# **Certificate of Analysis**

## pF4K CMV Flexi® Vector:

**Part No. Size (units)** C849A 20µg

**Description:** The pF4K CMV Flexi® Vector(a.b.d) is designed for use with the Flexi® System, Entry/Transfer (Cat.# C8640), and the Flexi® System, Transfer (Cat.# C8820). The vector contains a CMV immediate early enhancer/promoter plus a chimeric intron for mammalian expression of the protein-coding region and a T7 promoter for in vitro protein expression of the 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 pF4K CMV 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	pF4A CMV Flexi® Vector	Protein expression	Mammalian (CMV promoter)	Ampicillin
C8491	pF4K CMV Flexi® Vector		and in vitro (T7 promoter)	Kanamycin

## **Usage Information**

Concentration: 100ng/µl.

GenBank® Accession Number: AY753581.

Storage Buffer: The pF4K CMV 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 pF4K CMV 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:**  $A_{260}/A_{280} > 1.80$ .

**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 vector for 2 hours with 10 units of Pmel, Sgfl and Bgl II.

(D)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)The CMV promoter and its use are covered under U.S. Pat. Nos. 5,168,062 and 5,385,839 owned by the University of Iowa Research Foundation, Iowa City, Iowa, and licensed FOR RESEARCH USE ONLY. Commercial users must obtain a license to these patents directly from the University of Iowa Research Foundation.

Signed by:

R Wheeler Quality Assurance

# Part# 9PIC849 Revised 4/18



AF9PIC849 0418C849



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.

© 2004–2018 Promega Corporation. All Rights Reserved.

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

Flexi and HaloTag are registered trademarks 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# 9PIC849

Printed in USA. Revised 4/18.

<sup>(</sup>a)Patent Pending



# **Usage Information**

## pF4K CMV Flexi® Vector Features and Circle Map

The following features are present in the vector based on nucleotide sequence. CMV immediate early enhancer/promoter 1-742 chimeric intron 857-989 1033-1052 T7 RNA polymerase promoter (-17 to +3) 1056-1063 Sgfl site barnase coding region 1087-1422 Pmel site 1424-1431 SV40 late poly(A) signal 1583-1804 kanamycin resistance coding region 2112-2906 ColE1-derived plasmid origin of replication 3075-3111 cer site (site for E. coli XerCD recombinase) 3782-4067

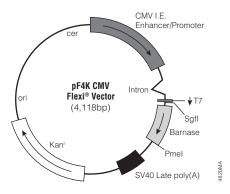


Figure 1. pF4K CMV Flexi® Vector circle map and sequence reference points.

### **Related Products**

Product	Size	Cat.#
Flexi® System, Entry/Transfer 5 ent	ry 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)	25µl	R1851
	100µl	R1852
Carboxy Flexi Enzyme Blend (Sgfl & EcolCRI)	) 50µl	R1901
HaloTag® Flexi® Vectors—CMV Dilution Series Sample Pack 9 × 2μς		G3780