## pF4A CMV Flexi® Vector:

Part No.	Size (units)
C848A	20µg

**Description:** The pF4A CMV Flexi® Vector<sup>(a,b,c)</sup> 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, 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 pF4A 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 <sup>®</sup> Vector	Utility	Expression	Drug Selection
C8441	pF1A T7 Flexi <sup>®</sup> Vector	Protein expression	E. coli and in vitro (T7 promoter)	Ampicillin
C8451	pF1K T7 Flexi <sup>®</sup> Vector			Kanamycin
C8461	pFN2A (GST) Flexi <sup>®</sup> Vector	Protein expression	E. coli and in vitro (T7 promoter)	Ampicillin
C8471	pFN2K (GST) Flexi <sup>®</sup> Vector	and purification		Kanamycin
L5671	pF3A WG (BYDV) Flexi® Vector	Protein expression	Wheat Germ in vitro (T7, SP6)	Ampicillin
L5681	pF3K WG (BYDV) Flexi <sup>®</sup> 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: AY753580.

Storage Buffer: The pF4A CMV Flexi® Vector is supplied in 10mM Tris-HCI (pH 8.0), 1mM EDTA.

**Storage Conditions:** Store the vector at  $-20^{\circ}$ 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 pF4A CMV Flexi<sup>®</sup> 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<sub>260</sub>/A<sub>280</sub> >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 Bgl II.

(a)Patent Pending.

<sup>(b)</sup>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.

<sup>(c)</sup>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.

Fin Wheeler

Signed by:

R. Wheeler, Quality Assurance

# Part# 9PIC848 Revised 4/18



AF9PIC848 0418C848



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.

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 WAR-RANTIES 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, MER-CHANTABILITY, CONDITION, OR ANY OTHER MAT-TER 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 profil), whether based upon warranty, contract, tot (including but not limited to loss of use, revenue or profil, whether based upon warranty, contract, tot 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# 9PIC848 Printed in USA. Revised 4/18.



# **Usage Information**

#### pF4A 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			
T7 RNA polymerase promoter (-17 to +3)	1033-1052			
Sgfl site	1056-1063			
barnase coding region	1087-1422			
Pmel site	1424-1431			
SV40 late poly(A) signal	1583-1804			
$\beta$ -lactamase (Amp <sup>r</sup> ) coding region	2065-2925			
ColE1-derived plasmid origin of replication	3080-3116			
cer site (site for E. coli XerCD recombinase)	3787-4072			



Product		Size	Cat.#
Flexi <sup>®</sup> System, Entry/Transfer	5 entry and 20 transfe	r reactions	C8640
Flexi <sup>®</sup> System, Transfer	100 transfe	r reactions	C8820
Carboxy Flexi <sup>®</sup> System, Transfer	50 transfe	r reactions	C9320
10X Flexi® Enzyme Blend (Sgfl & Pme	l)	25µl	R1851
		100µl	R1852
Carboxy Flexi Enzyme Blend (Sgfl & Ec	colCRI)	50µl	R1901
HaloTag <sup>®</sup> Flexi <sup>®</sup> Vectors–CMV Dilution	n Series Sample Pack	9 × 2µg	G3780

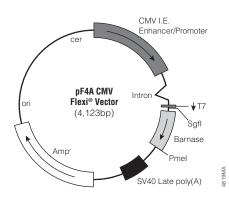


Figure 1. pF4A CMV Flexi® Vector circle map and sequence reference points.

Part# 9PIC848 Printed in USA. Revised 4/18.