Certificate of Analysis

pF1A T7 Flexi® Vector:

Part No. Size (units) C844A 20µg

Description: The pF1A T7 Flexi® Vector(a-c) is designed for use with the Flexi® System, Entry/Transfer (Cat.# C8640), and the Flexi® System, Transfer (Cat.# C8820). The vector contains a T7 promoter for bacterial or in vitro protein expression of a protein-coding region. The vector also contains the lethal barnase gene for positive selection of the insert, an ampicillin-resistance gene for selection of the plasmid and unique Sgfl and Pmel sites that allow easy insertion or transfer of the sequence of interest. Inserts containing a protein-coding region can easily be transferred from the pF1A T7 Flexi® Vector to other Flexi® Vectors with different expression options. For more information, see the *Flexi® Vector Systems Technical Manual* #TM254 available online at: **www.promega.com/protocols**

Concentration: 100ng/µl.

GenBank® Accession Number: AY753576.

Storage Buffer: The pF1A T7 Flexi® Vector is supplied in 10mM Tris-HCI (pH 8.0), 1mM EDTA.

 $\textbf{Storage Conditions:} Store \ the \ vector \ at \ -20 ^{\circ}\text{C}. \ A void \ multiple \ freeze-thaw \ cycles \ and \ exposure \ to \ frequent \ temperature$

changes. These fluctuations can greatly alter product stability.

Usage Notes: Concentration gradients may form in frozen products and should be dispersed upon thawing. Mix well prior to

use.

Quality Control Assays

Nuclease Assay: Following incubation of 1µg of pF1A T7 Flexi® Vector in Restriction Enzyme Buffer B at 37°C for 16 hours, no evidence of nuclease activity is detected by agarose gel electrophoresis.

Physical Purity: A₂₆₀/A₂₈₀ > 1.80.

Restriction Digestion: The presence of unique restriction sites for Pmel and Sgfl is confirmed by showing that the vector yields the expected fragment sizes after digesting 1µg of vector for 2 hours with 10 units of Pmel, Sgfl and Bglll.

(a)Patent Pending

(e)For research use only. Persons wishing to use this product or its derivatives in other fields of use, including without limitation, commercial sale, diagnostics or therapeutics, should contact Promega Corporation for licensing information.

(c)U.S. Pat. Nos. 8,293,503 and 8,367,403, European Pat. No. 1685247 and other patents and patents pending

Signed by:

R Wheeler Quality Assurance

Part# 9PIC844 Revised 10/16



AF9PIC844 1016C844



Promega Corporation	on
2800 Woods Hollow Road	
Madison, WI 53711-5399	USA
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	www.promega.com

PRODUCT USE LIMITATIONS, WARRANTY, DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be armful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

© 2004–2014, 2016 Promega Corporation. All Rights Reserved.

Flexi is a registered trademark of Promega Corporation.

GenBank is a registered trademark of US Dept of Health and Human Services.

All specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PIC844 Printed in USA. Revised 10/16.



Usage Information

pF1A T7 Flexi® Vector Features and Circle Map

The following features are present in the vector based on nucleotide sequence.

T7 RNA polymerase promoter (-17 to +3)	21-40
Sgfl site	100-107
barnase coding region	109-444
Pmel site	446-453
T7 terminator	573-620
β-lactamase (Amp ^r) coding region	954-1814
ColE1-derived plasmid origin of replication	1969-2005
cer site (site for <i>E. coli</i> XerCD recombinase)	2676-2961
rrnB transcription terminator	3012-3413

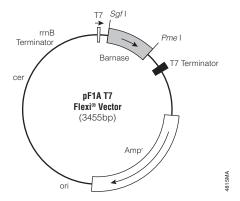


Figure 1. pF1A T7 Flexi® Vector circle map and sequence reference points.

Note: Maps of all the Flexi® Vectors are available at: **www.promega.com/vectors/cloning_vectors.htm**

Related Products

Product	Size	Cat.#
Flexi® System, Entry/Transfer	5 entry and 20 transfer reactions	C8640
Flexi® System, Transfer	100 transfer reactions	C8820
Carboxy Flexi® System, Transfer	50 transfer reactions	C9320
10X Flexi® Enzyme Blend (Sgfl & Pmel) <u>25</u> µl	R1851
	100µl	R1852
Carboxy Flexi Enzyme Blend (Sgfl & Ec	olCRI) 50µl	R1901
HaloTag® Flexi® Vectors—CMV Dilution	on Series Sample Pack 9 × 2μg	G3780
Single Step (KRX) Competent Cells	20 × 50µl	L3002

There are Flexi® Vectors available for many different applications. Visit: **www.promega.com/applications/cloning** to find out more.

Summary of Changes

The following changes were made to the 12/14 revision of this document: 1. Expired patent or license statements were removed.