

pSELECT-zeo-HSV1-tk::Sh

An expression plasmid coding for a CpG-free cytotoxic/resistance fusion gene

Catalog code: psetz-hsv1tksh

For research use only

Version 20L01-MM

PRODUCT INFORMATION

Contents:

- 20 µg of pSELECT-zeo-HSV1-tk::Sh provided as lyophilized DNA
- 1 ml of Zeocin™ (100 mg/ml)

Storage and stability:

- Product is shipped at room temperature.
- Lyophilized DNA should be stored at -20 °C.
- Resuspended DNA should be stored at -20 °C and is stable for 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 full-length ORF sequencing.
- Plasmid DNA was purified by ion exchange chromatography.

GENERAL PRODUCT USE

pSelect-zeo plasmids contain genes that have been chemically synthesized. The DNA sequence of these genes was modified by optimizing the codon usage, reducing or eliminating the CpG motifs and avoiding secondary DNA structures without changing the amino acid sequence of the wild type proteins.

pSELECT-zeo plasmids may be used:

To subclone the synthetic gene into another vector. To facilitate subcloning, the HSV1-tk::Sh gene is flanked by two unique restriction sites: Nco I at the 5' end that encompasses the Start codon, and Nhe I at the 3' end.

As a gene reporter plasmid. pSELECT-zeo is a mammalian expression plasmid selectable in *E. coli* and mammalian cells with Zeocin™, as the *Sh ble* gene in the second expression cassette is driven by the eukaryote CMV enhancer/promoter in tandem with the bacterial EM7 promoter.

PLASMID FEATURES

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

First expression cassette

- hEF1-HTLV prom is a composite promoter comprising the Elongation Factor-1alpha (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.
- HSV1-tk::Sh CpG-free: InvivoGen has engineered a fusion gene between the Herpes Simplex Virus 1 (HSV1) thymidine kinase gene and the *Sh ble* gene conferring Zeocin™ resistance. Both genes have been modified and contain no CpG motifs.
- SV40 pAn: the Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA³.

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*.
- Zeo: Resistance to Zeocin™ is conferred by the *Sh ble* gene from *Streptoalloteichus hindustanus*. 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*.
- βGlo pAn: The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription⁴.

1. Kim D. *et al.*, 1990. Use of the human elongation factor 1α promoter as a versatile and efficient expression system. *Gene* 91(2):217-23.
2. Takebe, Y. *et al.*, 1988. R 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.* 1:466-72.
3. Carswell S. & Alwine J., 1989. Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. *Mol. Cell Biol.* 9(10):4248-58.
4. Yu J. & Russell J. 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.

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.

TECHNICAL SUPPORT

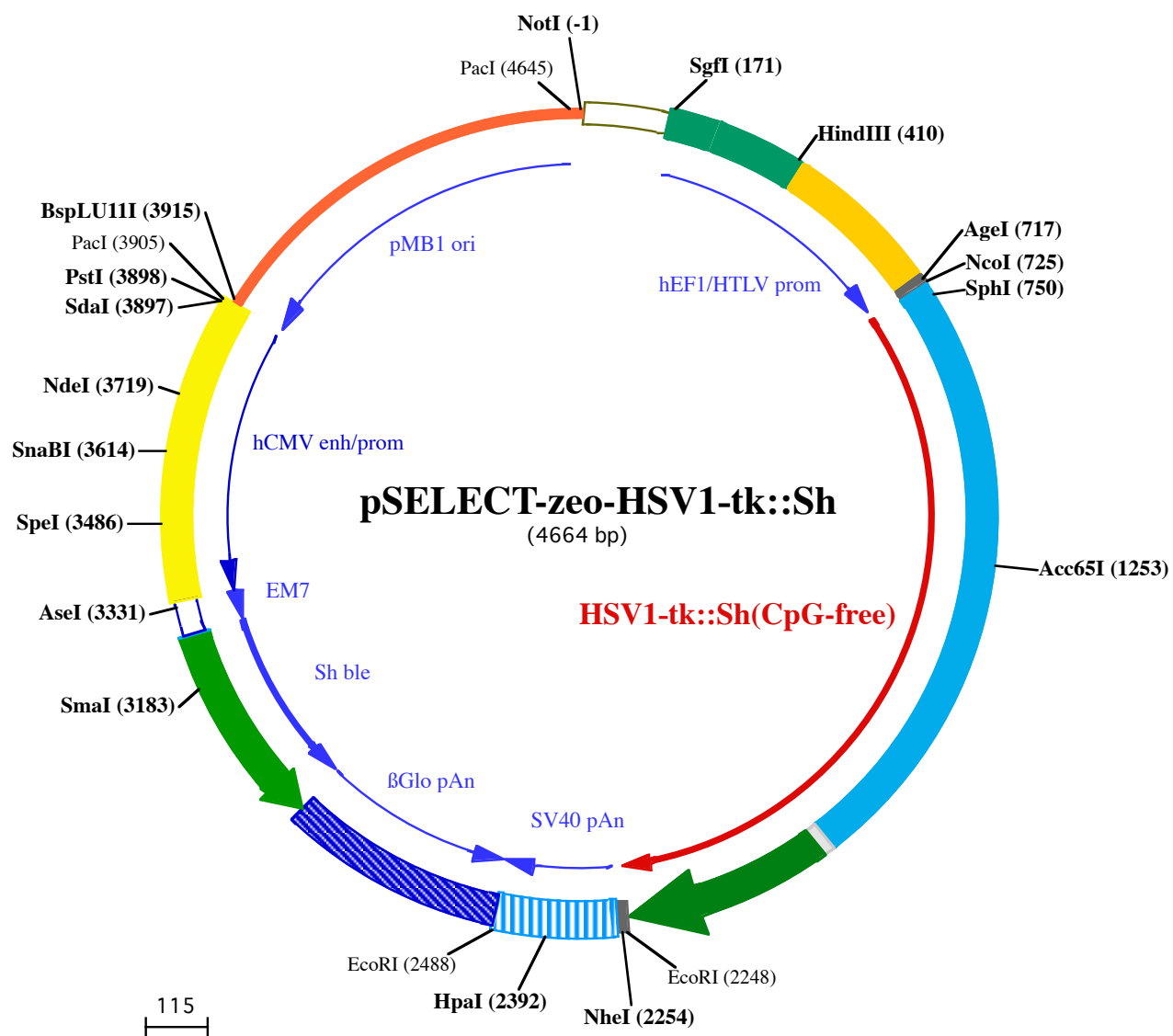
InvivoGen USA (Toll-Free): 888-457-5873

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



NotI (-1)

1 **CGGCCGCA**ATAAAAATCTTTATTTTCATTACATCTGTGTGTTGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAAACAAACGAAACA

SgfI (171)

101 AAACAAACTAGCAAAATAGGCTGTCCCCAGTGCAGTGCAGGTGCCAGAACATTTCTCTATCGAA**GGATCTGCGATCGCTCCGGTCCCGCTCAGTGGGCA**

201 **GAGCGCACATCGCCACAGTCCCGAGAAGTTGGGGGAGGGGTCCGCAATTGAACGGGTGCCTAGAGAAGTGGCCGGGGTAAACTGGGAAAGTATG**

301 **TCGTGACTGGCTCCGCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCCGTGAACGTTCTTTTCGCAACGGGTTTGCCGCCAG**

HindIII (410)

401 **AACACAGCTGAAGTTCGAGGGCTCGCATCTCTCTTCCAGCGCCGCCCTACCTGAGGCCGCCATCCACGCCGTTGAGTCGCGTTCTGCCGCT**

501 **CCCGCTGTGGTGCCTCTGAAGTGCCTCCGCCCTAGGTAAGTTAAAGCTCAGTGCAGACCGGGCTTTGTCCGGGCTCCCTTGAGCCTACTA**

601 **GACTCAGCCGGCTCTCCACGCTTTCCTGACCTGCTTGTCAACTCTACGCTTTGTCTGTTTTCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACC**

AgeI (717)

AgeI (717)

SphI (750)

701 **GGCGCTAC**CTGAGATCACCGTCCACCATGGCTTCTTACCCTGGACACCGATGCTTCTGCCTTTGACCAGGCTGCCAGATCCAGGGGCCACTCCAACA



1 M A S Y P G H Q H A S A F D Q A A R S R G H S N

801 **GGAGAAGTCCCTAAGACCCAGAAGACAGCAGGAAGCCACTGAGGTGAGGCTGAGCAGAAGATGCCAACCTGTGAGGGTGTACATTGATGGACCTCA**

25 R R T A L R P R R Q Q E A T E V R P E Q K M P T L L R V Y I D G P H

901 **TGGCATGGCAAGACCACCCTCAACTGCTGGTGGCACTGGCTCCAGGGATGACATTGTGTATGTGCTGAGCCAATGACCTACTGGAGAGTGTCTA**

58 G M G K T T T T Q L L V A L G S R D D I V Y V P E P M T Y R V L

1001 **GGAGCTCTGAGACATTGCCAACATCTACACCACCCAGCAGGCTGGACAGGGAGAATCTCTGCTGGAGATGCTGCTGGTGTGACCTCTGCC**

92 G A S E T I A N I Y T T Q H R L D Q G E I S A G D A A V V M T S A

1101 **AGATCACAATGGGAATGCCATGTCTGTGACTGATGCTTCTGGCTCCTCACATTGGAGGAGGGTGGCTTCTCATGCCCTCCACTGCCCTGAC**

125 Q I T M G M P Y A V T D A V L A P H I G G E A G S S H A P P P A L T

Acc65I (1253)

1201 **CCTGATCTTTGACAGACACCCATTGCAGCCCTGCTGTGCTACCAGCAGCAAGGTACCTCATGGGCTCCATGACCCACAGGCTGTGCTGGCTTTTGTG**

158 L I F D R H P I A A L L C Y P A A R Y L M G S M T P Q A V L A F V

1301 **GCCCTGATCCCTCAACCTCCCTGGCACCACATTGTTCTGGGAGCACTGCCTGAAGACAGACACATTGACAGGCTGGCAAAGAGGCAGAGACCTGGAG**

192 A L I P P T L P G T N I V L G A L A L P E D R H I D R L A K R Q R P G

1401 **AGAGACTGGACTGGCCATGCTGGCTGCAATCAGAAGGGTGTATGGACTGCTGGCAAACACTGTGAGATACCTCCAGTGTGGAGGCTCTGGAGAGGA**

225 E R L D L A M L A A I R R V Y G L L A N T V R Y L Q C G G S W R E D

1501 **CTGGGACAGCTCTGGAACAGCAGTGCCTCCTAAGGAGCTGAGCCCAAGTCCAATGCTGGTCCAAGACCCACATTGGGGACACCTGTTCCACCTG**

258 W G Q L S G T A V P P Q G A E P Q S N A G P R P H I G D T L F T L

1601 **TTGAGAGCCCTGAGCTGCTGGCTCCCAATGGAGCTGTACAATGTTTGCCTGGGCTGATGTTCTAGCCAAGAGGCTGAGTCCATGCTGATG**

292 F R A P E L L A P N G D L Y N V F A W A L D V L A K R L R S M H V

1701 **TCATCTGGACTATGACCACTCCCTGCTGGATGACAGATGCTCTGCTGCAACTAACCTCTGGCATGGTGCAGACCCATGTGACACCCCTGGCAGCAT**

325 F I L D Y D Q S P A G C R D A L L Q L T S G M V Q T H V T T P G S I

1801 **CCCCACCATCTGTGACCTAGCCAGAACCTTTGCCAGGAGATGGGAATCTCTGGAGCCAATGGAGTATGGCCAAGTTGACCAAGTGTGCTCCAGTGTCTC**

358 P T I C D L A R T F A R E M G I S G A N G V M A K L T S A V P V L

1901 **ACAGCCAGGATGTGGCTGGAGTGTGAGTCTGGACTGACAGGTTGGGTTCTCCAGAGATTTTGTGGAGGATGACTTTGCAGGTGTGGTGCAGAGATG**

392 T A R D V A G A V E F W T D R L G F S R D F V E D D F A G V V R D

2001 **ATGTACCTGTTTATCTCAGCAGTCCAGGACAGGTGGTGGCTGACAACACCTGGCTTGGTGTGGTGGAGGACTGGATGAGCTGTATGCTGAGTG**

425 D V T L F I S A V Q D Q V V P D N T L A W V W V R G L D E L Y A E W

2101 **GAGTGGTGTCTCCACCACTTCAGGATGCCAGTGGCCCTGCATGACAGAGATTGGAGAGCAGCCCTGGGGAGAGAGTTTGCCTGAGAGACCCA**

458 S E V V S T N F R D A S G P A M T E I G E Q P W G R E F A L R D P

NheI (2254)

EcoRI (2248)

2201 **GCAGCAACTGTGTGCACTTTGTGGCAGAGGAGCAGGACTGAGGATAAGAATTCGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAA**

492 A G N C V H F V A E E Q D •

HpaI (2392)

2301 **CCACAAGTAAATGACAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGTGAATAAACAGTTAAACA**

EcoRI (2488)

2401 **CAACAATTGCATTATTTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGTTTTTAAAGCAAGTAAAACCTCTACAATGTGGTATGGAAATCTAAAAT**



2501 **ACAGCATAGCAAACTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGCTGTTGCCAATGTGCA**

2601 **TTAGCTGTTTGACGCTCACCTCTTTCATGGAGTTAAGATATAGTGTATTTCCCAAGGTTTGAAGTACTAGCTTTCATTTCTTTATGTTTTAAATGCAC**

2701 **TGACCTCCACATCCCTTTTATGATAAATATTCAGAAATAATTAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAAATCCAGATGCT**

2801 **CAAGGCCCTTATAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAAGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTAT**

2901 **CCTCAGTCTGCTCTCTGCCACAAAGTGCACGAGTTGCCGGCCGGTCCGCGCAGGGGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTGATGG**

125 D 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

3001 **CCGGCCCGGAGGCTCCCGAAGTTCGTGGACACGACTCCGACCACTCGGCGTACAGCTCGTCCAGGCCCGCACCCACACCCAGGCCAGGGTGTGTTGTC**

92 P G 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

SmaI (3183)

3101 **CGGCACCCTGGTCTGGACCCGCTGATGAACAGGGTACGCTGCTCCGGACACACCCGCGAAGTCTGCTCCACGAAGTCCCGGGAGAACCCGAGC**

59 P V 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

3201 **CGGTGCTCAGAACTCGACCGCTCCGGCGAGCTGCGCGCGGTGAGCACCAGGAAACGGCACTGGTCAACTTGGCCATGATGGCCCTCTATAGTGAGTCG**

25 R D T W F E V A G A V D R A T L V P V A S T L K A M

AseI (3331)

3301 **TATTATACTATGCCATATACTATGCCGATGATTAATTGTCAAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCACTAAACGAGCTCTGCTTA**



3401 TATAGACCTCCCACCGTACACGCCTACCGCCATTTCGCTCAATGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTACTAGTCAAACAA **SpeI (3486)**

3501 ACTCCCATTGACGCTCAATGGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCACGCCATTGATGACTGCCAAAACCGCATCATCATGGTA

SnaBI (3614)

3601 ATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCATAAGGTCATGTAAGTGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGT

NdeI (3719)

3701 CAATAGGGGGCTACTTGGCATATGATACACTTGATGTAAGTGGCAGTTTACCCTAAATACTCCACCCATTGACGCTCAATGGAAAGTCCCTAT

PstI (3898)

SdaI (3897)

3801 TGGCGTTACTATGGGAACATACGTCATTATTGACGCTCAATGGGCGGGGTCGTTGGGCGGTCAGCCAGGCGGGCCATTTACCCTAAGTTATGTAACGCCT

PacI (3905) **BspLU11I (3915)**

3901 GCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGA

4001 CGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCTCGAAGCTCCCTCGTGCCTCT

4101 CCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGG

4201 TGTAGGTCGTTGCTCCAAGTGGGCTGTGTGCACGAACCCCGTTACGCCGACCCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGT

4301 AAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAAC

4401 TACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAACCA

4501 CCGCTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGCTGACGC

PacI (4645)

4601 TCAGTGAACGAAAACCTCACGTTAAGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA