Certificate of Analysis

pF1K T7 Flexi[®] Vector:

Part No. Size (units	
C845A	20µg

Description: The pF1K 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, a kanamycin-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 pF1K T7 Flexi® Vector to other Flexi® Vectors with different expression options (Table 1). For more information, see the *Flexi® Vector Systems Technical Manual* #TM254.

Table 1. Vectors Available for Use With the Flexi® Vector Systems.

Cat.#	Flexi® Vector	Utility	Expression	Drug Selection
C8441	pF1A T7 Flexi [®] Vector	Protein expression	E. coli and in vitro (T7 promoter)	Ampicillin
C8451	pF1K T7 Flexi [®] Vector			Kanamycin
C8461	pFN2A (GST) Flexi [®] Vector	Protein expression	E. coli and in vitro (T7 promoter)	Ampicillin
C8471	pFN2K (GST) Flexi [®] Vector	and purification		Kanamycin
L5671	pF3A WG (BYDV) Flexi® Vector	Protein expression	Wheat Germ in vitro (T7, SP6)	Ampicillin
L5681	pF3K WG (BYDV) Flexi® Vector	Protein expression	Wheat Germ in vitro (T7, SP6)	Kanamycin
C8481 C8491	pF4A CMV Flexi [®] Vector pF4K CMV Flexi [®] Vector	Protein expression	Mammalian (CMV promoter) and in vitro (T7 promoter)	Ampicillin Kanamycin

Usage Information

Concentration: 100ng/µl.

GenBank® Accession Number: AY753577.

Storage Buffer: The pF1K T7 Flexi® Vector is supplied in 10mM Tris-HCI (pH 8.0), 1mM EDTA.

Storage Conditions: Store the vector at –20°C. Avoid 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 pF1K 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 Pme I, Sgfl and BgIII.

^(a)Patent Pending.

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

Signed by:

for Wheeler

R. Wheeler, Quality Assurance

Part# 9PIC845 Revised 03/18



AF9PIC845 0318C845



Promega Corporation

2800 Woods Hollow Road	1
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 harmful 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 INDI-RECTLY, EXPRESS OR INPLIED, INCLUDING, WITH-OUT LIMITATION, AS TO THE SUITABILITY, PRODUC-TIVTY, 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 and HaloTag are registered trademarks of Promega Corporation.

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

Products may be covered by pending or issued patents or may have certain limitations. Please visit our web site for more information.

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# 9PIC845 Printed in USA. Revised 03/18.



Usage Information

pF1K 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			
kanamycin resistance coding region	1001-1795			
ColE1-derived plasmid origin of replication	1964-2000			
cer site (site for <i>E. coli</i> XerCD recombinase)	2671-2956			
rrnB transcription terminator	3007-3408			

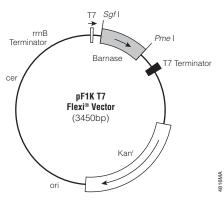


Figure 1. pF1K T7 Flexi® Vector circle map and sequence reference points.

Related Products

Size	Cat.#
entry and 20 transfer reactions	C8640
100 transfer reactions	C8820
50 transfer reactions	C9320
25µl	R1851
100µl	R1852
RI) 50µl	R1901
eries Sample Pack 9 × 2µg	G3780
20 × 50µl	L3002
	entry and 20 transfer reactions 100 transfer reactions 50 transfer reactions 25µl 100µl RI) 50µl rries Sample Pack 9 × 2µg

Summary of Changes

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

> Part# 9PIC845 Printed in USA. Revised 03/18.