

pUNO1-mSTING

Expression vector containing wild-type isoform mouse of STING

Catalog code: puno1-mstingwt
<https://www.invivogen.com/puno-sting>

For research use only

Version 20E18-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 AgeI which is compatible with BspEI, 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 as a 10 mg/ml colorless solution 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.

- **mouse STING-WT**

ORF size: 1137 bp

Cloning fragment size: 1148 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). Several variants of STING have been described in the human population, as well as various induced mutants of the human and mouse STING genes. Studies have revealed that STING variation can affect CDN recognition and signal transduction. Wild-type mouse STING (mSTING-WT) contains an arginine at position 231, similarly to human wild-type STING (hSTING-WT). Unlike hSTING-WT, mSTING-WT appears to have no preference for the cGAMP isomers¹ and efficiently binds DMXAA to produce type I IFNs².

- **EF-1α/HTLV hybrid promoter** is a composite promoter comprised of the Elongation Factor-1α (EF-1α) core promoter³ 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 Repeat⁴ 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 mRNA⁵.
- **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 interference⁶ and possible recombination events with the SV40 polyadenylation signal.

1. Gao P. *et al.*, 2013. Structure-function analysis of STING activation by c[G(2',5')pA(3',5')p] and targeting by antiviral DMXAA. *Cell*, 154(4):748-62. 2. Conlon J. *et al.*, 2013. Mouse, but not human STING, binds and signals in response to the vascular disrupting agent 5,6-dimethylxanthone-4-acetic acid. *J Immunol* 190(10):5216-25. 3. 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. 4. 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. 5. 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. 6. 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.

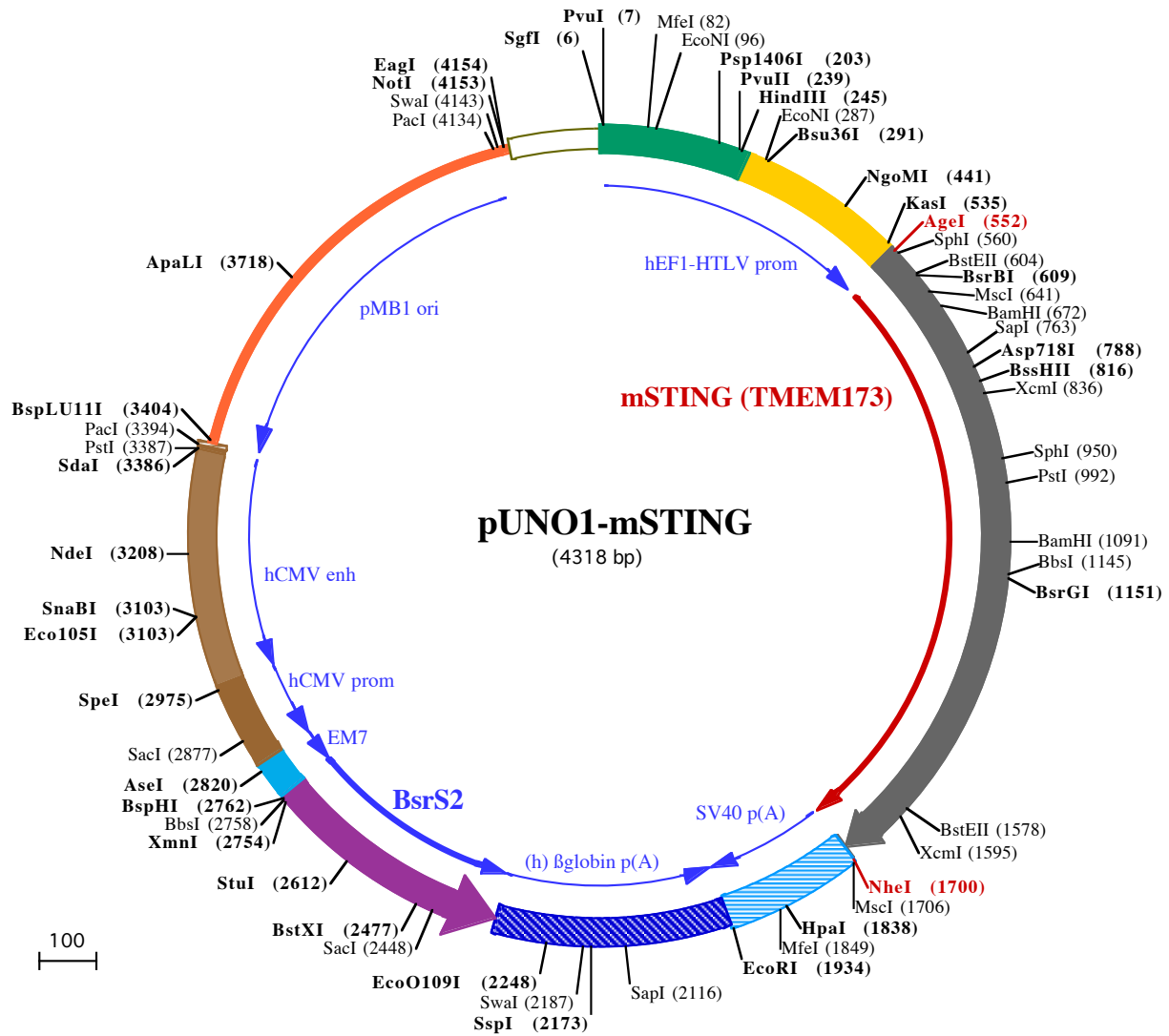
RELATED PRODUCTS

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) 1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA MfeI (82) EcoNI (96)

101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)
Psp1406I (203) 201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACAGCGCCCGCCGCTACCTGAGGCC EcoNI (287) **Bsu36I (291)**
PvuII (239)
EcoNI (287)

301 GCCATCCACGCGGTTGAGTCGCGTTCTGCCGCTCCCGCCTGTGGTGCTCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMI (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCGCTACCTAGACTCAGCGGCTCTCCACGCTTGGCTGACCCTGCTTGTCTAACTCTACGCTTTTGTTCGTTT

KasI (535) **AgeI (552)** **SphI (560)**
501 TCTGTTCTGCGCGGTTACAGATCCAAGCTGTGACCGCGGCTACCTGAGATCACCGGTCAGCATGCCATACTCCAACCTGCATCCAGCCATCCCACGGCC 1 M P Y S N L H P A I P R P

BsrBI (609)
601 CAGAGGTCACCGTCCAAATATGTAGCCCTCATCTTTCTGGTGGCCAGCCTGATGATCCTTTGGTGGCAAAGGATCCACCAAATCACACTCTGAAGTAC 13 R G H R S K Y V A L I F L V A S L M I L W V A K D P P N H T L K Y **BstEII (604)** **MscI (641)** **BamHI (672)**

701 CTAGCACTTACCTAGCCTGCACGAACCTGGACTACTGTTGAAAACTCTGCTGTCTGGCTGAAGAGCTGTGCCATGTCCAGTCCAGGTACCAGGGCA SapI (763) **Asp718I (788)**
47 L A L H L A S H E L G L L L K N L C C L A E E L C H V Q S R Y Q G

BssHII (816) **XcmI (836)**
801 GCTACTGGAAGGCTGTGCGCGCTGCCTGGGATGCCCATCCACTGTATGGCTATGATTCTACTATCGTCTTATTTCTATTTCTCCAAAACACTGCTGA 80 S Y W K A V R A C L G C P I H C M A M I L L S S Y F Y F L Q N T A D

901 CATATACCTCAGTTGGATGTTTGGCCTTCTGGTCTCTATAAGTCCCTAAGCATGCTCCTGGGCTTCAGAGCTTGACTCCAGCGGAAGTCTCTGCAGTC SphI (950) **PstI (992)**
113 I Y L S W M F G L L V L Y K S L S M L L G L Q S L T P A E V S A V

1001 TGTGAAGAAAAGAAGTTAAATGTTGCCACGGGCTGGCCTGGTACATACTACATTGGTACTTGGCGTTGATCTTACCAGGGCTCCAGGCCCGGATCCGAA BamHI (1091)
147 C E E K K L N V A H G L A W S Y Y I G Y L R L I L P G L Q A R I R

BsrGI (1151)
1101 TGTTCAATCAGTACATAACAACATGCTCAGTGGTGCAGGGAGCCGAAGACTGTACATCCTCTTTCCATTGGACTGTGGGGTGCCTGACAACCTGAGTGT BbsI (1145)
180 M F N Q L H N N M L S G A G S R R L Y I L F P L D C G V P D N L S V

1201 AGTTGACCCCAACATTTCGATTCCGAGATATGCTGCCCCAGCAAACATCGACCGTGTGGCATCAAGAATCGGGTTTATTCCAACAGCGTCTACGAGATT 213 V D P N I R F R D M L P Q Q N I D R A G I K N R V Y S N S V Y E I

1301 CTGGAGAACGGACAGCCAGCAGCGTCTGTATCCTGGAGTACGCCACCCCTTGCAGACCCTGTTGGCATGTACAGGATGCCAAAGCTGGCTTCAGTC 247 L E N G Q P A G V C I L E Y A T P L Q T L F A M S Q D A K A G F S

1401 GGGAGGATCGGCTTGAGCAGGCTAAACTCTTCTGCCGGACACTTGAGGAAATCCTGAAAGATGTCCCGAGTCTCGAAAATAACTGCCGCTCATTGTCTA 280 R E D R L E Q A K L F C R T L E E I L E D V P E S R N N C R L I V Y

1501 CCAAGAACCACAGACGGAAACAGTTTCTACTGTCTCAGGAGGTGCTCCGGCACATTCGTGAGGAAAGGAGGTTACCATGAATGCCCCCATG BstEII (1578) **XcmI (1595)**
313 Q E P T D G N S F S L S Q E V L R H I R Q E E K E E V T M N A P M

1601 ACCTCAGTGGCACCTCCTCCCTCGTACTGTCCCAAGAGCCTCCTCATCAGTGGTATGGATCAGCCTCTCCACTCCGCACTGACCTCATCTGAA MscI (1706) **NheI (1700)**
347 T S V A P P P S V L S Q E P R L L I S G M D Q P L P L R T D L I •

1701 GCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTA

HpaI (1838) **MfeI (1849)**
1801 TTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAACATTCATTTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTT

EcoRI (1934)
1901 TTAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAGCCTCTACTGAACTCTTTTCT

2001 GAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTTAAAGATATAGTGATTTTT

SapI (2116) **SspI (2173)** **SwaI (2187)**
2101 CCCAAGGTTTGAAGTCTCTTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTATAGTAAATATTCAGAATAATTTAAATACATC

EcoO109I (2248)
2201 ATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTGTAGTGTGGACTTAGGGAACAAGGA

2301 ACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTGGTGTACTTGGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTT 141 N R T Y K L P I L E E I T T K V L K **SacI (2448)** **BstXI (2477)**

2401 GCCATTCTCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGCACATGCCACAGGGGCTGACCCACTGATGGATCTGTCCACC 122 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 R I S R D V

2501 TCATCAGAGTAGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCGTTGCTCACAGCAGACCCAATGGCAATGGCTTCCAGCAGACAGTGA 88 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 A E A C V T V

StuI (2612)
 2601 CCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCACTTTGGTCTGATGGCCGCCCGACATGGTGCTTGTGCTCATAGAGCAT
 55 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 N D E Y L M

BspHI (2762)
BbsI (2758)
XmnI (2754)
 2701 GGTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCTGCTGAGAGATGTTGAAGTCTTCATGATGGCCCTCTATAGTGAGTCGTATTATACTAT
 22 T I K E T A V E V L E L D Q Q S I N F T K M

AseI (2820) **SacI (2877)**
 2801 GCCGATATACTATGCCGATGATTAATTGTCAAACACAGCGTGGATGGCGTCTCCAGCTATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCC

SpeI (2975)
 2901 CACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCCATG
 3000 ACGTCAATGGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGA

SnaBI (3103)
Eco105I (3103)
 3100 CTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCCATAAAGTCAATGTACTGGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGG

NdeI (3208)
 3200 CGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGAGTTTACCCTAAATACTCCACCATTGACGTCAATGGAAAGTCCCTATTGGCGTACT

PstI (3387)
SdaI (3386)**PacI (3394)**
 3300 ATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTACGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCC T G C A G G T T A

BspLU11I (3404)
 3398 A TTAAGAACATGTGAGCAAAAGGCCAGCAAAAGCCAGGAACCGTAAAAAGGCCGCTTGTGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATC
 3498 ACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCTGGAAGCTCCCTCGTGCGCTCTCTGTTCC
 3598 GACCCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTC

ApaLI (3718)
 3698 GTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGTAAGACAGC
 3798 ACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTA
 3898 CACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAAACAAACCACCGCTGGT
 3998 AGCGGTGGTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAA

EagI (4154)
PacI (4134) **SwaI (4143)** **NotI (4153)**
 4098 ACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC AGCGGCCGCAATAAAATATCTTTATTTTTCATTACATCTGTGTGT
 4198 GGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACAACTAGCAAAATAGGCTGTCCCCAGTGCAAGTGCAGGTG
 4298 CCAGAACATTTCTATCGAA