

Intramuscular inoculation

Plasmid DNA solution

- Prepare the vaccine plasmid solution by resuspending 10 μg of the vaccine plasmid DNA in 50 μl saline solution.
- Prepare the pBOOST2 solution by mixing 10 μg of pBOOST2-sahIRF73 and 90 μg of the mock plasmid pBOOST2-null in 50 μl saline solution for low dose, or 100 μg of pBOOST2-sahIRF73 in 50 μl saline solution for high dose.
- Combine both solutions to obtain a total of 110 μg DNA in 100 μl saline solution.

Note: The quantities are per mouse.

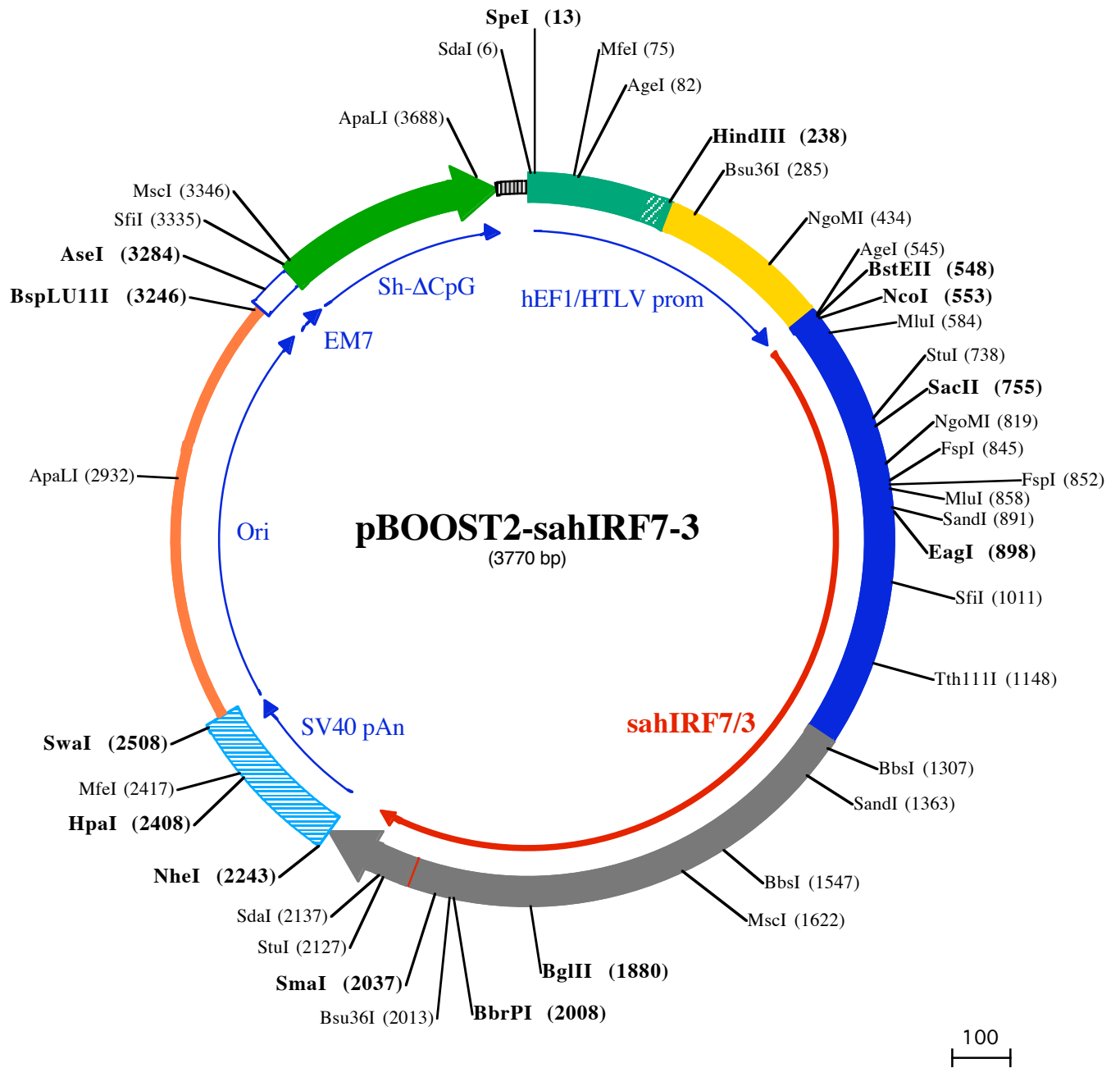
Intramuscular injections

- Inoculate 6 to 8-week old female BALB/c mice with 100 μl plasmid DNA solution (described above) into the quadriceps at 0 and 4 weeks.
- Collect sera and analyze for antibodies at 8 weeks.

Note: For more information see the article by Sasaki S. et al.¹

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



SdaI (6) SpeI (13) MfeI (75) AgeI (82)
1 CCTGCAGGGCCACTAGTCAGTGGGCAGAGCGACATCGCCACAGTCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACCGTGCCTAGAGAAGGT
101 GGC CGGGGTA AACTGGGAAAGTATGTCGTACTGGCTCCGCTTTTTCCCGAGGGTGGGGGAGAACC GTATATAAGTGCAGTAGTCGCCGTGAACGT
HindIII (238) Bsu36I (285)
201 TCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACGCGCCCGCCCTACCTGAGGCCGCATCCA
301 CGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAAGTCTCGCCGCTAGGTAAAGTTAAAGCTCAGGTCGAGACCGGGCCTTT
NgoMI (434)
401 GTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTGGCTGACCCTGCTTGTCAACTCTACGCTTTTGTTCGTTTTCTGTTCT
NcoI (553)
BstEII (548) MluI (584)
501 GCGCCGTTACAGATCCAAGCTGTGACCGCGCCCTACCTGAGATCACCGTACCATGGCCTTGGCTCCTGAGAGGGACGCCACGCTGCTTCCGGAG
1MetAl aLeuAl aProGl uArgAl aAl aProArgVal LeuPheGl yG
601 AGTGGCTCCTTGAGAGATCAGCAGCGGCTGCTATGAGGGGCTGCAGTGGCTGGACGAGGCCCGCACCTGTTCCGCGTGCCTGGAAGCACTTCGCGCG
16I uTrpLeuLeuGl yGl uI eSer Ser Gl yCysTyrGl uGl yLeuGl nTrpLeuAspGl uAl aArgThr CysPheArgVal P roT rpLysHi sPheAl aAr
StuI (738) SacII (755)
701 CAAGGACCTGAGCGAGGCCGACGCGCATCTTCAAGGCTGGCTGTGGCCGCGGAGGTGGCCGCTAGCAGCAGGGGAGGTGGCCCGCCCGGAG
49I gLysAspLeuSer Gl uAl aAspAl aArgI l ePheLysAl aTrpAl aVal Al aArgGl yArgT rpP roP roSer Ser ArgGl yGl yI yProP roP roGl u
EagI (898)
NgoMI (819) FspI (845) FspI (852) MluI (858) SandI (891)
801 GCTGAGACTGCGGAGCGCGCGGCTGGAAAACCACTTCCGCTGCGCACTGCGCAGCAGCGCTCGCTTCGTGATGCTGCGA GATAACTCGGGGACCCGG
83I Al aGl uThr Al aGl uArgAl aGl yTrpLysThr AsnPheArgCysAl aLeuArgSer Thr ArgArgPheVal MetLeuArgAspAsnSer Gl yAspProA
901 CCGACCCGACAAGGTGTACGCGCTCAGCCGGAGCTGTGCTGGCGAGAAGGCCAGGCAGGCAGACTGAGGCAGAGGCCCGCAGCTGTCCACC
116I l aAspP roHi sLysVal I TyrAl aLeuSer ArgGl uLeuCysT rpArgGl uGl yProGl yThr AspGl nThr Gl uAl aGl uAl aProAl aAl aVal P roP r
SfiI (1011)
1001 ACCACAGGGTGGGCCCCAGGGCCATTCTGGCACACACATGCTGGACTCCAAGCCCCAGGCCCTCCCTGCCCCAGCTGGTGACAAGGGGACCTC
149I oP roGl nGl yGl yProP roGl yProPheLeuAl aHi sThr Hi sAl aGl yLeuGl nAl aP roGl yProLeuP roAl aP roAl aGl yAspLysGl yAspLeu
Tth111I (1148)
1101 CTGCTCAGGCAGTGAACAGAGCTGCCTGGCAGCCATCTGCTGACAGCGTCATGGGGGAGATCCAGTCCCAACCAAGGCTCTGGAGAGGGACAAG
183I LeuLeuGl nAl aVal Gl nGl nSer CysLeuAl aAspHi sLeuLeuThr Al aSer TrpGl yAl aAspP roVal P roThr LysAl aP roGl yGl uGl yGl nG
1201 AAGGCTTCCCTGACTGGGCTGTGCTGGAGGCCAGGGCTCCCTGCTGGGAGCTGTACGGTGGCAGTAGAGACGCCCGCAGCCCACTTCTGA
216I l uGl yLeuP roLeuThr Gl yAl aCysAl aGl yGl yProGl yLeuP roAl aGl yGl uLeuTyrGl yTrpAl aVal I Gl uThr Thr P roSer P roThr SerAs
BbsI (1307) SandI (1363)
1301 TACCCAGGAAGACATTCTGGATGAGTACTGGTAAATGGTGTGGCCCACTCCAGATCCGGGACCCCAAGCCCTGGCTGTAGCCCTGAGCCCTGC
249I pThr Gl nGl uAspI l eLeuAspGl uLeuLeuGl yAsnMeTVal LeuAl aP roLeuP roAspP roGl yProP roSer LeuAl aVal Al aP roGl uP roCys
1401 CCTCAGCCCTTGGGAGCCCAAGCTTGGACAATCCCACTCCCTTCCAAACCTGGGGCCCTCTGAGAACCCTGAAGCGGCTGTTGGTGGCGGGGAAG
283I P roGl nP roLeuArgSer P roSer LeuAspAsnP roThr P roPheP roAsnLeuGl yProSer Gl uAsnP roLeuLysArgLeuLeuVal P roGl yGl uG
BbsI (1547)
1501 AGTGGGAGTTCAGGTGACAGCCTTCTACCGGGGCCCAAGTCTCCAGCAGACCATCTCTGCCCGGAGGGCTGCGGCTGGTGGGGTCCGAAGTGGG
316I l uTrpGl uPheGl uVal I Thr Al aPheTyrArgGl yArgGl nVal I LeuGl nGl nThr I l eSer CysP roGl uGl yLeuArgLeuVal Gl ySer Gl uVal I Gl
MscI (1622)
1601 AGACAGGACGCTGCCTGGATGGCCAGTCACTGCCAGACCCTGGCATGTCCCTGACAGACAGGGGAGTGATGAGCTACGTGAGGCATGTGCTGAGCTGC
349I yAspArgThr LeuP roGl yTrpP roVal I Thr LeuP roAspP roGl yMetSer LeuThr AspArgGl yVal MetSer TyrVal I r gHi sVal I LeuSer Cys
1701 CTGGTGGGGGACTGGCTCTCTGGCGGGCCGGCAGTGGCTCTGGGCCAGCGGCTGGGGCACTGCCACATACTGGCAGTGAGCGAGGAGCTGCTCC
383I LeuGl yGl yGl yLeuAl aLeuT rpArgAl aGl yGl nTrpLeuT rpAl aGl nArgLeuGl yHi sCysHi sThr TyrT rpAl aVal I Ser Gl uGl uLeuLeuP
BgIII (1880)
1801 CCAACAGCGGCATGGCCCTGATGGCGAGTCCCAAGGACAAGGAAGGAGCGTGTGGACCTGGGGCCCTCATTGTAGATCTGATTACCTTACCGGA
416I r oAsnSer Gl yHi sGl yProAspGl yGl uVal P roLysAspLysGl uGl yGl yVal I PheAspLeuGl yProPheI l eVal I AspLeuI l eThr PheThr Gl
1901 AGGAAGCGGACGCTACCCAGCTATGCCCTCTGGTCTGTGGGGAGTCAATGGCCAGGACAGCCGTGGACCAAGAGGCTCGTGATGGTCAAGGTT
449I uGl ySer Gl yArgSer P roArgTyrAl aLeuT rpPheCysVal Gl yGl uSer TrpP roGl nAspGl nP roT rpThr LysArgLeuVal I MetVal I LysVal
Bsu36I (2013)
BbrPI (2008) SmaI (2037)
2001 GTGCCACGTGCTCAGGCCTTGGTAGAAATGGCCCGGTAGGGGTCCTCCTCCCTGGAGAATACTGTGGACCTGCACATTGACAACGCCACCCAC
483I Val P roThr CysLeuArgAl aLeuVal Gl uMe tAl aArgVal I Gl yGl yAl aSer Ser LeuGl uAsnThr Val AspLeuHi s I l eAspAsnSer Hi sP roL
SdaI (2137)
StuI (2127)
2101 TCTCCCTCACCTCCGACCAGTACAAGGCCTACCTGCAGGACTTGGTGGAGGGCATGGATTTCCAGGGCCCTGGGGAGACTGAGCCCTCGCTCCTCATGG
516I euSer LeuThr SerAspGl nTyrLysAl aTyrLeuGl nAspLeuVal I Gl uGl yMetAspPheGl nGl yProGl yGl uThr ●●●
NheI (2243)
2201 TGTGCCCAACCCCTGTTCCCAACCTCAACCAATAAGCTAGCTCGACATGATAAGATACATTGATGAGTTTGACAAACCACAACCTAGAATGCA
2301 GTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACATTATAAGCTGCAATA
HpaI (2408) MfeI (2417)
2401 AACAAGTTAACAACAACAATTGCATTATTATGTTTCAGGTTGAGGGGAGGTGGGAGTTTTTAAAGCAAGTAAACCTTACAATGTGGTGA
SwaI (2508)
2501 ATCCATTTAAATGTTAATTAAGTACCATGACCAAAATCCCTTAACGTGAGTTTTCGTTCCACTGAGCGTCAGACCCCGTAGAAAAGATCAAAGGATCTT
2601 CTTGAGATCCTTTTTTCTGCGGTAATCTGCTGCTTGAACAAAAAACCACCGCTACAGCGGTGGTTTTGTTGCCGGATCAAGAGCTACCAACTCT
2701 TTTTCCGAAGGTAAGTGGCTTACGACAGCGCAGATACCAATACTGTTCTTCTAGTGTAGCCGTAGTTAGGCCACCACTTCAAGAAGCTGTAGCACCG

2801 CCTACATACCTCGCTCTGCTAATCCTGTTACCAGTGGCTGCTGCCAGTGGCGATAAGTCGTGTCTTACCGGGTTGGACTCAAGACGATAGTTACCGGATA

2901 AGGCGCAGCGGTTCGGGCTGAACGGGGGTTCTGTGCACACAGCCAGCTTGGAGCGAACGACCTACACCGAACTGAGATACCTACAGCGTGAGCTATGAGA
 ApaLI (2932)

3001 AAGCGCCACGCTTCCGAAGGGAGAAAGCGGACAGGTATCCGGTAAGCGGCAGGGTCGGAACAGGAGAGCGCACGAGGGAGCTTCCAGGGGAAACGCC

3101 TGGTATCTTTATAGTCTGTGCGGTTTCGCCACCTCTGACTTGAGCGTCGATTTTTGTGATGCTCGTCAGGGGGCGGAGCCTATGAAAAACGCCAGCA

3201 ACGCGGCCTTTTTACGGTTCCTGGCCTTTTGTGGCCTTTTGTCTCACATGTTCTTAATTAATTTTTCAAAGTAGTTGACAATTAATCATCGGCATAGT
 BspLU11I (3246) AseI (3284)

3301 ATATCGGCATAGTATAATACGACTCACTATAAGGAGGGCCATCATGGCCAAGTTGACCAAGTGGAGTGGCTGGAGCTG
 SfiI (3335) MseI (3346)

3401 TTGAGTTCTGGACTGACAGGTTGGGTTCTCCAGAGATTTGTGGAGGATGACTTTGCAGGTGTGGTCAGAGATGATGTCACCTGTTTCATCTCAGCAGT
 1▶MetAl aLysLeuThr SerAl aVal ProVal LeuThrAl aArgAspValAl aGl yAl aV

3501 CCAGGACCAGGTGGTGCCTGACAACACCCTGGCTTGGGTGGGTGAGAGGACTGGATGAGCTGTATGCTGAGTGGAGTGAGGTGGTCTCCACCACTTC
 20▶a l Gl uPheTrpThrAspArgLeuGl yPheSerArgAspPheVal Gl uAspAspPheAl aGl yVal ValA r gAspAspVal Thr LeuPhe l eSerAl aVa
 53▶l Gl nAspGl nVal Val ProAspAsnThr LeuAl aTrpVal l TrpVal A r gGl yLeuAspGl uLeuTyrAl aGl uTrpSer Gl uVal Val Ser ThrAsnPhe

3601 AGGGATGCCAGTGGCCCTGCCATGACAGAGATTGGAGAGCAGCCCTGGGGGAGAGAGTTGCCCTGAGAGACCCAGCAGGCAACTGTGTGCACCTTTGTGG
 87▶A r gAspAl aSer Gl yProAl aMe tThr Gl ul eGl yGl uGl nP roTrpGl yA r gGl uPheAl aLeuA r gAspP roAl aGl yAsnCysVal l Hi sPheVal A

3701 CAGAGGAGCAGGACTGAGGATAAGAATTGTAACAAAAACCCCGCCCGCGGGTTTTTTGTTAATTAA
 120▶l aGl uGl uGl nAsp●●●
 ApaLI (3688)