

# pBOOST2-mcs

Negative control plasmid for pBOOST2 DNA vaccine adjuvant plasmids

Catalog # pbst2-mcs

For research use only

Version 20K16-MM

## PRODUCT INFORMATION

### Content:

- 20 µg of lyophilized pBOOST2-mcs plasmid
- 1 ml of Zeocin™ (100 mg/ml)

### Shipping and storage:

Products are shipped at room temperature.

Lyophilized DNA is stable for 12 months when stored at -20°C.

Resuspended DNA is stable for 6 months when stored at -20°C.

Store Zeocin™ 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.

Plasmid DNA was purified by ion exchange chromatography and lyophilized.

## GENERAL PRODUCT USE

pBOOST2 plasmids were developed as genetic adjuvants for DNA vaccines to potentiate the immune response to a specific antigen. They feature different genes from the interferon regulatory factor family (IRF). IRFs are transcriptional activators for IFN- $\alpha$ , IFN- $\beta$  and IFN-stimulated genes. In particular IRF-1, IRF-3 and IRF-7 act as direct transducers of virus-mediated signaling pathways activating IFN- $\alpha$  and IFN- $\beta$  in infected cells. Recently, IRF-1, IRF-3 and IRF-7 were shown to be able to bias T cells towards type 1 or type 2 immune responses, leading to the activation of cytotoxic T cells and/or the production of antibodies. The method of plasmid DNA vaccine delivery is known to bias the immune response to a specific antigen towards a type 1 (T-cell) or type 2 (antibody) response<sup>1</sup>. These biases can be further enhanced by the codelivery of IRFs to increase the efficacy of the vaccination<sup>2,3</sup>.

## PLASMID FEATURES

- **hEF1 / HTLV prom** is a composite promoter comprising the Elongation Factor-1 $\alpha$  (EF-1 $\alpha$ ) core promoter<sup>4</sup> and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat<sup>5</sup>. The EF-1 $\alpha$  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 $\alpha$  core promoter to enhance stability of RNA.
- **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.
- **EM7** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*.

• **Sh- $\Delta$ CpG (Synthetic Zeocin' gene):** The *Sh ble* gene from *Streptoalloteichus hindustanus* encodes a small protein that confers resistance to Zeocin™ by binding to the antibiotic. To reduce the amount of CpG motifs that may skew the raised antigen-specific immune response, pBOOST2 contains a CpG-free allele of the Zeo<sup>s</sup> gene. All CpGs from the wild-type gene (50) were removed by synthesizing a new allele that contains no CpGs but encodes the exact same protein sequence.

### References:

1. Robinson HL., 1999. DNA vaccines: basic mechanism and immune responses (Review). *Int J Mol Med.* 4(5):549-55.
2. Sasaki S. *et al.*, 2002. Regulation of DNA-raised immune responses by cotransfected interferon regulatory factors. *J Virol.* 76(13):6652-9.
3. Bramson JL. *et al.*, 2003. Super-activated interferon-regulatory factors can enhance plasmid immunization. *Vaccine.* 21(13-14):1363-70.
4. Kim, D.W. *et al.*, 1990. Use of the human elongation factor 1 alpha promoter as a versatile and efficient expression system. *Gene* 2: 217-223.
5. Takebe, Y. *et al.*, 1988. R 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.

## 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 H<sub>2</sub>O. 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 $\alpha$ .

### **Zeocin™ usage**

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

---

## TECHNICAL SUPPORT

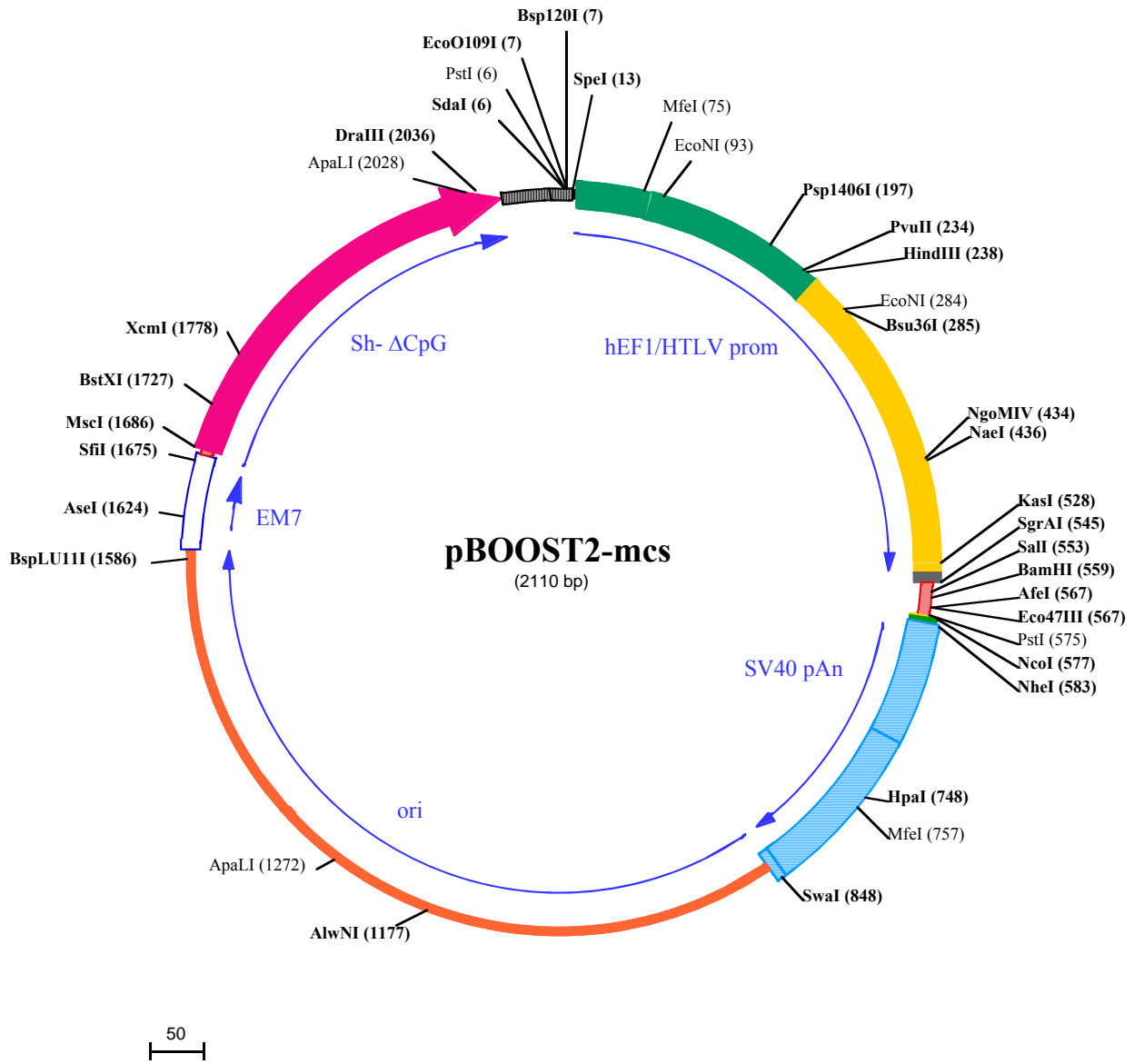
InvivoGen USA (Toll-Free): 888-457-5873

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)



**Bsp120I (7)**  
**EcoO109I (7)**  
PstI (6)  
**SdaI (6)**      **SpeI (13)**      MfeI (75)      EcoNI (93)

1 CCTGCAGGGCCCACTAGTCAGTGGGCGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTAGAGAAGGT

---

**Psp1406I**

101 GGC GCGGGTAAACTGGGAAAGTGATGTCGTGTACTGGCTCCGCTTTTTTCCCGAGGGTGGGGGAGAACCGTATATAAGTGCAGTAGTCCCGTGAACGT

---

**HindIII (238)**      **Bsu36I (285)**  
**PvuII (234)**      EcoNI (284)

201 TCTTTTTTCGCAAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCCTTCAAGCGCCCGCCCTACCTGAGGCCCATCCA

---

301 CGCCGGTTGAGTCGCGTTTCTGCCCTCCCGCTGTGGTGCCTCCTGAACTGCGTCCGCGCTTAGGTAAGTTTAAAGCTCAGGTCGAGACCGGGCTTT

---

**NgoMIV (434)**  
**NaeI (436)**

401 GTCCGGCGCTCCCTTGGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTGCCTGACCTGCTTGCCTCAACTCTACGCTCTTTGTTTCGTTTCTGTTCT

---

**Eco47III (567)**  
**KasI (528)**      **SgrAI (545)**      **SalI (553)**      **AfeI (567)**      **NcoI (577)**  
**BamHI (559)**      PstI (575)      **NheI (583)**

501 GCGCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCAccggcgtgtcgaccggatccagcgtctgagcCATGGCTAGCTCGACATGATAA

---

601 GATACATTGATGAGTTTGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTTATTTGTAAATTTGTGATGCTATTGCTTTATTTGTGAAATTTGT

---

**HpaI (748)**      MfeI (757)

701 GATGCTATTGCTTTATTTGTAAACCATTATAAGCTGCAATAAACCAAGTTAACCAACAATTGCATTCATTTTATGTTCAGGTTTCAGGGGAGGTGTGG

---

**SwaI (848)**

801 AGGTTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTAGATCCATTTAAATGTTAATTAAGTCCATGACCAAAATCCCTTAACGTGAGTTTTCGTTC

---

901 CACTGAGCGTCAGACCCCGTAGAAAAGATCAAAGGATCTTCTTGAGATCCTTTTTTCTGCGCGTAATCTGCTGCTTGCAAACAAAAAACCCCGCTAC

---

1001 CAGCGGTGGTTTGTTCGCCGATCAAGAGCTACCAACTCTTTTTTCCGAAGGTAAGTGGCTTCAGCAGAGCGCAGATACCAAATACTGTTCTTCTAGTGTA

---

**AlwNI (1177)**

1101 GCCGTAGTTAGGCCACCCTCAAGAACTCTGTAGCACCGCTACATACCTCGCTCTGCTAATCCTGTACACAGTGGCTGCTGCCAGTGGCGATAAGTCCG

---

**ApaLI (1272)**

1201 TGTCTTACCGGTTGGACTCAAGACGATAGTTACCGGATAAGGCGCAGCGGTGGGCTGAACGGGGGTTCTGTGCACACAGCCAGCTTGGAGCGAACGA

---

1301 CCTACCCGAACTGAGATACCTACAGCGTGAGCTATGAGAAAGCGCCACGCTTCCGAAAGGAGAAAGCGGACAGGTATCCGTAAGCGGCAGGGTCCGG

---

1401 AACAGGAGAGCGCAGAGGGAGCTTCCAGGGGAAACGCCCTGGTATCTTTATAGTCTGTGCGGGTTTCGCCACCTCTGACTTGAGCGTCGATTTTTGTGA

---

**BspLU11I (1586)**

1501 TGCTGTCAGGGGGCGGAGCCTATGGAAAAACGCCAACCGCGCCTTTTTACGGTTCCTGGCCTTTTGTGTCCTTTTGTGTCATGTTCTTAATTA

---

**AseI (1624)**      **MscI (1686)**  
**SfiI (1675)**

1601 AATTTTTCAAAGTAGTTGACAATTAATCATCGGCATAGTATATCGGCATAGTATAATACGACTCACTATAGGAGGGCCATCATGGCCAAGTTGACCACT

---

**BstXI (1727)**      **XcmI (1778)**

1701 GCTGTCCAGTGTCAAGCCAGGGATGTGGCTGGAGCTGTTGAGTTCTGGACTGACAGGTTGGGGTTCTCCAGAGATTTTGTGGAGGATGACTTTGCAG

7▶ A V P V L T A R D V A G A V E F W T D R L G F S R D F V E D D F A

1801 GTGTGGTCAGAGATGATGTCACCCTGTTTCATCTCAGCAGTCCAGGACAGGTGGTGCCTGACAAACCCCTGGCTTGGGTGTGGGTGAGAGGACTGGATGA

40▶ G V V R D D V T L F I S A V Q D Q V V P D N T L A W V W V R G L D E

1901 GCTGTATGCTGAGTGGAGTGAGTGGTCTCCACCAACTTCAGGGATGCCAGTGGCCTGCCATGACAGAGATTGGAGAGCAGCCCTGGGGGAGAGATTT

73▶ L Y A E W S E V V S T N F R D A S G P A M T E I G E Q P W G R E F

---

**DraIII (2036)**  
**ApaLI (2028)**

2001 GCCCTGAGAGACCAGCAGGCAACTGTGTGCACTTTGTGGCAGAGGAGCAGGACTGAGGATAAGAATTGTAACAAAAACCCCGCCCCGGGGTTTTTT

107▶ A L R D P A G N C V H F V A E E Q D •

2101 TGTTAATTAA