

pUNO1-hSTING-MRP

Expression vector containing an isoform of human STING lacking exon 7

Catalog code: puno1-hsting-mrp

<https://www.invivogen.com/hsting-mrp>

For research use only

Version 19K10-MM

PRODUCT INFORMATION

Contents

- 20 µg of lyophilized plasmid DNA
- 2 x 1 ml blasticidin at 10 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 at least for 1 year.
- 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 full-length open reading frame (ORF) sequencing.
- Plasmid DNA was purified by ion exchange chromatography.

GENERAL PRODUCT USE

- **Subclone gene into another vector.** Two unique restriction sites flank the gene, allowing convenient excision. The 5' site is BspEI which is compatible with AgeI, XmaI, NgoMIV and SgrAI. The 3' site is NheI which is compatible with XbaI, SpeI, and AvrII.
- **Stable gene expression in mammalian cells.** pUNO1 plasmids can be used directly in transfection experiments both *in vitro* and *in vivo*. pUNO1 plasmids contain the blasticidin-resistance gene (*bsr*) driven by the CMV promoter/enhancer in tandem with the bacterial EM7 promoter. This allows the amplification of the plasmid in *E. coli*, as well as the selection of stable clones in mammalian cells using the same selective antibiotic. pUNO1 allows high levels of expression and secretion of the gene product.

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 water. 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.

PLASMID FEATURES

- **Bsr (blasticidin resistance gene):** The *bsr* gene from *Bacillus cereus* encodes a deaminase that confers resistance to the antibiotic blasticidin. The *bsr* gene is driven by the CMV promoter/enhancer in tandem with the bacterial EM7 promoter. Therefore, blasticidin can be used to select stable mammalian cells transfectants and *E. coli* transformants.
- **CMV promoter & enhancer** drives the expression of the blasticidin resistance in mammalian cells.

• Human STING-MRP

ORF size: 852 bp

Cloning fragment size: 961 bp

STING (stimulator of interferon genes; also known as TMEM173, MITA, MPYS, and ERIS) is essential for the IFN response to microbial or self-DNA, and acts as a direct sensor of cyclic dinucleotides (CDNs). CDNs are important messengers in bacteria, affecting numerous responses of the prokaryotic cell, but also in mammalian cells, acting as agonists of the innate immune response. hSTING-MRP (MITA-related protein), discovered and identified in HEK293T cells¹, is an alternatively spliced isoform of hSTING lacking exon 7 that acts as a dominant negative mutant of STING. It was recently reported to block STING-mediated IFN response while retaining the ability to activate NF-κB¹.

• **EF-1α/HTLV hybrid promoter** is a composite promoter comprised of the Elongation Factor-1α (EF-1α) core promoter² and the 5' untranslated region of the Human T-Cell Leukemia Virus (HTLV). EF-1α utilizes a type 2 promoter that encodes for a «house keeping» gene. It is expressed at high levels in all cell cycles and lower levels during G0 phase. The promoter is also non-tissue specific; it is highly expressed in all cell types. The R segment and part of the U5 sequence (R-U5') of the HTLV Type 1 Long Terminal Repeat³ has been coupled to the EF-1α promoter to enhance stability of DNA and RNA. This modification not only increases steady state transcription, but also significantly increases translation efficiency possibly through mRNA stabilization.

• **SV40 pAn:** The Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions, resulting in high levels of steady-state mRNA⁴.

• **pMB1 ori** is a minimal *E. coli* origin of replication to limit vector size, but with the same activity as the longer Ori.

• **Human beta-Globin polyA** is a strong polyadenylation (pAn) signal placed downstream of *bsr*. The use of beta-globin pAn minimizes interference⁵ and possible recombination events with the SV40 polyadenylation signal.

1. Chen H. *et al.*, 2014. An alternative splicing isoform of MITA antagonizes MITA-mediated induction of type I IFNs. *J Immunol* 192(3):1162-70. 2. 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. 3. 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*. 8(1):466-72. 4. 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. 5. Yu J. & Russell J., 2001. Structural and functional analysis of an mRNP complex that mediates the high stability of human β-globin mRNA. *Mol Cell Biol*. 21(17):5879-88.

RELATED PRODUCTS

Product	Description	Cat. Code
Blasticidin	Selection antibiotic	ant-bl-1
ChemiComp GT116	Competent <i>E. coli</i>	gt116-11

TECHNICAL SUPPORT

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

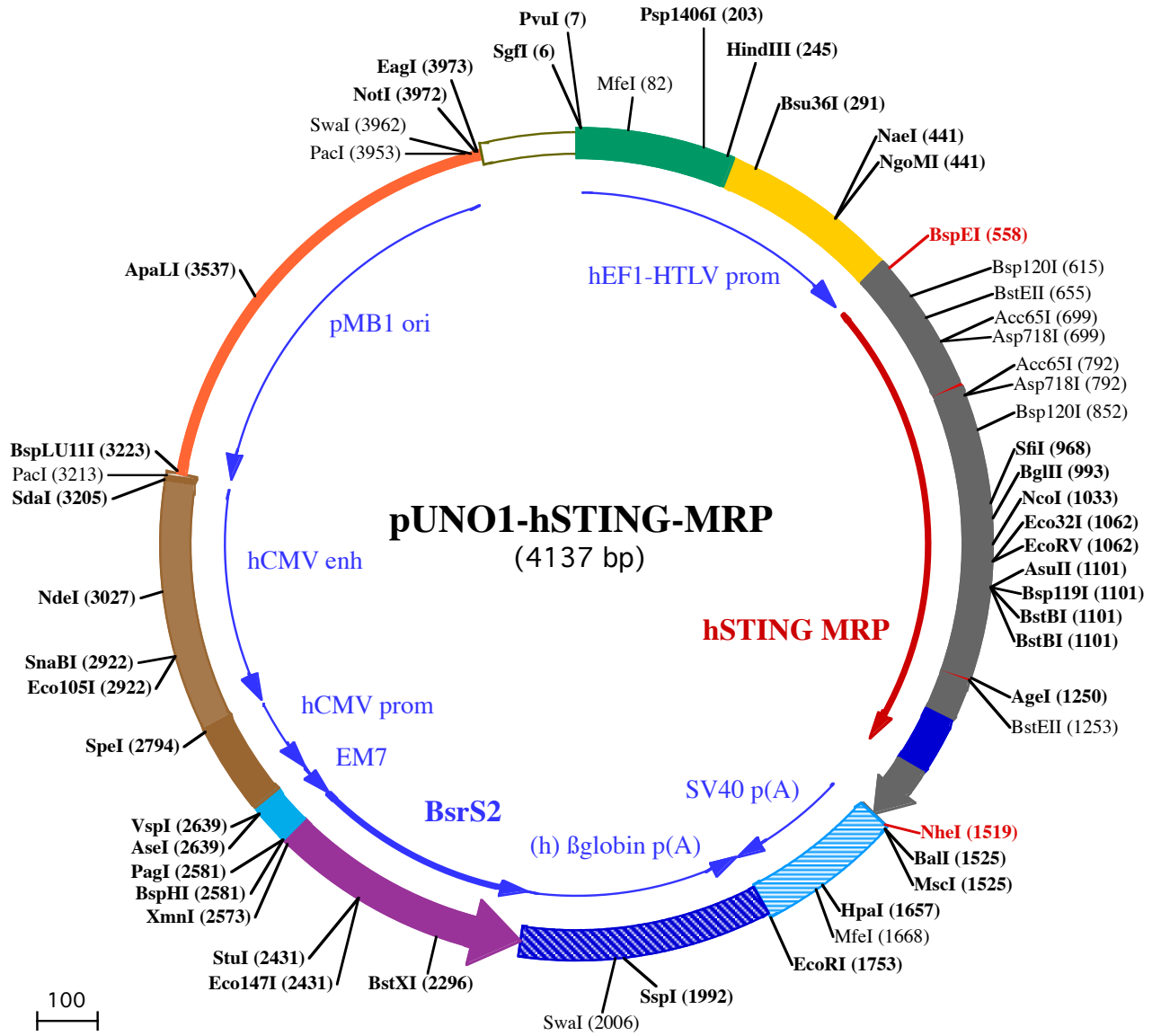
InvivoGen USA (International): +1 (858) 457-5873

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

InvivoGen Hong Kong: +852 3622-3480

E-mail: info@invivogen.com

 **InvivoGen**
www.invivogen.com



PvuI (7)
SgfI (6)
MfeI (82)

1 GGATCTGCATCGCTCCCGTGCCCGTCAGTGGGCAGAGCGACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCCGGCAATTGAACGGGTGCCTA

101 GAGAAGGTGGCGCGGGTAAACTGGGAAAGTATGTCGTGTACTGGCTCCGCCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203)
HindIII (245)
Bsu36I (291)

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCCGCCCTACCTGAGGCC

301 GCCATCCACGCGGGTTGAGTCGCGTCTGCGCCCTCCCGCTGTGGTGCCTCCTGAAGTCCGCTCCGCGCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMI (441)
NaeI (441)

401 GGGCCTTTGTCGGCGCTCCCTTGAGGCTACCTAGACTCAGCCGGCTCTCCAGCTTTGCTGACCCTGCTTGTCAACTCTACGCTTTTGTTTCGTTT

BspEI (558)

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGCCCTACCTGAGATCACCGGCTCCGGAAGATGCCCACTCCAGCTGCATCCATCCATCCCGT

1 ▶ M P H S S L H P S I P

Asp718I (699)
Acc65I (699)

601 GTCCAGGGGTACGGGCCAGAAAGCAGCCTTGGTTCTGTGAGTGCCTGCCTGGTGACCTTTGGGGGCTAGGAGAGCCACAGACACTCTCCG

12▶ C P R G H G A Q K A A L V L L S A C L V T L W G L G E P P E H T L R

Asp718I (792)
Acc65I (792)

701 GTACCTGGTCTCCACCTAGCCTCCCTGCAGCTGGGACTGCTGTTAAACGGGTCTGCAGCCTGGCTGAGGAGTGCACCACATCCACTCCAGGTACCCG

45▶ Y L V L H L A S L Q L G L L L N G V C S L A E E L R H I H S R Y R

Bsp120I (852)

801 GGCAGTACTGGAGACTGTGCGGGCTGCCTGGGCTGCCCCCTCCGCGTGGGGCCTGTTGCTGCTGCATCTATTTCTACTACTCCCTCCCAAATG

79▶ G S Y W R T V R A C L G C P L R R G A L L L L S I Y F Y Y S L P N

SfiI (968)
BglII (993)

901 CGGTGCGCCCGCCTTCACTTGGATGCTTGCCTCCTGGCCTCTCGAGGACTGAACATCCTCCTGGCCTCAAGGGCCTGGCCCAAGCTGAGATCTC

112▶ A V G P P F T W M L A L L G L S Q A L N I L L G L K G L A P A E I S

NcoI (1033)
EcoRV (1062)
Eco32I (1062)

1001 TGCAGTGTGAAAAAGGAATTTCAACGTGGCCATGGGCTGGCATGGTCAATTACATCGGATATCTGCGGCTGATCTGCCAGAGCTCCAGGCCCGG

145▶ A V C E K G N F N V A H G L A W S Y Y I G Y L R L I L P E L Q A R

BstBI (1101)
BstBI (1101)
Bsp119I (1101)
AsuII (1101)

1101 ATTCGAACCTTACAATCAGCATTACAACACCTGCTACGGGGTGCAGTGAAGCAGCGGCTGTATTTCTCTCCATTGGACTGTGGGTGCCTGATAACC

179▶ I R T Y N Q H Y N N L L R G A V S Q R L Y I L L P L D C G V P D N

BstEII (1253)

1201 TGAGTATGGCTGACCCCAACATTCGCTTCTCGGATAAACTGCCCCAGCAGACCGTGACCGTGTCTGGCATCAAGGATCGGGTTTACAGCAACAGCATCTA

212▶ L S M A D P N I R F L D K L P Q Q T G D R A G I K D R V Y S N S I Y

1301 TGAGCTTCTGGAGAACGGGACGCGAACCTGACAGTACAGCAGCTTCTCGCTGCTCCAGGAGGTTCTCCGACCTCGCGCAGGAGAAAAGGAAAGG

245▶ E L L E N G Q R N L Q M T A A S R C P R R F S G T C G R R K R K R

1401 TTACTGTGGGACGTTGAGACCTCAGCGGTGCCAGTACCTCCAGATGTCCAAGAGCCTGAGTCTCTCAGTGAATGAAAAGCCCTCCCTCT

279▶ L L W A A •

MscI (1525)
BalI (1525)
NheI (1519)

1501 CCGCACGGATTCTCTTGAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTGGACAACCACTAGAATGCAGTGAAAAAATGCTTTATT

HpaI (1657)
MfeI (1668)

1601 TGTGAAATTTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACATTGCATTCATTTTATGTTTCAGGTCAGG

EcoRI (1753)

1701 GGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTCATAAATACAGCATAGCAAACTTTAACCTCCAAATCAAGCC

1801 TCTACTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTCTTTTCATGGAT

SspI (1992)

1901 TTAAGATATAGTGTATTTTCCCAAGGTTTGAAGTCTCTTCAATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAATATTTCA

SwaI (2006)

2001 GAAATAATTTAAATACATCATTGCAATGAAAATAATGTTTTTATTAGGCAGAACCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTT

2101 GGACTTAGGGAACAAGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTACTTGAGGGGGATGAGTTCCTCAA

141 ◀ • N R T Y K L P I L E E I

BstXI (2296)

2201 TGGTGGTTTTGACCAGCTGCCATTCTCATATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGCACATGCCACAGGGCTGACCAC

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

2301 CCTGATGGATCTGCCACCTCATCAGAGTAGGGTGCCTGACAGCCAAATGGTGTCAAAGTCTTCTGCCGTTGCTCACAGCAGACCAATGGCAATG

95▶ R I S R D V E D S Y P H R V A V I T D F D K Q G N S V A S G I A I

StuI (2431)
Eco147I (2431)

2401 GCTTCAGCAGACAGTACCTGCCAATGTAGGCTCAATGTGGACAGCAGAGATGATCTCCCACTTGGTCTGATGGCCGCCCGACATGGTGTCT

61▶ A E A C V T V R G I Y A E I H V A S I I E G T K T R I A A G V H H K

PagI (2581)

BspHI (2581)

XmnI (2573)

2501 TGTGTCTCATAGAGCATGGTATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGGTCTTCATGATGGCCCTCTATA
28 N D E Y L M T I K E T A V E V L E L D Q Q S I N F T K M

VspI (2639)

AseI (2639)

2601 GTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTCACTAAACGAGC

SpeI (2794)

2701 TCTGCTTATATAGACCTCCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGT

2801 CAAAACAACTCCCATTGACGTCAATGGGTGGAGACTTGGAAATCCCGTGGTCAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCA

SnaBI (2922)

Eco105I (2922)

2901 TCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTAAGTGGCATAATGCCAGGCGGGCCATTTACCGTC

NdeI (3027)

3001 ATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGTACTGCAAGTGGGCGAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAA

3101 GTCCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGCGGTGAGCCAGGCGGGCCATTTACCGTAAGTTATG

PacI (3213)

SdaI (3205)

BspLU11I (3223)

3201 TAACGCTGCAGGTTAATTAAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCGTTGCTGCGTTTTTCCATAGGCTCCCG

3301 CCCCCTGACGAGCATCACAATAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCG

3401 TGGCCTCTCTGTTCCGACCTGCCGCTTACCGGATACCTGTCCGCTTTTCTCCCTCGGGAAGCGTGGCGTTTTCTCATAGCTCACGCTGTAGGTATCT

ApaLI (3537)

3501 CAGTTCGGTGTAGGTCGTTCCGCTCAAGTGGGCTGTGTGCACGAACCCCGTTCCAGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCC

3601 AACCCCGTAAGACACGACTTATGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGT

3701 GGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAA

3801 ACAAAACCACCGCTGGTAGCGGTGTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGG

EagI (3973)

PacI (3953) SwaI (3962) NotI (3972)

3901 TCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGAATAAAATATCTTTATTTTC

4001 ATTACATCTGTGTGTTGTTTTTTGTGTGAATCGTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCCC

4101 AGTGCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA