

# pUNO1-SpikeV8

Expression vector encoding the SARS-CoV-2 Delta variant (B.1.617.2 lineage) Spike gene

Catalog code: p1-spike-v8

<https://www.invivogen.com/delta-b1617-spike-expression-vectors>

For research use only

Version 21G07-ED

## PRODUCT INFORMATION

### Contents

- 20 µg of lyophilized pUNO1-SpikeV8 (plasmid DNA)
- 2 x 1 ml of **Blasticidin** (10 mg/ml)

### Storage and Stability

- Product is shipped at room temperature.
- Lyophilized DNA should be stored at -20°C.
- Resuspended DNA is stable for 1 year at -20°C.
- Store Blasticidin at 4°C or -20°C. The expiry date is specified on the product label.

### Quality control

- Plasmid construct is confirmed by restriction analysis and full-length open reading frame (ORF) sequencing.
- After purification by ion exchange chromatography, predominant supercoiled conformation is verified by electrophoresis.

## PLASMID FEATURES

### Delta Variant SARS-CoV-2 Spike cassette

• **EF-1 $\alpha$ /HTLV hybrid promoter** is a composite promoter comprised of the Elongation Factor-1 $\alpha$  (EF-1 $\alpha$ ) core promoter<sup>1</sup> and the 5' untranslated region of the Human T-Cell Leukemia Virus (HTLV). EF-1 $\alpha$  utilizes a type 2 promoter that encodes a "house-keeping" gene. It is expressed at high levels in all cell cycles and lower levels during the G0 phase. Additionally, since the promoter is not 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>2</sup> has been coupled to the EF-1 $\alpha$  promoter to enhance stability of DNA and RNA. This modification not only increases steady state transcription, but also significantly increases translation efficiency.

- **Codon-optimized Spike ORF**

pUNO1-SpikeV8 contains the Spike coding sequence from the Delta SARS-CoV-2 variant (B.1.617.2 lineage), first identified in India. This variant is characterized by a number of mutations and deletions within the Spike coding sequence (see below)<sup>3</sup>. Additionally, to improve expression of the S protein in cell lines, the gene is codon-optimized and the last 19 amino acids, which contain an endoplasmic reticulum (ER)-retention motif (KxHxx), have been removed<sup>4,5</sup>.

pUNO1-SpikeV8 includes the following sequence features:

- **S1 domain:** T19R, T95I, G142D, E156G,  $\Delta$ F157-R158, D614G, P681R
- **RBD:** L452R, T478K
- **S1/S2 boundary:** Functional furin cleavage site
- **S2 domain:** D950N

Spike (S) is a structural glycoprotein expressed on the surface of SARS-CoV-2. It mediates membrane fusion and viral entry into target cells upon binding to the host receptor ACE2 and the proteolytic activity of host proteases such as furin and TMPRSS2<sup>6</sup>.

For more information visit: <https://www.invivogen.com/sars2-spike>

- **SV40 pAn** is the Simian Virus 40 late polyadenylation (pAn) signal and it enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA<sup>7</sup>.

### Antibiotic selection cassette

- **hCMV (human cytomegalovirus) enhancer & promoter** drive the expression of the blasticidin resistance gene (*bsr*) in mammalian cells.
- **EM7** is a bacterial promoter that enables the constitutive expression of the blasticidin resistance gene (*bsr*) in *E. coli*.
- ***bsr* (blasticidin resistance gene)** encodes a deaminase from *Bacillus cereus* that confers resistance to the antibiotic blasticidin. The expression of the *bsr* gene is driven by the CMV promoter/enhancer and the bacterial EM7 promoter. Therefore, **Blasticidin** can be used to select stable clones in mammalian cells and *E. coli* transformants.
- **Human  $\beta$ -Globin pAn** is a strong polyadenylation (pAn) signal placed downstream of *bsr*. The use of  $\beta$ -globin pAn minimizes interference and possible recombination events with the SV40 pAn signal<sup>8</sup>.

### General features of pUNO1-SpikeV8

- **pMB1 ori** is a minimal *E. coli* origin of replication.

## APPLICATIONS

### Stable gene expression in mammalian cells.

pUNO1 plasmids are designed for both transient and stable transfection in mammalian cell lines by selection with **Blasticidin**. Furthermore, they facilitate high levels of expression of the gene of interest.

### Cell fusion assays

pUNO1-SpikeV8 has been specifically designed for mammalian cell expression of the SARS-CoV-2 S protein. This plasmid features a functional furin cleavage site, and therefore when expressed by a host cell (e.g. 293T cells) it facilitates cell-cell fusion assays with ACE2-expressing cells (e.g. **HEK-Blue™ hACE2 Cells**)<sup>9</sup>.

## 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 the 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 DH5 $\alpha$ .

### 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.

## 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 Asia: +852 3622-3480

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

## REFERENCES

1. Kim D. *et al.*, 1990. Use of the human elongation factor 1 $\alpha$  promoter as a versatile and efficient expression system. *Gene* 91(2):217-23

2. 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.

3. Davis, C. *et al.* 2021. Reduced neutralisation of the Delta (B.1.617.2) SARS-CoV-2 variant of concern following vaccination. medRxiv doi:10.1101/2021.06.23.21259327.

4. Johnson, M.C. *et al.* 2020. Optimized Pseudotyping Conditions for the SARS-COV-2 Spike Glycoprotein. *J Virol* 94.

5. Ou, X. *et al.* 2020. Characterization of spike glycoprotein of SARS-CoV-2 on virus entry and its immune cross-reactivity with SARS-CoV. *Nat Commun* 11, 1620.

6. Hoffmann M. *et al.*, 2020. SARS-CoV-2 cell entry depends on ACE2 and TMPRSS2 and is blocked by a clinically proven protease inhibitor. *Cell.* 181:1-16.

7. 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.

8. 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.

9. Papa, G. *et al.* 2021. Furin cleavage of SARS-CoV-2 Spike promotes but is not essential for infection and cell-cell fusion. *PLoS Pathog* 17, e1009246.

## RELATED PRODUCTS

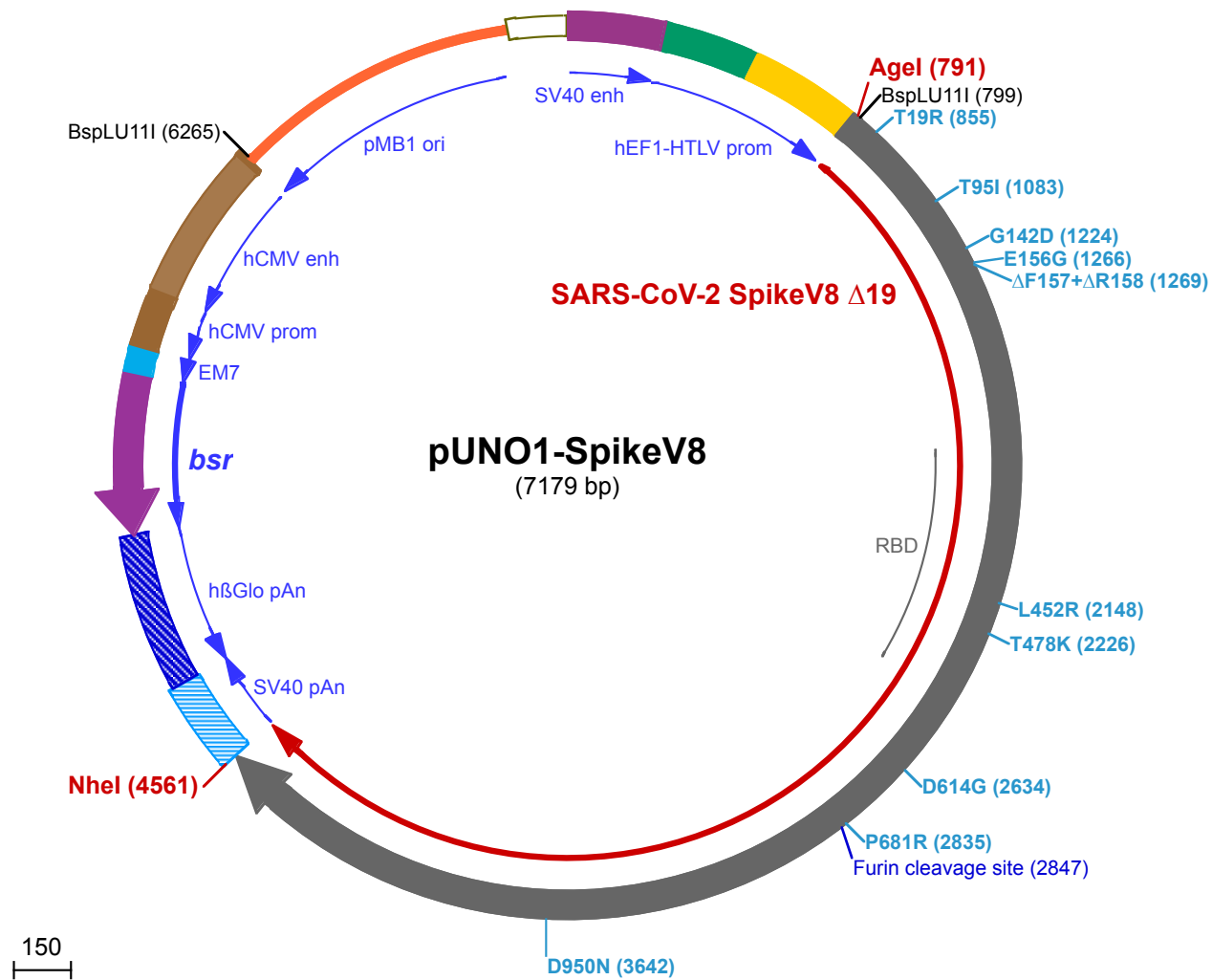
Product	Description	Cat. Code
Blasticidin	Selection antibiotic	ant-bl-1
ChemiComp GT116	Competent <i>E. coli</i>	gt116-11
<b>COVID-19 Product Range</b>		
293-hMyD88 Cells	Cell line	293-hmyd
HEK-Blue™ hACE2 Cells	Cell line	hkb-hace2
A549-hACE2-TMPRSS2 Cells	Cell Line	a549-hace2-tpsa
pUNO1-hACE2	Expression vector	puno1-hace2
pUNO1-hTMPRSS2a	Expression vector	puno1-htp2a
Anti-CoV2RBD-c1-hlgG1	Recombinant Antibody	cov2rbdc1-mab1

For a complete list of InvivoGen's COVID-19 related products visit:  
<https://www.invivogen.com/covid-19>

### 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 Asia: +852 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)





Agel (791)

1 GGACCTGCAGGGCTGAATAACCTCTGAAAGAGGAAGCTTGGTTAGGTACCTTCTGAGGCGGAAAGAACAGCTGTGGAATGTGTGCAGTTAGGGTGTG  
 101 GAAAGTCCCAGGCTCCCAGCAGGCAGAAGTATGCAAAGCATGCATCTCAATTAGTCAGCAACCAGGTGTGGAAAGTCCCAGGCTCCCAGCAGGCAG  
 201 AAGTATGCAAAGCATGCATCTCAATTAGTCAGCAACCATAGTCCCACTAGTCTCCGGTGGCCGTGAGTGGGAGAGCGCACATCGCCACAGTCCCCGA  
 301 GAAAGTGGGGGGAGGGGTGCGCAATTGAACGGGTGCCTAGAGAAGGTGGCGGGGTAACCTGGGAAAGTGTGCTGTACTGGCTCCGCTTTTTCCC  
 401 GAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCCGTGAACGTTCTTTTTTCGCAACGGGTTTCCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTC  
 501 GCATCTCTCTTCACGCGCCCGCCGCTACCTGAGGCCGCCATCCACGCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAACTGC  
 601 GTCCGCGCTTAGGTAAGTTTAAAGCTCAGGTCGAGACCGGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGC

701 CTGACCTGCTTGCTCAACTCTACGTCTTTGTTTCTGTTTCTGTTTCTGCGCAGTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTCAA

T19R (855)

801 CATGTTTGTGTTCTTGGTGTGCTTCCACTGGTCAGTCCCAATGCGTTAATCTCAGAACCCGAACTCAACTCCCACCCGCATATACAAATCCTTCCAC  
 1 M F V F L V L L P L V S S Q C V N L R T R T Q L P P A Y T N S F T  
 901 AGAGGAGTGTACTATCCTGACAAAGTGTTCGGTCAAGTGTCTCCACTCTACTCAGGACCTCTTCTGCCTTCTTTTCAACGTTACATGGTTTCATG  
 34 R G V Y Y P D K V F R S S V L H S T Q D L F L P F F S N V T W F H

T95I (1083)

1001 CAATCCATGTGTCTGGGACAAACGGCACAAACGCTTCGACAACCCTGTATTGCCATTCAATGATGGGGTGTACTTTGCTCCATTGAGAAATCCAACAT  
 67 A I H V S G T N G T K R F D N P V L P F N D G V Y F A S I E K S N I  
 1101 CATTGAGGATGGATTTTCGGGACTACTCTGGACTCAAAGACACAGAGCCTGCTGATCGTTAAACAACGCCACAAACGTTGTCATCAAAGTGTGCGAATTC  
 100 I R G W I F G T T L D S K T Q S L L I V N N A T N V V I K V C E F

ΔF157+ΔR158 (1269)

G142D (1224)

E156G (1266)

1201 CAGTTTTGCAATGATCCCTTCTGGATGTGACTATCACAGAATAACAAGTCTGGATGGAGAGCGGAGTCTACAGCAGCGCAAACAACCTGCACCTTCG  
 134 Q F C N D P F L D V Y Y H K N N K S W M E S G V Y S S A N N C T F  
 1301 AGTACGTGAGTCAACCTTTCTGATGGACCTGGAAGGGAAACAGGGAACTTCAAGAACCTGAGAGAGTTTGTCTTTAAGAACATCGACGGCTATTTTAA  
 167 E Y V S Q P F L M D L E G K Q G N F K N L R E F V F K N I D G Y F K  
 1401 GATCTATAGTAAGCATACGCCTATCAACTGGAAGGGATCTTCCCAGGGCTTTTACGCCCTGGAACCTTTGGTTGACTTGCCTATTGGTATCAATATC  
 200 I Y S K H T P I N L V R D L P Q G F S A L E P L V D L P I G I N I  
 1501 ACCAGATTTAGACCTTCTGGCATTGCAICGGTCTTACTCTCAGGATGATTCCTCCTCCGGGTGGACTGCCGGCGCCGCTGCCTACTATGTCCGGT  
 234 T R F Q T L L A L H R S Y L T P G D S S S G W T A G A A A Y Y V G  
 1601 ATCTGCAACCAAGAAGCTTCTGCTCAAGTACAACGAAACGGCAGTACTATTACGGATGCTGTTGATTGTGCCCTGGACCTCTGTCTGAGACTAAATGCAC  
 267 Y L Q P R T F L L K Y N E N G T I T D A V D C A L D P L S E T K C T  
 1701 CCTCAAGAGCTTACCCTTGAGAAGGGGATTTACCAAACAGTAATTTCCGGGTCCAACCCACCGAAAGCATTGTGCGGTTCCCAAATATACCAATCTG  
 300 L K S F T V E K G I Y Q T S N F R V Q P T E S I V R F P N I T N L

1801 TGTCCCTTTGGCGAAGTGTCAATGCTACAAGTTTGTCTTCTGTGTACGCATGGAATAGGAAACGCATCTCCAATTGTGTCGCTGATTACTCCGTGCTGT  
 334 C P F G E V F N A T R F A S V Y A W N R K R I S N C V A D Y S V L

1901 ACAATCCGCCTCTTCTCAACCTTCAAGTGTATGGCGTTTACCTACCAAACCTTAAACGACCTGTGCTTCACTAATGTGTATGCCACTCTTTTGTGAT  
 367 Y N S A S F S T F K C Y G V S P T K L N D L C F T N V Y A D S F V I

2001 ACGAGGCGATGAAGTGAACAGATTGCACCAGGGCAGACCGGCAAAATTGCCGACTACAACCTACAAGCTTCCAGATGACTTTACCGGATGTGTTATTGCA  
 400 R G D E V R Q I A P G Q T G K I A D Y N Y K L P D D F T G C V I A

L452R (2148)

2101 TGGAACTCAAACAATCTGGATTCCAAGTGGGTGGCAACTATAACTACCGCTATAGACTGTTCAAGAAATCCAACCTGAAACCATTGAGCGAGATATAA  
 434 W N S N N L D S K V G G N Y N Y R Y R L F R K S N L K P F E R D I

T478K (2226)

2201 GCACAGAAATCTACCAGGCTGGAAGTAAAACCTGCAACGGCGTGAAGGGTTCAACTGCTACTTCCCATTGACAGATTACGGATTCCAGCCTACAAACGG  
 467 S T E I Y Q A G S K P C N G V E G F N C Y F P L Q S Y G F Q P T N G

2301 GGTGGGTTACCAACCCTATCGTGTCTGATGCTCCTGAGTTTTGAGCTCCTCCATGCCCCAGCCACAGTCTGTGGCCCCAAGAAAAGCACCAATCTGGTGAAG  
 500 V G Y Q P Y R V V V L S F E L L H A P A T V C G P K K S T N L V K

2401 AACAAATGCGTGAACCTTAACTTAAACGGACTCACAGGAACCGGCGTATTGACGGAGAGTAACAAGAAGTTCTGCCATTCCAGCAGTTCCGGTTCGCGATA  
 534 N K C V N F N F N G L T G T G V L T E S N K K F L P F Q Q F G R D

2501 TTGCCGACACTACCGACGCTGTCCGAGATCCCCAGACATTGGAGATTCTTGATATCACACCCTGTAGTTTCCGGCGAGTGAGCGTGATTACGCCCGGAAC  
 567 I A D T T D A V R D P Q T L E I L D I T P C S F G G V S V I T P G T

D614G (2634)

2601 CAATACCAGCAATCAGGTTGCCGTCCTGTATCAGGGCGTGAATTGCACCGAGGTACCTGTGCGCCATCCACGCTGACCAACTTACCCACATGGCGAGTA
600 N T S N Q V A V L Y Q G V N C T E V P V A I H A D Q L T P T W R V
2701 TATTCCACCGGCTCCAACGTCTTTCAGACACGTGCTGGATGTCTGTATCGGTGCAGAACACGTTAATAATAGCTACGAGTGTGATATCCCCATCGGTGCTG
634 Y S T G S N V F Q T R A G C L I G A E H V N N S Y E C D I P I G A

P681R (2835) Furin cleavage site (2847)

2801 GAATATGCGCCTCTTATCAAACCTCAAACCAACTCTCGTAGGGGGCACGTAGTGTAGCATCCCAAAGTATCATTGCCTACACAATGAGCCTCGGTGCTGA
667 G I C A S Y Q T Q T N S R R R R A R S V A S Q S I I A Y T M S L G A E
2901 GAATTCTGTGCGCTACAGCAACAACTCCATTGTATCCCTACTAATCTCACAATCAGTGTGACAACTGAAATCTGCCGATCTATGACCAAAAACAAGC
700 N S V A Y S N N S I A I P T N F T I S V T T E I L P V S M T K T S
3001 GTTGACTGCACCATGTACATCTGTGGCATTCTACCGAATGTAGCAATCTCCTCGCAATACGGATCATTCTGCATCAGTGAATCGCCCTCACAG
734 V D C T M Y I C G D S T E C S N L L L Q Y G S F C T Q L N R A L T
3101 GTATTGCAAGTTGAGCAGGACAAGAATACGCAGGAAGTGTGGCCAGGTGAAGCAAATCTACAAAACCTCCACCATAAAAAGACTTTGGCGGATTCAATTT
767 G I A V E Q D K N T Q E V F A Q V K Q I Y K T P P I K D F G G F N F
3201 CTCACAGATCCTGCCGATCCCTCAAACCTCAAGCGTAGCTTTATCGAGGATCTGCTCTTCAACAAGGTAACCCCTCGCAGATGCCGGTTTCATCAAG
800 S Q I L P D P S K P S K R S F I E D L L F N K V T L A D A G F I K
3301 CAGTATGGCGATTGTCTGGGAGACATCGCCGCTCGGGACCTGATCTGTGCACAGAAGTTCAATGGACTGACCGTGTGCTCCCTTGTGACCGACGAGA
834 Q Y G D C L G D I A A R D L I C A Q K F N G L T V L P P L L T D E
3401 TGATAGCCAATACACTAGCGCCCTGTGGCCGACCATCTTCTGGGTGACATTCGGAGTTCGGAGTTCGGGCTGCCCTCAGATTCCTTTTGTATGCAGAT
867 M I A Q Y T S A L L A G T I T S G W T F G A G A A L Q I P F A M Q M
3501 GGCCTACCGCTTAAACGGCATCGGTGTGACACAAAACGTTCTGTATGAAAACAGAACTCATCGCCAACAGTTCAACAGTGTATCGGTAAGATACAG
900 A Y R F N G I G V T Q N V L Y E N Q K L I A N Q F N S A I G K I Q

D950N (3642)

3601 GATAGCCTGTCATCCACTGCCAGCGCATTGGGAAAGTTGCAGAAATGTAGTGAACCCAGGCAATGCCAGGCACTTAACACCCTGGTGAACAGCTCTCTTCAA
934 D S L S S T A S A L G K L Q N V N Q N A Q A L N T L V K Q L S S
3701 ATTTTGGTGCCATTTCTAGCGTCTGAATGACATACTAGCCGGTGGCAAGGTGGAGGCTGAAGTGCAGATTGATAGGCTGATAACTGGGCGCCTTCA
967 N F G A I S S V L N D I L S R L D K V E A E V Q I D R L I T G R L Q
3801 GTCTCTCAGACCTATGTGACCCAGCAGCTCATCCGCGCTGCTGAAATTCGCGCATCCGCTAACCTGGCAGCAACCAAAATGTCGAGTGTGTGCTGGGT
1000 S L Q T Y V T Q Q L I R A A E I R A S A N L A A T K M S E C V L G
3901 CAGTCTAAGAGAGTGGACTTTTGCGGGAAGGGGTATCACTGATGTCTTTTCTCAGTCTGCACCCCATGGTGTGGTCTTTCTGCACGTGACTTATGTCC
1034 Q S K R V D F C G K G Y H L M S F P Q S A P H G V V F L H V T Y V
4001 CAGCTCAGGAAAAGAACTTCACTACAGCCCATCTGCCACGATGGGAAAGCCCACTTCCAGGGAAGGCGTATTCGTGTCCAATGGTACTCATTG
1067 P A Q E K N F T T A P A I C H D G K A H F P R E G V F V S N G T H W
4101 GTTCGCTACTCAGAGAAATTTCTACGAGCCCCAGATTATAACCACTGACAATACATTTGTATCCGGCAATTGTGATGTGGTTATCGGGATTGTGAATAAT
1100 F V T Q R N F Y E P Q I I T T D N T F V S G N C D V V I G I V N N
4201 ACTGTTTACGATCCTTTGCGCCAGAGCTGGACTCCTTCAAGGAGGAGCTTGACAAAATATTTAAGAATCACACATCACTGACGTGACCTCGGAGATA
1134 T V Y D P L Q P E L D S F K E E L D K Y F K N H T S P D V D L G D
4301 TTTCAGGAATCAATGCTTCCGTGGTCAATATTCAGAAGGAGATAGACAGGCTGAATGAGGTTGCCAAGAACCTCAACGAGTCTCTGATCGATCTGCAGGA
1167 I S G I N A S V V N I Q K E I D R L N E V A K N L N E S L I D L Q E
4401 GTTGGGCAAGTACGAACAGTATATCAAATGGCCATGGTACATTTGGCTTGGGTTTCATTGCTGGGCTGATAGCTATCGTCATGGTGACAATTATGTTGTG
1200 L G K Y E Q Y I K W P W Y I W L G F I A G L I A I V M V T I M L C

NheI (4561)

4501 TGCATGACATCCTGCTGTAGTTGTCTGAAGGGCTGCTGCTCATGCGGCAGCTGTTGCTAAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTT
1234 C M T S C C S C L K G C C S C G S C C •
4601 GGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAG
TTAACAACAACAATTGCATTCAATTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATT
CTAAAATACAGCATAGCAAAAACCTTAACTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCA
ATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCCAAGGTTTGAAGTACTCTTCATTTCTTTATGTTTTA
AATGCACTGACCTCCACATTCCCTTTTATGTAATAATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCC
AGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTGTGGACTTAGGGAACAAAGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCT
AGCTTTAGTTCCTGGTGTACTTGAGGGGATGAGTTTCTCAATGGTGGTTTTGACCAGCTTGCCATTCATCTCAATGAGCACAAAAGCAGTCAGGAGCATA
141 • N R T Y K 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
5301 GTCAGAGATGAGCTCTCTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGCTGCAGCCACAATGGTGTCA
109 D S I L E R C M G C P S V V R I S R D V E D S Y P H R V A V I T D
5401 AAGTCTTCTGCCGTTGCTCACAGCAGACCAATGGCAATGGCTTACGACAGACAGTGCACCTGCAATGTAGGCCTCAATGTGGACAGCAGAGATGA
75 F D K Q 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 V A S I I
5501 TCTCCCAGTCTTGGTCTGTAGGGCCGCGGACATGGTGTGTTGCTCCTCATAGAGCATGGTGTATCTCAGTGGCGACCTCCACCAGCTCCAGATC
42 E G T K T R I A A G V H K N D E Y L M T I K E T A V E V L E L D
5601 CTGCTGAGAGATGTTGAAGGCTTTCATGATGGCCCTCTATAGTGAGTCTGATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCG
9 Q Q S I N F T K M
5701 TGGATGGCGTCTCCAGCTTATCTGACGGTTCATAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCTACCGCCATTTGCGTCAATGGGGCG

5801 GAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAACAAACTCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCGTGAGTCAA  
5901 ACCGCTATCCACGCCATTGATGTAAGTCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTAAGTCCCAAGTAGGAAAGTCCATAA  
6001 GGTCATGTAAGTGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGTACTGCAAGTGG  
6101 GCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCAATTATTGACGTCAATGGGCGGGGT  
6201 CGTTGGGCGGTGAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAAC  
6301 CGTAAAAAGGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCTGACGAGCATCAGAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAG  
6401 GACTATAAAGATACCAGGCGTTTCCCTGGAAGTCCCTCGTGCCTCTCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTC  
6501 GGAAGCGTGGCGTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCCGTTCAG  
6601 CCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGTAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCA  
6701 GAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCC  
6801 AGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGTATCCGCAAAACAAACCACCGCTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGA  
6901 AAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAAACGAAACTCACGTTAAGGGATTTTGGTCATGGCTAGTTAAT  
7001 TAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAA  
7101 AACAAAACGAAACAAAACAAACTAGCAAATAGGCTGTCCCAAGTGAAGTGCAGGTGCCAGAACATTTCTCTATCGAA