

Product usage

Before using this product, please read the Limited Use statement below

Important Limited Use information for pTiGer2-eGFP

The purchase of the pTiGer2-eGFP vector conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes.

The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes.

Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic, or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research.

If the purchaser is unwilling to accept the limitations of this limited use statement, InvivoGen is willing to accept return of the product with a full refund. The product must be returned in resaleable condition. For information on purchasing a license to this product for purposes other than research, contact us at outlicensing@invivogen.com.

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 Asia: +852 3622-3480

E-mail: info@invivogen.com



pTiGer2-eGFP

A multigenic plasmid for inducible enhanced green fluorescent protein expression, selectable with Zeocin®

Catalog code: ptg2-gfp

<https://www.invivogen.com/tet-on-ptiger-reporter>

For research use only

Version 24A16-NJ

PRODUCT INFORMATION

Contents:

- 20 µg of pTiGer2-eGFP provided as lyophilized DNA
- 1 ml of Zeocin® (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 for up to 1 year.
- Store Zeocin® at 4 °C or at -20 °C. The expiry date is specified on the product label.

Quality control

- Plasmid construct is confirmed by restriction analysis and full-length open reading frame (ORF) sequencing.
- After purification by ion exchange chromatography, predominant supercoiled conformation is verified by electrophoresis.

PRODUCT DESCRIPTION

InvivoGen provides a family of plasmids featuring a tetracycline-inducible reporter gene. The pTiGer2-eGFP plasmid encodes the enhanced green fluorescent protein (eGFP) and the Zeocin® resistance marker for selection in both mammalian cells and bacteria. This plasmid can be used as a transfection control for plasmids of the pTiGer-mcs family.

The eGFP expression is only possible upon transfection of cells featuring the tetracycline repressor (TetR) protein¹, such as InvivoGen's HEK-RepTor™ or A549-RepTor™ cells. These cells express TetR constitutively in the nucleus, where it binds to tetracycline operator (tetO) sequences and represses gene transcription. Upon incubation with doxycycline (a synthetic tetracycline derivative), TetR is released from the tetO sequences and the eGFP-encoding gene is transcribed.

pTiGer-SEAP and pTiGer-Lucia plasmids are also available.

PLASMID FEATURES

SEAP expression cassette

- **hCMV enh/ hEF1 prom:** This composite promoter combines the human cytomegalovirus (HCMV) enhancer and the core promoter of the human elongation factor-1 α (EF-1 α)².
- **tetOtetO:** This sequence is also known as TRE (Tetracycline Response Element). It is a repeat of the 19-nucleotide sequence of the tetracycline operator (tetO)¹.
- **eGFP (enhanced green fluorescent protein)** is a synthetic variant of the wild-type GFP. It features mutations that allow for better folding efficiency at 37°C and higher-intensity emission, compared to GFP. Moreover its codon sequence is optimized for expression in mammalian cells. The eGFP excitation peak is at ~ 490 nm, and its emission peak at ~ 509 nm.

- **SV40 pAn:** The Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA³.

Zeocin® antibiotic selection cassette

- **pMB1 Ori:** This minimal *E. coli* origin of replication with the same activity as the longer Ori.
- **hAldA enh/ hFerH prom:** This composite promoter combines the human aldehyde dehydrogenase (aldA) enhancer and the core promoter of the human ferritin heavy chain gene (FerH).
- **EM7:** This bacterial promoter enables the constitutive expression of the *Sh ble* gene in *E. coli*.
- **Sh ble:** The resistance to Zeocin® is conferred by the *Sh ble* gene from *Streptoalloteichus hindustanus*. The *Sh ble* gene is driven by the human AldA/FerH promoter in tandem with the bacterial EM7 promoter allowing selection in both mammalian cells and *E. coli*.
- **mEF1 5'UTR:** The 5'UTR (untranslated region) of the murine EF-1 α enhances *Sh ble*-encoding mRNA stability and protein translation.
- **hβGlo pAn:** The human β -Globin pAn is a strong polyadenylation signal placed downstream of *Sh ble*⁴.

1. Hillen, W., Wissmann, A. (1989). Tet repressor-tet operator interaction. Protein-Nucleic Acid Interaction. DOI: 10.1007/978-1-349-09871-2_7. 2. Kim DW, et al., 1990. Use of the human elongation factor 1 α promoter as a versatile and efficient expression system. Gene 91(2):217-23. 3. Carswell S. & Alwine JC., 1989. Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. Mol Cell Biol. 9(10):4248-58. 4. Yu J. & Russell JE., 2001. Structural and functional analysis of an mRNP complex that mediates the high stability of human β -globin mRNA. Mol Cell Biol. 21(17):5879-88.

METHODS

Plasmid resuspension

- Quickly spin the tube to pellet the DNA.
- To obtain a plasmid solution at 1 µg/µl, resuspend the DNA in 20 µl of sterile water.
- Store the resuspended plasmid at -20°C.

Plasmid amplification and cloning

Plasmid amplification and cloning can be performed in *E. coli* GT115 or other commonly used laboratory *E. coli* strains, such as DH5 α .

Zeocin® usage

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

Generation of Tet-inducible expression cells

For a general procedure using InvivoGen's RepTor™ cell lines, please visit: <https://www.invivogen.com/reptor-cells>.

RELATED PRODUCTS

Product	Description	Cat. Code
Zeocin®	Selection antibiotic	ant-zn-1
HEK-RepTor™ cells	TetR-expressing cells	hk-rtor

TECHNICAL SUPPORT

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

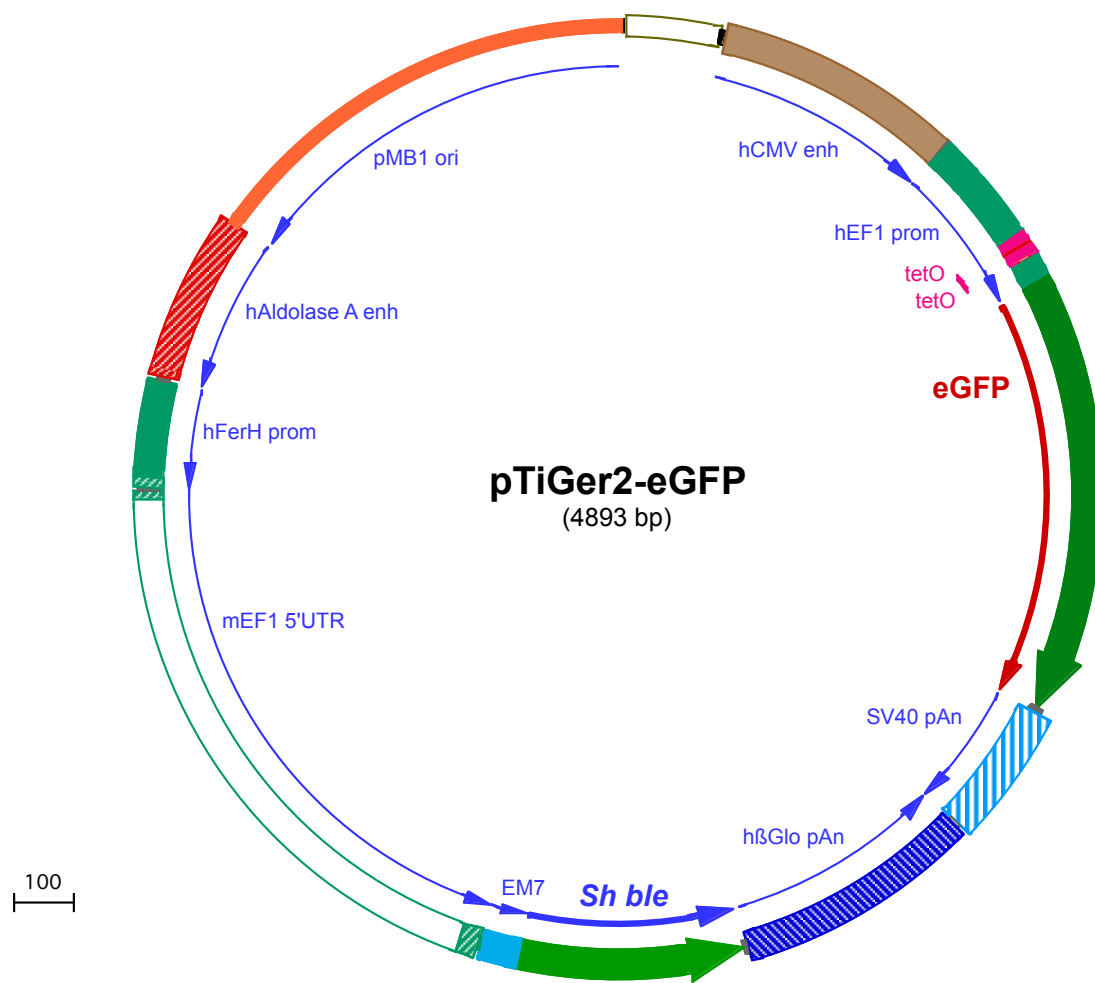
InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: info@invivogen.com

 **InvivoGen**
www.invivogen.com



1 CTCGAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACG
101 AAACAAAACAACTAGCAAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGACCTGCAGGCGTTACATAACTTACGGTA
201 AATGGCCCCGCTGGTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCATTGACGTC
301 AATGGGTGGAGTATTTACGGTAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC
401 CGCCTGGCATTATGCCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGATGATGCGGTTTTGG
501 CAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGTTTACTAGTCAGTGGCC
601 AGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGATCCGGTGCCTAGAGAAGTGGCGCGGGTAACTGGGAAAGTGAT
701 GTCGTGTACTGGCTCCGCTTTTCCGAGGTTGGGGGAGAACCCTATATAAGTGCAGTAGTTGCCGTGAACGTTTCCCTATCAGTGATAGAGATCTCCC
801 TATCAGTGATAGAGATCTTTCGAACGGGTTTCCGCCAGAACACAGCTGAAGCTTACCAGGTCACCATGGTGAAGCAAGGGCGAGGAGCTGTTACCAGGG
125 M V S K G E E L F T G
901 GTGGTGCCATCTGGTTCGAGCTGGACGGCGACGTAACGGCCACAAGTTACGCGTGTCCGGCAGGGCGAGGGCGATGCCACCTACGGCAAGCTGACCC
125 V V P I L V E L D G D V N G H K F S V S G E G E G D A T Y G K L T
1001 TGAAGTTCATCTGCACCACCGCAAGCTGCCGTGCCCTGGCCACCCTCGTACCACCTGACCTACGGCGTGCAGTGTTCAGCCGCTACCCCGACCA
45 L K F I C T T G K L P V P W P T L V T T L T Y G V Q C F S R Y P D H
1101 CATGAAGCAGCAGACTTCTTCAAGTCCGCCATGCCGGAAGGTACGTCCAGGAGCGACCATCTTCTTCAAGGACGACGGCAACTACAAGCCCGCGCC
78 M K Q H D F F K S A M P E G Y V Q E R T I F F K D D G N Y K T R A
1201 GAGGTGAAGTTCGAGGGCGACACCCTGGTGAACCGCATCGAGCTGAAGGGCATCGACTTCAAGGAGGACGGCAACATCCTGGGCGACAAGCTGGAGTACA
112 E V K F E G D T L V N R I E L K G I D F K E D G N I L G H K L E Y
1301 ACTACAACAGCCACAACGCTATATCATGCGCCACAAGCAGAAGAAGCGCATCAAGTGAACCTCAAGATCCGCCACAACATCGAGGACGGCAGCGTGCA
145 N Y N S H N V Y I M A D K Q K N G I K V N F K I R H N I E D G S V Q
1401 GCTCGCCGACCACTACCAGCAGAACACCCCATCGGCGACGGCCCGTGTGTGCCGACAACCACTACCTGAGCACCCAGTCCGCCCTGAGCAAAGAC
178 L A D H Y Q Q N T P I G D G P V L L P D N H Y L S T Q S A L S K D
1501 CCCAACGAGAAGCGGATCACATGGTCTGCTGGAGTTCGTGACCGCCCGGGATCACTCTCGGCATGGACGAGCTGTACAAGTAAAGCTAGCTGGCCA
212 P N E K R D H M V L L E F V T A A G I T L G M D E L Y K
1601 GACATGATAAGATACATTGATGAGTTTGACAAACCACAAC TAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTG
1701 TAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGTTTTTAAAGCAAGTA
1801 AAACCTCTACAAATGTGGTATGGAATCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATA
1901 AGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGACGCTCACCTCTTTTATGAGTTTAAAGATATAGTGTATTTTCCAAGGTTTGA
2001 ACTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTATGAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAA
2101 ATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAGGAACCTTTAATAGA
2201 AATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTCTGCCACAAAGTGCACGCAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCC
125 D Q E E A V F H V C N G A P D R L A F E
2301 CGCCCCACGGTGTCTCGCCGATCTCGGTATGCGCCGGCCGGAGGCGTCCCGGAAGTTCTGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCA
103 R G W P Q E G I E T M A P G S A D R F N T S V V E S W E A Y L E D L
2401 GGCCGCGACCCACCCAGGCGGTTGTCGGCACCACTGGTCTGGACCGCGCTGATGAACAGGGTACGTCGTCCCGGACCAACCCGGCGAA
70 G R V W V W A L T N D P V V Q D Q V A S I F L T V D D R V V G A F
2501 GTCGTCTCCACGAAGTCCCGGAGAACCAGCGGTCGGTCCAGAACTCGACCGTCCGGCGACGTCGCGCGGGTGAACCCGGAACGGCACTGGTC
37 D D E V F D R S F G L R D T W F E V A G A V D R A T L V P V A S T
2601 AACTTGGCCATGATGGCCCTCTATAGTGAAGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCATCCGGTTGCTTTGAATTAGCGGT
3 L K A M
2701 GGTTCCTACAACCTAAAAAGGGTTAAAAGATACCTTTGAACCGTAAGAAGCCCGAGAATTAGCTCCGCTCAAAACTCAAGGGGACAAATTCAAA
2801 AATGACTTCCAGCGCCAGGCTGGCCTGACTAGTCTCCACCCACCAATGTGAACAACTCCAACGCCATTACATCCCTCCCCCGCCGACTAGCCGT
2901 GCTCAAAAGCCGAGGTGACTATTGCGGCCGATAGGACCACGGGTACAGGAAGCAGCAGCCGGTGAAGGACCAGGCCCTCTTCTTTGTGGTGGTAC
3001 TCACCCGCCCCTCCACCGGCTGCCGCTCCTCATTGAGCTCCTTGAACAGGGCCCGGAGCGCCATCTTCCACGCACGCAACTGGTGCCGGA
3101 CGGGATGGCTCACCTAGTTAGGAGGAGGCAACGCGGCCCAAGCCAGATCGTCCGGTGTGGGACACATGGCTCGGCACGCTAACCC
3201 AGCCTGTTGCTTCGGGAAAAACCCAGGCTCGCCCATCCAGTGGCGTGGACATGTGCTCCGAAGGCGGGGCGCCAGCCGCACTCCTGTCCC

3301 TCCATTCTCCCAACCATGACCTCTCCGGGCTCCGGGCGAGCAAGCCCCGACCCCTCCCTTTGTTAGCCCTATTGCTGAACGGCAATCGAAGGCAGCA
3401 GGGCAACAACAACAAAAAAAAAAGACCAGAGTGCGGCCGGAGTAGCACGCGGGCGGGCGGACACCACGCTAGGCCTCAAGCCGGACACGAGGCGA
3501 GGCTACGGGGTTGCCGCTAGGCCTCGCACTCTGCCTCCCGCGCCGCCCAACTCGAAGCGGGAATGCTCGCAGCTAATCCCCGCCGACGACAGCGGGGC
3601 CCGGCCGCTCGGAGCAGGACCTCCAGCTCGGCGGCCCGGGAAGCCACCCGCCCTCACCTGCGTTCTGACGGCAAGCTTCGGCGAAGAAGCTCTGG
3701 CCCTGCGGGTCGCTTGTGGTCTTTATAGCCGCTCGGCGTCAGGCCCGCCCGCCAATCAGCGCCCGCCCGAGCCCGCTCCTTCGGTGGGCGC
3801 GGGACCCCGCCCTGCTGTGGGGAGGGGCGGCCGCTGGAGGCCCTCGCGCTCTGGCGAACTAGT**CGACGGACTGGGCTACGGGCCGCCCGAGAG**
3901 **CGCAGCCAGAGGCCGCTAGGAAGGGGCGGGCGCCGAGAACACGATCCCTCCCCACCCCTCGGACGTGACTCGGACCACATCCCGGGTTCGCTA**
4001 **GGGCCCTCCCTTCTGCTCCTTTCCCCAGCCTGGCGGCTCTGGGGCGCCGTGACTCAGCCAGAATGTTGGCAATGGGGAGGGCGGAACGGGAAGTGG**
4101 **AGGACCGGATGAAAAAGTCGGAACGAAGGAAGCTGAGTTTCGCCTGCAGGTTAATTAAGAACATGTGAGCAAAAGCCAGCAAAAGCCAGGAACCGT**
4201 **AAAAAGCCCGTGTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGAC**
4301 **TATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCCTTCTCCCTTCGGG**
4401 **AAGCGTGGCGTTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTAGCCC**
4501 **GACCGCTGCGCCTTATCCGTAACCTATCGTCTTGAGTCCAACCCGTAAGACACGACTTATGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAG**
4601 **CGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGT**
4701 **TACCTTCGAAAAAGAGTTGGTAGCTTTGATCCGGCAAAACAAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGCAAGCAGCAGATTACGCGCAGAAAA**
4801 **AAAGGATCTCAAGAAGATCCTTTGATCTTTTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGT**