

Product Contents

pFC14K HaloTag[®] CMV Flexi[®] Vector:

Part No.	Size
G966A	20µg

Part# 9PIG966

Revised 10/16

Description: The pFC14K HaloTag[®] CMV Flexi[®] Vector^(a-9) is configured to append the HaloTag[®] tag to the carboxy-terminus of the protein fusion partner. The vector provides constitutive protein expression in mammalian cells using the human cytomegalovirus (CMV) intermediate early enhancer/promoter. The vector can be used for both stable and transient gene expression; for stable expression, cotransfection with a vector containing a selectable marker is required.

The pFC14K HaloTag[®] CMV Flexi[®] Vector contains the following features:

- A **CMV intermediate/early enhancer/promoter** for constitutive expression in mammalian cells.
- A **T7 RNA polymerase promoter** for in vitro HaloTag[®] fusion protein expression in cell-free systems (e.g., TnT[®] lysate reaction).
- The **C-terminal HaloTag[®] region**, which rapidly forms covalent bonds with HaloTag[®] ligands, enabling labeling or immobilization of expressed proteins.
- A **TEV protease site** for cleavage of the expressed protein from HaloTag[®] using ProTEV Protease (Cat.# V6051).
- The lethal **barnase gene** for positive selection of the insert. **Note: The pFC14K HaloTag[®] CMV Flexi[®] Vector can be propagated only in *E. coli* once the barnase gene is replaced with the protein-coding sequence of interest.**
- A **kanamycin-resistance gene** for selection of the plasmid.
- Unique **SgfI and EcoRI sites**, which allow easy insertion of the sequence of interest. These sites create a readthrough sequence that can be joined to a protein-coding region flanked by SgfI and PmeI sites, enabling easy transfer to the pFC14K HaloTag[®] CMV Flexi[®] Vector from other Flexi[®] Vectors with different expression options. **Once inserted in this vector, the sequence is no longer available for transfer.** For more information, see the *Flexi[®] Vector Systems Technical Manual #TM254*, available online at www.promega.com/protocols

Concentration: 100ng/µl.

GenBank[®] Accession Number: EU113047.

Storage Buffer: The pFC14K HaloTag[®] CMV Flexi[®] Vector is supplied in 10mM Tris-HCl (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. See label for expiration date.

Usage Notes:

1. Use this vector in conjunction with pFC15, pFC16 and pFC17 Flexi[®] Vectors to determine which vector provides the appropriate protein expression level for your particular application. The pFC14 Flexi[®] Vector carries the full-length CMV promoter while pFC15, pFC16 and pFC17 Flexi[®] Vectors contain various deletions of the CMV promoter. Since the full-length CMV promoter expresses highly in many cell types, it may be inappropriate for applications where high concentrations of fusion protein may affect physiological function.
2. This vector was designed to be used with the Flexi[®] Vector System, a directional cloning method to shuttle protein-coding sequences between compatible vectors. In this system, carboxy-terminal tag fusions cannot shuttle the insert to other expression vectors. To retain the capacity to transfer a protein-coding sequence to multiple vectors, first clone the protein-coding sequence into an ampicillin-resistant Flexi[®] Vector with no tag or an amino-terminal tag [e.g., pF4A CMV Flexi[®] Vector (Cat.# C8481) or pFN21A HaloTag[®] CMV Flexi[®] Vector (Cat.# G2821)] prior to transferring the insert to the pFC14K HaloTag[®] CMV Flexi[®] Vector. For more information, see the *Flexi[®] Vector Systems Technical Manual #TM254*, available online at: www.promega.com/protocols
3. Concentration gradients may form in frozen products and should be dispersed upon thawing. Mix well prior to use.

Quality Control Assays

Contaminant Assays

Contaminating Nucleic Acids: RNA, single-stranded DNA and chromosomal DNA are not evident in specified quantities of the vector as determined by agarose gel electrophoresis.

Nuclease Assay: Following incubation of 1µg of the vector in restriction enzyme buffer at 37°C for 16–24 hours, no evidence of nuclease activity is detected by agarose gel electrophoresis.

Physical Purity: $A_{260}/A_{280} \geq 1.80$, $A_{260}/A_{250} \geq 1.05$.

Functional Assays

Identity Assay: The vector has been sequenced completely and has 100% identity with the published sequence available at: www.promega.com/vectors/

Restriction Digestion: The functional purity of the vector DNA is verified by successful digestion with restriction enzymes at the optimal temperature for one hour. Samples are examined by agarose gel electrophoresis, comparing cut and uncut vector DNA with marker DNA.



AF9PIG966 1016G966



Promega

Promega Corporation

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 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 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.

© 2007–2009, 2013, 2016 Promega Corporation. All Rights Reserved.

Flexi, HaloTag, TnT and Wizard 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# 9PIG966
Printed in USA. Revised 10/16.

Signed by:

R. Wheeler, Quality Assurance

pFC14K HaloTag[®] CMV Flexi[®] Vector Features and Circle Map

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

CMV intermediate 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
EcoICRI site	1442–1447
TEV site	1462–1482
HaloTag [®] coding region	1492–2382
SV40 late polyadenylation signal	2516–2737
Kanamycin resistance (Kan ^r) coding region	3045–3839
Col/E1-derived plasmid origin of replication	4008–4044
cer site (site for <i>E. coli</i> XerCD recombinase)	4715–5000

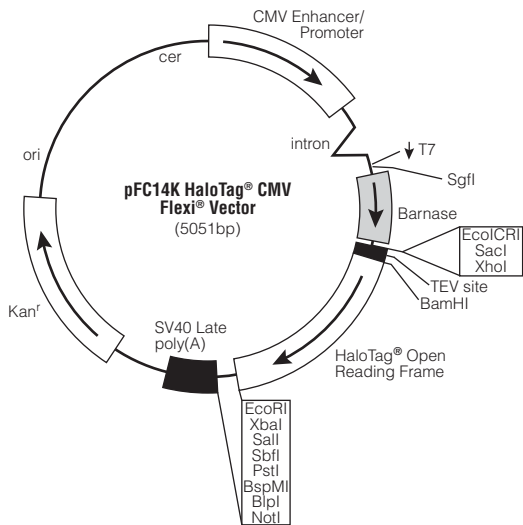


Figure 1. pFC14K HaloTag[®] CMV Flexi[®] Vector circle map and sequence reference points.

Note: Maps of all the Flexi[®] Vectors are available at: www.promega.com/vectors

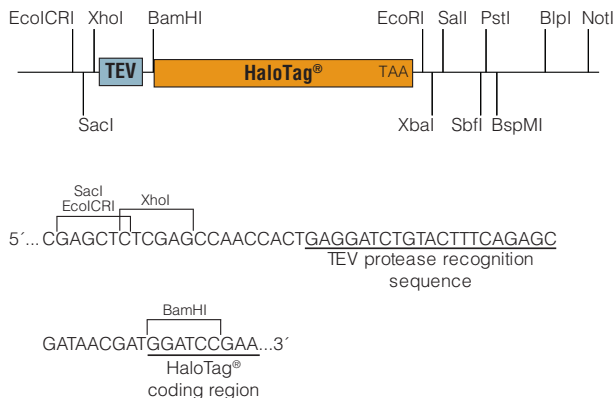


Figure 2. pFC14K HaloTag[®] CMV Flexi[®] Vector sequence upstream and downstream of the HaloTag[®] gene.

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)	25µl	R1851
	100µl	R1852
Carboxy Flexi Enzyme Blend (Sgfl & EcoICRI)	50µl	R1901
HaloTag [®] Flexi [®] Vectors—CMV Dilution Series	9 × 2µg	G3780
Single Step (KRX) Competent Cells	5 × 200µl	L3001

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

©READ THIS FIRST BEFORE OPENING PRODUCT

The sale of this product and its use are subject to the terms of a limited use label license, the full text of which is available at: www.promega.com/LULL. That text must be read by the purchaser prior to opening this product to determine whether the purchaser agrees that all use of the product shall be in accordance with the license terms. If the purchaser is not willing to accept the terms of the limited use label license, Promega is willing to accept the return of the unused product and provide the purchaser with a full refund. However, if the product is opened for any reason, then the purchaser agrees to be bound by the terms of the limited use label license.

©U.S. Pat. Nos. 7,425,436 and 7,935,803 and other patents pending.

©European Pat. No. 1685247 and other patents pending.