

pNiFty2-IFB-SEAP

Inducible reporter plasmid feature mouse IFN β promoter

Catalog code: pnf2-ifbsp

For research use only

Version 20L03-MM

PRODUCT INFORMATION

Content:

- 20 μ g of pNiFty2-IFB-SEAP provided as lyophilized DNA.
- 1 ml of ZeocinTM (100 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 up to 1 year.
- Store ZeocinTM at 4 °C or at -20 °C. The expiry date is specified on the product label.

Quality control:

Plasmid construct has been confirmed by restriction analysis and sequencing.

GENERAL PRODUCT USE

Interferons are key modulators of the immune response. Their pleiotropic activities are mediated by the induction of many IFN-stimulated genes (ISGs). To help study the transcriptional regulation and signal transduction of type I IFNs, InvivoGen provides several reporter systems, called **pNiFty2**, based on the inducible expression of the secreted embryonic alkaline phosphatase (SEAP) gene. The SEAP gene is cloned under the control of three different promoters that are activated by various transcription factors, such as IRF3, IRF5, IRF7 and NF- κ B.

pNiFty2-IFB-SEAP features the mouse IFN β promoter.

PLASMID FEATURES

- **mIFN β prom:** The mouse interferon beta minimal promoter comprises several positive regulatory domains (PRDs) that bind different cooperating transcription factors such as NF- κ B, IRF3 and IRF7¹. Co-expression of IFN β -SEAP with constitutively activated IRF3 (saIRF3) or IRF7 (saIRF7) in HEK293 cells led to a strong increase in SEAP expression.
- **5U-140** is a synthetic 5'UTR containing an intron.
- **SEAP** is a secreted form of human embryonic alkaline phosphatase. Unlike endogenous alkaline phosphatases, SEAP is extremely heat stable and resistant to the inhibitor L-homoarginine. It catalyses the hydrolysis of pNitrophenyl phosphate (pNpp) producing a yellow end product. SEAP expression can be readily quantified by collecting samples of culture medium and measuring the hydrolysis of pNpp with a spectrophotometer at 405 nm.
- **SV40 pAn:** The Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA.
- **Ori** is a minimal *E. coli* origin of replication with the same activity as the longer Ori
- **hEF1/HTLV prom** is a composite promoter comprising the Elongation Factor-1 α (EF-1 α) core promoter² and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat³. The EF-1 α promoter exhibits a strong activity and yields long lasting expression of a transgene *in vivo*. The R-U5' has been coupled to the EF-1 α core promoter to enhance stability of RNA.
- **EM7** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*.

- **Zeo:** Resistance to the antibiotic ZeocinTM is conferred by the *Sh ble* gene from *Streptoalloteichus hindustanus*. The *Sh ble* gene is driven by the hEF1-HTLV promoter in tandem with the bacterial EM7 promoter allowing selection in both mammalian cells and *E. coli*.
- **β Glo pAn:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription⁴.

1. Vodjdani G. et al. 1998. Structure and characterization of a murine chromosomal fragment containing the interferon beta gene. *J Mol Biol.* 204(2):221-31. **2. Kim DW. et al., 1990.** Use of the human elongation factor 1 alpha 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.* 1: 466-472. **4. Yu J & Russell JE. 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.

METHODS

Plasmid resuspension

Quickly spin the tube containing the lyophilized plasmid to pellet the DNA. To obtain a plasmid solution at 1 μ g/ μ l, resuspend the DNA in 20 μ l of sterile H₂O. Store resuspended plasmid at -20 °C.

Plasmid amplification and cloning

Plasmid amplification and cloning can be performed in *E. coli* GT116 or in other commonly used laboratory *E. coli* strains, such as DH5 α .

ZeocinTM usage

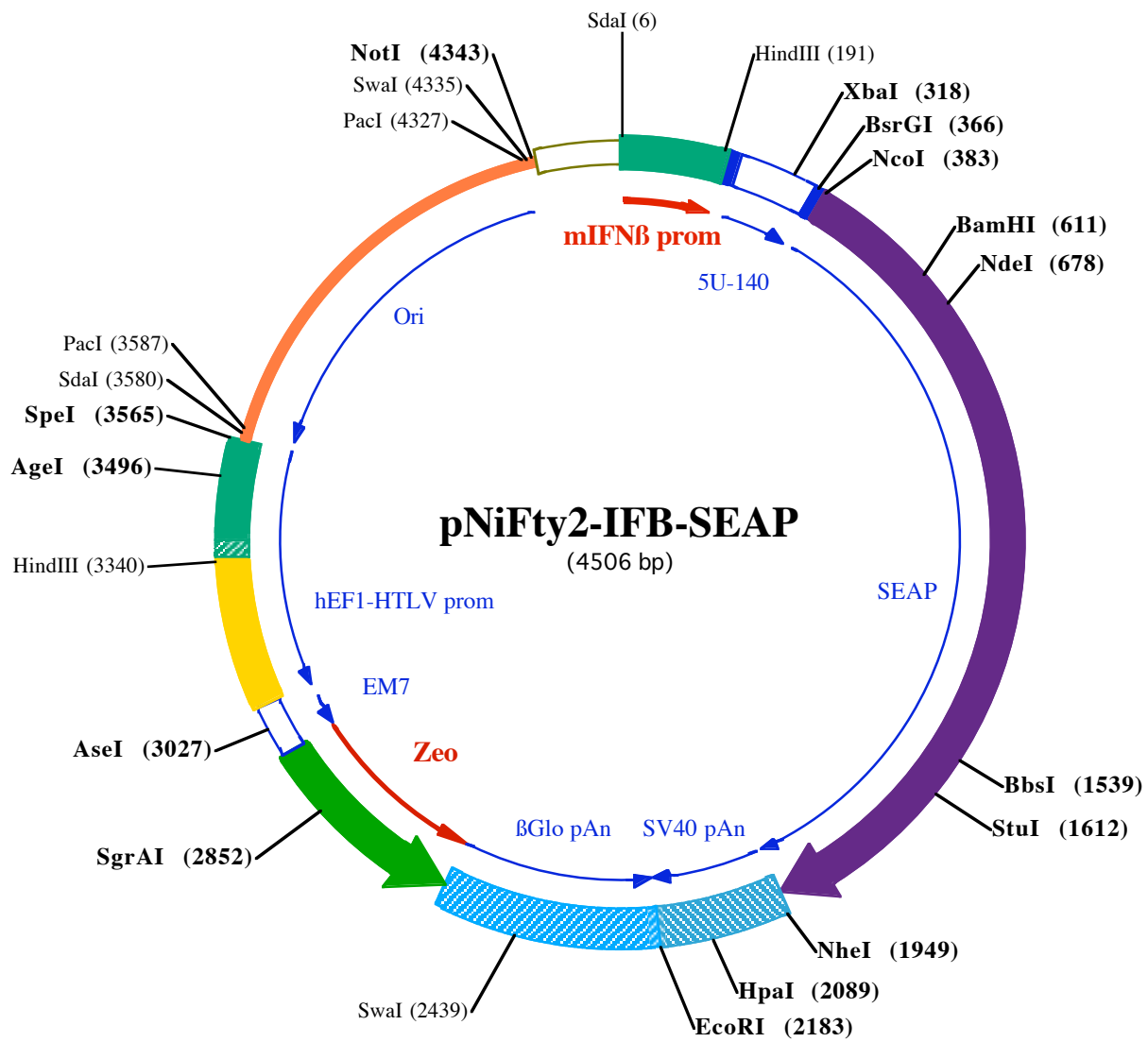
This antibiotic can be used for *E. coli* at 25 μ g/ml in liquid or solid media and at 50-200 μ g/ml to select ZeocinTM-resistant mammalian cells.

RELATED PRODUCTS

Product	Catalog Code
pUNO-sahIRF3	puno-sahirf3
pUNO-sahIRF7 Δ	puno-sahirf7d
pNiFty2-IFA-SEAP	pnf2-ifasp
pNiFty2-56K-SEAP	pnf2-56ksp
Zeocin TM	ant-zn

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 3-622-34-80
E-mail: info@invivogen.com



SdaI (6)
1 CCTGCAGGAGCTTGAATAAAATGAATATTAGAAGCTGTTAGAATAAGAGAAAATGACAGAGGAAAACCTGAAAGGGAGAACTGAAAGTGGGAAATTCCTCT

HindIII (191)
101 GAGGCAGAAAGGCATCCCTTATAATAGCACAGCCATGAAGGAAGATCATTCTACTGCAGCCTTTGACAGCCTTTGCCTCATCTTGAAGCTTCTGC

201 CTTCTCCCTCTGTGAGTTTgtaagtcactgactgtctatgctgtggaaaggggtggcaggagatgggacagtcaggaaaagtgccactatgaaccT

XbaI (318) BsrGI (366) NcoI (383)
301 GCAGCCCTAGGAATGCATAGAcactgtactaaccttctctcttctctctctctctgacagGTTGTTGTACAGTAGCTTCCACATGGTTCTGGGGCCCT
401 GCATGCTGCTGCTGCTGCTGCTGGCCTGAGGCTACAGCTCTCCCTGGGCATCATCCAGTTGAGGAGGAGAACCAGGACTTCTGGAACCCGCGAGGC
61>ysMe tLeuLeuLeuLeuLeuLeuLeuLeuLeuGl yLeuArgLeuGl nLeuSer LeuGl y l e l l eP roVal Gl uGl uAsnProAspPheTrpAsnArgGl uAl
501 AGCCGAGGCCCTGGGTGCCCAAGAAGCTGCAGCCTGCACAGACAGCCGCAAGAACCTCATCATCTTCTGGCGATGGATGGGGTGTCTACGGTG
39>aAl aGl uAl aLeuGl yAl aAl aLysLysLeuGl nP roAl aGl nThr Al aAl aLysAsnLeu l e l l ePheLeuGl yAspGl yMetGl yVal Ser Thr Val

BamHI (611) NdeI (678)
601 ACAGCTGCCAGGATCCTAAAAGGGCAGAAGAAGGACAACTGGGCCTGAGATACCCCTGGCTATGGACCCTCCCATATGTGGCTCTGTCCAAGACAT
73>Thr Al aAl aArg l l eLeuLysGl yGl nLysLysAspLysLeuGl yProGl u l eP roLeuAl aMetAspArgPheP roTyrVal l Al aLeuSer LysThr T
701 ACAATGTAGACAAACATGTGCCAGACAGTGGAGCCACAGCCACGGCCTACCTGTGCGGGTCAAGGGCACTTCCAGACCATTGGCTTGAGTGAGCCGC
106>yrAsnVal l AspLysH i sVal P roAspSer Gl yAl aThr Al aThr Al aTyrLeuCysGl yVal LysGl yAsnPheGl nThr l l eGl yLeuSer Al aAl aAl
801 CCGCTTTAACAGTGAACACGACACGCGGCAACGAGGTCATCTCCGTGATGAATCGGGCCAAAGAAAGCAGGGAAGTCAGTGGGAGTGGTAACACCACA
139>aArgPheAsnGl nCysAsnThr Thr ArgGl yAsnGl uVal l l eSer Val MetAsnArgAl aLysLysAl aGl yLysSer Val Gl yVal Val Thr Thr
901 CGAGTGCAGCAGCCTGCCAGCCGACCTACGCCACAGGTAACCGAACTGGTACTCGGACGCGACGTGCTGCTCGGCCCGCAGGAGGGGT
173>ArgVal Gl nHi sAl aSer P roAl aGl yThr TyrAl aHi sThr Val l AsnArgAsnTrpTyrSerAspAl aAspVal l P roAl aSer Al aArgGl nGl uGl yC
1001 GCCAGGACATCGCTACGACGCTCATCTCCAACATGGACATTTGATGTGATCTGGGTGGAGGCGGAAAGTACATGTTTCCGATGGGAACCCAGACCCTGA
206>ysGl nAsp l l eAl aThr Gl nLeu l l eSerAsnMetAsp l l eAspVal l l eLeuGl yGl yGl yA rgLysTyrMetPheArgMetGl yThr P roAspP roGl
1101 GTACCCAGATGACTACAGCCAAGGTGGGACCAGGCTGGACGGGAGAATCTGGTGCAGGAATGGCTGGCGAAGCGCCAGGGTCCCGGATGTGTGGAAC
239>uTyrP roAspAspTyrSer Gl nGl yGl yThr ArgLeuAspGl yLysAsnLeuVal Gl nGl uTrpLeuAl aLysArgGl nGl yAl aArgTyrVal l TrpAsn
1201 CGCACTGAGCTCATGCAGGCTTCCCTGGACCCGTCTGTGACCATCTCATGGTCTCTTTGAGCCTGGAGACATGAAATACGAGATCCACCAGACTCCA
273>ArgThr Gl uLeuMetGl nAl aSer LeuAspP roSer Val Thr Hi sLeuMetGl yLeuPheGl uP roGl yAspMetLysTyrGl u l l eHi sArgAspSer T
1301 CACTGGACCCCTCCTGATGGAGATGACAGAGGCTGCCCTGCTGAGCAGGAACCCCGCGCTTCTTCTTCTTCTGCGAGGGTGGTGCATCGA
306>hrLeuAspP roSerLeuMetGl uMetThr Gl uAl aAl aLeuArgLeuLeuSerArgAsnP roArgGl yPhePheLeuPheVal Gl uGl yGl yA rg l l eAs
1401 CCACGGTTCATCAGAAAGCAGGGCTTACCGGCCTGACTGAGACGATCATGTTCTGACGACGCCATTGAGAGGGCGGGCCAGCTCACAGCGAGGAGGAC
339>phi sGl yHi sHi sGl uSerArgAl aTyrArgAl aLeuThr Gl uThr l l eMetPheAspAspAl a l l eGl uArgAl aGl yGl nLeuThr Ser Gl uGl uAsp

BbsI (1539)
1501 ACGCTGAGCCTCGTCACTGCCGACCACTCCACGCTTCTTCTTCCGGAGGCTACCCCTGCGAGGGAGCTCCATCTTCCGGCTGGCCCTGGCAAGGCC
373>ThrLeuSerLeuVal Thr Al aAspHi sSer Hi sVal l PheSer PheGl yGl yTyrP roLeuArgGl ySer Ser l l ePheGl yLeuAl aP roGl yLysAl aAl

StuI (1612)
1601 GGGACAGGAAGGCTACACGGTCTCTTATACGAAACGGTCCAGGCTATGTGCTCAAGGACGGCGCCGGCCGGATGTTACCAGAGAGCGAGAGCGGGAG
406>rgAspArgLysAl aTyrThr Val l LeuLeuTyrGl yAsnGl yProGl yTyrVal l LeuLysAspGl yAl aArgP roAspVal Thr Gl uSer Gl uSer Gl ySe
1701 CCCGAGTATCGGCAGCAGTCAAGCAGTGGCCCTGGACGAAGAGACCCACGAGCGAGGACGTGGCGGTGTTCCGCGCGGCCCGCCAGGCGCACTGGTT
439>rP roGl uTyrArgGl nGl nSer Al aVal P roLeuAspGl uGl uThr Hi sAl aGl yGl uAspVal l Al aVal l PheAl aArgGl yP roGl nAl aHi sLeuVal
1801 CACGGCTGACAGGAGCAGACTTCATAGCGCAGCTGATGGCTTCCCGCCTGCTGGAGCCCTACACCCTGCGACTGGCGCCCCCGCGGCACCA
473>Hi sGl yVal Gl nGl uGl nThr Phe l l eAl aHi sVal l MetAl aPheAl aAl aCysLeuGl uP roTyrThr Al aCysAspLeuAl aP roP roAl aGl yThr T

NheI (1949)
1901 CCGACCGCCGACCCGGGGCGTCCCGTCAAGCGTGGATTGAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAA
506>hrAspAl aAl aHi sP roGl yArgSerArgSer LysArgLeuAsp•••

HpaI (2089)
2001 CTAGAATCGAGTGAATAAAATGCTTTATTGTGAAATTTGTGATGCTATTGCTTTATTGTAAACATTATAAGCTGCAATAAACAAGTTAAACAACAACA

EcoRI (2183)
2101 TTGCATTCATTTATGTTTCAGGTTCAAGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTTCAAAATACAGCA
2201 TAGCAAACTTTAACTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGATGAATAAGGCATAGGCATCAGGGCTGTGGCAATGTGCATTAGCT
2301 GTTTGCAGCCTCACCTTCTTATGAGTAAAGATATAGTATTTTCCCAAGGTTTGAAGTACTCTTCTTCTTATGTTTTAAATGCAGTACCT

Swal (2439)
2401 CCCACATTCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGC
2501 CCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAG
1274•••Gl y•••A

SgrAI (2852)
2601 TCCTGCTCCTTGCACAAAGTGCACGAGTTGCCGCGGGTGCAGAGGGGCAACTCCCGCCCCAGGCTGCTCGCCGATCTCGGTCATGGCCGGCC
123>spGl nGl uGl uAl aVal l PheHi sVal l CysAsnGl yAl aP roAspArgLeuAl aPheGl uArgGl yTrpP roGl nGl uGl y l l eGl uThrMetAl aP roGl
2701 CGGAGCGTCCCGAAGTTCGTGGACACGACTCCGACCACTCGGCTACAGCTGCTCAGGCGCGCACCCACCCAGGCGAGGTTGTCTCGGCAC
904>ySer Al aAspArgPheAsnThr Ser Val Val Gl uSer TrpGl uAl aTyrLeuGl uAspLeuGl yA rgVal l TrpVal l TrpAl aLeuThrAsnAspP roVal l

AseI (3027)
2801 CACTGGTCTTGACCGCTGATGAACAGGGTCACTGCTCCCGGACCACCCGCGAAGTCTGCTCCACGAAGTCCCGGAGAACCAGCCGCTCG
57>Val Gl nAspGl nVal l Al aSer l l ePheLeuThr Val l AspAspArgVal l Val Gl yAl aPheAspAspGl uVal l PheAspArgSer PheGl yLeuArgAspT
2901 GTCCAGAAGCTCGACCGCTCCGGCGACGTCGCGCGGGTGGACCGGAAAGCAGCTGGTCAACTTGGCCATGATGGCCCTCTATAGTGAAGTCTATTAT
234>hrTrpPheGl uVal l Al aGl yAl aVal l AspArgAl aThrLeuVal l P roVal l Al aSer ThrLeuLysAl aMet

HindIII (3340)
3001 ACTATGCCATATACTATGCCGATGATTAATTGTCAACTACTGTTTGTAGGCCCGGTCACAGCTTGGATCTGTAACGGCGCAGAACAGAAAACGAAACA
3101 AAGACGTAGAGTTGAGCAAGCAGGTTGAGCAAGCGTGGAGAGCGGCTGAGTCTAGGTAGGCTCCAAGGGAGCGCCGGACAAAGGCCCGCTCTGCACC
3201 TGAGCTTTAACTTACCTAGACGGCGGACGAGTTCAGGAGGCACACAGCGGGAGGCGGAGCGGACTCAACCGCGCTGGATGGCGGCTCAGGT
3301 AGGGCGCGGGCGGTGAAGGAGAGATGCGAGCCCTCGAAGCTTCAGCTGTGTTCTGGCGCAAAACCCGTTGCGAAAAGAACGTTACGGCGACTACT

3401 GCACTTATATACGGTTCTCCCCACCTCGGGAAAAAGGCGGAGCCAGTACACGACATCACTTTCCCAAGTTTACCCCGGCCACCTTCTCTAGGCACCGG **AgeI (3496)**

3501 TTCAATTGCGGACCCCTCCCCCACTTCTCGGGGACTGTGGCGATGTGGCTCTGCCCACTGACTAGTGGGCCCTGCAGGTTAATTAAGAACATGTGA **SpeI (3565)** SdaI (3580) PacI (3587)

3601 GCAAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCTGACGAGCATCACAAAAATCGACGCT

3701 CAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAGATACCAGGCTTTCCCTGGAAGTCCTCGTGGCTCTCTGTTCCGACCTGCGGCTTAC

3801 CGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGTTTTTCATAGCTCACGCTGTAGGTATCTCAGTTCGGGTAGGTCGTTCCGCTCAAGCTG

3901 GGCTGTGTGCACGAACCCCGTTTCAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCGGTAAGACACGACTTATCGCCACTGG

4001 CAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGT

4101 ATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTT

4201 GTTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGCTGACGCTCAGTGGAACGAAACTCACGTT

4301 AAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGAA PacI (4327) SwaI (4335) **NotI (4343)**

4401 TCGTAACATAACATACGCTCTCCATCAAAACAAACGAAACAAACAAACTAGCAAATAGGCTGTCCCAGTGCAAGTGCAAGGTGCCAGAACATTTCTCT

4501 ATCGAA