

# pBROAD3-mcs

An optimized vector for mouse and rat transgenesis

Catalog # pbroad3

For research use only

Version # 05A12-SV

## PRODUCT INFORMATION

### Content:

- 20 µg of pBROAD3-mcs provided as lyophilized DNA

### Storage and Stability:

- Products are shipped at room temperature.
- Lyophilized DNA is stable for 1 year when stored at -20°C
- Resuspended DNA is stable for 6 months at -20°C.

### Quality control:

- Plasmid construct has been confirmed by restriction analysis and sequencing.
- Plasmid DNA were purified by ion exchange chromatography and lyophilized.

## GENERAL PRODUCT USE

The pBROAD3-mcs plasmid was designed for expression of a transgene in virtually all tissues of transgenic mice and rats. This feature is brought by the ROSA promoter. The murine ROSA26 promoter was initially identified by random retroviral gene trapping in mouse embryonic stem cells<sup>1</sup>. This high CpG content promoter was shown to drive ubiquitous expression of the human placental alkaline phosphatase and enhanced green fluorescent protein during embryonic and postnatal development in mouse and rat<sup>2</sup>.

A multiple cloning site (MCS) has been added downstream of the ROSA promoter for convenient cloning of your gene of interest. The MCS contains several restriction sites that are compatible with many other enzymes, thus facilitating cloning. Furthermore, the *E. coli* region is flanked on either side by the well cutting 8 bp-recognizing restriction enzyme *Pac* I that enables linearization and easy excision of the *E. coli* region.

## PLASMID FEATURES

- **mROSA prom:** This TATA-less promoter was found to be very effective *in vitro* in a very broad range of mammalian cell lines. The strength of the murine ROSA promoter is ascribed to the 10 potential Sp1 sites found within the CpG island extending from the core promoter to the first half of 5'untranslated region (5'UTR), the highest number of Sp1 sites ever recorded in any natural promoter. The 5'UTR contains an engineered intron of 350 bp which increases the transcription of the transgene<sup>3</sup>.
- **βGlo pAn:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription<sup>4</sup>.
- **pMB1 ori:** a minimal *E. coli* origin of replication to limit vector size but with the same activity as the longer Ori.
- **Amp:** The ampicillin resistance gene allows the selection of transformed *E. coli* carrying a pBROAD plasmid.

- **MCS:** The multiple cloning site contains the following restriction sites:

*Age* I, *Bsp* LU11I, *Bgl* II, *Hind* III, *Eco* RI, and *Msc* I

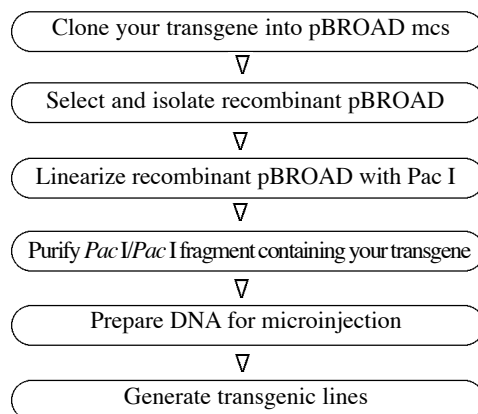
*Age* I is compatible with *Bsp* EI and *Sgr* AI.

*Bsp* LU11I is compatible with *Bsp* HI and *Nco* I.

*Bgl* II is compatible with *Bam* HI, *Bst* YI and *Bcl* I.

*Eco* RI is compatible with *Apo* I and *Mfe* I

## EXPERIMENTAL OUTLINE



## METHODS

### Plasmid resuspension:

Quickly spin the tube containing the lyophilized plasmid to pellet the DNA. To obtain a plasmid solution at 1µg/µl, resuspend the DNA in 20µl of sterile H<sub>2</sub>O. Store resuspended plasmid at -20°C.

### Pac I linearization of recombinant pBROAD:

1- Digest 10 µg recombinant pBROAD3 plasmid with 1 to 5 units of *Pac* I restriction enzyme.

**Note:** *Pac* I may be purchased from New England Biolabs and used at 0.1-0.5 unit per µg plasmid DNA.

2- Incubate at 37°C for 1-2 hours.

3- Purify the fragment containing the ROSA26 prom-transgene-βGlo pAn cassette by agarose gel following your usual protocol.

### References:

1. Zambrowicz BP, et al. 1997. Disruption of overlapping transcripts in the ROSA beta geo 26 gene trap strain leads to widespread expression of beta-galactosidase in mouse embryos and hematopoietic cells. *Proc Natl Acad Sci USA*. 94:3789-94.
2. Kisseberth WC, et al. 1999. Ubiquitous expression of marker transgenes in mice and rats. *Dev Biol*. 214:128-38.
3. Brinster RL, et al. 1988. Introns increase transcriptional efficiency in transgenic mice. *Proc Natl Acad Sci USA* 85(3):836-40
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.

## TECHNICAL SUPPORT

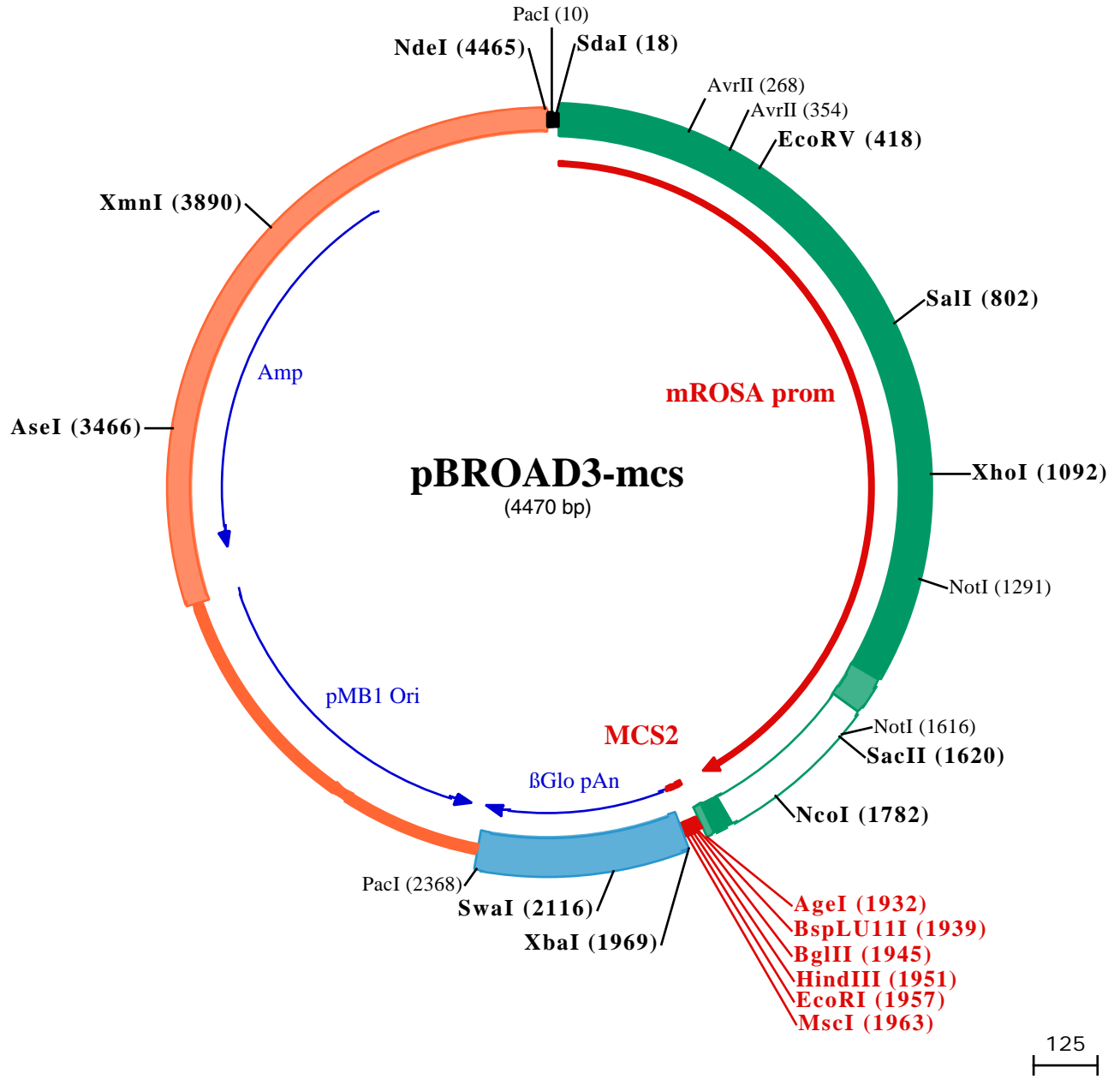
InvivoGen USA (Toll-Free): 888-457-5873

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



PacI (10) SdaI (18)  
1 GATCTCGACTTAATTAACCTGCAGGTGAAGACGTTACACAAGTAACATGAGAAAGCAGAAAAATGCAGGTCATCCACGCACCCTGACCCAGGCCAGCAGG  
101 GCGGGCTGCAGCATCAGTACACAGGAGAAAGATCCTTATTCTAAGAATGAGAAAGGCAAAGGCGCCGATAGAATAAATTAGCATAGAAGGGCTTTCC  
AvrII (268)  
201 CAGGAGTAAAACCTTCTTCTGAGCGATTACCTACTAAAACAGGGCTTTTGCCACTACCATTACCTAGGATCTTGCTTGCACGGATTCATAGGGG  
AvrII (354)  
301 CATATCCCTCCCCTCTTCTTAGAGTCGTTCTTAAAAGATCGCTCTCCACGCCCTAGGCAGGGAAAAACGACAAAATCTGGCTCAATTCAGGCTAGAAC  
EcoRV (418)  
401 CCTACAAATTCACAGGGATATCGCAAGGATACTGGGGCATACGCCACAGGGAGTCCAAGAATGTGAGGTGGGGGTGGCGAAGTAATGTCTTTGGTGTG  
501 GGAAGCAGCAGCCATCTGAGATAGGAAGTGGAAAACAGAGGAGGCGTTCAGGAAGATTATGGAGGGGAGGACTGGCCCCACGAGCGACCAGAG  
601 TTGTCACAAGGCCGAAGAACAGGGAGGTGGGGGCTCAGGGACAGAAAAAAGTATGTGTATTTGAGAGCAGGGTTGGGAGGCTCTCTGAAAAG  
701 GGTATAAACGTGGAGTAGGCAATACCAGGCAAAAAGGGAGACCAGAGTAGGGGGAGGGGAAGAGTCTGACCCAGGGAAGACATTA AAAAGGTAGTGG  
SalI (802)  
801 GGTTCGACTAGATGAAGGAGAGCCTTCTCTCTGGCAAGAGCGGTGCAATGCTGTGTAAGGTAGCTGAGAAGACGAAAAGGCAAGCATCTTCTGCTA  
901 CCAGGCTGGGGAGGCCAGGCCACGACCCCGAGGAGAGGGAACGCAGGAGACTGAGGTGACCCTTCTTTCCCGGGGGCCGGTCTGTGTGTTGCTG  
XhoI (1092)  
1001 TCTCTTTTCTGTTGGACCCTTACCTTGACCCAGGCGCTGCCGGGCTGGGCCCGGGTGGCGGCACGGCACTCCCGGAGGCAGCGAGACTCGAGTTA  
1101 GGCCCAACGCGGCGCCACGGCGTTTCTGCGCGGAATGGCCGTACCCGTGAGGTGGGGTGGGGGCAGAAAAGCGGAGCGAGCCGAGCGGGGAG  
NotI (1291)  
1201 GGGGAGGGCCAGGGCGGAGGGGCCGCGACTACTGTGTTGGCGGACTGGCGGACTAGGGTGCCTGAGTCTCTGAGCGAGCGGGCGCGCCGCC  
1301 CTCCCCGGCGGCGGACGCGGCGGACGCGGCGGAGCTACTCAGCCCGTGCCTGAGCGGAAACGCCACTGACCGACGGGGATTCCCAGTGCCGGCCG  
1401 CAGGGGCAGCGGGACAGCCCCCTCCGCGCGCATTGGCCTCTCCGCCACCGCCACACTTATTGGCCGGTGCGCCGCAATCAGCGGAGGCTGC  
1501 CGGGGCCGCTAAAGAAGAGGCTGTGCTTTGGGGCTCCGGCTCCTCAGAGAGCTCGGCTAGgtagggatcgggactctggcgggagggcggttggtg  
SacII (1620)  
NotI (1616)  
1601 cgtttgccccgatggcgcccgccgagccctccgagcgtggtggagccgttctgtgagacagccgggtacgagtcgtgacgctggaaggggcaagcgg  
NcoI (1782)  
1701 gtggtgggcaggaatgcggtccgcccctgcagcaaccggagggggagggagaaggagcggaaaagtctccaccggacgcggccatggctcgggggggggg  
1801 gggcagcggaggagcgttccggccgacgtctcgtcgtgatggcttcttttctccgcccgtgtgtgaaaacacaattgtactaaccttctctcttt  
HindIII (1951)  
BspLU11I (1939) MscI (1963)  
AgeI (1932) BglII (1945) EcoRI (1957) XbaI (1969)  
1901 cctctcctgacagGTGTGAAACAGGAAGAGAACCGGTGACATGTAGATCTAAGCTTGAATTC TGCCATCTAGAAGCTCGCTTCTTGCTGTCCAATTC  
2001 TATTAAGGTTCTTTGTTCCCTAAGTCCAACACTAACTGAGGGATATTATGAAGGCCCTTGAGCATCTGGATTCTGCCTAATAAAAAACATTTATT  
SwaI (2116)  
2101 TCATTGCAATGATGTATTTAAATTATTTCTGAATATTTTACTAAAAAGGGAATGTGGGAGGTCAGTGCATTTAAAACATAAAGAAATGAAGAGCTAGTTC  
2201 AAACCTTGGGAAAATACACTATATCTTAACTCCATGAAAAGGTTGAGGCTGCAAACAGCTAATGCACATTGGCAACAGCCCCTGATGCCTATGCCTTA  
PacI (2368)  
2301 TTCATCCCTCAGAAAAGGATTCAAGTAGAGGCTTGATTGGAGGTTAAAGTTTTGCTATGCTGTATTTAATTA AAAACCCGCTTCGGCGGTTTTTTTA  
2401 TGCATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCAAAAA  
2501 ATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTGCGCTCTCTGTTCCGACCCT  
2601 GCCGCTTACCGGATACCTGTCCGCTTCTCCCTTCGGGAAGCGTGGCGTTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGTCTGTTCCG  
2701 TCCAAGCTGGGCTGTGTGCACGAACCCCGTTCCAGCCGACCGTGGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGTAAGACAGCACTTAT  
2801 CGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGTGGCCTAACTACGGCTACACTAG  
2901 AAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAACAAACCACCGTGGTAGCGGT  
3001 GGTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGCTGACGCTCAGTGGAAACGAAA

3101 ACTCACGTTAAGGGATTTTGGTCATGAGATTATCAAAAAGGATCTTCACCTAGATCCTTTTAAATTA AAAATGAAGTTTAAATCAATCTAAAGTATATA

3201 TGAGTAAACTTGGTCTGACAGTTACCAATGCTTAATCAGTGAGGCACCTATCTCAGCGATCTGTCTATTTTCGTTTCATCCATAGTTGCCTGACTCCCCGTC  
287 ••• TrpHisLysI leLeuSerAlaGlyI leGluAlaI leGlnArgAsnArgGluAspMetThrAlaGlnSerGlyThrT

3301 GTGTAGATAACTACGATACGGGAGGGCTTACCATCTGGCCCCAGTGCTGCAATGATACCGCGAGACCCACGCTCACCGGCTCCAGATTTATCAGCAATAA  
260 hrTyrI leValVal I leArgSerProLysGlyAspProGlyLeuAlaAlaI leI leGlyArgSerGlyArgGluGlyAlaGlySerLysAspAlaI lePh  
AseI (3466)

3401 ACCAGCCAGCCGGAAGGGCCGAGCGCAGAAGTGGTCTGCAACTTTATCCGCCTCCATCCAGTCTATTAATTGTTGCCGGAAGCTAGAGTAAGTAGTTC  
227 eTrpGlyAlaProLeuAlaSerArgLeuLeuProGlyAlaValLysAspAlaGluMetTrpAspI leLeuGlnGlnArgSerAlaLeuThrLeuLeuGlu

3501 GCCAGTTAATAGTTTGGCAACGTTGTTGCCATTGCTACAGGCATCGTGGTGTACGCTCGTCTGTTGGTATGGCTTCATTCAGCTCCGGTCCCAACGA  
194 GlyThrLeuLeuLysArgLeuThrThrAlaMetAlaValProMetThrThrAspArgGluAspAsnProI leAlaGluAsnLeuGluProGluTrpArgA

3601 TCAAGCGAGTTACATGATCCCCATGTTGTGCAAAAAAGCGGTTAGCTCCTTCGGTCTCCGATCGTTGTGAGAAGTAAGTTGCCCGCAGTGTATCAC  
160 spLeuArgThrValHisAspGlyMetAsnHisLeuPheAlaThrLeuGluLysProGlyGlyI leThrThrLeuLeuLeuAsnAlaAlaThrAsnAspSe

3701 TCATGGTTATGGCAGCACTGCATAATTCTCTTACTGTCATGCCATCCGTAAGATGCTTTCTGTGACTGGTGAGTACTCAACCAAGTCATTCTGAGAATA  
127 rMetThrl leAlaAlaSerCysLeuGluArgValThrMetGlyAspThrLeuHisLysGluThrValProSerTyrGluValLeuAspAsnGlnSerTyr  
XmnI (3890)

3801 GTGTATGCGGCGACCGAGTTGCTCTTGGCCGGCGTCAATACGGGATAATACCGCGCCACATAGCAGAACTTTAAAAGTGCTCATCATTGGAAAACGTTCT  
94 HisI leArgArgGlyLeuGlnGluGlnGlyAlaAspI leArgSerLeuValAlaGlyCysLeuLeuValLysPheThrSerMetMetProPheArgGluG

3901 TCGGGCGAAAACCTCTCAAGGATCTTACCGCTGTTGAGATCCAGTTCGATGTAACCCACTCGTGCACCCAACCTGATCTTCAGCATCTTTTACTTTCACCA  
60 luProArgPheSerGluLeul leLysGlySerAsnLeuAspLeuGluI leTyrGlyValArgAlaGlyLeuGlnAspGluAlaAspLysValLysValLe

4001 CCGTTTCTGGGTGAGCAAAAACAGGAAGGCAAAATGCCGCAAAAAGGGAATAAGGGCGACACGGAAATGTTGAATACTCATACTCTTCCTTTTCAATA  
27 uThrGluProHisAlaPheValProLeuCysPheAlaAlaPhePheProI leLeuAlaValArgPheHisGlnI leSerMet

4101 TTATTGAAGCATTATCAGGGTTATTGTCTCATGAGCGGATACATATTGAATGTATTTAGAAAAATAAACAAATAGGGGTTCCGCGCACATTTCCCGGA  
4201 AAAGTGCCACCTGACGTCTAAGAAACCATTATTATCATGACATTAACCTATAAAAAATAGGCGTATCACGAGGCCCTTTCGTCTCGCGGTTTCGGTGATG  
4301 ACGGTGAAAACCTCTGACACATGCAGCTCCCGGAGACGGTACAGCTTGTCTGTAAGCGGATGCCGGAGCAGACAAGCCGTCAGGGCGGTCAGCGGG  
NdeI (4465)

4401 TGTTGGCGGGTGTCCGGGCTGGCTTAACTATGCGGCATCAGAGCAGATTGTACTGAGAGTGCACCATATG