pFN21A HaloTag® CMV Flexi® Vector:

Part No.	
G282A	

Description: The pFN21A HaloTag[®] CMV Flexi[®] Vector^(a,b,c) is configured to append the HaloTag[®] tag to the aminoterminus of the protein fusion partner and 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 pFN21A HaloTag® CMV Flexi® Vector contains the following features:

Size 20µg

- · A CMV intermediate/early enhancer/promoter for constitutive expression in mammalian cells.
- A T7 RNA polymerase promoter for in vitro HaloTag® fusion protein expression.
- The N-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 pFN21A HaloTag[®] CMV Flexi[®] Vector can be
 propagated only in *E. coli* once the barnase gene is replaced with the protein-coding sequence of interest.
- An ampicillin-resistance gene for selection of the plasmid.
- Unique Sgfl and Pmel 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 Sgfl and Pmel sites, enabling easy transfer to the pFN21A HaloTag® CMV Flexi® Vector from other Flexi® Vectors with different expression options. Once inserted in this vector, the sequence is available for transfer to other Flexi® Vectors. For more information, see the Flexi® Vector Systems Technical Manual #TM254, available online at: www.promega.com/protocols/

Concentration: 100ng/µl

GenBank® Accession Number: EU621374.

Storage Buffer: The pFN21A 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:

- Use this vector in conjunction with pFN22, pFN23 and pFN24 Flexi[®] Vectors to determine which vector provides the appropriate protein expression level for your particular application. The pFN21 Flexi[®] Vector carries the full-length CMV promoter while pFN22, pFN23 and pFN24 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. See Table 1 on reverse side for additional information.
- This vector was designed to be used with the Flexi® Vector System, a directional cloning method to shuttle protein-coding sequences between compatible vectors. To prepare the HaloTag® fusion protein, the protein coding region is cloned into the pFN21A HaloTag® CMV Flexi® Vector using the Flexi® System, Entry/Transfer (Cat.# C8640). For more information, see the *Flexi*® *Vector Systems Technical Manual* #TM254.
- 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: $A_{260}/A_{280} \ge 1.80$, $A_{260}/A_{250} \ge 1.05$.

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.

Signed by:

Ren Wheeler

R. Wheeler, Quality Assurance

Part# 9PIG282 Revised 10/16



AF9PIG282 1016G282



Promega Corporation

2800 Woods Hollow Road	ł
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 WAR-RANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, DIRCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, FONDUCTIVITY, 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 any other damages, whether direct, incidental, forese-able, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tor (including the the failure of Promega products to perform in accordance with the stated specifications.

 \circledast 2011, 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# 9PIG282 Printed in USA. Revised 10/16.



Usage Information

pFN21A 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
HaloTag [®] protein coding region	1067-1957
TEV site	1970-1990
Sgfl site	1997-2004
barnase coding region	2028-2363
Pmel site	2365-2372
SV40 late polyadenylation signal	2524-2745
β -lactamase (Amp ^r) coding region	3006-3866
Co/E1-derived plasmid origin of replication	4021-4057
cer site (site for <i>E. coli</i> XerCD recombinase)	4728-5013

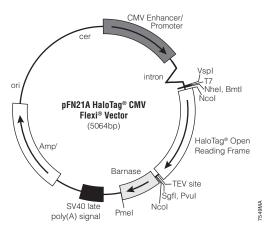


Figure 1. pFN21A HaloTag® CMV Flexi® Vector circle map and sequence reference points.

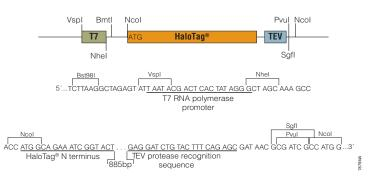


Figure 2. pFN21A HaloTag $^{\otimes}$ CMV Flexi $^{\otimes}$ Vector sequence upstream and downstream of the HaloTag $^{\otimes}$ gene.

Table 1. Relative Mammalian Expression Levels for HaloTag® Flexi® Vectors.

Vector Name	Cat.#	Expression Level*	
pFC14A HaloTag® CMV Flexi® Vector	G9651	- High	
pFC14K HaloTag [®] CMV Flexi [®] Vector	G9661		
pFC15A HaloTag® CMV <i>d1</i> Flexi® Vector	G1611	- Medium	
pFC15K HaloTag® CMV <i>d1</i> Flexi® Vector	G1601		
pFC16A HaloTag [®] CMV <i>d2</i> Flexi [®] Vector	G1591	- Low	
pFC16K HaloTag® CMV <i>d2</i> Flexi® Vector	G1571		
pFC17A HaloTag® CMV <i>d3</i> Flexi® Vector	G1551	- Ultralow	
pFC17K HaloTag [®] CMV <i>d3</i> Flexi [®] Vector	G1321		
pFN21A HaloTag® CMV Flexi® Vector	G2821	- High	
pFN21K HaloTag® CMV Flexi® Vector	G2831		
pFN22A HaloTag [®] CMV <i>d1</i> Flexi [®] Vector	G2841	Medium	
pFN22K HaloTag® CMV <i>d1</i> Flexi® Vector	G2851		
pFN23A HaloTag® CMV <i>d2</i> Flexi® Vector	G2861	- Low	
pFN23K HaloTag [®] CMV <i>d2</i> Flexi [®] Vector	G2871		
pFN24A HaloTag® CMV <i>d3</i> Flexi® Vector	G2881	Ultralow	
pFN24K HaloTag® CMV <i>d3</i> Flexi® Vector	G2981		

*Expression level depends on the cell type and the protein fused to HaloTag® coding region.

Related Products

Product	Size	Cat.#
HaloTag [®] Cloning Starter System		G6050
HaloTag [®] Flexi [®] Vectors-CMV Dilution S	eries 9 × 2µg	G3780
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 & Ecol	CRI) 50µl	R1901

(a)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 terms of the purchaser agrees to be bound by the terms of the limited use label license.

^(b)U.S. Pat. Nos. 7,425,436 and 7,935,803 and other patents pending.

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

Part# 9PIG282 Printed in USA. Revised 10/16.