pF5A CMV-neo Flexi[®] Vector:

Size (units) Part No. C940A

20ua

Description: The pF5A CMV-neo Flexi® Vector(a.b.c) is designed for use with the Flexi® System, Entry/Transfer (Cat.# C8640) and the Flexi® System, Transfer (Cat.# C8820). The pF5A CMV-neo Flexi® Vector contains the CMV immediate early enhancer/ promoter region plus a chimeric intron for mammalian expression and a T7 promoter for in vitro expression of the proteincoding region. The vector also contains the barnase gene for positive selection of the insert, an ampicillin resistance gene for selection of the plasmid, and unique Sqfl and Pmel sites that allow easy insertion and transfer of the insert of interest. The pF5A CMV-neo Flexi® Vector also contains a neomycin resistance gene under the control of the SV40 promoter for long-term selection of transfected mammalian cells. Transfection of a pF5A CMV-neo Flexi® Vector containing a cloned protein-coding region into mammalian cells provides resistance to the antibiotic G-418 (Cat.# V7981, V8091), allowing selection of transfectants. Do not use the pF5A CMV-neo Flexi® Vector without an insert as a negative control because the barnase gene decreases the viability of the transfected cells. Cloned protein-coding regions can be transferred from the pF5A CMV-neo Flexi® Vector to other Flexi® Vectors with different expression options using the Flexi® Entry/Transfer Systems (Cat.# C8640 and C8820). For more information, see the Flexi® Vector Systems Technical Manual #TM254.

Usage Information

Concentration: 100ng/µl.

GenBank® Accession Number: DQ487155

Storage Buffer: The pF5A CMV-neo Flexi® Vector is supplied in 10mM Tris-HCI (pH 8.0), 1mM EDTA.

Storage Conditions: See the Product Information Label for storage recommendations. 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 pF5A CMV-neo 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: $A_{260}/A_{280} \ge 1.80$

Restriction Digestion: The presence of unique restriction sites for Pmel and Sgfl is confirmed by showing that the vector is linearized and 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

(b)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)European Pat. No. 1685247 and other patents pending

Kin Wheeler

Part# 9PIC940 Revised 10/16



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 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 replace-ment 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, foresee-able, consequential, or special (including but not lim-ited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including ne gligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

© 2006–2013, 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# 9PIC940 Printed in USA. Revised 10/16.

Signed by:



Usage Information

pF5A CMV-neo 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
SV40 enhancer and early promoter	1903-2320
neomycin phosphotransferase coding region	2366-3160
synthetic polyadenylation signal	3224-3272
β-lactamase coding region	3533-4393
ColE1-derived plasmid replication of origin	4548-4584
cer site (site for <i>E. coli</i> XerCD recombinase)	5255-5540

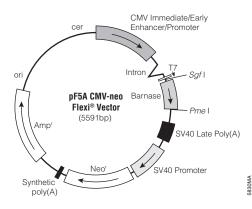


Figure 1. pF5A CMV-neo Flexi® Vector circle map.

Related Products

Product	S	ize	Cat.#
Flexi® System, Entry/Transfer	5 entry and 20 transfer reaction	ons	C8640
Flexi [®] System, Transfer	100 transfer reaction	ons	C8820
Carboxy Flexi [®] System, Transfer	50 transfer reaction	ons	C9320
10X Flexi® Enzyme Blend (Sgfl & Pm	el)2	5µl	R1851
	10	0µI	R1852
Carboxy Flexi Enzyme Blend (Sgfl & E	EcolCRI) 5	0µI	R1901
HaloTag® Flexi® Vectors-CMV Diluti	on Series Sample Pack 9 × 2	2µg	G3780
Single Step (KRX) Competent Cells	5 x 20	0µl	L3001

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

> Part# 9PIC940 Printed in USA. Revised 10/16