

# pUNO1-hSTING-M155

Expression vector containing M155 isoform human STING (V155M) open reading frame

Catalog code: puno1-hsting-m155  
<https://www.invivogen.com/hsting-m155>

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-M155**

**ORF size:** 1140 bp

**Cloning fragment size:** 1181 bp

STING (stimulator of interferon genes; also known as TMEM173, MITA, MPYS, and ERIS) is essential for the interferon (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. The M155 (V155M) variant results in a gain-of-function mutation with the constitutive activation of STING and an upregulation of IFN production<sup>1</sup>. This mutation is associated with a chronic autoinflammatory disease, known as STING-associated vasculopathy with onset in infancy (SAVI)<sup>2</sup>.

- **EF-1α/HTLV hybrid promoter** is a composite promoter comprised of the Elongation Factor-1α (EF-1α) core promoter<sup>3</sup> 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 Repeat<sup>4</sup> 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 mRNA<sup>5</sup>.

- **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 interference<sup>6</sup> and possible recombination events with the SV40 polyadenylation signal.

1. **Jeremiah N. et al., 2013.** Inherited STING-activating mutation underlies a familial inflammatory syndrome with lupus-like manifestations. *J Clin Invest.* 124(12):5516-20.
2. **Liu Y. et al., 2014.** Activated STING in a vascular and pulmonary syndrome. *N Engl J Med.* 371(6):507-18.
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

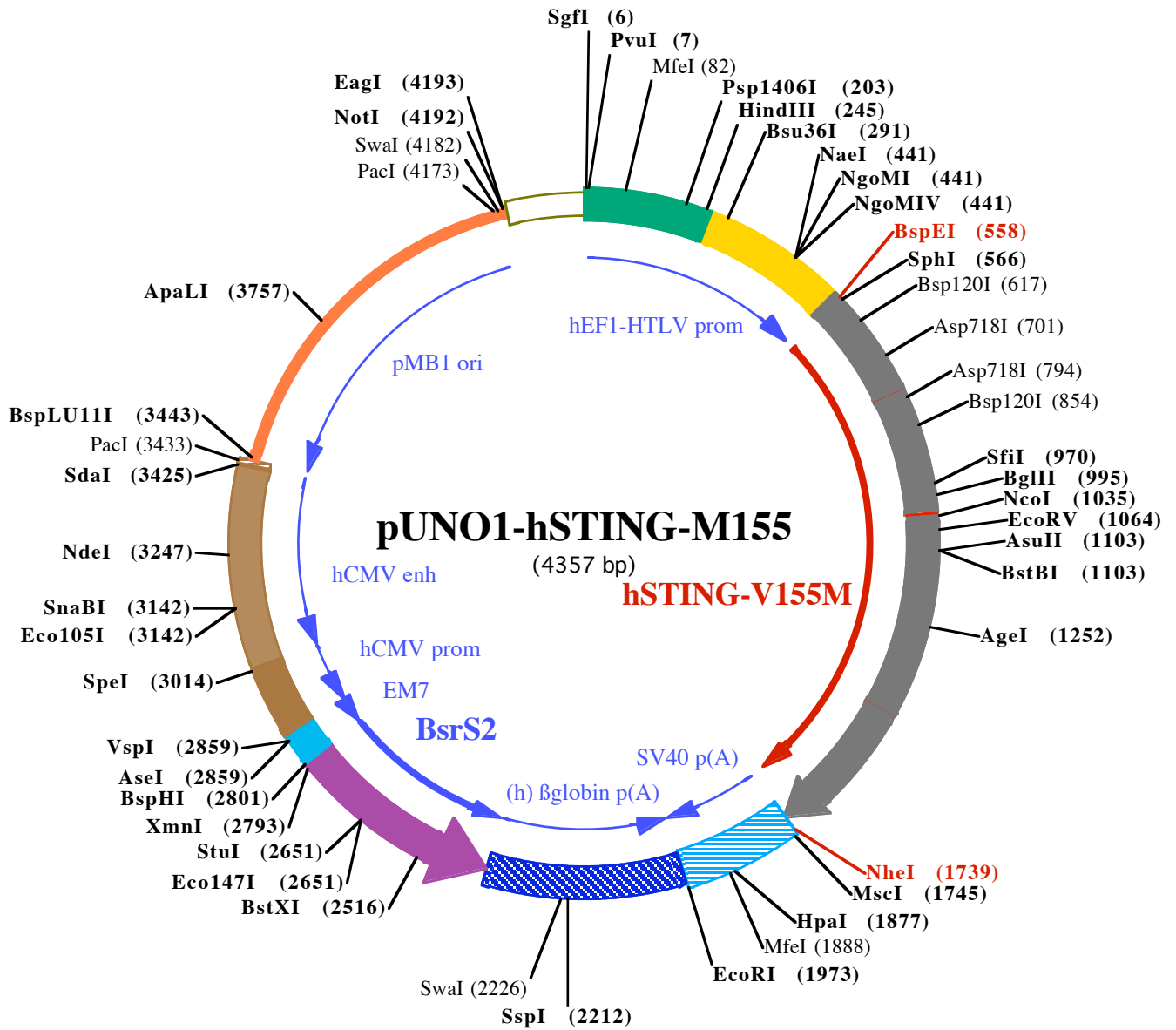
InvivoGen USA (International): +1 (858) 457-5873

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

InvivoGen Hong Kong : +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)





**PvuI (7)**  
**SgfI (6)**  
1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGA

---

**MfeI (82)**  
75 GGGGTCGGCAATTGAACGGGTGCCTAGAGAAGGTGGCGGGGTAAACTGGGAAAGTGATGTCGTGACTGGCT

---

**Psp1406I (203)**  
149 CCGCCTTTTTCCCGAGGGTGGGGGAGAACCGTATATAAGTGCAGTAGTCGCCGTGAACGTTCTTTTTCGCAACG

---

**HindIII (245)** **Bsu36I (291)**  
223 GGTGGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCCTTCACGCGCCCGCCGCTACCTGA

---

297 GGCCGCCATCCACGCGGTTGAGTCGCGTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCGCT

---

**NgoMIV (441)**  
**NgoMI (441)**  
**NaeI (441)**  
371 CTAGGTAAGTTTAAAGCTCAGGTCGAGACCGGGCCTTTGTCCGGCGCTCCCTTGGAGCCTACCTAGACTCAGCC

---

445 GGCTCTCCACGCTTTGCTGACCCTGCTTGCTCAACTCTACGTCTTTGTTTCGTTTTCTGTTCTGCGCCGTTAC

---

**SphI (566)**  
**BspEI (558)**  
519 AGATCCAAGCTGTGACCGGCGCCTACCTGAGATCACCGGCTCCGGACAGCATGCCCCACTCCAGCCTGCATCCA

---

1 M P H S S L H P

**Bsp120I (617)**  
593 TCCATCCCGTGTCCAGGGGTCACGGGGCCAGAAGGCAGCCTTGGTTCTGCTGAGTGCCCTGCCTGGTGACCCT

---

9 S I P C P R G H G A Q K A A L V L L S A C L V T L

**Asp718I (701)**  
667 TTGGGGCTAGGAGAGCCACCAGAGCACACTCTCCGGTACCTGGTCTCCACCTAGCCTCCCTGCAGCTGGGAC

---

33 W G L G E P P E H T L R Y L V L H L A S L Q L G

**Asp718I (794)**  
741 TGCTGTAAACGGGGTCTGCAGCCTGGCTGAGGAGCTGCCACATCCACTCCAGGTACCGGGGCAGCTACTGG

---

58 L L L N G V C S L A E E L R H I H S R Y R G S Y W

**Bsp120I (854)**  
815 AGGACTGTGCGGGCCTGCCTGGGCTGCCCCCTCCGCCGTGGGGCCCTGTTGCTGCTGTCCATCTATTTCTACTA

---

83 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

889 CTCCTCCCAAATGCGGTGCGCCCGCCCTTCACTTGGATGCTTGCCTCCTGGGCCTCTCGAGGCACTGAACA

---

107 S L P N A V G P P F T W M L A L L G L S Q A L N

**SfiI (970)** **BglIII (995)** **NcoI (1035)**  
963 TCCTCCTGGGCCTCAAGGGCCTGGCCCCAGCTGAGATCTCTGCAGTGTGTGAAAAAGGGAATTTCAACATGGCC

---

132 I L L G L K G L A P A E I S A V C E K G N F N M A

**BstBI (1103)**  
**AsuII (1103)**  
1037 CATGGGCTGGCATGGTCATATTACATCGGATATCTGCGGCTGATCCTGCCAGAGCTCCAGGCCCGGATTTCGAAC

---

157 H G L A W S Y Y I G Y L R L I L P E L Q A R I R T

1111 TTACAATCAGCATTACAACAACCTGCTACGGGGTGCAGTGAGCCAGCGGCTGTATATTCTCTCCCATTGGACT

---

181 Y N Q H Y N N L L R G A V S Q R L Y I L L P L D

**AgeI (1252)**  
1185 GTGGGGTGCCTGATAACCTGAGTATGGCTGACCCCAACATTGCTTCTGGATAAACTGCCCCAGCAGACCGGT

---

206 C G V P D N L S M A D P N I R F L D K L P Q Q T G

1259 GACCGTGTGGCATCAAGGATCGGGTTTACAGCAACAGCATCTATGAGCTTCTGGAGAACGGGCAGCGGGCGGG

---

231 D R A G I K D R V Y S N S I Y E L L E N G Q R A G

1333 CACCTGTGCTGGAGTACGCCACCCCTTGCAGACTTTGTTTGCATGTCAATAACAGTCAAGCTGGCTTTA

---

255 T C V L E Y A T P L Q T L F A M S Q Y S Q A G F

1407 GCCGGGAGGATAGGCTTGGAGCAGGCCAACTCTTCTGCCGACACTTGAGGACATCCTGGCAGATGCCCTGAG

---

280 S R E D R L E Q A K L F C R T L E D I L A D A P E

1481 TCTCAGAACAACCTGCCGCTCATTGCCTACCAGGAACCTGCAGATGACAGCAGCTTCTCGCTGTCCCAGGAGGT

---

305 S Q N N C R L I A Y Q E P A D D S S F S L S Q E V

1555 TCTCCGGCACCTGCGGCAGGAGGAAAAGGAAGAGGTTACTGTGGGCAGCTTGAAGACCTCAGCGGTGCCAGTA

---

329 L R H L R Q E E K E E V T V G S L K T S A V P S

1629 CCTCCACGATGTCCCAAGAGCCTGAGCTCCTCATCAGTGAATGGAAAAGCCCTCCTCCTCCGCACGGATTTTC

---

354 T S T M S Q E P E L L I S G M E K P L P L R T D F

MscI (1745)  
NheI (1739)

1703 TCTTGAGACCCAGGGTCACCAGGCCAGAGCCTCCAGTGCTAGCTGGCCAGACATGATAAGATACATTGATGAGT  
379 S •  
1777 TTGGACAAACCACAACACTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTT

HpaI (1877) MfeI (1888)

1851 GTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCATTTTATGTTTTCAGGTTCCAGGGGGA

EcoRI (1973)

1925 GGTGTGGGAGGTTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCAAA  
1999 ACTTTAACCTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTG  
2073 TTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTTAAGATATAGTGTATTTTCCCAAGG

SspI (2212)

2147 TTTGAACTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAAAATATTCA

SwaI (2226)

2221 GAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCT  
2295 TCATAATATCCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAA

2369 GCGAGCTTCTAGCTTTAGTTCCTGGTGTACTTGAGGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCC  
141 • N R T Y K L P I L E E I T T K V L K G

2443 ATTCATCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTGCACATGCCACAGGGGCTGA  
121 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

BstXI (2516)

2517 CCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTTCTGC  
96 V R I S R D V E D S Y P H R V A V I T D F D K Q

StuI (2651)

Eco147I (2651)

2591 CCGTTGCTCACAGCAGACCCAATGGCAATGGCTTCAGCACAGACAGTGACCCTGCCAATGTAGGCCCTCAATGTG  
71 G N S V A S G I A I A E A C V T V R G I Y A E I H

2665 GACAGCAGAGATGATCTCCCCAGTCTTGGTCTGATGGCCGCCCCGACATGGTGCTTGTTCCTCATAGAGCA  
47 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 (2801)

XmnI (2793)

2739 TGGTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGGTCTTCATGATGGCC  
22 T I K E T A V E V L E L D Q Q S I N F T K M

VspI (2859)

AseI (2859)

2813 CTCCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACACAGCGTGGATGGC

2887 GTCTCCAGCTATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCCACCGTACACGCCTACCGCCC

SpeI (3014)

2961 ATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAAACTCC

3034 CATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGCTATCCACGCCCAATTGATGTACTGC

SnaBI (3142)

Eco105I (3142)

3108 CAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTAAGTCCCAAGTAGGAAAGTCCCATAAGGT

NdeI (3247)

3182 CATGTACTGGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGAT

3256 ACACTTGATGTACTGCCAAGTGGGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATT

3330 GCGTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTGCTTGGGCGGTGAGCCAGGCGGGCC

**SdaI (3425)** PacI (3433) **BspLU11I (3443)**

3404 ATTTACCGTAAGTTATGTAACGCCTGCAGGTTAA TTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGG



3476 AACCGTAAAAAGGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACG

3550 CTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCCTGGAAGCTCCCTCGTGC

3624 GCTCTCCTGTTCCGACCCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCT

**ApaLI (3757)**

3698 CATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCC

3772 CGTTCAGCCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGC

3846 CACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGG

3920 TGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAA

3994 AAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTTGTTTGAAGCAGCAGA

4068 TTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGGAACGAA

**EagI (4193)**

PacI (4173) SwaI (4182) **NotI (4192)**

4142 AACTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC AGCGGCCGCAATAAAATATCTTTA

4216 TTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAA

4290 ACAAACAAACTAGCAAATAGGCTGTCCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA