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGCTGGCCAGACATGAT
 185 P Q R C T T F A S K I Q G R V D K I K G L A G D R •
 1201 AAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAGAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAAACCAT

1301 ATAAGCTCAATAAACAAGTTAACAACAACAATGCATTCATTTATGTTTCAGGTTGAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAACCTCT

EcoRI (1414)

1401 ACAAATGTGGTATGGAATCAGTCAATATGTTACCCCAAAAAAGCTGTTTGTAACTGCCAACCTCATTCTAAAATGTATATAGAAGCCAAAAAGACA

1501 ATAACAAAAATATTCTGTAGAACAAAATGGGAAAGAATGTTCCACTAAATATCAAGATTTAGAGCAAAAGCATGAGATGTGTGGGATAGACAGTGAAGCC

SacI (1615)

1601 TGATAAAATAGAGTAGAGCTCAGAAACAGACCCATTGATATATGTAAGTGACCTATGAAAAAATATGGCATTTTACAATGGGAAAATGATGGTCTTTTT

1701 CTTTTTAGAAAAACAGGAAATATATTTATATGTAAAAAATAAAGGGAACCCATATGTCATACCATACACAAAAAATCCAGTGAATTATAAGTC

1801 TAAATGGAGAAGGCAAACTTTAAATCTTTAGAAAATAATAGAAGCATGCCATCAAGACTTCAGTGTAGAGAAAAATTTCTATGACTCAAGTCTCT

1901 AACCAAAAGAAAAGATTGTTAATTAGATTGCATGAATATTAAGACTATTTTTAAATTAATAAACCATTAAGAAAAGTCAAGCCATAGAAATGACAGAA

2001 AATATTTGCAACACCCAGTAAAGAGAATTGTAATATGCAGATTATAAAAAAGAGTCTTACAATCAGTAAAAAATAAACTAGACAAAAATTTGAACAG

2101 ATGAAAGAGAACTCTAATAATCATTACACATGAGAACTCAATCTCAGAAATCAGAGAACTATCATTGCATATACACTAAATAGAGAAATATTAATA

PacI (2220)

2201 GGCTAAGTAACATCTGTGGCTTAATTA~~AAATCAGCAGTTCAACCTGTTGATAGTATGTAAGCTCTCATGTTAATGTAAGCTCTCATGTTAAT~~

AseI (2395)

2301 GAACTAAACCCTCATGGCTAATGTAAGCTCTCATGGCTAATGTAAGCTCTCATGTTTATGTAAGCTCTCATGTTTGAACATAAAATTA

2401 TATAAATCAGCAACTTAAATAGCCTCTAAGGTTTTAAGTTTTATAAGAAAAAAGAATATATAAGGCTTTTAAAGGTTTTAAGGTTTCTAGCTTTAGT
 125 • D

2501 CCTGTTCTCAGTACAAAATGGACACAATTTCCAGCAGGGTCTCTGAGGGCAAATCCCTTCCCAAGGTTGTTACCAATTTCTGTCATGGCTGGGCC
 123 Q E E A V F H V C N G A P D R L A F E R G W P Q E G I E T M A P G
 2601 AGAGGCATCCCTGAAATTTGTGCTGACTACTTCTGACCACTCTGCATAAAGCTCATCTAGGCCTCTGACCCAGACCCAAGCAAGGGTGTGTGAGGACCA
 90 S A D R F N T S V V E S W E A Y L E D L G R V W V W A L T N D P V
 2701 ACTTGGTCTGAATGCTGAGATGAAGAGGGTGACATCATCTGCAACACCAAGCAAAATCATCTTCAACAAAGTCTCTGGAGAAATCCTAATCTGTGACG
 56 V Q D Q V A S I F L T V D D R V V G A F D D E V F D R S F G L R D T

SfiI (2864)

2801 TCCAGAACTCTACAGCCCTGCAACATCCCTTGTGAGGACTGGGACTGCAGAAGTGAAGTTTGGCCATGATGGCCCTCCTATAGTGAAGTTGATTATA
 23 W F E V A G A V D R A T L V P V A S T L K A M
 2901 CTATGCAGATATACTATGCCAATGTTAATTGTCAACTACCTGTT