

pBLAST42-hAngio

An anti-angiogenic plasmid expressing the human Angiostatin gene

Catalog # pbla-hangio

For research use only

Version # 05L21-SV

PRODUCT INFORMATION

Content:

- 20 µg of lyophilized pBLAST42-hAngio plasmid DNA.
- 4 pouches of *E. coli* Fast-Media® Blas (2 for agar media, 2 for liquid media).

Storage and stability:

- Products are shipped at room temperature.
- Upon receipt, resuspend lyophilized DNA and store at -20°C. Avoid repeated freeze-thaw cycles.
- Store *E. coli* Fast-Media® Blas at room temperature. Fast-Media® pouches are stable 18 months when stored properly.

Quality control:

- Plasmid DNA was prepared using affinity column and lyophilized.
- Plasmid construct has been confirmed by restriction analysis sequencing.

GENERAL PRODUCT USE

Anti-angiogenic proteins (also called angiostatic) are utilized in cancer therapy to inhibit tumor angiogenesis, i.e. the formation of blood vessels that feed tumor cells. However, angiostatic proteins are difficult to obtain and purify. Therefore, the use of **pBLAST**, a family of plasmids expressing human or murine angiostatic genes should simplify the production of angiostatic proteins *in vitro* and *in vivo*.

pBLASTs may be used for:

- convenient production of angiostatic proteins *in vitro*
- anti-angiogenic gene therapy

pBLAST allows a high level of expression and secretion of the gene product: genes coding for secreted proteins are cloned with their native signal sequence. However, proteins that lack a natural signal sequence (e.g. angiostatin) are engineered for secretion by addition of the hIL-2 signal sequence.

pBLAST carries a single antibiotic resistance gene, blasticidin, which allows very rapid and convenient selection of both bacteria and mammalian cell transformants.

PLASMID FEATURES

- **EF-1α / HTLV hybrid promoter** is a composite promoter comprised of the Elongation Factor-1α (EF-1α) promoter¹ and 5' untranslated region of the Human T-Cell Leukemia Virus (HTLV). EF-1α is an 'housekeeping' gene ubiquitously expressed in eukaryotic cells. The EF-1α promoter exhibits a strong activity, higher than viral promoters and, on the contrary to the CMV promoter, yields persistent expression of the transgene *in vivo*. The R segment and part of the U5 sequence (R-U5') of the HTLV Type 1 Long Terminal Repeat² has been coupled to the EF-1α promoter to enhance stability of DNA and RNA. This modification not only increases steady state transcription, but also significantly increases translation efficiency possibly through mRNA stabilization.
- **III7** is a bacterial promoter that is spliced out as an intron in mammalian cells. Expression of the transgene in *E. coli* can be further increased by the addition of IPTG when working with bacteria that constitutively express *Lacl*
- **SV40 polyA**: The Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA. The efficiency of this signal was first described by Carswell et al.³
- **SpAn**: A synthetic polyadenylation site and a strong pause site are placed downstream of the pMB1 Ori to limit transcriptional interference between both transcriptional units. The synthetic polyA site is based on the highly efficient polyA signal of the rabbit β-globin gene⁴.

• **Angiostatic gene hAngiostatin**

Size: 1155 bp.

- **pMB1 Ori** is a minimal *E. coli* origin of replication with the same activity as the longer Ori.
- **CMV prom:** The CMV promoter allows the expression of the blasticidin resistance gene in mammalian cells.
- **Bsr (blasticidin resistance gene):** The *bsr* gene from *Bacillus cereus* encodes a deaminase that confers resistance to the antibiotic Blasticidin S. The *bsr* gene is driven by the SV40 promoter in tandem with the bacterial EM7 promoter. Therefore each pBLAST plasmid can be used to select stable mammalian cells transfectants and *E. coli* transformants.
- **bGh pAn:** The bovine growth hormone (bGh) polyadenylation (pAn) signal and a transcriptional pause are placed 3' of the blasticidin gene. The bGh pAn has been shown to be as efficient as SV40 and HSV1tk polyadenylation signals in many different cell types⁵. The use of bGh pAn minimizes interference and possible recombination events with the SV40 polyadenylation signal. The pause site prevents transcriptional interference and read-through.

References

- 1- Kim et al (1990). Gene 2: 217-223.
- 2- Takebe et al (1988). Mol. Cell Biol. 1: 466-472.
- 3- Carswell et al (1989). Mol. Cell Biol. 10: 4248-4258.
- 4- Levitt et al. (1989). Genes Dev. 7: 1019-1025.
- 5- Goodwin et al. (1992). J. Biol. Chem. 23: 16330-16334.

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 H₂O. Store resuspended plasmid at -20°C.

Selection of bacteria with *E. coli* Fast-Media Blas:

E. coli Fast-Media® Blas is a **new, fast and convenient** way to prepare liquid and solid media for bacterial culture by using only a microwave. *E. coli* Fast-Media® Blas is a TB (liquid) or LB (solid) based medium with blasticidin, and contains stabilizers. *E. coli* Fast-Media® Blas can be ordered separately (catalog code # fas-bl-l, fas-bl-s).

Method:

- 1- Pour the contents of a pouch into a clean borosilicate glass bottle or flask.
- 2- Add 200 ml of distilled water to the flask
- 3- Heat in a microwave on MEDIUM power setting (about 400Watts), until bubbles start appearing (approximately 3 minutes). **Do not heat a closed container. Do not autoclave Fast-Media®.**
- 4- Swirl gently to mix the preparation. **Be careful, the bottle and media are hot, use heatproof pads or gloves and care when handling.**
- 5- Reheat the media for 30 seconds and gently swirl again. Repeat as necessary to completely dissolve the powder into solution. But be careful to avoid overboiling and volume loss.
- 6- Let agar medium cool to 45°C before pouring plates. Let liquid media cool to 37°C before seeding bacteria.

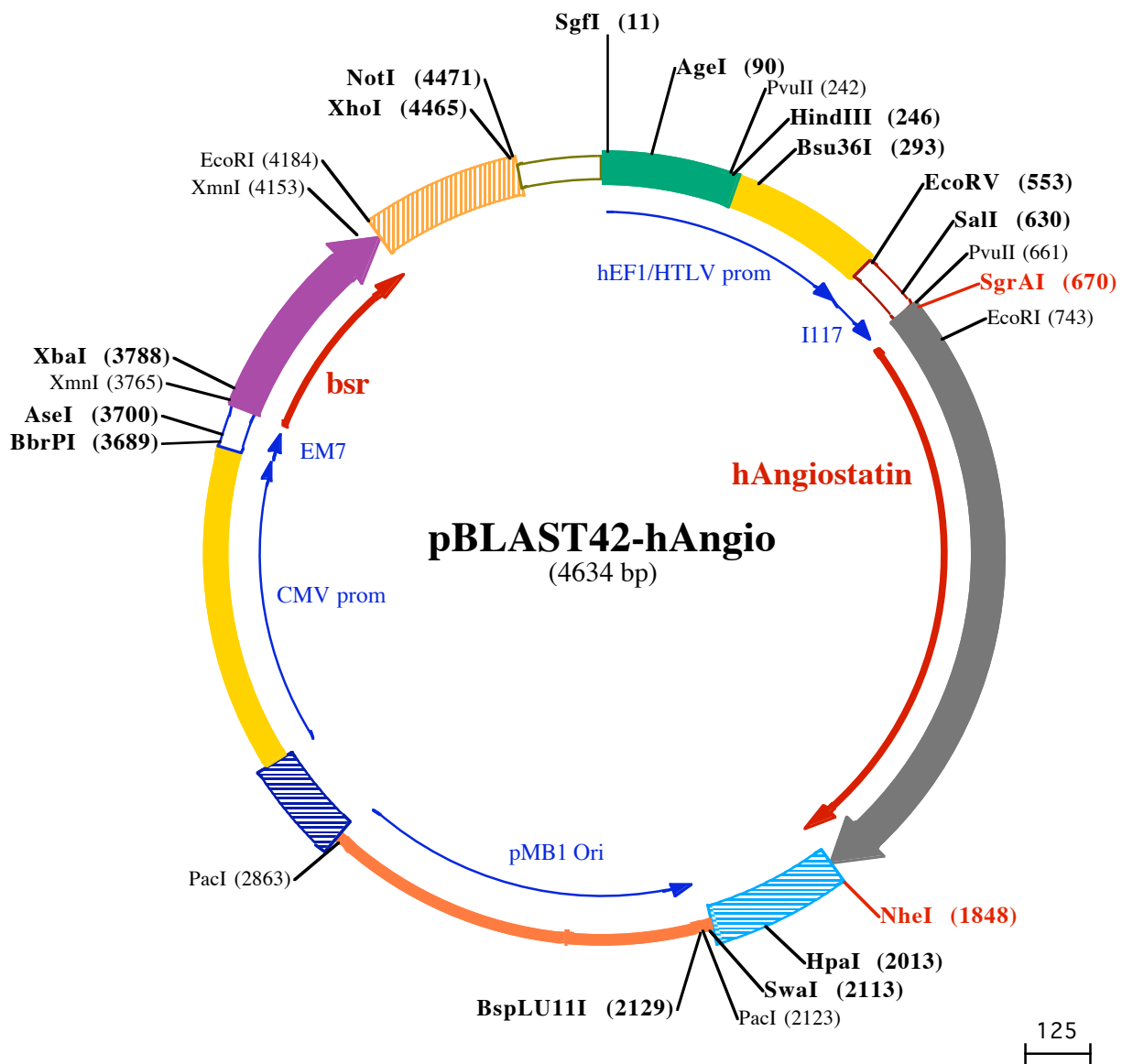
Note: Do not reheat solidified Fast-Media® as the antibiotic will be permanently destroyed by the procedure.

TECHNICAL SUPPORT

Toll free (US): 888-457-5873
Outside US: (+1) 858-457-5873
E-mail: info@invivogen.com
Website: www.invivogen.com

InvivoGen

3950 Sorrento Valley Blvd. Suite A
San Diego, CA 92121 - USA



SgfI (11) AgeI (90)
1 GGATCTGCGATCGCTCCCGTCCCGTCCAGTGGCGAGAGCGCACATGCCCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGCGCAATTGAACCGGTGCCTA
101 GAGAAGGTGGCGCGGGTAAACTGGGAAAGTGATGTCGTGACTGGCTCCGCTTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (246) PvuII (242) Bsu36I (293)
201 GTGAACGTTCTTTTTTCAACGGGTTTGGCCGCAAGACACAGCTGAAGCTTCAGAGGGCTCGCATCTCTCCTTACGCGCCCGCCCTACCTGAGGCC
301 GCCATCCACGCGGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCTGAACTGCGTCCGCGCTAGGTAAGTTTAAAGCTCAGGTCGAGACC
401 GGGCTTTGTCCGGCTCCCTTGAGCTACCTAGACTCAGCGGCTCTCCACGCTTTCCTGACCCTGCTTGTCAACTCTACGTTCTTTGTTTCGTTT

EcoRV (553)
501 TCTGTTCTGCGCGTTACAGATCAAGCTGTGACCGCGCTACGtaagtgatactactagatttatcaaaagagtgttgacttgtgagcgctcaca

SalI (630) PvuII (661) SgrAI (670)
601 ttgatacttagattcatcgagaggacgctcgactactaaccttcttcttcttctctacagCTGAGATCACCGGCGAAGGAGGGCCACCATGTACAGGAT
1MetTyrArgMe

EcoRI (743)
701 GCAACTCTGTCTTGCATTGCACATAAGTCTTGCACCTTGTCAAGTTCGGTGTATCTCTCAGAGTGCAGACTGGGAATGGAAGAACTACAGAGGGACG
4tGlnLeuLeuSerCysIleAlaLeuSerLeuAlaLeuValThrAsnSerValTyrLeuSerGluCysLysThrGlyAsnGlyLysAsnTyrArgGlyThr
801 ATGTCAAAAACAAAAATGGCATCACCTGTCAAAAATGGAGTTCACCTTCTCCACAGACCTAGATTCTCACCTGCTACACCCCTCAGAGGGACTGG
38MetSerLysThrLysAsnGlyIleThrCysGlnLysTrpSerSerThrSerProHisArgProArgPheSerProAlaThrHisProSerGluGlyLeuG
901 AGGAGAAGTACTGCAGGAATCCAGACAACGATCCGAGGGGCGCTGGTGTATATACTACTGATCCAGAAAAGAGATATGACTACTGCGACATCTTGTAGTG
71luGluAsnTyrCysArgAsnProAspAsnAspProGlnGlyProTrpCysTyrThrThrAspProGluLysArgTyrAspTyrCysAspIleLeuGluuCy
1001 TGAAGAGGAATGTATGCATTGCAGTGGAGAAAATATGACGGCAAAATTTCCAAGACCATGTCTGGACTGGAATGCCAGGCTGGGACTCTCAGAGCCCA
104sGluGluGluCysMetHisCysSerGlyGluAsnTyrAspGlyLysIleSerLysThrMetSerGlyLeuGluCysGlnAlaTrpAspSerGlnSerPro
1101 CACGCTCATGGATACATTCCTTCAAATTTCCAACAAGAACCTGAAGAAGAAATTAAGTGTCTGTAACCCGATAGGAGCTGCGGCTTGGTGTTCACCA
138HisAlaHisGlyTyrIleProSerLysPheProAsnLysAsnLeuLysLysAsnTyrCysArgAsnProAspArgGluLeuArgProTrpCysPheThrT
1201 CCGACCCCAACAGCGCTGGAACTTTGTGACATCCCGCTGCACAACACCTCCACCATCTTGTGTCACCTACCAGTGTCTGAAGGGAACAGGTGA
171hrAspProAsnLysArgTrpGluLeuCysAspIleProArgCysThrThrProProProSerSerGlyProThrTyrGlnCysLeuLysGlyThrGlyY
1301 AAACATATCGGGGAATGTGGCTGTACCGTGTCCGGGCACACCTGTGACAGCTGGAGTGCACAGACCCCTCACACACATAACAGGACCCAGAAAATTC
204uAsnTyrArgGlyAsnValAlaValThrValSerGlyHisThrCysGlnHisTrpSerAlaGlnThrProHisThrHisAsnArgThrProGluAsnPhe
1401 CCCTGCAAAAATTTGGATGAAAATCTACTGCCGCAATCTGACGAAAAAGGGCCCATGGTGCATACAACCAACAGCCAAAGTGGGTTGGGAGTACTGTA
238ProCysLysAsnLeuAspGluAsnTyrCysArgAsnProAspGlyLysArgAlaProTrpCysHisThrThrAsnSerGlnValArgTrpGluTrpCysL
1501 AGATACCGTCTGTGACTCTCCAGTATCCACGGAACAATTTGGCTCCACAGCACCTGAGCTAACCCCTGTGGTCCAGGACTGCTACCATGGTGA
271ysIleProSerCysAspSerSerProValSerThrGluGlnLeuAlaProThrAlaProProGluLeuThrProValValGlnAspCysTyrHisGlyYAs
1601 TGGACAGAGTACCGAGGCACATCTCCACCACCACAGGAAAGAGTGTAGTCTTGGTCATCTATGACACCACCCGACCCAGAAAGACCCAGAA
304pGlyGlnSerTyrArgGlyThrSerSerThrThrThrGlyLysLysCysGlnSerTrpSerSerMetThrProHisArgHisGlnLysThrProGlu
1701 AACTACCCAAATGCTGGCCTGACAATGAACTACTGCAAGATCCAGATGCCGATAAAGGCCCTGGTGTTTTACCACAGACCCAGCGTCAAGTGGGAGT
338AsnTyrProAsnAlaGlyLeuThrMetAsnTyrCysArgAsnProAspAlaAspLysGlyProTrpCysPheThrThrAspProSerValArgTrpGluT
NheI (1848)
1801 ACTGCAACCTGAAAAATGCTCAGGAACGAGAAGCGAGTGTGTATAGCTAGCTCGCATGATAAGATACATTGATGAGTTTGGACAAACACAACCTAGA
371yrCysAsnLeuLysLysCysSerGlyThrGluAlaSerValVal●●●
1901 ATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCTATAAGCTG
HpaI (2013)
2001 CAATAAACAAGTTAACAACAATTCATTCTTTATGTTTCAGGTTTCAGGGGAGGTGGGGAGTTTTTAAAGCAAGTAAACCTCTACAATGT
PaeI (2123)
2101 GGTAGATCCATTTAAATGTTAATTAAGAATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCTAAAAAGCCGCTTGTGGCTTTTTCCATAGGC
2201 TCCGCCCCCTGACGAGCATCAAAAATCGACGCTCAAGTCAGAGGTGGGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTTGGAAAGCTC
2301 CCTCGTGCCTCTCTGTTCCGACCTTCCGCTTACCGGATACCTGTCGCTTCTCCCTTGGGAAGCGTGGCGTTTCTCATAGCTCACGCTGAGG
2401 TATCTCAGTTCGGTGTAGTCTGTTCCGCTCAAGCTGGGCTGTGTGCAGAACCCCGCTTACGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTG
2501 AGTCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCTGTTAAGGATTAGCAGAGCGAGGTATGAGGCGGTGCTACAGAGTTCTTGAA
2601 GTGGTGGCTAACTACGGCTACTAGAAGAACGATTTGGTATCTCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTTTGATCC
2701 GGCAACAACCCCGCTGGTAGCGGTTTTTTTTGTTTGAAGCAGCAGATTACGGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTA
PaeI (2863)
2801 CGGGTCTGACGCTCAGTGGAACTCACGTTAAGGGATTTGGTCTAGTAAATTAAGCTGCAATAAACAATCATTATTTTTCATTGGATCT
2901 GTGTGTTGTTTTTTGTGTGGCTTGGGGAGGGGAGCCAGAATGACTCAAGAGCTACAGGAAGCCAGGTGAGAGACCCACTGGACAAACAGTGGC
3001 TGGACTCTGCACATAACACAATCAACAGGGGAGTGAAGTGGATCGAGCTAGAGTCCGTTACATAACTTACGGTAAATGGCCGCTGGCTGACCGCC
3101 CAACGACCCCGCCATTGACGTCATAATGACGTATGTTCCATAGTAACGCAATAGGGACTTTCCATTGACGTCATGGTGGAGTATTTACGGTAA
3201 ACTGCCACTTGGCAGTACATCAAGGTATCATATGCCAAGTACGCCCTTATTGACGTCATGACGGTAAATGGCCGCTGGCATTATGCCAGTACA
3301 TGACCTTATGGGACTTCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGGTGTGCGGTTTTGGCAGTACATCAATGGGCGTGGATA
3401 GCGGTTTGTACTCACGGGATTTCCAAGTCTCCACCCATTGACGTCATGGGAGTTTGTTTGGACCAAAATCAACGGGACTTTCAAAAATGCTGTAAC
3501 AACTCCGCCCCATTGACGCAAAATGGCGGTAGCGGTGACGGTGGGAGGTCTATATAAGCAGAGCTGTTTGTGAAACCGTCAAGTCCGCTGGAGACGCC
BbrPI (3689) AseI (3700)
3601 ATCCACGCTTTTTGACCTCCATAGAAGACCCGGACCGATCCAGCTCCGCGGCGGGAACGGTGCATTGGAACCGACCTGCAGCAGCTGTTGACAA

3701 TAATCATCGGCATAGTATATCGGCATAGTATAATACGACTCACTATAGGAGGGCCACCATGAAGACCTTCAACATCTCCAGCAGGATCTAGAATTAGTA
XmnI (3765)
XbaI (3788)
1▶MetLysThr PheAsnI l eSer Gl nGl nAspLeuGl uLeuVal
 3801 GAAGTAGCGACAGAGAAGATTACAATGCTTTATGAGGATAATAAACATCATGTGGGAGCGGCAATTCGTACGAAAACAGGAGAAATCATTTCGGCAGTAC
15▶Gl uValAl aThr Gl uLysI l eThr Me tLeuTyrGl uAspAsnLysHi sHi sVal Gl yAl aAl al l eArgThr LysThr Gl yGl u l l e l l eSer Al aVal H
 3901 ATATTGAAGCGTATATAGGACGAGTAACTGTTTGTGAGAAGCCATTGCGATTGGTAGTGCAGTTTCGAATGGACAAAAGGATTTTGACACGATTGTAGC
48▶i s l l eGl uAl aTyr l l eGl yArgVal Thr Val CysAl aGl uAl al l eAl al l eGl ySer Al aVal SerAsnGl yGl nLysAspPheAspThr l l eValAl
 4001 TGTTAGACACCCTTATTCTGACGAAGTAGATAGAAGTATTCGAGTGGTAAGTCCTTGTGGTATGTGTAGGGAGTTGATTTAGACTATGCACCAGATTGT
81▶aValArgHi sProTyrSerAspGl uValAspArgSer l l eArgVal Val Ser ProCysGl yMetCysArgGl uLeu l l eSerAspTyrAl aProAspCys
XmnI (4153)
EcoRI (4184)
 4101 TTTGTGTTAATAGAAATGAATGGCAAGTTAGTCAAACTACGATTGAAGAAGTCTATTCCACTCAAATATACCCGAAATTAAGAAATTCGCTAGAGGGCC
115▶PheVal l Leul l eGl uMe tAsnGl yLysLeuVal l LysThr Thr l l eGl uGl uLeu l l eProLeuLysTyrThr ArgAsn••••
 4201 CTATTCTATAGTGTACCTAAATGCTAGAGCTCGCTGATCAGCCTCGACTGTGCCTTCTAGTTGCCAGCCATCTGTTGTTTGCCCTCCCCCGTGCCTTC
 4301 CTTGACCCTGGAAGGTGCCACTCCACTGTCTTTCTAATAAAATGAGGAAATTGCATCGCATTGTCTGAGTAGGTGCATTCTATTCTGGGGGTGGG
NotI (4471)
XhoI (4465)
 4401 GTGGGGCAGGACAGCAAGGGGGAGGATTGGGAAGACAATAGCAGGCATGCGCAGGGCCCAATTGCTCGAGCGGCCCGCAATAAAATATCTTTATTTTCATT
 4501 ACATCTGTGTGTTGTTTTTTGTGTGAATCGTAACATAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAATAGGGTGTCCCCAGT
 4601 GCAAGTGCAGGTGCCAGAACATTCTCTATCGAA