

pUNO1-hSTING-A230

Expression vector containing A230 isoform human STING (G230A) open reading frame

Catalog code: puno1-hsting-a230

<https://www.invivogen.com/hsting-a230>

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-A230

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 IFN response to microbial or self-DNA, and acts as a direct sensor of cyclic dinucleotides (CDNs), such as c-di-GMP. 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. Several non-synonymous variants of STING have been described in the human population. Studies have shown that STING variation can affect CDN recognition and signal transduction. A230 contains a single amino acid substitution G232A. G230 is located in the flexible loop that forms a lid above the c-di-GMP binding pocket. The G230A variant is able to respond to lower concentrations of CDNs due to a different binding to the ligand¹.

• **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. Yi G. *et al.*, 2013. Single nucleotide polymorphisms of human STING can affect Innate immune response to cyclic dinucleotides. PLoS One 8(10):e77846. 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. 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. 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

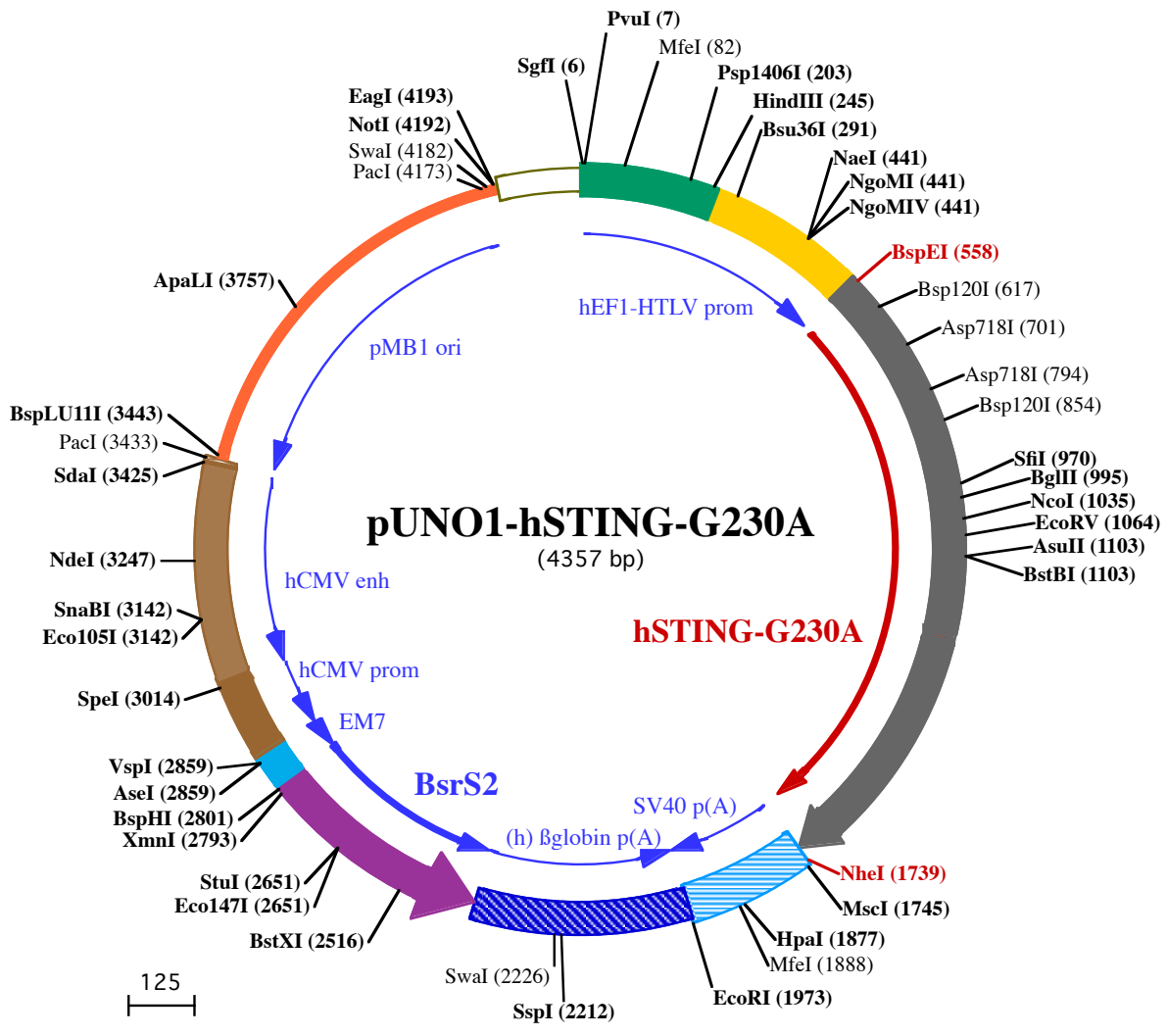
InvivoGen USA (Toll-Free): 888-457-5873

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



PvuI (7)
SgfI (6) **MfeI (82)**
1 GGATCTGCGATCGCTCCGGTCCCGTCACTGGGCGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
101 GAGAAGGTGGCGGGGTAAGTGGGAAAGTGTCTGTACTGGCTCCGCCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203) **HindIII (245)** **Bsu36I (291)**
201 GTGAACGTTCTTTTTTCGCAACGGGTTTCCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGGCCCGCCGCTACCTGAGGCC
301 GCCATCCACGGCGTTGAGTCGGCTTCTGCGCCCTCCCGCTGTGGTGCCTCCTGAAGTCCGCTCCGCTTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
NgoMI (441)
NaeI (441)
401 GGGCCTTTGTCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTTCCACGCTTTCCTGACCTGCTTGCTCAACTCTACGCTTTGTTTCGTTT

BspEI (558)
501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCTACCTGAGATCACGGCTCCGACAGCATGCCCACTCCAGCCTGCATCCATCCATCCC
1► M P H S S L H P S I P

Bsp120I (617)
601 GTGTCCAGGGTACGGGCCAGAAAGCAGCCTTGGTCTGCTGAGTGCCTGCCTGTTGACCTTTGGGGCTAGGAGAGCCACCAGAGCACACTCTC
11► 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

Asp718I (701) **Asp718I (794)**
701 CGGTACCTGGTCTCCACCTAGCTCCCTGCAGTGGGACTGCTGTTAAACGGGGTCTGCAGCCTGGCTGAGGAGCTGCGCCACATCCACTCCAGGTACC
45► R 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

Bsp120I (854)
801 GGGCAGCTACTGGAGGACTGTGCGGGCTGCCTGGGCTGCCCTCCGCGTGGGGCTGTGCTGCTGCCATCTATTCTACTACTCCCTCCAAA
78► R 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 (970) **BglII (995)**
901 TCGGTCGCGCCGCTTCACTTGGATGCTTCCCTCCTGGCCTCTCGCAGGCACTGAACATCCTCCTGGCCTCAAGGGCTGGCCCAAGCTGAGATC
111► 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

NcoI (1035) **EcoRV (1064)**
1001 TCTGCAGTGTGTGAAAAAGGAATTTCAACGTGGCCATGGGCTGGCATGGTCAATTACATCGGATATCTCGGGCTGATCCTGCCAGAGCTCCAGGCC
145► S 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

BstBI (1103)
AsuII (1103)
1101 GGATTCGAACCTACAATCAGCATTACAACAACCTGTACTCGGGGTGCGATGAGCCAGCGGCTGTATATTCTCCTCCATTGGACTGTGGGGTGCCTGATAA
178► R 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
1201 CCTGAGTATGGCTGACCCCAACATTCTGCTTCTGGATAAACTGCCCGCAGCAGCCGCTGACCGTGGTGGATCAAGGATCGGGTTTACAGCAACAGCAGC
211► L S M A D P N I R F L D K L P Q Q T A D R A G I K D R V Y S N S I
1301 TATGAGTCTTGGAGAACGGGAGCGGGGCGGACCTGTGTCTGGATAGCCACCCCTTGCAGACTTTGTTGCCATGTCACAATACAGTCAAGCTG
245► Y E L L E N G Q R A G T C V L E Y A T P L Q T L F A M S Q Y S Q A
1401 GCTTTAGCCGGGAGGATAGGCTTGAGCAGGCCAACTCTTCTGCCGACACTTGAGGACATCCTGGCAGATGCCCTGAGTCTCAGAACAACCTGCCCT
278► G F 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 S Q N N C R L
1501 CATTGCCTACCAGGAACCTGCAGATGACAGCAGCTTCTCGTGTCCAGGAGTTCTCCGGCACCTGCGGCAGGAGGAAAAGGAGGTTACTGTGGC
311► I A Y Q E P A D D S S F S L S Q E V L R H L R Q E E K E E V T V G
1601 AGCTTGAAGACTCAGCGGTGCCAGTACCTCCACGATGTCCAAAGACCTGAGCTCCTCATCAGTGAATGAAAAGCCCTCCCTCTCCGACGGATT
345► S L K T S A V P S T S T M S Q E P E L L I S G M E K P L P L R T D

MscI (1745)
NheI (1739)
1701 TCTCTGAGACCCAGGGTACCAGGCCAGAGCTCCAGTGTAGCTGGCCAGCATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGC
378► F S •

HpaI (1877) **MfeI (1888)**
1801 AGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAACCTGATTC

EcoRI (1973)
1901 TTTTATGTTTCAGTTACAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATTTAAAATACAGCATAGCAAAA
2001 TTTAACCTCAAATCAAGCCTCTACTTGAATCTTTTCTGAGGATGAATAAGGCATAGGCATCAGGGGCTGTGCCAATGTGCATTAGCTGTTGACG
2101 CTCACCTCTTTTATGAGTTAAGATATAGTGTATTTTCCAAAGTTTGAAGTCTCTCATTCTTTATGTTTAAATGCACTGACCTCCACATTC

SspI (2212) **SwaI (2226)**
2201 CCTTTTTAGTAAAATATTCAGAAAATAATTAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAA
2301 TATCCCCAGTTTAGTGTGGACTTAGGGAACAAAGAACCTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTGTACTT
141► • N R T Y K

BstXI (2516)
2401 GAGGGGATGAGTCTCAATGGTGGTTTTGACCAGCTTGCATTCATCTCAATGAGCACAAGCAGTCAAGGAGCATAGTCAAGATGAGCTCTCTGCAC
135► L P I L E E I 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

StuI (2651)
Eco147I (2651)
2601 CAGCAGCCCAATGGCAATGGCTCAGCACAGACAGTACCCTGCCAATGTAGCCCTCAATGTGGACAGCAGAGATGATCCTCCAGCTTGGTCTGAT
68► A S G I A I 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

XmnI (2793)
2701 GGCCGCCCCGACATGGTCTTGTGCTCATAGAGCATGGTGTCTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGGTC
35► A A G V H H K 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

2801 **BspHI (2801)** **VspI (2859)**
AseI (2859)
TTCATGATGGCCCTCCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATC
1 K M
2901 TGACGGTTCACATAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCTACCGCCATTGCGTCAATGGGCGGAGTTGTTACGACATTTGGAA

3001 **SpeI (3014)**
AGTCCCCTTGATTTACTAGTCAAAACAACTCCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGAT

3101 **SnaBI (3142)**
Eco105I (3142)
GTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGTGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGGCATAATGC

3201 **NdeI (3247)**
CAGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGTACTGCCAAGTGGGCAGTTTACCGTAAATACTCCA
3301 CCCATTGACGTCAATGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTATTATTGACGTCAATGGGCGGGGTCGTTGGCGGTCAGCCAGCGCG

3401 **PacI (3433)**
SdaI (3425) **BspLU11I (3443)**
GCCATTTACCGTAAGTTATGTAACGCCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGAAACCGTAAAAGGCCGCTTGTCTGG
3501 CGTTTTCCATAGGCTCCGCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGCGCTT
3601 TCCCCCTGGAAGCTCCCTCGTGCCTCTCTGTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCAT

3701 **ApaLI (3757)**
AGCTCACGCTGTAGTATCTCAGTTCGGTGTAGGTGTTGCTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCCGACCGCTGCGCCTTATCCG
3801 GTAACATCGTCTTGAGTCCAACCCGTAAGACACGACTTATGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGC
3901 TACAGAGTCTTGAAGTGGTGCCTAACTACGGTCACTAGAAAGCAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAGAGTT
4001 GGTAGCTTTGATCCGGCAAAACAAACACCGCTGGTAGCGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATC

4101 **EagI (4193)**
PacI (4173) **SwaI (4182)** **NotI (4192)**
CTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGGAAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGC
4201 AATAAAATATCTTTATTTTATTACATCTGTGTGGTTTTTTGTGTGAATCGTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAAAC
4301 TAGCAAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAATTTCTATCGAA