

# STOP

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

## **Important Limited Use License information for pSELECT-NLucia-blasti**

The purchase of the pSELECT-NLucia-blasti 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](mailto:info@invivogen.com)



# pSELECT-NLucia-blasti

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

Catalog code: psetb-nlucia

<https://www.invivogen.com/pselect-lucia-tag>

For research use only

Version 20117-MM

## PRODUCT INFORMATION

### Contents

- 20 µg of pSELECT-NLucia-blasti plasmid provided as lyophilized DNA
- 2 x 1 ml blasticidin at 10 mg/ml

### Storage and stability

- Product is shipped at room temperature.
- Upon receipt, store lyophilized DNA at -20°C.
- Resuspended DNA should be stored at -20°C.
- Store blasticidin at 4°C or -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 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-1α (EF-1α) core promoter<sup>1</sup> and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat<sup>2</sup>. 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' - BamHI, Eco47III, NcoI, NheI - 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 mRNA<sup>3</sup>.
- **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*.
- **Bsr:** Resistance to Blasticidin is conferred by the *bsr* gene from *Bacillus cereus*. The *bsr* gene is driven by the CMV enhancer/promoter in tandem with the bacterial EM7 promoter allowing selection in both mammalian cells and *E. coli*.
- **βGlo pAn:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription<sup>4</sup>.

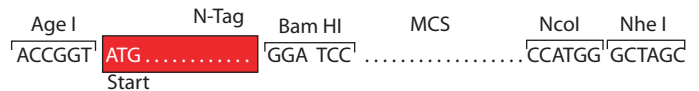
## CLONING STRATEGY

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 BamHI/NheI 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 BamHI restriction site is compatible with BglII.



If it is not possible to use the NheI restriction site, it is possible to use another restriction site such as NcoI.



Alternatively, blunt end cloning can be achieved using the Eco47III site, which is downstream of the BamHI 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](mailto: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 H<sub>2</sub>O. Store resuspended plasmid at -20°C.

### Plasmid amplification and cloning

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

### Blasticidin usage

Blasticidin should be used at 25-100 µg/ml in bacteria and 1-30 µg/ml in mammalian cells. Blasticidin is supplied at 10 mg/ml in HEPES buffer.

### References:

1. Kim, D.W. et al., 1990. Use of the human elongation factor 1 alpha promoter as a versatile and efficient expression system. *Gene* 2: 217-223. 2. Takebe, Y. et al., 1988. SR alpha promoter: an efficient and versatile mammalian cDNA expression system composed of the simian virus 40 early promoter and the R-U5 segment of human T-cell leukemia virus type 1 long terminal repeat. *Mol. Cell Biol.* 1: 466-472. 3. Carswell, S., & Alwine, J.C., 1989. Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. *Mol. Cell Biol.* 10: 4248-4258. 4. Yu J & Russell JE., 2001. Structural and functional analysis of an mRNP complex that mediates the high stability of human beta-globin mRNA. *Mol Cell Biol.* 21(17):5879-88.

## RELATED PRODUCTS

Product	Catalog Code
ChemiComp GT116	gt116-11
QUANTI-Luc™	rep-qlc1
QUANTI-Luc™ Gold	rep-qlcg1

---

### TECHNICAL SUPPORT

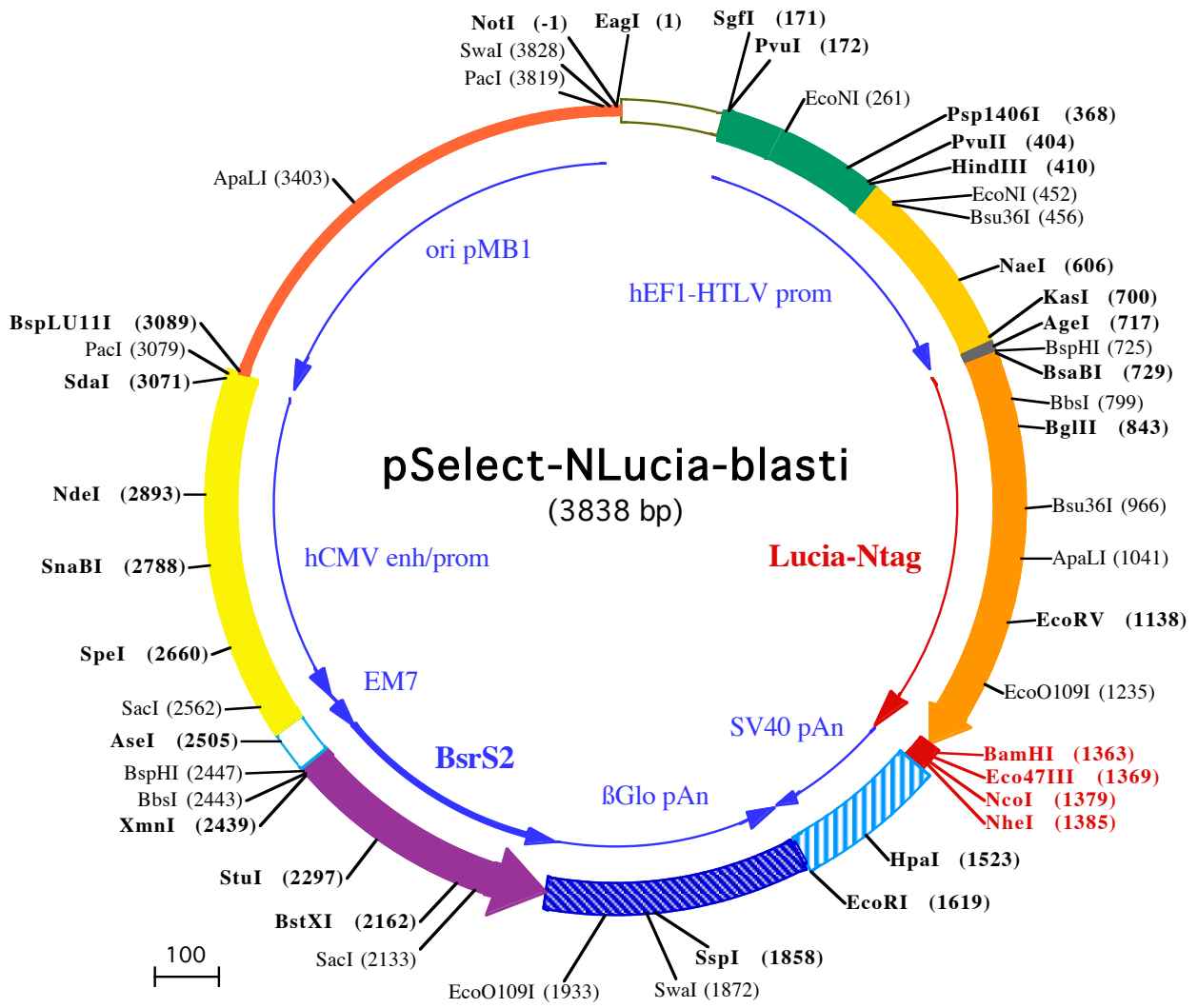
InvivoGen USA (Toll-Free): 888-457-5873

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](mailto:info@invivogen.com)



EagI (1)  
NotI (-1)  
1 GCGGCCGCAATAAAATATCTTTATTTTTCATTACATCTGTGTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACA

PvuI (172)  
SgfI (171)  
101 AAACAACTAGCAAAATAGGCTGTCCCGAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGATCTGCGATCGCTCCGGTCCCGTCAGTGGCA

EcoNI (261)  
201 GAGCGCACATCGCCACAGTCCCGAGAAGTTGGGGGAGGGTCCGCAATTGAACGGTGCCTAGAGAAGTGGCGGGGTAACGGGAAAGTGTATG

Psp1406I (368)  
301 TCGTGTACTGGTCCGCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCCGCTGAACGTTCTTTTTCCGCAACGGGTTTCCGCCAG

HindIII (410)  
PvuII (404)  
401 AACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACCGCGCCCGCCCTACCTGAGGCCGCCATCCACGCCGTTGAGTCGCGTTCTGCCGCCCT

Bsu36I (456)  
EcoNI (452)  
501 CCCCCTGTGGTCCCTCTGAACTGCGTCCGCCGTCTAGTAAAGTTAAAGTCAAGTGCAGACCGGGCCTTTGTCGGCGCTCCCTTGAGGCTACCTA

NaeI (606)  
601 GACTCAGCCGCTCTCCACGCTTTGCCTGACCCTGCTTCTCAACTCTACGCTTTTGTTCGTTTTCTGTTCTGCCCGTTACAGATCCAAGCTGTGACC

KasI (700)  
AgeI (717)  
BspHI (725)  
BsaBI (729)  
701 GGGCCTACCTGAGATCACCGGTTCATCATGATGGAAATCAAGGTGCTGTTTGCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCCACTGAAATCAATG

BbsI\*(799)  
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

BglIII (843)  
801 AAGACCTCAATATAGCTGCTGTGGCCTCAACTTTGCCACCACAGATCTTGAGACTGACCTGTTACCAACTGGGAGACCATGAATGTGATTAGCACTGA

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 CACAGAGCAGGTGAACACAGATGTGACAGGGGAAGTGCCTGGCAAAAACCTCCCGCAGATGCTCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGCT

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

ApaLI (1041)  
1001 GTTGCAACAAGAGGCTGCTCATTGCTCTCCACATTAAGTGCACCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAA

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 AGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCTGGCTCAAGGATAAGGAGCCACTGGACCAGTTTATTGCTCAAGT

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 GGACCTCTGTGCTGATTGCCACTGGCTGTCTGAAGGGCCTTGCCAAATGTCAGTGTCTGACCTCCTGAAGAAGTGGCTTCCCAGAGGTGTACCCT

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 TTTGCCAGCAAGATTGAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGAGGGTGGATCCAGCGCTGCAGCCATGGGCTAGCTGGCCAGAC

192 F A S K I Q G R V D K I K G L A G D R G G G S S A A A M G •

1401 ATGATAAGATACATTGATGAGTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTGTAA

HpaI (1523)  
1501 CCATTATAAGCTGCAATAAACAAGTTAACAAACAATTCGATTCATTTTATGTTTCAGGTTCCAGGGGAGGTGGGAGTTTTTTAAAGCAAGTAAAA

EcoRI (1619)  
1601 CCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGG

1701 CATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCCAAGGTTTGAAC

SspI (1858)  
SwaI (1872)  
1801 AGCTTTCATTTCTTTATGTTTTAAATGCACTGACCTCCCACATTCCTTTTTAGTAAATATTCAGAAATATTTAAATACATCATTGCAATGAAAATA

EcoO109I (1933)  
1901 AATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAAT

2001 TGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTATCTCAAT

141 • N R T Y K L P I L E E I T T K V L K G N M E I

SacI (2133)  
BstXI (2162)  
2101 GAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGG

117 L V F C D P A Y D S I L E R C M G C P S V V R I S R D V E D S Y P

StuI (2297)  
2201 TGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGCTCACAGCAGACCCAATGGCAATGGCTTCAGCAGACAGTACCCTGCCAATGTAGG

83 H R V A V I T D F D K Q G N S V A S G I A I A E A C V T V R G I Y A

2301 CCTCAATGTGGACAGCAGAGATGATCTCCAGTCTGGTCTGATGGCCGCCGACATGGTCTGTTGCTCCTCATAGAGCATGGTATCTTCTCAGT

50 E I H V A S I I E G T K T R I A A G V H H K N D E Y L M T I K E T

BspHI (2447)  
BbsI (2443)  
XmnI (2439)  
2401 GCGACCTCCACAGCTCCAGATCCTGCTGAGAGATGTTGAAGTCTTCATGATGGCCCTCCTATAGTGAAGTCTGATTATACTATGCCGATATACTATGC

17 A V E V L E L D Q Q S I N F T K M

AseI (2505)  
SacI (2562)  
2501 CGATGATTAATTGTCAAACACGCTGGATGGCGTCTCCAGCTTATCTGACGGTTCCTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCTT

SpeI (2660)  
2600 ACCGCCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGAAAGTCCCGTTGATTTACTAGTCAAAACAAACTCCCATTGACGTCAATGGGGT

SnaBI (2788)  
2699 GAGACTTGAAATCCCGTGAAGTCAAACCGCTATCCACGCCATTGATGACTGCCAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATG

NdeI (285)  
2799 TACTGCCAAGTAGGAAAGTCCATAAGGTGATGACTGGGCATAATGCCAGGCGGGCCATTACCCTGATTGACGTCAATAGGGGCGTACTTGGCATAT

2899 GATACACTTGTACTGCCAAGTGGGCGAGTTTACCCTAAATACTCCACCCATTGACGTCAATGAAAGTCCCTATTGGCGTTACTATGGGAACATACGT

SdaI (3071)  
PacI (3079)  
BspLU11I (3089)  
2999 CATTATTGACGTCAATGGGCGGGGCTGTTGGGCGGTACGCCAGCGGGCCATTTACCCTAAGTTATGTAACCGCTGCAGGTTAATAAGAACAATGTG

3097 AGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGC

3197 TCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAGATACCAGGCGTTTCCCCCTGGAAGCTCCCTCGTGCGCTCTCCTGTTCCGACCCTGCCGTTA

3297 CCGGATACCTGTCCGCCTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTTCGCTCCAAGCT

3397 GGGCTGTGTGCACGAACCCCCGTTACGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTG

3497 GCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAG

3597 TATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGATTGGTAGCTCTTGATCCGGCAAACAACCACCGCTGGTAGCGGTGTTTTTT

3697 TGTTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACCTCACGT

3797 TAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC A

    ApaLI (3403)  
    PacI (3819)  Swal (3828)