

STOP

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

Important Limited Use License information for pSELECT-NLucia-zeo

The purchase of the pSELECT-NLucia-zeo 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 3622-3480

E-mail: info@invivogen.com



pSELECT-NLucia-zeo

Plasmid for the expression of Lucia-N-terminal tagged proteins

Catalog # psetz-nlucia

For research use only

Version 20L01-MM

PRODUCT INFORMATION

Content:

- 20 µg of pSELECT-NLucia-zeo plasmid provided as lyophilized DNA
- 1 ml of Zeocin™ (100 mg/ml)

Storage and Stability:

Product is shipped at room temperature. Lyophilized DNA should be resuspended upon receipt and stored at -20°C. Lyophilized DNA is stable for 3 months at -20°C. Resuspended DNA is stable more than one year at -20°C.

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.

GENERAL PRODUCT USE

pSELECT plasmids are specifically designed for strong and constitutive expression of a gene of interest in a wide variety of cell lines. They allow the selection of stable transfectants and offer a variety of selectable markers. pSELECT plasmids contain two expression cassettes: the first drives the expression of the gene of interest and the second drives the expression of a large choice of dominant selectable markers for both *E. coli* and mammalian cells. They are both terminating with a strong polyadenylation signal (polyA) that separates the two expression cassettes thus preventing any transcription interference. The late SV40 polyA terminates the transcription of the gene of interest while the human β-globin polyA terminates the transcription of the selectable marker.

pSELECT Lucia-Tag is a family of expression plasmids designed to generate tagged proteins in mammalian cells with Lucia luciferase in order to detect and quantify the tagged protein of interest by bioluminescence assay. Lucia luciferase is a novel secreted luciferase with strong bioluminescent activity. The luciferase activity of Lucia luciferase can be detected and quantified directly in the culture medium of transfected cells using InvivoGen's QUANTI-Luc™ detection reagent.

N-terminal tag: the tag encompasses the Start codon and is followed by a multiple cloning site (MCS).

PLASMID FEATURES

First expression cassette

• **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1a (EF-1α) core promoter¹ and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat². The EF-1α promoter exhibits a strong activity and yields long lasting expression of a transgene *in vivo*. The R-U5' has been coupled to the EF-1α core promoter to enhance stability of RNA.

• **MCS:** The multiple cloning site contains the following restriction sites: 5' - BamH I, Eco47III, Nco I, Nhe I - 3'

Each restriction site is compatible with many other enzymes.

• **SV40 pAn:** the Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA³.

• **ori:** a minimal *E. coli* origin of replication to limit vector size, but with the same activity as the longer Ori.

Second expression cassette

• **CMV enh/prom:** The human cytomegalovirus immediate-early gene 1 promoter/enhancer was originally isolated from the Towne strain and was found to be stronger than any other viral promoters.

• **EM7** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*.

• **Sh ble gene** confers Zeocin™ resistance. The *Sh ble* gene is driven by the CMV enhancer/promoter in tandem with the bacterial EM7 promoter allowing selection in both mammalian cells and *E. coli*.

• **BGlo pAn:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription⁴.

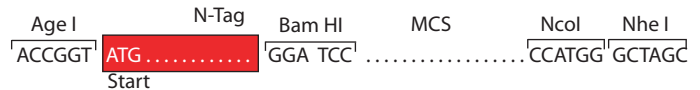
CLONING STRATEGY

For expression of a tagged protein, it is important to ensure to clone your gene-of-interest into the correct reading frame. In general it is recommended to use Bam HI/Nhe I restriction site combination for cloning into plasmids with the N-Tag. For the plasmids with N-Tag, check whether your gene of interest and the start codon of the N-Tag are in the correct reading frame.

Note: The Bam HI restriction site is compatible with Bgl II.



If it is not possible to use the Nhe I restriction site, it is possible to use another restriction site such as Nco I.



Alternatively, blunt end cloning can be achieved using the Eco 47 III site, which is downstream of the Bam HI site. Blunt end PCR fragments can be obtained using T4 DNA polymerase or Klenow.



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 3622-3480

E-mail: info@invivogen.com

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.

Plasmid amplification and cloning

Plasmid amplification and cloning can be performed in *E. coli* GT116 or in other commonly used laboratory *E. coli* strains, such as DH5α.

Zeocin™ usage

This antibiotic can be used for *E. coli* at 25 µg/ml in liquid or solid media and at 50-200 µg/ml to select Zeocin™-resistant mammalian cells.

References:

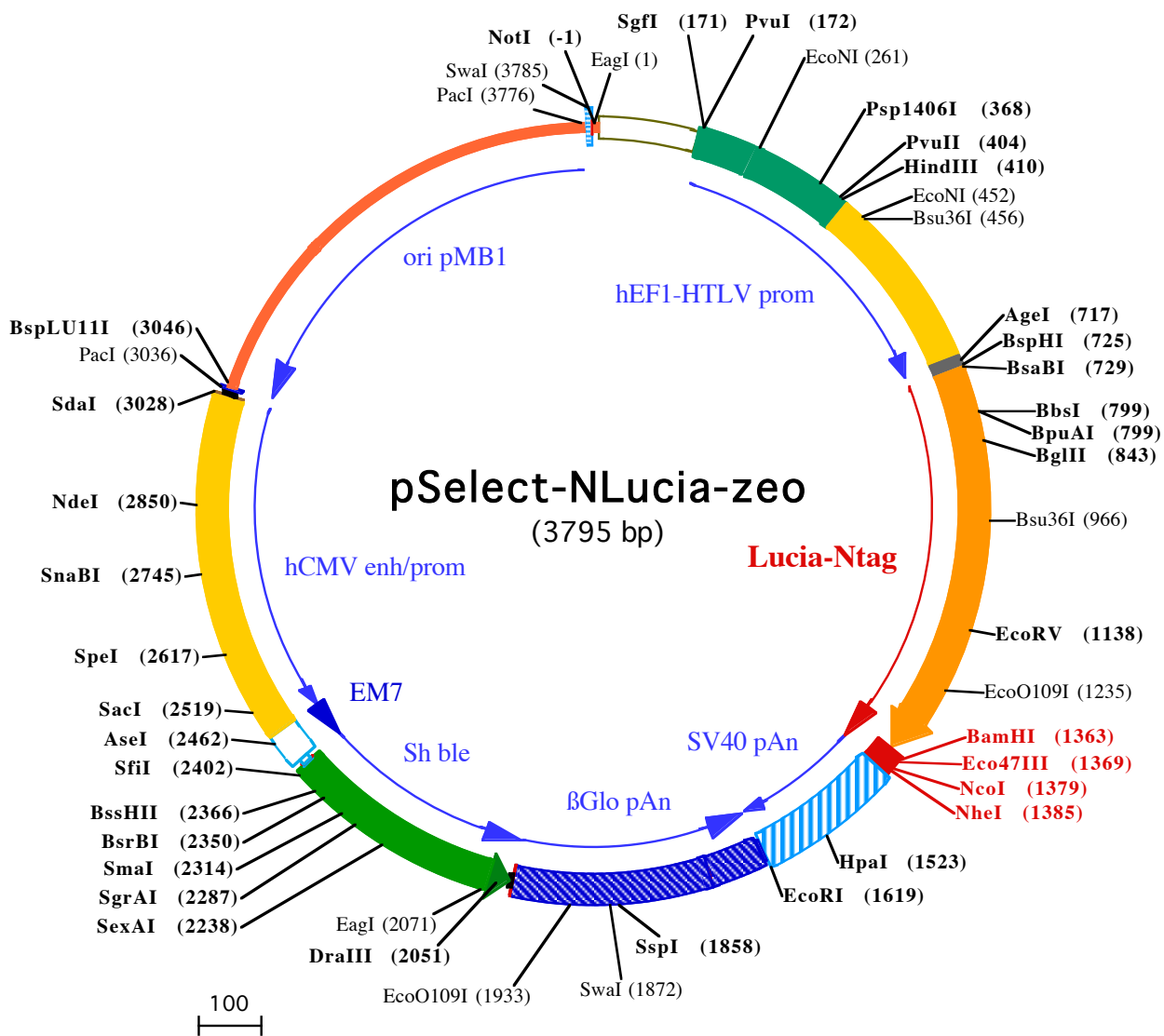
1. Kim, D.W. *et al.* (1990). *Gene* 2: 217-223.
2. Takebe, Y. *et al.* (1988). *Mol. Cell Biol.* 1: 466-472.
3. Carswell, S., and Alwine, J.C. (1989). *Mol. Cell Biol.* 10: 4248-4258.
4. Yu J & Russell JE. (2001). *Mol Cell Biol*, 21(17):5879-88.

RELATED PRODUCTS

Product	Catalog Code
pSELECT-NLucia-blasti	psetb-nlucia
QUANTI-Luc™	rep-qlc1
ChemiComp GT116	gt116-11

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 3622-3480
E-mail: info@invivogen.com



EagI (1)
NotI (-1)
1 CCGGCCCAATAAAATATCTTTATTTTATTACATCTGTGTGTTGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACA
PvuI (172)
SgfI (171)
101 AAACAACTAGCAAATAGGCTGTCCCCAGTGAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGATCTGCGATCGCTCCGGTGCCCGTCAGTGGCA
EcoNI (261)
201 GAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTCCGCAATTGAACGGTGCCTAGAGAAGTGGCGGGGTAACCTGGGAAAGTGATG
Psp1406I (368)
301 TCGTGTAAGTGGCTCCGCTTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCCGTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAG
HindIII (410) Bsu36I (456)
PvuII (404) EcoNI (452)
401 AACACAGCTGAAGCTTCAGGGGCTCGCATCTCTCCTTACCGCCCGCCCTACCTGAGGCCCATCCACGCCGGTTGAGTCGCTTCTGCCGCT
501 CCGCCTGTGGTGCCTCCTGAACGCTCCGCCGCTAGGTAAGTTAAAGTCCAGTGCAGACCGGGCCTTTGTCGGCGCTCCCTGGAGCTACCTA
601 GACTCAGCCGGCTCTCCACGCTTGGCTGACCCTGCTCAACTCTACGTCTTTGTTTCGTTTTCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACC
BspHI (725) BpuAI
AgeI (717) BsaBI (729) BbsI
701 GCGCCTACCTGAGATCACCGGTATCATGATGGAATCAAGGTGCTGTTGGCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCACTGAAATCAATG
1 M M E I K V L F A L I C I A V A E A K P T E I N
BgIII (843)
801 AAGACCTCAATATAGCTGTGTGGCCTCCAACCTTTGCCACCACAGATCTTGAGACTGACCTGTTCACTGGGAGACCATGAATGTGATTAGCACTGA
25 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 N V I S T D
Bsu36I (966)
901 CACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACTCCCCCAGATGCTCGAGGGAGCTGGAGGCAATGCCAGAAGGGCT
58 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 A N A R R A
1001 GGTTCACAAGAGGCTGCCTCATTGCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAA
92 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 H T Y E G E
EcoRV (1138)
1101 AGGAGTCTGCTCAGGGAGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCTGGCTCAAGGATAAGGAGCCACTGGACCAGTTTATTGCTCAAGT
125 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 Q F I A Q V
EcoO109I (1235)
1201 GGACCTCTGTGCTGATTGCACCCTGGCTGTGAAGGGCCTTCCAATGTCCAGTGTCTGACCTCTGAAGAAGTGGCTTCCCAGAGGTGTACCACT
158 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 P Q R C T T
Eco47III (1369) NheI (1385)
BamHI (1363) NcoI (1379)
1301 TTTGCCAGCAAGATTGAGGGTGGGACAAATCAAGGGTCTGGCTGGGACAGAGGAGGTGGATCCAGCGCTGCAGCCATGGGCTAGCTGGCCAGAC
192 F A S K I Q G R V D K I K G L A G D R G G S S A A A M G
1401 ATGATAAGATACATTGATGAGTTTGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTTGTGAATTTGTGATGCTATTGCTTTATTTGTAA
HpaI (1523)
1501 CCATTATAAGCTGCAATAAACAAGTAAACAACAACAAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAA
EcoRI (1619)
1601 CCTCTACAAATGTGGTATGGAATTCATAAATACAGCATAGCAAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGG
1701 CATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCCAAGTTTGAAC
SspI (1858) SmaI (1872)
1801 AGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTATGAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATA
EcoO109I (1933)
1901 AATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAAT
DraIII (2051) EagI (2071)
2001 TGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCCTGCTCCTCTGCCACAAGTGCACGAGTTGCCGGCGGGTCCGCGAGGGCGAACTCCCGC
127 G • D Q E E A V F H V C N G A P D R L A F E R
2101 CCCCACGGCTGCTCGCCGATCTCGGTGATGGCCGGCCCGAGGCGTCCCGAAAGTTCGTGGACACGACCTCCGACCCTCGCGGTACAGCTCGTCCAGG
102 G W P Q E G I E T M A P G S A D R F N T S V V E S W E A Y L E D L G
SexAI (2238) SgrAI (2287)
2201 CGCGACCCACACCAGGCCAGGGTGTGTCGGCACCACTGGTCTGGACCGCGTGTGAACAGGGTACGTCGTCGCGGACCACCCGCGAAGTC
69 R V W V W A L T N D P V V Q D Q V A S I F L T V D D R V V G A F D
SmaI (2314) BsrBI (2350) BssHIII (2366)
2301 GTCCTCCACGAAGTCCCGGAGAACCAGCCGGTCCGACTCGACCGCTCCGGCAGCTCGCGCGGTGAGCACCAGGACGGCACTGGTCAAC
36 D E V F D R S F G L R D T W F E V A G A V D R A T L V P V A S T L
SfiI (2402) AseI (2462)
2401 TTGGCCATGATGGCCCTCTATAGTGAGTCGATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTT
2 K A M
SacI (2519)
2501 ATCTGACGGTTCACATAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTG

2601 **SpeI (2617)**
GAAAGTCCC GTT GATTTAC TAGTCAA AACAAACTCCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCCGTGAGTCAAACCGCTATCCACGCCATT

2701 **SnaBI (2745)**
GATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGGCATAA

2801 **NdeI (2850)**
TGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGCAGTTTACCGTAAATACT

2901
CCACCCATTGACGTCAATGGAAAGTCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTCAGCCAGG

3001 **SdaI (3028)** **PacI (3036)** **BspLU11I (3046)**
CGGGCCATTTACCGTAAGTTATGTAACGCCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAGGCCGCGTTGC

3101
TGGCGTTTTTCCATAGGCTCCGCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGC

3201
GTTTCCCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCTGCCGCTTACCGGATACCTGTCCGCCTTCTCCCTTCGGAAGCGTGGCGCTTCT

3301
CATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCGCTGCGCTTAT

3401
CCGGTAACTATCGTCTT GAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGG

3501
TGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGA

3601
GTTGGTAGCTCTTGATCCGGCAAACAACACCGCTGGTAGCGGTGGTTTTTTTTGTTTGCAAGCAGCAGATTACGGCAGAAAAAAGGATCTCAAGAAG

3701 **PacI (3776)** **SwaI (3785)**
ATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA